
Selected CF® Dye References, by Dye Color and Application

Click on the blue link to see journal article

Contents:

The use of CF® Dyes have been published in hundreds of journal articles. This list of publications using CF® Dyes is organized below by the dye color and by application that the CF® Dyes were used with.

CF® Dyes by Color

- [CF®350](#) p. 2
- [CF®405M](#) p. 2
- [CF®405S](#) p. 4
- [CF®405L](#) p. 6
- [CF®488A](#) p. 6
- [CF®514](#) p. 22
- [CF®532](#) p. 22
- [CF®543](#) p. 22
- [CF®555](#) p. 24
- [CF®568](#) p. 30
- [CF®583R](#) p. 40
- [CF®594](#) p. 40
- [CF®597R](#) p. 45
- [CF®620R](#) p. 45
- [CF®633](#) p. 45
- [CF®640R](#) p. 50
- [CF®647](#) p. 55
- [CF®660C](#) p. 60
- [CF®660R](#) p. 61
- [CF®680](#) p. 62
- [CF®680R](#) p. 66
- [CF®750](#) p. 67
- [CF®770](#) p. 69
- [CF®790](#) p. 74

CF® Dyes by Application

- [Mix-n-Stain™](#) p. 74
- [Flow Cytometry](#) p. 79
- [Western Blotting](#) p. 80
- [In-Cell Western®](#) p. 83
- [In Vivo Imaging](#) p. 84
- [Super-Resolution Microscopy](#) p. 86
- [TIRF Microscopy](#) p. 92
- [Fluorescence Polarization Assay](#) p. 93
- [2-Photon Microscopy](#) p. 93

Abbreviations: FC: flow cytometry; FIONA: fluorescence imaging with one-nanometer accuracy; FISH: fluorescence in situ hybridization; FLImP: fluorophore localization imaging with photobleaching; FRAP: fluorescence recovery after photobleaching; gSHRImP: generalized single molecule high-resolution imaging with photobleaching; IF: immunofluorescence; ICW: in-cell western®; IHF: immuno-histofluorescence; SIM: structured illumination microscopy; SMLM: single-molecule localization microscopy; STED: stimulated emission depletion; STORM: stochastic optical reconstruction microscopy; TIRF: total internal reflection microscopy; WB: Western blotting.

LI-COR, Odyssey, and In-Cell Western are registered trademarks of LI-COR Biosciences in the United States and other countries. IVIS is a registered trademark of Caliper Life Sciences.

CF® Dye References by Dye Color

CF®350

Ahi, Y. S., et al. [Adenoviral E2 IVa2 Protein Interacts with L4 33K Protein and E2 DNA-Binding Protein](#). J Gen Virol 94(6), 1325, (2013), DOI: DOI:10.1099/vir.0.049346-0 (Mix-n-Stain CF350, Mix-n-Stain CF488A, CF568 goat anti-mouse IgG; ICC)

Bojarczuk, A., et al. [Cryptococcus neoformans Intracellular Proliferation and Capsule Size Determines Early Macrophage Control of Infection](#). Sci Rep 6, 21489, (2016), DOI: 10.1038/srep21489 (CF350 goat anti-mouse fragment)

Fan, J. X., et al. [Targeting epithelial-mesenchymal transition: Metal organic network nano-complexes for preventing tumor metastasis](#). Biomaterials 139, 116-126, (2017), DOI: 10.1016/j.biomaterials.2017.06.007 (CF350 succinimidyl ester)

Jin, S., et al. [Tetherin Suppresses Type I Interferon Signaling by Targeting MAVS for NDP52-Mediated Selective Autophagic Degradation in Human Cells](#). Mol Cell 68(2), 308-322 e304, (2017), DOI: 10.1016/j.molcel.2017.09.005 (CF350 donkey anti-rabbit)

Lee, W., et al. [Synergy between Piezo1 and Piezo2 channels confers high-strain mechanosensitivity to articular cartilage](#). Proc Natl Acad Sci U S A 111(47), E5114-5122, (2014), DOI: 10.1073/pnas.1414298111 (CF350 phalloidin)

Pu, Q., et al. [Identifying drug resistant cancer cells using microbubble well arrays](#). Biomed Microdevices 19(3), 17, (2017), DOI: 10.1007/s10544-017-0160-9 (CF350 goat anti-mouse IgG1)

Shanks, H. R., et al. [Including fluorescent nanoparticle probes within injectable gels for remote strain measurements and discrimination between compression and tension](#). Soft Matter 17(4), 1048-1055, (2021), DOI: 10.1039/d0sm01635g (CF350 amine)

Takenaka, N., et al. [Involvement of the protein kinase Akt2 in insulin-stimulated Rac1 activation leading to glucose uptake in mouse skeletal muscle](#). PLoS One 14(2), e0212219, (2019), DOI: 10.1371/journal.pone.0212219, PONE-D-18-28417 (CF350 CF543 CF647 secondaries)

Takenaka, N., et al. [Role of the guanine nucleotide exchange factor in Akt2-mediated plasma membrane translocation of GLUT4 in insulin-stimulated skeletal muscle](#). Cell Signal 26(11), 2460-2469, (2014), DOI: S0898-6568(14)00221-6, 10.1016/j.cellsig.2014.07.002 (CF350, CF543 & CF647; ICC)

Tschumper, R. C., et al. [Molecular analysis of immunoglobulin genes reveals frequent clonal relatedness in double monoclonal gammopathies](#). Blood Cancer J 3, e112, (2013), DOI: DOI:10.1038/bcj.2013.12 (Mix-n-Stain CF350; ICC)

Zhang, M., et al. [A high M1/M2 ratio of tumor-associated macrophages is associated with extended survival in ovarian cancer patients](#). J Ovarian Res 7(1), 19, (2014), DOI: 10.1186/1757-2215-7-19, 10.1186/1757-2215-7-19 (CF350, CF488A labeled primaries; IHC)

CF®405M

Barbirou, M., et al. [Single Circulating-Tumor-Cell-Targeted Sequencing to Identify Somatic Variants in Liquid Biopsies in Non-Small-Cell Lung Cancer Patients](#). Current Issues in Molecular Biology 44(2), 750-763, (2022), DOI: 10.3390/cimb44020052 (CF405 succinimidyl ester & CF647 succinimidyl ester)

Bolduc, J. F., et al. [Epigenetic Metabolite Acetate Inhibits Class I/II Histone Deacetylases, Promotes Histone Acetylation, and Increases HIV-1 Integration in CD4\(+\) T Cells](#). J Virol 91(16), (2017), DOI: 10.1128/JVI.01943-16 (CF4505M Mix-n-Stain flow)

Chiva-Blanch, G., et al. [Microparticle Shedding from Neural Progenitor Cells and Vascular Compartment Cells Is Increased in Ischemic Stroke](#). PLoS One 11(1), e0148176, (2016), DOI: 10.1371/journal.pone.0148176 (CF405M Annexin V)

Comrie, W. A., et al. [F-actin flow drives affinity maturation and spatial organization of LFA-1 at the immunological synapse](#). J Cell Biol 208(4), 475-491, (2015), DOI: jcb.201406121, 10.1083/jcb.201406121 (CF405M phalloidin)

Demmerle, J., et al. [Strategic and practical guidelines for successful structured illumination microscopy](#). Nat Protoc 12(5), 988-1010, (2017), DOI: nprot.2017.019, 10.1038/nprot.2017.019 (CF405S and CF405M SIM)

Gielen, P. R., et al. [Connexin43 confers Temozolomide resistance in human glioma cells by modulating the mitochondrial apoptosis pathway](#). Neuropharmacology, (2013), DOI: DOI:10.1016/j.neuropharm.2013.05.002 (CF405M Annexin V; FC)

Hadassa Y. Holzapfel, and Marc R. Birtwistle. [Creating Complex Fluorophore spectra on Antibodies Through Combinatorial Labeling](#). Transl Sci, 1-10, (2016). (CF405M, CF488A, CF568, CF640R, FRET, MnS)

Kim, Y. R., et al. [Neutrophils Return to Bloodstream Through the Brain Blood Vessel After Crosstalk With Microglia During LPS-Induced Neuroinflammation](#). Front Cell Dev Biol 8, 613733, (2020), DOI: 10.3389/fcell.2020.613733 (CF405M wheat-germ agglutinin, 2-photon)

Kraus, F., et al. [Quantitative 3D structured illumination microscopy of nuclear structures](#). Nat Protoc 12(5), 1011-1028, (2017), DOI: nprot.2017.020, 10.1038/nprot.2017.020 (CF568 azide, CF405M azide, SIM)

Lee, A. T., et al. [A class of GABAergic neurons in the prefrontal cortex sends long-range projections to the nucleus accumbens and elicits acute avoidance behavior](#). J Neurosci 34(35), 11519-11525, (2014), DOI: 34/35/11519, 10.1523/JNEUROSCI.1157-14.2014 (CF405 mouse anti-biotin; IHC)

Lee, K. S., et al. [Functional Synaptic Architecture of Callosal Inputs in Mouse Primary Visual Cortex](#). Neuron 101(3), 421-428 e425, (2019), DOI: 10.1016/j.neuron.2018.12.005 (CF405M CF647 expansion microscopy)

Lindstrom, A., et al. [Fusion between M2-macrophages and cancer cells results in a subpopulation of radioresistant cells with enhanced DNA-repair capacity](#). Oncotarget 8(31), 51370-51386, (2017), DOI: 10.18632/oncotarget.17986 (CF405M succinimidyl ester)

Loh, L. N., et al. [Dissecting Bacterial Cell Wall Entry and Signaling in Eukaryotic Cells: an Actin-Dependent Pathway Parallels Platelet-Activating Factor Receptor-Mediated Endocytosis](#). MBio 8(1), (2017), DOI: mBio.02030-16, 10.1128/mBio.02030-16 (CF405M wheat germ agglutinin, CF488A succinimidyl ester, CF640R succinimidyl ester; CF640R SIM)

Markaki, Y., et al. [Fluorescence In Situ Hybridization Applications for Super-Resolution 3D Structured Illumination Microscopy](#). Methods Mol Biol 950, 43-64, (2013), DOI: DOI:10.1007/978-1-62703-137-0_4 (CF405M secondary antibodies; super resolution microscopy (3D-SIM); 3D-FISH)

Martinko, A. J., et al. [Switchable assembly and function of antibody complexes in vivo using a small molecule](#). Proc Natl Acad Sci U S A, (2022), DOI: 119. 10.1073/pnas.2117402119 (CF405M succinimidyl ester)

Miron, E., et al. [In Vivo and In Situ Replication Labeling Methods for Super-resolution Structured Illumination Microscopy of Chromosome Territories and Chromatin Domains](#). Methods Mol Biol 1431, 127-140, (2016), DOI: 10.1007/978-1-4939-3631-1_10 (CF405M SIM)

Nam, J. H., et al. [TRPV1 on astrocytes rescues nigral dopamine neurons in Parkinson's disease via CNTF](#). Brain 138(Pt 12), 3610-3622, (2015), DOI: 10.1093/brain/awv297 (CF405M secondaries)

Namburi, P., et al. [A circuit mechanism for differentiating positive and negative associations](#). Nature 520(7549), 675-678, (2015), DOI: nature14366, 10.1038/nature14366 (CF405 Streptavidin)

Nolke, T., et al. [Septins guide microtubule protrusions induced by actin-depolymerizing toxins like Clostridium difficile transferase \(CDT\)](#). Proc Natl Acad Sci U S A 113(28), 7870-7875, (2016), DOI: 10.1073/pnas.1522717113 (CF405M secondary)

Ohgomori, T., et al. [Differential involvement of vesicular and glial glutamate transporters around spinal alpha-motoneurons in the pathogenesis of SOD1\(G93A\) mouse model of amyotrophic lateral sclerosis](#). Neuroscience 356, 114-124, (2017), DOI: 10.1016/j.neuroscience.2017.05.014 (CF405M streptavidin)

Seo, J., et al. [PICASSO allows ultra-multiplexed fluorescence imaging of spatially overlapping proteins without reference spectra measurements](#). Nat Commun 13, 2475, (2022), DOI: 10.1038/s41467-022-30168-z (CF405S, CF405M, CF405L, CF488A, CF568, CF633, CF660R, CF680R)

Sutherland, D. R., et al. [High-sensitivity 5-, 6-, and 7-color PNH WBC assays for both Canto II and Navios platforms](#). Cytometry B Clin Cytom 94(4), 637-651, (2018), DOI: 10.1002/cyto.b.21626 (CF405M succinimidyl ester, flow)

Trofimenko, E., et al. [The endocytic pathway taken by cationic substances requires Rab14 but not Rab5 and Rab7](#). Cell Rep 37, 109945, (2021), DOI: 10.1016/j.celrep.2021.109945 (CF405M, CF488A, CF594, CF647, Mix-n-Stain)

Trofimenko, E., et al. [Identification of a new, Rab14-dependent, endo-lysosomal pathway](#). bioRxiv, (2021), DOI: 10.1101/2021.03.25.436964 (CF405M, CF594, Mix-n-Stain)

Yamada, J., et al. [Perineuronal nets affect parvalbumin expression in GABAergic neurons of the mouse hippocampus](#). Eur J Neurosci 41(3), 368-378, (2015), DOI: 10.1111/ejn.12792 (CF405 streptavidin)

Zhang, R., et al. [The mechanisms of dynamin-actin interaction](#). bioRxiv, (2019), DOI: 10.1101/586461 (CF405M goat anti-rabbit, STED)

CF®405S

Demmerle, J., et al. [Strategic and practical guidelines for successful structured illumination microscopy](#). Nat Protoc 12(5), 988-1010, (2017), DOI: nprot.2017.019 [pii] 10.1038/nprot.2017.019 (CF405S and CF405M SIM)

Essig, K., et al. [Roquin Suppresses the PI3K-mTOR Signaling Pathway to Inhibit T Helper Cell Differentiation and Conversion of Treg to Tfr Cells](#). Immunity, (2017), DOI: 10.1016/j.immuni.2017.11.008 (CF405S SIM)

Gothe, H. J., et al. [Spatial Chromosome Folding and Active Transcription Drive DNA Fragility and Formation of Oncogenic MLL Translocations](#). Mol Cell 75(2), 267-283 e212, (2019), DOI: 10.1016/j.molcel.2019.05.015 (CF405S dUTP nick translation FISH)

He, X., et al. [Targeting the Endocannabinoid/CB1 Receptor System For Treating Major Depression Through Antidepressant Activities of Curcumin and Dexanabinol-Loaded Solid Lipid Nanoparticles](#). Cell Physiol Biochem 42(6), 2281-2294, (2017), DOI: 10.1159/000480001 (CF405S succinimidyl ester, RedDot)

Irie, T. [Essential Role of Somatic Kv2 Channels in High-Frequency Firing in Cartwheel Cells of the Dorsal Cochlear Nucleus](#). eNeuro, (2021), DOI: 10.1523/ENEURO.0515-20.2021 (CF405S IgG2a CF568 IgG2b)

Hwang, J., et al. [Aptamer-conjugated live human immune cell based biosensors for the accurate detection of C-reactive protein](#). Sci Rep 6, 34778, (2016), DOI: 10.1038/srep34778 (CF594 Mix-n-Stain)

Jankowska, K. I., et al. [Integrins Modulate T Cell Receptor Signaling by Constraining Actin Flow at the Immunological Synapse](#). Front Immunol 9, 25, (2018), DOI: 10.3389/fimmu.2018.00025 (CF488A Mix-n-Stain, CF405 phalloidin)

Kohara, K., et al. [Cell type-specific genetic and optogenetic tools reveal hippocampal CA2 circuits](#). Nat Neurosci 17(2), 269-279, (2014), DOI: nn.3614 [pii] 10.1038/nn.3614 (CF633 goat anti-mouse, CF633 goat anti-rabbit, CF633 SA, CF405S goat anti-mouse; IHC)

MacDonald, A. B. [Lyme ELISA seronegative patients with borrelia antibody coated borrelia microbes in their blood](#). Med Clin Res 7, (2022), ISSN: 2577 - 8005 (CF405S, CF660R, Mix-n-Stain)

Nazari, B., et al. [Altered Dermal Fibroblasts in Systemic Sclerosis Display Podoplanin and CD90](#). Am J Pathol 186(10), 2650-2664, (2016), DOI: 10.1016/j.ajpath.2016.06.020 (CF405S tyramide)

Ruger, M., et al. [Species-specific viability analysis of Pseudomonas aeruginosa, Burkholderia cepacia and Staphylococcus aureus in mixed culture by flow cytometry](#). BMC Microbiol 14, 56, (2014), DOI: 1471-2180-14-56 [pii] 10.1186/1471-2180-14-56 (CF405S Mix-n-Stain; Flow Cytometry)

Ruger, M., et al. [A flow cytometric method for viability assessment of Staphylococcus aureus and Burkholderia cepacia in mixed culture](#). Cytometry A 81(12), 1055, (2012), DOI: 10.1002/cyto.a.22219 (CF405S Mix-n-Stain labeling of wheat germ agglutinin; Flow Cytometry)

Schneider, K., et al. [Dissection of cell cycle-dependent dynamics of Dnmt1 by FRAP and diffusion-coupled modeling](#). Nucleic Acids Res 41(9), 4860-4876, (2013), DOI: gkt191 [pii] 10.1093/nar/gkt191 (CF405S goat anti-mouse; FRAP)

Seo, J., et al. [PICASSO allows ultra-multiplexed fluorescence imaging of spatially overlapping proteins without reference spectra measurements](#). Nat Commun 13, 2475, (2022), DOI: 10.1038/s41467-022-30168-z (CF405S, CF405M, CF405L, CF488A, CF568, CF633, CF660R, CF680R)

Torkzadeh, H., et al. [Quantification and modeling of the response of surface biofilm growth to continuous low intensity UVC irradiation](#). Water Res 193, 116895, (2021), DOI: 10.1016/j.watres.2021.116895 (CF405S ConA Biofilm)

Ugolini, I., et al. [Chromatin localization of nucleophosmin organizes ribosome biogenesis](#). Mol Cell 82, 4443-4457 e4449, (2022), DOI: 10.1016/j.molcel.2022.10.033 (CF405S maleimide)

Van Dyke, J. M., et al. [Macrophage-mediated inflammation and glial response in the skeletal muscle of a rat model of familial amyotrophic lateral sclerosis \(ALS\)](#). *Exp Neurol* 277, 275-282, (2016), DOI: 10.1016/j.expneurol.2016.01.008 (CF405S bungarotoxin)

Vasilyeva, N. A., et al. [Influence of Membrane Receptor Lateral Diffusion on the Short-Term Depression of Acetylcholine-Induced Current in Helix Neurons](#). *Cell Mol Neurobiol* 37(8), 1443-1455, (2017), DOI: 10.1007/s10571-017-0475-3 (CF405S bungarotoxin FRAP)

Yamada, J., et al. [Alterations in expression of Cat-315 epitope of perineuronal nets during normal ageing, and its modulation by an open-channel NMDA receptor blocker, memantine](#). *J Comp Neurol* 525(9), 2035-2049, (2017), DOI: 10.1002/cne.24198 (CF405 streptavidin)

Yu, C. H., et al. [Integrin-beta3 clusters recruit clathrin-mediated endocytic machinery in the absence of traction force](#). *Nat Commun* 6, 8672, (2015), DOI: 10.1038/ncomms9672 (CF680R maleimide and CF405 phalloidin)

CF®405L

Seo, J., et al. [PICASSO allows ultra-multiplexed fluorescence imaging of spatially overlapping proteins without reference spectra measurements](#). *Nat Commun* 13, 2475, (2022), DOI: 10.1038/s41467-022-30168-z (CF405S, CF405M, CF405L, CF488A, CF568, CF633, CF660R, CF680R)

CF®488A

Agrawal, A., et al. [Manipulation of coacervate droplets with an electric field](#). *Proc Natl Acad Sci U S A* 119, e2203483119 (2022), DOI: 10.1073/pnas.2203483119 (CF488A BSA, CF640R BSA, fluorescence microscopy)

Ahi, Y. S., et al. [Adenoviral E2 IVa2 Protein Interacts with L4 33K Protein and E2 DNA-Binding Protein](#). *J Gen Virol* 94(6), 1325, (2013), DOI: DOI:10.1099/vir.0.049346-0 (Mix-n-Stain CF350, Mix-n-Stain CF488A, CF568 goat anti-mouse IgG; ICC)

Ahi, Y. S., et al. [Functional Interplay Between Murine Leukemia Virus Glycag, Serinc5, and Surface Glycoprotein Governs Virus Entry, with Opposite Effects on Gammaretroviral and Ebolavirus Glycoproteins](#). *mBio* 7(6), (2016), DOI: 10.1128/mBio.01985-16 (CF488A goat anti-rabbit & CF594 donkey anti-mouse)

Akhmetshina, A., et al. [Activation of canonical Wnt signalling is required for TGF-beta-mediated fibrosis](#). *Nat Commun* 3, 735, (2012), DOI: DOI:10.1038/ncomms1734 (CF488A and CF594 streptavidin; IHC, ICC)

Al Tanoury, Z., et al. [Genes involved in cell adhesion and signaling: a new repertoire of retinoic acid receptor target genes in mouse embryonic fibroblasts](#). *J Cell Sci* 127(Pt 3), 521-533, (2014), DOI: jcs.131946 [pii] 10.1242/jcs.131946 (CF488A phalloidin; ICC)

Alissafi, T., et al. [De novo-induced self-antigen-specific Foxp3+ regulatory T cells impair the accumulation of inflammatory dendritic cells in draining lymph nodes](#). *J Immunol* 194(12), 5812-5824, (2015), DOI: 10.4049/jimmunol.1500111 (CF488A goat anti-rabbit)

Amorosi, C. J., et al. [Massively parallel characterization of CYP2C9 variant enzyme activity and abundance](#). *Am J Hum Genet* 108, 1735-1751, (2021), DOI: 10.1016/j.ajhg.2021.07.001 (CF488A picolyl azide)

Anderson, M. S., et al. [Maturation and vesicle-mediated egress of primate gammaherpesvirus rhesus monkey rhadinovirus require inner tegument protein ORF52](#). J Virol 88(16), 9111-9128, (2014), DOI: JVI.01502-14 [pii] 10.1128/JVI.01502-14 (CF488A Mix n Stain; ICC)

Anestopoulos, I., et al. [A Novel Role of Silibinin as a Putative Epigenetic Modulator in Human Prostate Carcinoma](#). Molecules 22(1), (2016), DOI: 10.3390/molecules22010062 (GaR CF405S, goat anti-rabbit CF488A)

Angelov, B. and Angelova, A. [Nanoscale clustering of the neurotrophin receptor TrkB revealed by super-resolution STED microscopy](#). Nanoscale 9(28), 9797-9804, (2017), DOI: 10.1039/c7nr03454g (CF488A STED)

Antolín-Urbaneja, J. C., et al. [Compact and cost effective instrument for detecting drug precursors in different environments based on fluorescence polarization](#). Paper presented at: Optical Sensors 2013, (Prague, Czech Republic, Proc SPIE), (2013), DOI: DOI:10.1117/12.2017257 (CF488A succinimidyl ester; fluorescence polarization assay)

Aravindan, S., et al. [Polyphenols from marine brown algae target radiotherapy-coordinated EMT and stemness-maintenance in residual pancreatic cancer](#). Stem Cell Res Ther 6, 182, (2015), DOI: 10.1186/s13287-015-0173-3 (Mix-n-Stain CF488A)

Arul, M., et al. [Culture of low passage colorectal cancer cells and demonstration of variation in selected tumour marker expression](#). Cytotechnology 66(3), 481, (2013), DOI: 10.1007/s10616-013-9600-4 (CF488A Mix-n-Stain, CF594 Mix-n-Stain, CF488A goat anti-mouse; ICC)

Bargul, J. L., et al. [Species-Specific Adaptations of Trypanosome Morphology and Motility to the Mammalian Host](#). PLoS Pathog 12(2), e1005448, (2016), DOI: 10.1371/journal.ppat.1005448 (anti-Biotin CF488A)

Bekiaris, C., et al. [Neurogenesis in the septal and temporal part of the adult rat dentate gyrus](#). Hippocampus 25(4), 511-523, (2015), DOI: 10.1002/hipo.22388 (CF488A goat anti-mouse & CF555 goat anti-rabbit)

Bohara, M., et al. [C-type natriuretic peptide modulates permeability of the blood-brain barrier](#). J Cereb Blood Flow Metab 34(4), 589-596, (2014), DOI: jcbfm2013234 [pii] 10.1038/jcbfm.2013.234 (CF488A goat anti-mouse; ICC)

Bonifacio, G., et al. [Protonation controls ASIC1a activity via coordinated movements in multiple domains](#). J Gen Physiol 143(1), 105-118, (2014), DOI: jgp.201311053 [pii] 10.1085/jgp.201311053 (CF488A maleimide; voltage clamp)

Briceno, M. P., et al. [Toxoplasma gondii Infection Promotes Epithelial Barrier Dysfunction of Caco-2 Cells](#). J Histochem Cytochem 64(8), 459-469, (2016), DOI: 10.1369/0022155416656349 (CF488A goat anti-rabbit)

Burgueno, J. F., et al. [TLR2 and TLR9 modulate enteric nervous system inflammatory responses to lipopolysaccharide](#). J Neuroinflammation 13(1), 187, (2016), DOI: 10.1186/s12974-016-0653-0 (CF488A donkey anti-chicken & CF488A donkey anti-mouse)

Cantuti-Castelvetri, L., et al. [Neuropilin-1 facilitates SARS-CoV-2 cell entry and infectivity](#). Science 370(6518), 856-860, (2020), DOI: 10.1126/science.abd2985 (CF488A, CF555, CF647)

Cao, F. Y., et al. [Evaluating the effects of charged oligopeptide motifs coupled with RGD on osteogenic differentiation of mesenchymal stem cells](#). ACS Appl Mater Interfaces 7(12), 6698-6705, (2015), DOI: 10.1021/acsami.5b00064 (CF488A phalloidin)

Cao, F. Y., et al. [A novel function of BMHP1 and cBMHP1 peptides to induce the osteogenic differentiation of mesenchymal stem cells](#). Biomater Sci 3(2), 345-351, (2015), DOI: 10.1039/c4bm00300d (CF488A phalloidin)

Chen, Z. H., et al. [Characterization of interstitial Cajal progenitors cells and their changes in Hirschsprung's disease](#). PLoS One 9(1), e86100, (2014), DOI: 10.1371/journal.pone.0086100 PONE-D-13-26924 [pii] (CF488A donkey anti-rabbit, CF543 BAG, CF633 donkey anti-mouse; IHC)

Daly, D. T. and Ariel, M. [A novel cerebellar commissure and other myelinated axons in the Purkinje cell layer of a pond turtle \(Trachemys scripta elegans\)](#). J Comp Neurol 526(17), 2802-2823, (2018), DOI: 10.1002/cne.24528 (CF488A and CF594 10K dextran)

Darvish-Damavandi, M., et al. [Towards the development of an enzyme replacement therapy for the metabolic disorder propionic acidemia](#). Mol Genet Metab Rep 8, 51-60, (2016), DOI: 10.1016/j.ymgmr.2016.06.009, S2214-4269(16)30048-9 [pii] (CF488A Mix-n-Stain)

Dasgupta, P., et al. [Rapid assessment of G protein signaling of four opioid receptors using a real-time fluorescence-based membrane potential assay](#). Eur J Pharmacol 890, 173640, (2021), DOI: 10.1016/j.ejphar.2020.173640 (CF488A Anti-6X His CF640R Anti-FLAG)

Decembre, E., et al. [Sensing of immature particles produced by dengue virus infected cells induces an antiviral response by plasmacytoid dendritic cells](#). PLoS Pathog 10(10), e1004434, (2014), DOI: 10.1371/journal.ppat.1004434 PPATHOGENS-D-14-00241 [pii] (CF488A phalloidin)

Di Giovanni, S., et al. [Determination of benzyl methyl ketone - a commonly used precursor in amphetamine manufacture](#). Analytical Methods 4(11), 3558-3564, (2012), DOI: DOI:10.1039/c2ay25772f (CF488A succinimidyl ester; fluorescence polarization assay)

Diamantopoulou, A., et al. [Denial or receipt of expected reward through maternal contact during the neonatal period differentially affect the development of the rat amygdala and program its function in adulthood in a sex-dimorphic way](#). Psychoneuroendocrinology 38(9), 1757, (2013), DOI: DOI:10.1016/j.psyneuen.2013.02.012 (Mix-n-Stain CF488A, Mix-N-Stain CF555; IHC)

Donskow-Lyonewska, K., et al. [Heligmosmoides polygyrus fourth stages induce protection against DSS-induced colitis and change opioid expression in the intestine](#). Parasite Immunol 34(11), 536, (2012), DOI: DOI:10.1111/pim.12003 (Mix-n-Stain CF488A; IHC)

Droste, D., et al. [Ca\(2+\)-permeable AMPA receptors in mouse olfactory bulb astrocytes](#). Sci Rep 7, 44817, (2017), DOI: 10.1038/srep44817 (CF488A donkey anti-guinea pig)

Egashira, A., et al. [Kid depletion in mouse oocytes associated with multinucleated blastomere formation and inferior embryo development](#). Anim Sci J 87(8), 1048-1054, (2016), DOI: 10.1111/asj.12530 (CF488A goat anti-rabbit)

Endo, H., et al. [Ibuprofen enhances the anticancer activity of cisplatin in lung cancer cells by inhibiting the heat shock protein 70](#). Cell Death Dis 5, e1027, (2014), DOI: cddis2013550 [pii] 10.1038/cddis.2013.550 (CF488A Annexin V; FC)

Ersek, B., et al. [CD3zeta-Chain Expression of Human T Lymphocytes Is Regulated by TNF via Src-like Adaptor Protein-Dependent Proteasomal Degradation](#). J Immunol 189(4), 1602, (2012), DOI: DOI:10.4049/jimmunol.1102365 (Mix-n-Stain CF488A; ICC)

Falibene, A., et al. [Long-term avoidance memory formation is associated with a transient increase in mushroom body synaptic complexes in leaf-cutting ants](#). Front Behav Neurosci 9, 84, (2015), DOI: 10.3389/fnbeh.2015.00084 (CF488A goat anti-mouse)

Falibene, A., et al. [Daily Thermal Fluctuations Experienced by Pupae via Rhythmic Nursing Behavior Increase Numbers of Mushroom Body Microglomeruli in the Adult Ant Brain](#). Front Behav Neurosci 10, 73, (2016), DOI: 10.3389/fnbeh.2016.00073 (CF488A goat anti-mouse)

Falter, C. and Voigt, C. A. [The Eurotiomycete Apinisia graminicola as the causal agent of a leaf spot disease on the energy crop Miscanthus x giganteus in Northern Germany](#). European Journal of Plant Pathology 149(4), 797-806, (2017), DOI: 10.1007/s10658-017-1226-5 (CF488A wheat germ agglutinin)

Fang, S., et al. [Synthesis, antitumor activity, and mechanism of action of 6-acrylic phenethyl ester-2-pyranone derivatives](#). Org Biomol Chem 13(16), 4714-4726, (2015), DOI: 10.1039/c5ob00007f (CF488A goat anti-rabbit IF)

Farias, G. G., et al. [Imaging the Polarized Sorting of Proteins from the Golgi Complex in Live Neurons](#). Methods Mol Biol 1496, 13-30, (2016), DOI: 10.1007/978-1-4939-6463-5_2 (CF488A, CF555 & CF640R Mix-n-Stain IF)

Francois, A., et al. [The Low-Threshold Calcium Channel Cav3.2 Determines Low-Threshold Mechanoreceptor Function](#). Cell Rep 10(3), 370-382, (2015), DOI: 10.1016/j.celrep.2014.12.042 (CF488A donkey anti-chicken)

Frangeul, L., et al. [A cross-modal genetic framework for the development and plasticity of sensory pathways](#). Nature 538(7623), 96-98, (2016), DOI: nature19770 [pii] 10.1038/nature19770 CF488A hydrazide

Fu, X., et al. [In Vivo Single-Molecule Detection of Nanoparticles for Multiphoton Fluorescence Correlation Spectroscopy to Quantify Cerebral Blood Flow](#). Nano Lett 20(8), 6135-6141, (2020), DOI: 10.1021/acs.nanolett.0c02280 (CF488A dextran 10K 70K 250K CF568 dextran 10K)

Gan, E. S., et al. [Jumonji demethylases moderate precocious flowering at elevated temperature via regulation of FLC in Arabidopsis](#). Nat Commun 5, 5098, (2014), DOI: ncomms6098 [pii] 10.1038/ncomms6098 (CF488A goat anti-rabbit, CF555 goat anti-mouse; ICC)

Geherin, S. A., et al. [IL-10+ Innate-like B Cells Are Part of the Skin Immune System and Require alpha4beta1 Integrin To Migrate between the Peritoneum and Inflamed Skin](#). J Immunol 196(6), 2514-2525, (2016), DOI: 10.4049/jimmunol.1403246 (CF488A CF594 Mix-n-Stain)

Giuntini, S., et al. [Combined roles of human IgG subclass, alternative complement pathway activation, and epitope density in the bactericidal activity of antibodies to meningococcal factor h binding protein](#). Infect Immun 80(1), 187-194, (2012), DOI: DOI:10.1128/IAI.05956-11 (CF488A goat anti-human IgG; FC)

Hadjiantoniou, S. V., et al. [Physical confinement signals regulate the organization of stem cells in three dimensions](#). J R Soc Interface 13(123), (2016), DOI: 10.1098/rsif.2016.0613 (CF488A ARt)

Halstead, J. M., et al. [TRICK: A Single-Molecule Method for Imaging the First Round of Translation in Living Cells and Animals](#). Methods Enzymol 572, 123-157, (2016), DOI: 10.1016/bs.mie.2016.02.027 (CF488A anti-GFP)

Hamazaki, N., et al. [Gene activation-associated long noncoding RNAs function in mouse preimplantation development](#). Development 142(5), 910-920, (2015), DOI: 10.1242/dev.116996 (CF488A donkey anti-mouse IF)

Hadassa Y. Holzapfel, and Marc R. Birtwistle. [Creating Complex Fluorophore spectra on Antibodies Through Combinatorial Labeling](#). Transl Sci, 1-10, (2016). (CF405M, CF488A, CF568, CF640R, FRET, MnS)

He, J., et al. [Cucurbitacin Ila induces caspase-3-dependent apoptosis and enhances autophagy in lipopolysaccharide-stimulated RAW 264.7 macrophages](#). Int Immunopharmacol, (2013), DOI: DOI:10.1016/j.intimp.2013.03.013 (CF488A goat anti-mouse IgG, CF568 goat anti-rabbit IgG; ICC)

Higa, M., et al. [TRF2 recruits ORC through TRFH domain dimerization](#). Biochim Biophys Acta Mol Cell Res 1864(1), 191-201, (2017), DOI: 10.1016/j.bbamcr.2016.11.004 (CF488A goat anti-rabbit, CF594 DRRt, CF594 goat anti-mouse)

Hirano, K. and Namihira, M. [LSD1 Mediates Neuronal Differentiation of Human Fetal Neural Stem Cells by Controlling the Expression of a Novel Target Gene, HEYL](#). Stem Cells 34(7), 1872-1882, (2016), DOI: 10.1002/stem.2362 (CF488A, CF568, CF647 donkey secondaries)

Hirtz, M., et al. [Click-Chemistry Immobilized 3D-Infused Microarrays in Nanoporous Polymer Substrates](#). Advanced Materials Interfaces, n/a-n/a, (2016), DOI: 10.1002/admi.201500469 (CF488A Mix-n-Stain)

Ho, A. K., et al. [The topology, structure and PE interaction of LITAF underpin a Charcot-Marie-Tooth disease type 1C](#). BMC Biol 14(1), 109, (2016), DOI: 10.1186/s12915-016-0332-8 (CF488A wheat germ agglutinin)

Hoepflinger, M. C., et al. [Molecular and biochemical analysis of the first ARA6 homologue, a RAB5 GTPase, from green algae](#). J Exp Bot 64(18), 5553-5568, (2013), DOI: ert322 [pii] 10.1093/jxb/ert322 (CF488A goat anti-rabbit; ICC)

Home, P., et al. [Altered subcellular localization of transcription factor TEAD4 regulates first mammalian cell lineage commitment](#). Proc Natl Acad Sci U S A 109(19), 7362-7367, (2012), DOI: DOI:10.1073/pnas.1201595109 (Mix-n-Stain CF488A, Mix-n-Stain CF568; ICC)

Hui, T. K., et al. [Ablation of Lrp4 in Schwann Cells Promotes Peripheral Nerve Regeneration in Mice](#). Biology (Basel) 10(6), (2021), DOI: 10.3390/biology10060452 (CF568 CF488A alpha-bungarotoxin)

Inutsuka, A., et al. [Concurrent and robust regulation of feeding behaviors and metabolism by orexin neurons](#). Neuropharmacology 85, 451-460, (2014), DOI: S0028-3908(14)00237-8 [pii] 10.1016/j.neuropharm.2014.06.015 (CF488 CF594 CF647; IHC)

Inutsuka, A., et al. [The integrative role of orexin/hypocretin neurons in nociceptive perception and analgesic regulation](#). Sci Rep 6, 29480, (2016), DOI: 10.1038/srep29480 (CF488A, CF594 & CF647 secondaries)

Ise, H., et al. [Engulfment and clearance of apoptotic cells based on a GlcNAc-binding lectin-like property of surface vimentin](#). Glycobiology 22(6), 788, (2012), DOI: DOI:10.1093/glycob/cws052 (CF488A goat anti-mouse IgG, CF488A donkey anti-rat IgG, CF555 donkey anti-rabbit IgG; ICC, IHC)

Ishizuka, Y. and Bramham, C. R. [A simple DMSO-based method for cryopreservation of primary hippocampal and cortical neurons](#). J Neurosci Methods 333, 108578, (2020), DOI: 10.1016/j.jneumeth.2019.108578 (AM1-43, CF488A donkey anti-chicken)

Ito, H., et al. [Expression analyses of Phactr1 \(phosphatase and actin regulator 1\) during mouse brain development](#). Neurosci Res 128, 50-57, (2018), DOI: 10.1016/j.neures.2017.08.002 (CF488A goat anti-rabbit & CF543 goat anti-mouse)

Jankowska, K. I., et al. [Integrins Modulate T Cell Receptor Signaling by Constraining Actin Flow at the Immunological Synapse](#). Front Immunol 9, 25, (2018), DOI: 10.3389/fimmu.2018.00025 (CF488A Mix-n-Stain, CF405 phalloidin)

Jones, R. A., et al. [NMJ-morph reveals principal components of synaptic morphology influencing structure-function relationships at the neuromuscular junction](#). Open Biol 6(12), (2016), DOI: 10.1098/rsob.160240 (CF488A bungarotoxin)

Kaneda, M., et al. [Synaptic localisation of SRF coactivators, MKL1 and MKL2, and their role in dendritic spine morphology](#). Sci Rep 8(1), 727, (2018), DOI: 10.1038/s41598-017-18905-7 (CF488A goat anti-rabbit, CF488A goat anti-mouse and CF594 GAGp)

Kanno, T., et al. [A novel small molecule, N-\(4-\(2-pyridyl\)\(1,3-thiazol-2-yl\)\)-2-\(2,4,6-trimethylphenoxy\)acetamide, selectively protects against oxidative stress-induced cell death by activating the Nrf2-ARE pathway: Therapeutic implications for ALS](#). Free Radic Biol Med, (2012), DOI: DOI:10.1016/j.freeradbiomed.2012.09.010 (CF488A goat anti-mouse IgG; ICC)

Karagounis, I. V., et al. [Repression of the autophagic response sensitises lung cancer cells to radiation and chemotherapy](#). Br J Cancer 115(3), 312-321, (2016), DOI: 10.1038/bjc.2016.202 (CF488A & CF568 secondaries)

Kasagi, Y., et al. [Peritoneal Dissemination Requires an Sp1-Dependent CXCR4/CXCL12 Signaling Axis and Extracellular Matrix-Directed Spheroid Formation](#). Cancer Res 76(2), 347-357, (2016), DOI: 10.1158/0008-5472.CAN-15-1563 (CF488A donkey anti-goat)

Kawagishi, T., et al. [Reverse Genetics for Fusogenic Bat-Borne Orthoreovirus Associated with Acute Respiratory Tract Infections in Humans: Role of Outer Capsid Protein sigmaC in Viral Replication and Pathogenesis](#). PLoS Pathog 12(2), e1005455, (2016), DOI: 10.1371/journal.ppat.1005455 (CF488A goat anti-mouse)

Kawase, A., et al. [Decreased radixin function for ATP-binding cassette transporters in liver in adjuvant-induced arthritis rats](#). J Pharm Sci 103(12), 4058-4065, (2014), DOI: 10.1002/jps.24210, S0022-3549(15)30300-2 [pii] (CF488A goat anti-rat & CF568 goat anti-mouse)

Kern, F., et al. [Nogo-A couples with Apg-1 through interaction and co-ordinate expression under hypoxic and oxidative stress](#). Biochem J 455(2), 217-227, (2013), DOI: BJ20130579 [pii] 10.1042/BJ20130579 (CF488A GAC; IHC)

Kida, A., et al. [Cell surface-fluorescence immunosorbent assay for real-time detection of hybridomas with efficient antibody secretion at the single-cell level](#). Anal Chem 85(3), 1753-1759, (2013), DOI: DOI:10.1021/ac303067k (CF594 amine; ICC)

Kijani, S., et al. [Filter-Dense Multicolor Microscopy](#). PLoS One 10(3), e0119499, (2015), DOI: 10.1371/journal.pone.0119499 PONE-D-14-20661 [pii] (CF488A, CF543, CF594)

Kikuchi, K., et al. [Rho signaling inhibitor, CCG-1423, inhibits axonal elongation and dendritic complexity of rat cortical neurons](#). Biochem Biophys Res Commun 492(3), 474-479, (2017), DOI: 10.1016/j.bbrc.2017.08.071 (CF488A goat anti-rabbit)

Kim, H. J. and Selvin, P. R. [Fluorescence Imaging with One Nanometer Accuracy](#). in: SpringerReference, Springer, (2013), (CF633; Fluorescence Imaging with One Nanometer Accuracy (FIONA))

Kiratitanaporn, W., et al. [3D printing a biocompatible elastomer for modeling muscle regeneration after volumetric muscle loss](#). Biomater Adv 142, 213171, (2022), DOI: 10.1016/j.bioadv.2022.213171 (CF555 GAMs, CF555 GARb, CF488A phalloidin, fluorescence microscopy)

Kishi, J. Y., et al. [Light-Seq: light-directed in situ barcoding of biomolecules in fixed cells and tissues for spatially indexed sequencing](#). Nat Methods 19, 1393-1402, (2022), DOI: 10.1038/s41592-022-01604-1 (CF488A WGA, fluorescence microscopy)

Kiyono, M., et al. [Expression of the bacterial heavy metal transporter MerC fused with a plant SNARE, SYP121, in Arabidopsis thaliana increases cadmium accumulation and tolerance](#). Planta 235(4), 841-850, (2012), DOI: 10.1007/s00425-011-1543-4 (CF488A goat anti-rabbit IgG; IHC)

Klimas, A., et al. [Nanoscopic Imaging of Human Tissue Sections via Physical and Isotropic Expansion](#). J Vis Exp (151), (2019), DOI: 10.3791/60195 (CF488A CF633 goat anti-chicken)

Knighton, N. J., et al. [Toward cardiac tissue characterization using machine learning and light-scattering spectroscopy](#). J Biomed Opt 26, (2021), DOI: 10.1117/1.JBO.26.11.116001 (CF488A WGA, fluorescence microscopy)

Koman, V. B., et al. [Atomically Thin 2D Interfaces as Sensors for Molecular Permeability through Cellular Layers and Thin Tissues](#). Advanced Functional Materials, (2022), DOI: 10.1002/adfm.202109598 (CF488A wheat germ agglutinin)

Komura, K., et al. [Dynamic behaviors of vimentin induced by interaction with GlcNAc molecules](#). Glycobiology 22(12), 1741, (2012), DOI: DOI:10.1093/glycob/cws118 (CF488A goat anti-mouse IgG; ICC)

Kubickova, J., et al. [Haloperidol Affects Plasticity of Differentiated NG-108 Cells Through sigma1R/IP3R1 Complex](#). Cell Mol Neurobiol 38(1), 181-194, (2018), DOI: 10.1007/s10571-017-0524-y (CF488A goat anti-rabbit)

Kumaran, G. K. and Hanukoglu, I. [Identification and classification of epithelial cells in nephron segments by actin cytoskeleton patterns](#). FEBS J 287(6), 1176-1194, (2019), DOI: <https://doi.org/10.1111/febs.15088> (CF488A phalloidin & CF555 wheat germ agglutinin)

Kume, K., et al. [alpha-Amanitin Restrains Cancer Relapse from Drug-Tolerant Cell Subpopulations via TAF15](#). Sci Rep 6, 25895, (2016), DOI: 10.1038/srep25895 (CF488A goat anti-mouse)

Kume, K. and Nishizuka, S. S. [Colony Lysate Arrays for Proteomic Profiling of Drug-Tolerant Persisters of Cancer Cell](#). Anal Chem 89(17), 8626-8631, (2017), DOI: 10.1021/acs.analchem.7b01215 (CF488A goat anti-mouse & CF594 goat anti-rabbit)

Lasher, R. A., et al. [Electrical stimulation directs engineered cardiac tissue to an age-matched native phenotype](#). J Tissue Eng 3(1), 2041731412455354, (2012), DOI: DOI:10.1177/2041731412455354 (CF488A wheat germ agglutinin; IHC)

Lencesova, L., et al. [Changes and role of adrenoceptors in PC12 cells after phenylephrine administration and apoptosis induction](#). Neurochem Int 57(8), 884-892, (2010), DOI: DOI:10.1016/j.neuint.2010.09.007 (CF488A goat anti-rabbit IgG; ICC)

Li, G., et al. [Patching of Lipid Rafts by Molecular Self-Assembled Nanofibrils Suppresses Cancer Cell Migration](#). Chem 2(2), 283-298, (2017), DOI: 10.1016/j.chempr.2017.01.002 (CF488A goat anti-mouse)

Li, W., et al. [The Effect of Layer-by-Layer Assembly Coating on the Proliferation and Differentiation of Neural Stem Cells](#). ACS Appl Mater Interfaces 7(5), 3018-3029, (2015), DOI: 10.1021/am504456t (CF488A Phalloidin)

Li, W., et al. [Elevation of brain magnesium prevents and reverses cognitive deficits and synaptic loss in Alzheimer's disease mouse model](#). J Neurosci 33(19), 8423-8441, (2013), DOI: DOI:10.1523/JNEUROSCI.4610-12.2013 (CF488A goat anti-mouse IgG, CF488A goat anti-rabbit IgG, and CF488A rabbit anti-goat IgG; IHC)

Liedmann, S., et al. [Localization of a TORC1-eIF4F translation complex during CD8\(+\) T cell activation drives divergent cell fate](#). Mol Cell 82, 2401-2414 e2409, (2022), DOI: 10.1016/j.molcel.2022.04.016 (CF488A DARt, CF568 DAGt, CF568 DARt, super-resolution, STORM)

Lim, A. K., et al. [The nuage mediates retrotransposon silencing in mouse primordial ovarian follicles](#). Development 140(18), 3819-3825, (2013), DOI: dev.099184 [pii] 10.1242/dev.099184 (CF488A Mix-n-Stain; IHC)

Lindberg, O. R., et al. [Epidermal growth factor treatment of the adult brain subventricular zone leads to focal microglia/macrophage accumulation and angiogenesis](#). Stem Cell Reports 2(4), 440-448, (2014), DOI: 10.1016/j.stemcr.2014.02.003, S2213-6711(14)00055-1 [pii] (CF488A donkey anti-mouse CF555 donkey anti-rabbit, donkey anti-goat CF633 donkey anti-goat; IHC)

Lindberg, O. R., et al. [EGF-Induced Expansion of Migratory Cells in the Rostral Migratory Stream](#). PLoS One 7(9), e46380, (2012), DOI: DOI:10.1371/journal.pone.0046380 (CF488A donkey anti-mouse IgG; CF555 donkey anti-goat IgG; CF555 donkey anti-rabbit IgG; CF633 donkey anti-goat IgG; IHC)

Linhoff, M. W., et al. [A high-resolution imaging approach to investigate chromatin architecture in complex tissues](#). Cell 163(1), 246-255, (2015), DOI: S0092-8674(15)01122-8 [pii] 10.1016/j.cell.2015.09.002 (CF488A goat anti-mouse CF488A goat anti-chicken CF647 goat anti-mouse)

Liu, K. P., et al. [LC3B-II deacetylation by histone deacetylase 6 is involved in serum-starvation-induced autophagic degradation](#). Biochem Biophys Res Commun 441(4), 970-975, (2013), DOI: S0006-291X(13)01880-9 [pii] 10.1016/j.bbrc.2013.11.007 (CF488A goat anti-mouse CF568 goat anti-rabbit; ICC)

Locatelli, M., et al. [DNA damage reduces heterogeneity and coherence of chromatin motions](#). Proc Natl Acad Sci U S A 119, e2205166119, (2022), DOI: 10.1073/pnas.2205166119 (CF488A-dUTP)

Loh, L. N., et al. [Dissecting Bacterial Cell Wall Entry and Signaling in Eukaryotic Cells: an Actin-Dependent Pathway Parallels Platelet-Activating Factor Receptor-Mediated Endocytosis](#). MBio 8(1), (2017), DOI: mBio.02030-16 [pii] 10.1128/mBio.02030-16 (CF405M wheat germ agglutinin, CF488A succinimidyl ester, CF640R succinimidyl ester; CF640R SIM)

Loftus, M. S., et al. [A Conserved Leucine Zipper Motif in Gammaherpesvirus ORF52 Is Critical for Distinct Microtubule Rearrangements](#). J Virol 91(17), (2017), DOI: 10.1128/JVI.00304-17 (CF488A Mix-n-Stain)

Lourbopoulos, A., et al. [Changes in thyroid hormone receptors after permanent cerebral ischemia in male rats](#). J Mol Neurosci 54(1), 78-91, (2014), DOI: 10.1007/s12031-014-0253-3 (CF488A goat anti-rabbit, CF488A donkey anti-rabbit, CF555 goat anti-mouse, CF555 donkey anti-goat, EverBrite; IHC)

Maenohara, S., et al. [Role of UHRF1 in de novo DNA methylation in oocytes and maintenance methylation in preimplantation embryos](#). PLoS Genet 13(10), e1007042, (2017), DOI: 10.1371/journal.pgen.1007042 (CF488A donkey anti-rabbit, CF488A donkey anti-rat and CF594 donkey anti-rabbit)

Mahassni, S. H. and Al-Reemi, R. M. [Apoptosis and necrosis of human breast cancer cells by an aqueous extract of garden cress \(Lepidium sativum\) seeds](#). Saudi Journal of Biological Sciences 20(2), 131-139, (2013), DOI: DOI: http://dx.doi.org/10.1016/j.sjbs.2012.12.002 (CF488A Annexin V; ICC, FC)

Martineau, L., et al. [Lymphoblastoids cell lines - Derived iPSC line from a 26-year-old myotonic dystrophy type 1 patient carrying \(CTG\)200 expansion in the DMPK gene: CHUQi01-A](#). Stem Cell Res 26, 103-106, (2018), DOI: 10.1016/j.scr.2017.12.010 (CF488A goat anti-mouse)

Matsuda, K., et al. [Granule cells control recovery from classical conditioned fear responses in the zebrafish cerebellum](#). Sci Rep 7(1), 11865, (2017), DOI: 10.1038/s41598-017-10794-0 (CF488A goat anti-rabbit)

Meadows, M. K., et al. [Boronic Acid Mediated Coupling of Catechols and N-Hydroxylamines: A Bioorthogonal Reaction to Label Peptides](#). Org Lett 19(12), 3179-3182, (2017), DOI: 10.1021/acs.orglett.7b01198 (CF488A aminoxy)

Mercier, L., et al. [In vivo imaging of skeletal muscle in mice highlights muscle defects in a model of myotubular myopathy](#). Intravital 5(1), e1168553, (2016), DOI: 10.1080/21659087.2016.1168553 (CF488A alpha-bungarotoxin, 2-photon)

Miyagishima, S. Y., et al. [DipM is required for peptidoglycan hydrolysis during chloroplast division](#). BMC Plant Biol 14, 57, (2014), DOI: 1471-2229-14-57 [pii] 10.1186/1471-2229-14-57 (CF488A Mix-n-Stain, CF555 Mix-n-Stain; ICC)

Model, M. A. [Methods for cell volume measurement](#). Cytometry A 93(3), 281-296, (2018), DOI: 10.1002/cyto.a.23152 (CF488A 10K Dextran)

Mohammadpour, R., et al. [Comparative Endocytosis Mechanisms and Anticancer Effect of HPMA Copolymer- and PAMAM Dendrimer-MTCP Conjugates for Photodynamic Therapy](#). Macromol Biosci 17(4), (2017), DOI: 10.1002/mabi.201600333 (CF488A wheat germ agglutinin)

Montesinos-Rongen, M., et al. [Primary Central Nervous System \(CNS\) Lymphoma B Cell Receptors Recognize CNS Proteins](#). J Immunol 195(3), 1312-1319, (2015), DOI: jimmunol.1402341 [pii] 10.4049/jimmunol.1402341 (Mix-n-Stain CF488A)

Morita, S., et al. [Targeted DNA demethylation in vivo using dCas9-peptide repeat and scFv-TET1 catalytic domain fusions](#). Nat Biotechnol 34(10), 1060-1065, (2016), DOI: 10.1038/nbt.3658 (CF488A donkey anti-chicken & CF543 donkey anti-rabbit)

Myler, L. R., et al. [Single-molecule imaging reveals the mechanism of Exo1 regulation by single-stranded DNA binding proteins](#). Proc Natl Acad Sci U S A 113(9), E1170-1179, (2016), DOI: 10.1073/pnas.1516674113 (CF488A anti-biotin)

Nabeshima, Y., et al. [Calpain 1 inhibitor BDA-410 ameliorates alpha-klotho-deficiency phenotypes resembling human aging-related syndromes](#). Sci Rep 4, 5847, (2014), DOI: srep05847 [pii] 10.1038/srep05847 (CF594 goat anti-rat & CF488 goat anti-chicken; IHC)

Nagata, Y., et al. [Interleukin-1beta \(IL-1beta\)-induced Notch ligand Jagged1 suppresses mitogenic action of IL-1beta on human dystrophic myogenic cells](#). PLoS One 12(12), e0188821, (2017), DOI: 10.1371/journal.pone.0188821 (CF488A anti-Sheep)

Nahidazar, L. and Harkes, R. [Multicolor Localization-Based Super Resolution Microscopy](#). Methods Mol Biol 2350, 69-76, (2021), DOI: 10.1007/978-1-0716-1593-5_5 (CF488A, CF680, CF660C, CF514 SMLM)

Nakashima, H., et al. [Canonical TGF-beta Signaling Negatively Regulates Neuronal Morphogenesis through TGIF/Smad Complex-Mediated CRMP2 Suppression](#). J Neurosci 38(20), 4791-4810, (2018), DOI: 10.1523/JNEUROSCI.2423-17.2018 (CF488A, CF555, CF647 secondaries)

Nakashima, H., et al. [MeCP2 controls neural stem cell fate specification through miR-199a-mediated inhibition of BMP-Smad signaling](#). Cell Rep 35(7), 109124, (2021), DOI: 10.1016/j.celrep.2021.109124 (CF488A donkey anti-rabbit CF488A donkey anti-goat CF555 donkey anti-mouse CF555 donkey anti-rabbit CF555 donkey anti-guinea pig CF647 donkey anti-mouse and CF647 donkey anti-rat)

Niikura, K., et al. [Endosomal escape efficiency of fusogenic B18 and B55 peptides fused with anti-EGFR single chain Fv as estimated by nuclear translocation](#). J Biochem 159(1), 123-132, (2016), DOI: 10.1093/jb/mvv083 (CF488A goat anti-mouse)

Nishimura, T., et al. [DAAM1 stabilizes epithelial junctions by restraining WAVE complex-dependent lateral membrane motility](#). J Cell Biol 215(4), 559-573, (2016), DOI: 10.1083/jcb.201603107 (CF488A donkey anti-rat)

Noguchi, H., et al. [Expression of DNMT1 in neural stem/precursor cells is critical for survival of newly generated neurons in the adult hippocampus](#). Neurosci Res 95, 1-11, (2015), DOI: S0168-0102(15)00033-4 [pii] 10.1016/j.neures.2015.01.014 (CF488A donkey anti-mouse, CF543 donkey anti-rabbit, CF568 donkey anti-rat, CF647 donkey anti-goat, CF647 donkey anti-rabbit)

Nomura, R., et al. [Azidothymidine-triphosphate impairs mitochondrial dynamics by disrupting the quality control system](#). Redox Biol 13, 407-417, (2017), DOI: 10.1016/j.redox.2017.06.011 (CF488A goat anti-rabbit)

Oda, H., et al. [Plasma microRNAs Are Potential Biomarkers of Acute Rejection After Hindlimb Transplantation in Rats](#). Transplant Direct 2(11), e108, (2016), DOI: 10.1097/TXD.0000000000000620 (CF488A and CF543 secondaries)

Ohira, M., et al. [A novel anti-microtubule agent with carbazole and benzohydrazide structures suppresses tumor cell growth in vivo](#). Biochim Biophys Acta 1850(9), 1676-1684, (2015), DOI: 10.1016/j.bbagen.2015.04.013 (CF488A goat anti-rabbit, CF594 goat anti-mouse)

Okere, B., et al. [In vitro differentiation of human amniotic epithelial cells into insulin-producing 3D spheroids](#). Int J Immunopathol Pharmacol 28(3), 390-402, (2015), DOI: 10.1177/0394632015588439 (CF488A goat anti-mouse)

Orme, M. H., et al. [The unconventional myosin CRINKLED and its mammalian orthologue MYO7A regulate caspases in their signalling roles](#). Nat Commun 7, 10972, (2016), DOI: 10.1038/ncomms10972 (CF488A donkey anti-mouse, CF633 donkey anti-rabbit & CF633 donkey anti-goat)

Osteikoetxea, X., et al. [Improved characterization of EV preparations based on protein to lipid ratio and lipid properties](#). PLoS One 10(3), e0121184, (2015), DOI: 10.1371/journal.pone.0121184 PONE-D-14-38407 [pii] (CF488A Mix-n-Stain)

Pang, J., et al. [Step-by-step preparation of mouse eye sections for routine histology, immunofluorescence, and RNA in situ hybridization multiplexing](#). STAR Protoc 2, 100879, (2021), DOI: 10.1016/j.xpro.2021.100879 (CF488A DAMs, CF568 DARb, CF488A DAGt)

Parekh, N. J., et al. [Type I interferon-dependent CCL4 is induced by a cGAS/STING pathway that bypasses viral inhibition and protects infected tissue, independent of viral burden](#). PLoS Pathog 15(10), e1007778, (2019), DOI: 10.1371/journal.ppat.1007778 (CF488A wheat germ agglutinin)

Park, K. S., et al. [A crucial requirement for Hedgehog signaling in small cell lung cancer](#). Nat Med 17(11), 1504-1508, (2011), DOI: DOI:10.1038/nm.2473 (Mix-n-Stain CF488A, Mix-n-Stain CF568, and Mix-n-Stain CF647; ICC)

Petrelli, A., et al. [Efficient Conjugation of Oligosaccharides to Polymer Particles through Furan/Maleimide Diels-Alder Reaction: Application to the Capture of Carbohydrate-Binding Proteins](#). Chembiochem 18(2), 206-212, (2017), DOI: 10.1002/cbic.201600509 (CF488A, CF568 Mix-n-Stain)

Peixoto-Neves, D., et al. [KV7.1 channel blockade inhibits neonatal renal autoregulation triggered by a step decrease in arterial pressure](#). Am J Physiol Renal Physiol 322(2), F197-F207, (2022), DOI: 10.1152/ajprenal.00568.2020 (CF488A donkey anti-rabbit & CF555 donkey anti-goat)

Platt, J. L., et al. [C3d regulates immune checkpoint blockade and enhances antitumor immunity](#). JCI Insight 2(9), (2017), DOI: 10.1172/jci.insight.90201 (CF488A donkey anti-rat and CF555 donkey anti-goat)

Polidoro, G., et al. [Substance P and the neurokinin-1 receptor expression in dog ileum with and without inflammation](#). Res Vet Sci 114, 297-307, (2017), DOI: 10.1016/j.rvsc.2017.06.002 (CF488A donkey anti-mouse)

Pons, M., et al. [Phosphorylation of filamin A regulates chemokine receptor CCR2 recycling](#). J Cell Sci 130(2), 490-501, (2017), DOI: 10.1242/jcs.193821 (CF488A, CF568 Mix-n-Stain)

Prapa, E., et al. [Effect of Anti-Mullerian hormone \(AMH\) and bone morphogenetic protein 15 \(BMP-15\) on steroidogenesis in primary-cultured human luteinizing granulosa cells through Smad5 signalling](#). J Assist Reprod Genet 32(7), 1079-1088, (2015), DOI: 10.1007/s10815-015-0494-2 (CF488A donkey anti-rabbit & CF594 donkey anti-mouse)

Qian, M., et al. [Unraveling of a Neutralization Mechanism by Two Human Antibodies against Conserved Epitopes in the Globular Head of H5 Hemagglutinin](#). J Virol 87(6), 3571-3577, (2013), DOI: DOI:10.1128/JVI.01292-12 (CF488A goat anti-human IgG; IHC)

Rahman, A., et al. [GMAP is an Atg8a-interacting protein that regulates Golgi turnover in Drosophila](#). Cell Rep 39, 110903, (2022), DOI: 10.1016/j.celrep.2022.110903 (CF488A, CF568 secondaries, fluorescence microscopy)

Raynes, J. K., et al. [Coaggregation of kappa-Casein and beta-Lactoglobulin Produces Morphologically Distinct Amyloid Fibrils](#). Small 13(14), (2017), DOI: 10.1002/smll.201603591 (CF488A maleimide CF647 maleimide)

Ren, J., et al. [Sirt1 Protects Subventricular Zone-Derived Neural Stem Cells from DNA Double-Strand Breaks and Contributes to Olfactory Function Maintenance in Aging Mice](#). Stem Cells 40, 493-507, (2022), DOI: 10.1093/stmcls/sxac008 (CF488A, CF543, CF633 secondaries. fluorescence microscopy)

Ren, S., et al. [Anti-proliferative effect of 23,24-dihydrocucurbitacin F on human prostate cancer cells through induction of actin aggregation and cofilin-actin rod formation](#). Cancer Chemother Pharmacol 70(3), 415, (2012), DOI: DOI:10.1007/s00280-012-1921-z (CF488A goat anti-mouse, CF568 goat anti-rabbit; ICC)

Ritzenthaler, J. D., et al. [Nicotinic Acetylcholine Receptors are Sensors for Ethanol in Lung Fibroblasts](#). Alcohol Clin Exp Res 37(6), 914, (2013), DOI: DOI:10.1111/acer.12044 (CF488A a-Bungarotoxin; ICC)

Rizzi, F., et al. [Polyphenon E\(R\), a standardized green tea extract, induces endoplasmic reticulum stress, leading to death of immortalized PNT1a cells by anoikis and tumorigenic PC3 by necroptosis](#). Carcinogenesis 35(4), 828, (2014), DOI: bgt481 [pii] 10.1093/carcin/bgt481 (CF488A goat anti-mouse; IHC)

Ruiz, R., et al. [alpha-Synuclein A30P decreases neurodegeneration and increases synaptic vesicle release probability in CSPalpha-null mice](#). Neuropharmacology 76 Pt A, 106-117, (2014), DOI: S0028-3908(13)00399-7 [pii] 10.1016/j.neuropharm.2013.08.032 (CF488A donkey anti-rabbit & CF488A donkey anti-mouse; IHC)

Russell, D. F., et al. [Lectin binding and gel secretion within Lorenzinian electroreceptors of Polyodon](#). PLoS One 17, e0276854, (2022), DOI: 10.1371/journal.pone.0276854 (CF488A WGA, fluorescence microscopy)

Sadeghinezhad, J., et al. [Neurochemistry of myenteric plexus neurons of bank vole \(*Myodes glareolus*\) ileum](#). Res Vet Sci 95(3), 846-853, (2013), DOI: S0034-5288(13)00278-6 [pii] 10.1016/j.rvsc.2013.07.028 (CF488A goat anti-mouse, goat anti-rabbit, goat anti-rat; IHC)

Saito, M., et al. [Tctex-1 controls ciliary resorption by regulating branched actin polymerization and endocytosis](#). EMBO Rep 18(8), 1460-1472, (2017), DOI: 10.15252/embr.201744204 (CF488A anti-chicken)

Sawaguchi, S., et al. [O-GlcNAc on NOTCH1 EGF repeats regulates ligand-induced Notch signaling and vascular development in mammals](#). Elife 6, (2017), DOI: 10.7554/elife.24419 (CF488A donkey anti-goat, CF640R donkey anti-rat and CF488A streptavidin)

Schafer, N., et al. [Complement Regulator FHR-3 Is Elevated either Locally or Systemically in a Selection of Autoimmune Diseases](#). Front Immunol 7, 542, (2016), DOI: 10.3389/fimmu.2016.00542 (CF488A goat anti-mouse)

Schmidt, C., et al. [The Dual Role of a Polyvalent IgM/IgA-Enriched Immunoglobulin Preparation in Activating and Inhibiting the Complement System](#). Biomedicines 9(7), (2021), DOI: 10.3390/biomedicines9070817 (CF488A DAGt CF568 donkey anti-mouse, western blotting)

Seo, J., et al. [PICASSO allows ultra-multiplexed fluorescence imaging of spatially overlapping proteins without reference spectra measurements](#). Nat Commun 13, 2475, (2022), DOI: 10.1038/s41467-022-30168-z (CF405S, CF405M, CF405L, CF488A, CF568, CF633, CF660R, CF680R)

Sevenich, M., et al. [Development of an alpha-synuclein fibril and oligomer specific tracer for diagnosis of Parkinson's disease, dementia with Lewy bodies and multiple system atrophy](#). Neurochem Int 161, 105422, (2022), DOI: 10.1016/j.neuint.2022.105422 (CF488A SE, CF633 maleimide)

Sheth, V., et al. [Quantifying Intracellular Nanoparticle Distributions with Three-Dimensional Super-Resolution Microscopy](#). ACS Nano 17, 8376-8392, (2023), DOI: 10.1021/acsnano.2c12808 (CF488A WGA, fluorescence microscopy)

Sheu, J. R., et al. [A novel indication of platonin, a therapeutic immunomodulating medicine, on neuroprotection against ischemic stroke in mice](#). Sci Rep 7, 42277, (2017), DOI: 10.1038/srep42277 (CF488A donkey anti-mouse, CF488A donkey anti-rabbit, CF594 donkey anti-mouse and CF594 donkey anti-rabbit)

Shi, L., et al. [Non-thermal plasma induces a stress response in mesothelioma cells resulting in increased endocytosis, lysosome biogenesis and autophagy](#). Free Radic Biol Med 108, 904-917, (2017), DOI: 10.1016/j.freeradbiomed.2017.04.368 (CF488A & CF568 secondaries)

Shibata, T., et al. [Crosslinking of a Peritrophic Matrix Protein Protects Gut Epithelia from Bacterial Exotoxins](#). PLoS Pathog 11(10), e1005244, (2015), DOI: 10.1371/journal.ppat.1005244 (CF488A goat anti-mouse, CF568 goat anti-mouse and CF568 goat anti-rabbit)

Shiimori, M., et al. [A specific set of exon junction complex subunits is required for the nuclear retention of unspliced RNAs in *Caenorhabditis elegans*](#). Mol Cell Biol 33(2), 444-456, (2013), DOI: MCB.01298-12 [pii] 10.1128/MCB.01298-12 (CF488A phalloidin; ICC)

Shimizu, K., et al. [Three-Dimensional Culture Model of Skeletal Muscle Tissue with Atrophy Induced by Dexamethasone](#). Bioengineering (Basel) 4(2), (2017), DOI: 10.3390/bioengineering4020056 (CF488A goat anti-rabbit)

Shiozaki, H. M. and Kazama, H. [Parallel encoding of recent visual experience and self-motion during navigation in Drosophila](#). Nat Neurosci 20(10), 1395-1403, (2017), DOI: 10.1038/nn.4628 (CF488A anti-rat, CF633 anti-mouse and CF633 anti-rabbit)

Shuto, T., et al. [Increase in receptor activator of nuclear factor kappaB ligand/osteoprotegerin ratio in peri-implant gingiva exposed to Porphyromonas gingivalis lipopolysaccharide](#). J Dent Sci 11(1), 8-16, (2016), DOI: 10.1016/j.jds.2015.10.005 (CF488A goat anti-rabbit)

Smith, P. L., et al. [Neonatal peripheral immune challenge activates microglia and inhibits neurogenesis in the developing murine hippocampus](#). Dev Neurosci 36(2), 119-131, (2014), DOI: 000359950 [pii] 10.1159/000359950 (CF555 donkey anti-goat, CF488A donkey anti-chicken & CF647 donkey anti-mouse; IHC)

Somiya, M., et al. [Intracellular trafficking of bio-nanocapsule-liposome complex: Identification of fusogenic activity in the pre-S1 region of hepatitis B virus surface antigen L protein](#). J Control Release 212, 10-18, (2015), DOI: 10.1016/j.jconrel.2015.06.012 (CF488A succinimidyl ester)

Steijven, K., et al. [Learning performance and brain structure of artificially-reared honey bees fed with different quantities of food](#). PeerJ 5, e3858, (2017), DOI: 10.7717/peerj.3858 (CF488A goat anti-mouse)

Stein, E., et al. [In vitro and in vivo uptake study of Escherichia coli Nissle 1917 bacterial ghosts: cell-based delivery system to target ocular surface diseases](#). Invest Ophthalmol Vis Sci 54(9), 6326-6333, (2013), DOI: iovs.13-12044 [pii] 10.1167/iov.13-12044 (CF488A goat anti-rabbit; IHC)

Steinman, M. Q., et al. [Hypothalamic vasopressin systems are more sensitive to the long term effects of social defeat in males versus females](#). Psychoneuroendocrinology 51, 122-134, (2015), DOI: 10.1016/j.psyneuen.2014.09.009 (CF488A donkey anti-guinea pig)

Sundberg, S. C. and Granseth, B. [Cre-expressing neurons in the cortical white matter of Ntsr1-Cre GN220 mice](#). Neurosci Lett 675, 36-40, (2018), DOI: 10.1016/j.neulet.2018.03.053 (CF488A donkey anti-guinea pig)

Takenaka, N., et al. [A critical role of the small GTPase Rac1 in Akt2-mediated GLUT4 translocation in mouse skeletal muscle](#). FEBS J 281(5), 1493-1504, (2014), DOI: 10.1111/febs.12719 (CF488A, CF543, CF647 goat anti-mouse, goat anti-rabbit, goat anti-rat; ICC)

Takenaka, N., et al. [Rac1 Activation Caused by Membrane Translocation of a Guanine Nucleotide Exchange Factor in Akt2-Mediated Insulin Signaling in Mouse Skeletal Muscle](#). PLoS One 11(5), e0155292, (2016), DOI: 10.1371/journal.pone.0155292 (CF405, CF488A, CF543 & CF647 IHF)

Tanaka, K., et al. [A novel acylaminoimidazole derivative, WN1316, alleviates disease progression via suppression of glial inflammation in ALS mouse model](#). PLoS One 9(1), e87728, (2014), DOI: 10.1371/journal.pone.0087728, PONE-D-13-43248 [pii] (CF488A goat anti-mouse; ICC)

Tapia-Lopez, L. V., et al. [Plasma and chemical treatments of zirconia surface for cell response improvement: A comparative study](#). Materials Letters 303, (2021), DOI: 10.1016/j.matlet.2021.130518 (CF488A phalloidin)

Thiede, B. R. and Corwin, J. T. [Permeation of fluorophore-conjugated phalloidin into live hair cells of the inner ear is modulated by P2Y receptors](#). J Assoc Res Otolaryngol 15(1), 13-30, (2014), DOI: 10.1007/s10162-013-0425-9 (CF488A phalliodin, CF488A amine)

Toth, R., et al. [CpG distribution and methylation pattern in porcine parvovirus](#). PLoS One 8(12), e85986, (2013), DOI: 10.1371/journal.pone.0085986, PONE-D-13-31047 [pii] (CF488A goat anti-rabbit; ICC)

Trofimenko, E., et al. [The endocytic pathway taken by cationic substances requires Rab14 but not Rab5 and Rab7](#). Cell Rep 37, 109945, (2021), DOI: 10.1016/j.celrep.2021.109945 (CF405M, CF488A, CF594, CF647, Mix-n-Stain)

Troidl, K., et al. [Actin-binding rho activating protein \(Abra\) is essential for fluid shear stress-induced arteriogenesis](#). Arterioscler Thromb Vasc Biol 29(12), 2093-2101, (2009), DOI: DOI:10.1161/ATVBAHA.109.195305 (CF488A goat anti-rabbit IgG; ICC, IHC)

von Walden, F., et al. [Mechanical loading induces the expression of a Pol I regulon at the onset of skeletal muscle hypertrophy](#). Am J Physiol Cell Physiol 302(10), C1523-1530, (2012), DOI: DOI:10.1152/ajpcell.00460.2011 (CF488A goat anti-mouse IgG, CF633 goat anti-rabbit IgG; ICC, IHC)

Voulgaridou, G. P., et al. [Aldehyde dehydrogenase 3A1 promotes multi-modality resistance and alters gene expression profile in human breast adenocarcinoma MCF-7 cells](#). Int J Biochem Cell Biol 77(Pt A), 120-128, (2016), DOI: 10.1016/j.biocel.2016.06.004 (CF488A goat anti-mouse)

Vullo, S., et al. [Conformational dynamics and role of the acidic pocket in ASIC pH-dependent gating](#). Proc Natl Acad Sci U S A 114(14), 3768-3773, (2017), DOI: 10.1073/pnas.1620560114 (CF488A maleimide)

Wang, K., et al. [Ultralow-Fouling Zwitterionic Polyurethane-Modified Membranes for Rapid Separation of Plasma from Whole Blood](#). Langmuir 37(33), 10115-10125, (2021), DOI: 10.1021/acs.langmuir.1c01477 (CF488A wheat germ agglutinin)

Wang, X., et al. [Quantitative proteomic analysis of age-related subventricular zone proteins associated with neurodegenerative disease](#). Sci Rep 6, 37443, (2016), DOI: 10.1038/srep37443 (CF488A & CF543 secondaries)

Wang, Y., et al. [Cucurbitacin IIb exhibits anti-inflammatory activity through modulating multiple cellular behaviors of mouse lymphocytes](#). PLoS One 9(2), e89751, (2014), DOI: 10.1371/journal.pone.0089751, PONE-D-13-38697 [pii] (CF488A goat anti-mouse, CF568 goat anti-rabbit; IHC)

Watari, K., et al. [Tumor-derived interleukin-1 promotes lymphangiogenesis and lymph node metastasis through M2-type macrophages](#). PLoS One 9(6), e99568, (2014), DOI: 10.1371/journal.pone.0099568, PONE-D-13-45479 [pii] (CF488A donkey anti-rat & CF594 donkey anti-rabbit; IHC)

Wei, Y., et al. [Synergistic Entry of Individual Nanoparticles into Mammalian Cells Driven by Free Energy Decline and Regulated by Their Sizes](#). ACS Nano 16, 5885-5897, (2022), DOI: 10.1021/acsnano.1c11068 (CF488A, CF555, CF647 SE, nanoparticle labeling, fluorescence microscopy)

West, A. B., et al. [Differential LRRK2 expression in the cortex, striatum, and substantia nigra in transgenic and nontransgenic rodents](#). J Comp Neurol 522(11), 2465-2480, (2014), DOI: 10.1002/cne.23583 (CF555 donkey anti-rabbit & CF488A donkey anti-mouse; IHC)

Wieduwild, R., et al. [Minimal Peptide Motif for Non-covalent Peptide-Heparin Hydrogels](#). J Am Chem Soc 135(8), 2919-2922, (2013), DOI: DOI:10.1021/ja312022u (CF488A phalloidin; IHC)

Wienand, K. and Shires, K. [The use of MAGE C1 and flow cytometry to determine the malignant cell type in multiple myeloma](#). PLoS One 10(3), e0120734, (2015), DOI: 10.1371/journal.pone.0120734, PONE-D-14-44112 [pii] (CF488A flow)

Wu, S., et al. [AUF1 is recruited to the stress granules induced by coxsackievirus B3](#). Virus Res 192, 52-61, (2014), DOI: S0168-1702(14)00331-1 [pii] 10.1016/j.virusres.2014.08.003 (CF488A goat anti-rabbit, CF555 goat anti-mouse & CF555 goat anti-rabbit; ICC)

Xu, Y., et al. [Arabidopsis MRG domain proteins bridge two histone modifications to elevate expression of flowering genes](#). Nucleic Acids Res 42(17), 10960-10974, (2014), DOI: 10.1093/nar/gku781 (CF488A goat anti-rabbit & CF555 goat anti-mouse)

Xu, Z., et al. [Mucosal immunoglobulins at respiratory surfaces mark an ancient association that predates the emergence of tetrapods](#). Nat Commun 7, 10728, (2016), DOI: 10.1038/ncomms10728 (CF488A, CF555 & CF640R Mix-n-Stain)

Xu, Z., et al. [Teleost skin, an ancient mucosal surface that elicits gut-like immune responses](#). Proc Natl Acad Sci U S A 110(32), 13097-13102, (2013), DOI: 1304319110 [pii] 10.1073/pnas.1304319110 (CF488A Mix-n-Stain, CF555, CF640R; IHC)

Yamagami, K., et al. [Preimplantation embryo-secreted factors modulate maternal gene expression in rat uterus](#). Cell Tissue Res 364(2), 453-463, (2016), DOI: 10.1007/s00441-015-2336-z (CF488A goat anti-rabbit)

Yamamoto, S., et al. [Contribution of calumin to embryogenesis through participation in the endoplasmic reticulum-associated degradation activity](#). Dev Biol 393(1), 33-43, (2014), DOI: S0012-1606(14)00320-0 [pii] 10.1016/j.ydbio.2014.06.024 (CF488A Mix-n-Stain, CF568 Mix-n-Stain; IHC)

Yamashita, A., et al. [Increased metabolite levels of glycolysis and pentose phosphate pathway in rabbit atherosclerotic arteries and hypoxic macrophage](#). PLoS One 9(1), e86426, (2014), DOI: 10.1371/journal.pone.0086426, PONE-D-13-24660 [pii] (CF488A Mix-n-Stain, CF594 Mix-n-Stain; IHC)

Yang, J., et al. [Transcriptomic profiling of Alternaria longipes invasion in tobacco reveals pathogenesis regulated by AIHK1, a group III histidine kinase](#). Sci Rep 7(1), 16083, (2017), DOI: 10.1038/s41598-017-16401-6 (CF488A wheat germ agglutinin fungi in plants)

Yang, R., et al. [Cellular biophysical dynamics and ion channel activities detected by AFM-based nanorobotic manipulator in insulinoma beta-cells](#). Nanomedicine 9(5), 636, (2012), DOI: DOI:10.1016/j.nano.2012.10.011 (CF488A phalloidin; ICC)

Yang, W., et al. [Regulation of Meristem Morphogenesis by Cell Wall Synthases in Arabidopsis](#). Curr Biol 26(11), 1404-1415, (2016), DOI: 10.1016/j.cub.2016.04.026 (CF488A & CF568 secondaries)

Yang, Y., et al. [Regulation of Viral RNA Synthesis by the V Protein of Parainfluenza Virus 5](#). J Virol 89(23), 11845-11857, (2015), DOI: 10.1128/JVI.01832-15 (CF488A anti-FLAG)

Yasuda, M., et al. [In utero intraocular AAV injection for early gene expression in the developing rodent retina](#). STAR Protoc 2(3), 100742, (2021), DOI: 10.1016/j.xpro.2021.100742 (CF488A CTB)

Yasuga, H., et al. [Fluid interfacial energy drives the emergence of three-dimensional periodic structures in micropillar scaffolds](#). Nature Physics, (2021), DOI: 10.1038/s41567-021-01204-4 (CF488A GAMs1 and CF488A ConA)

Yi, W., et al. [CRISPR-assisted detection of RNA-protein interactions in living cells](#). Nat Methods 17(7), 685-688, (2020), DOI: 10.1038/s41592-020-0866-0 (CF488A secondary)

Yin, C. and Li, P. [Growth Suppression of Glioma Cells Using HDAC6 Inhibitor, Tubacin](#). Open Med (Wars) 13, 221-226, (2018), DOI: 10.1515/med-2018-0034 (CF488A goat anti-rabbit CF568 goat anti-mouse)

Yoshimura, T., et al. [Mouse GTSF1 is an essential factor for secondary piRNA biogenesis](#). EMBO Rep 19(4), (2018), DOI: 10.15252/embr.201642054 (CF488A succinimidyl ester)

Yuan, L., et al. [Crystal structures reveal catalytic and regulatory mechanisms of the dual-specificity ubiquitin/FAT10 E1 enzyme Uba6](#). Nat Commun 13, 4880, (2022), DOI: 10.1038/s41467-022-32613-5 (CF488A maleimide, enzime kenetics assay)

Zanetti-Domingues, L. C., et al. [Hydrophobic fluorescent probes introduce artifacts into single molecule tracking experiments due to non-specific binding](#). PLoS One 8(9), e74200, (2013), DOI: 10.1371/journal.pone.0074200, PONE-D-13-17594 [pii] (CF488A, CF568, CF633 & CF640R maleimide, TIRF)

Zha, Q. B., et al. [Cucurbitacin E Induces Autophagy via Downregulating mTORC1 Signaling and Upregulating AMPK Activity](#). PLoS One 10(5), e0124355, (2015), DOI: 10.1371/journal.pone.0124355 (CF488A goat anti-mouse & CF568 goat anti-rabbit)

Zhang, M., et al. [A high M1/M2 ratio of tumor-associated macrophages is associated with extended survival in ovarian cancer patients](#). J Ovarian Res 7(1), 19, (2014), DOI: 1757-2215-7-19 [pii], 10.1186/1757-2215-7-19 (CF350, CF488A labeled primaries; IHC)

Zhang, Y., et al. [Deep-brain magnetic stimulation promotes adult hippocampal neurogenesis and alleviates stress-related behaviors in mouse models for neuropsychiatric disorders](#). Mol Brain 7, 11, (2014), DOI: 1756-6606-7-11 [pii] 10.1186/1756-6606-7-11 (CF488A goat anti-mouse & CF555 donkey anti-mouse; IHC)

Zhang, Y. T., et al. [Formation of cofilin-actin rods following cucurbitacin-B-induced actin aggregation depends on slingshot homolog 1-mediated cofilin hyperactivation](#). J Cell Biochem 114(10), 2415, (2013), DOI: DOI:10.1002/jcb.24587 (CF488A goat anti-mouse, CF568 goat anti-rabbit; ICC)

Zhang, Y. T., et al. [VASP activation via the Galpha13/RhoA/PKA pathway mediates cucurbitacin-B-induced actin aggregation and cofilin-actin rod formation](#). PLoS One 9(4), e93547, (2014), DOI: 10.1371/journal.pone.0093547, PONE-D-13-34310 [pii] (CF488A goat anti-mouse, CF568 goat anti-rabbit; ICC)

Zhou, H. and Liu, G. [Regulation of density of functional presynaptic terminals by local energy supply](#). Mol Brain 8, 42, (2015), DOI: 10.1186/s13041-015-0132-z (CF488A, CF555 & CF640R secondaries, MitoView 633)

Zhu, J. S., et al. [Cucurbitacin B induces cell cycle arrest, apoptosis and autophagy associated with G actin reduction and persistent activation of cofilin in Jurkat cells](#). Pharmacology 89(5-6), 348-346, (2012), DOI: DOI:10.1159/000338757 (CF488A goat anti-mouse, CF568 goat anti-rabbit; ICC)

Zong, G., et al. [Total Synthesis and Biological Evaluation of Ipomoeassin F and Its Unnatural 11R-Epimer](#). J Org Chem 80(18), 9279-9291, (2015), DOI: 10.1021/acs.joc.5b01765 (CF488A Annexin V)

CF®514

Chen, K., et al. [Excitation spectral microscopy for highly multiplexed fluorescence imaging and quantitative biosensing](#). bioRxiv, (2021), DOI: 10.1101/2021.04.13.439576 (CF532 wheat germ agglutinin CF514SE and CF568SE)

Chen, K., et al. [Super-multiplexing excitation spectral microscopy with multiple fluorescence bands](#). Biomed Opt Express 13, 6048-6060, (2022), DOI: 10.1364/BOE.473241 (CF583R, CF620R, CF514, LipidSpot488, super-resolution, Excitation spectral microscopy)

Nahidazar, L. and Harkes, R. [Multicolor Localization-Based Super Resolution Microscopy](#). Methods Mol Biol 2350, 69-76, (2021), DOI: 10.1007/978-1-0716-1593-5_5 (CF488A, CF680, CF660C, CF514 SMLM)

Stoltzfus, C. R., et al. [CytoMAP: A Spatial Analysis Toolbox Reveals Features of Myeloid Cell Organization in Lymphoid Tissues](#). Cell Rep 31(3), 107523, (2020), DOI: 10.1016/j.celrep.2020.107523 (CF750, CF594, CF514, CF633 Mix-n-Stain)

CF®532

Chen, K., et al. [Excitation spectral microscopy for highly multiplexed fluorescence imaging and quantitative biosensing](#). bioRxiv, (2021), DOI: 10.1101/2021.04.13.439576 (CF532 wheat germ agglutinin CF514SE and CF568SE)

Sainski, A. M., et al. [Casp8p41 generated by HIV protease kills CD4 T cells through direct Bak activation](#). J Cell Biol 207(1), 159, (2014), DOI: jcb.20140505109252014c [pii] 10.1083/jcb.20140505109252014c (CF640R Mix-n-Stain; FC)

CF®543

Ara, A., et al. [Mechanism of Enhanced HIV Restriction by Virion Coencapsidated Cytidine Deaminases APOBEC3F and APOBEC3G](#). J Virol 91(3), (2017), DOI: 10.1128/JVI.02230-16 (CF543 Rabbit anti-V5 & CF640R rabbit anti-human)

Bonafede, R., et al. [ASC-Exosomes Ameliorate the Disease Progression in SOD1\(G93A\) Murine Model Underlining Their Potential Therapeutic Use in Human ALS](#). Int J Mol Sci 21(10), (2020), DOI: 10.3390/ijms21103651 (CF543 alpha-bungarotoxin)

Chen, Y., et al. [Turtle interacts with borderless in regulating glial extension and axon ensheathment](#). Mol Brain 10(1), 17, (2017), DOI: 10.1186/s13041-017-0299-6 (CF543 donkey anti-mouse)

Chen, Z. H., et al. [Characterization of interstitial Cajal progenitors cells and their changes in Hirschsprung's disease](#). PLoS One 9(1), e86100, (2014), DOI: 10.1371/journal.pone.0086100, PONE-D-13-26924 [pii] (CF488A donkey anti-rabbit, CF543 biotin anti-goat, CF633 donkey anti-mouse; IHC)

Held, R. G., et al. [Synapse and Active Zone Assembly in the Absence of Presynaptic Ca\(2+\) Channels and Ca\(2+\) Entry](#). Neuron 107(4), 667-683 e669, (2020), DOI: 10.1016/j.neuron.2020.05.032 (CF555 donkey anti-mouse 3D STORM)

Ito, H., et al. [Expression analyses of Phactr1 \(phosphatase and actin regulator 1\) during mouse brain development](#). Neurosci Res 128, 50-57, (2018), DOI: 10.1016/j.neures.2017.08.002 (CF488A goat anti-rabbit & CF543 goat anti-mouse)

Kijani, S., et al. [Filter-Dense Multicolor Microscopy](#). PLoS One 10(3), e0119499, (2015), DOI: 10.1371/journal.pone.0119499, PONE-D-14-20661 [pii] (CF488A, CF543, CF594)

Lannin, T. B., et al. [Comparison and optimization of machine learning methods for automated classification of circulating tumor cells](#). Cytometry A 89(10), 922-931, (2016), DOI: 10.1002/cyto.a.22993 (CF543 pan anti-cytokeratin & CF594 Mix-n-Stain)

Louzada, S., et al. [Fluorescence In Situ Hybridization onto DNA Fibres Generated Using Molecular Combing](#). Fluorescence In Situ Hybridization (FISH). pp. 275-293, (2017), DOI: 10.1007/978-3-662-52959-1_31 (CF543 succinimidyl ester)

Morita, S., et al. [Targeted DNA demethylation in vivo using dCas9-peptide repeat and scFv-TET1 catalytic domain fusions](#). Nat Biotechnol 34(10), 1060-1065, (2016), DOI: 10.1038/nbt.3658 (CF488A donkey anti-chicken & CF543 donkey anti-rabbit)

Noguchi, H., et al. [Expression of DNMT1 in neural stem/precursor cells is critical for survival of newly generated neurons in the adult hippocampus](#). Neurosci Res 95, 1-11, (2015), DOI: S0168-0102(15)00033-4 [pii] 10.1016/j.neures.2015.01.014 (CF488A donkey anti-mouse, CF543 donkey anti-rabbit, CF568 donkey anti-rat, CF647 donkey anti-goat, CF647 donkey anti-rabbit)

Oda, H., et al. [Plasma microRNAs Are Potential Biomarkers of Acute Rejection After Hindlimb Transplantation in Rats](#). Transplant Direct 2(11), e108, (2016), DOI: 10.1097/TXD.0000000000000620 (CF488A and CF543 secondaries)

Ren, J., et al. [Sirt1 Protects Subventricular Zone-Derived Neural Stem Cells from DNA Double-Strand Breaks and Contributes to Olfactory Function Maintenance in Aging Mice](#). Stem Cells 40, 493-507, (2022), DOI: 10.1093/stmcls/sxac008 (CF488A, CF543, CF633 secondaries, fluorescence microscopy)

Takenaka, N., et al. [A critical role of the small GTPase Rac1 in Akt2-mediated GLUT4 translocation in mouse skeletal muscle](#). FEBS J 281(5), 1493-1504, (2014), DOI: 10.1111/febs.12719 (CF488A, CF543, CF647 goat anti-mouse, goat anti-rabbit, goat anti-rat; ICC)

Takenaka, N., et al. [Rac1 Activation Caused by Membrane Translocation of a Guanine Nucleotide Exchange Factor in Akt2-Mediated Insulin Signaling in Mouse Skeletal Muscle](#). PLoS One 11(5), e0155292, (2016), DOI: 10.1371/journal.pone.0155292 (CF405, CF488A, CF543 & CF647 IHF)

Takenaka, N., et al. [Role of the guanine nucleotide exchange factor in Akt2-mediated plasma membrane translocation of GLUT4 in insulin-stimulated skeletal muscle](#). Cell Signal 26(11), 2460-2469, (2014), DOI: S0898-6568(14)00221-6 [pii] 10.1016/j.cellsig.2014.07.002 (CF350, CF543 & CF647; ICC)

Takeuchi, H., et al. [The efficacy of combining a vascularized biogenic conduit and a decellularized nerve graft in the treatment of peripheral nerve defects: An experimental study using the rat sciatic nerve defect model](#). Microsurgery 42, 254-264, (2022), DOI: 10.1002/micr.30853 (CF488A DARb, CF543 DAMs)

Thege, F. I., et al. [Microfluidic immunocapture of circulating pancreatic cells using parallel EpCAM and MUC1 capture: characterization, optimization and downstream analysis](#). Lab Chip 14(10), 1775-1784, (2014), DOI: 10.1039/c4lc00041b (CF543 succinimidyl ester; fluorescence microscopy)

Wang, X., et al. [Quantitative proteomic analysis of age-related subventricular zone proteins associated with neurodegenerative disease](#). Sci Rep 6, 37443, (2016), DOI: 10.1038/srep37443 (CF488A & CF543 secondaries)

CF®555

Accogli, G., et al. [A lectin-based cell microarray approach to analyze the mammalian granulosa cell surface glycosylation profile](#). Glycoconj J 33(5), 717-724, (2016), DOI: 10.1007/s10719-016-9666-2 10.1007/s10719-016-9666-2 [pii] (CF555 streptavidin)

Alves, A. A., et al. [Dynamics of the orphan myosin MyoF over Trypanosoma cruzi life cycle and along the endocytic pathway](#). Parasitol Int 86, 102444, (2022), DOI: 10.1016/j.parint.2021.102444 (CF555 BSA)

Avital-Shmilovici, M., et al. [A Mega-high-throughput Screening Platform for the Discovery of Biologically Relevant Sequence-defined Non-natural polymers](#). ChemRxiv (2020). DOI: 10.26434/chemrxiv.13106069.v1 (CF555 succinimidyl ester)

Bago, J. R., et al. [Therapeutically engineered induced neural stem cells are tumour-homing and inhibit progression of glioblastoma](#). Nat Commun 7, 10593, (2016), DOI: 10.1038/ncomms10593 (CF555 rabbit anti-goat)

Barboro, P., et al. [Androgen receptor activity is affected by both nuclear matrix localization and the phosphorylation status of the heterogeneous nuclear ribonucleoprotein K in anti-androgen-treated LNCaP cells](#). PLoS One 8(11), e79212, (2013), DOI: 10.1371/journal.pone.0079212 PONE-D-13-18847 [pii] (CF555 donkey anti-mouse, CF633 donkey anti-rabbit; ICC)

Barboro, P., et al. [The Role of Nuclear Matrix Proteins Binding to Matrix Attachment Regions \(MARs\) in Prostate Cancer Cell Differentiation](#). PLoS One 7(7), e40617, (2012), DOI: DOI:10.1371/journal.pone.0040617 (CF555 goat anti-mouse IgG, CF633 anti-rabbit IgG, CF555 streptavidin; ICC, FISH)

Bekiaris, C., et al. [Neurogenesis in the septal and temporal part of the adult rat dentate gyrus](#). Hippocampus 25(4), 511-523, (2015), DOI: 10.1002/hipo.22388 (CF488A goat anti-mouse & CF555 goat anti-rabbit)

Bertok, T., et al. [Carboxybetaine Modified Interface for Electrochemical Glycoprofiling of Antibodies Isolated from Human Serum](#). Langmuir 31, 7148-7157, (2015), DOI: 10.1021/acs.langmuir.5b00944 (CF555 streptavidin, microarray (STED))

Bostrom, M., et al. [The hippocampal neurovascular niche during normal development and after irradiation to the juvenile mouse brain](#). Int J Radiat Biol 90(9), 778-789, (2014), DOI: 10.3109/09553002.2014.931612 (CF555 donkey anti-goat; IHC)

Cantuti-Castelvetro, L., et al. [Neuropilin-1 facilitates SARS-CoV-2 cell entry and infectivity](#). Science 370(6518), 856-860, (2020), DOI: 10.1126/science.abd2985 (CF488A, CF555, CF647)

Cruz, A., and Joseph, S. [Interaction of the Influenza A Virus NS1 Protein with the 5'-m7G-mRNA.eIF4E.eIF4G1 Complex](#). Biochemistry 61, 1485-1494, (2022), DOI: 10.1021/acs.biochem.2c00019 (CF555 hydrazide (STED))

Dai, H. Q., et al. [TET-mediated DNA demethylation controls gastrulation by regulating Lefty-Nodal signalling](#). Nature 538(7626), 528-532, (2016), DOI: 10.1038/nature20095 (CF555 Streptavidin)

Deo, D. I., et al. [Biofunctionalization of PEGylated microcapsules for exclusive binding to protein substrates](#). Biomacromolecules 15(7), 2555-2562, (2014), DOI: 10.1021/bm500412d (CF555 Mix-n-Stain microcapsules; fluorescence microscopy)

Diamantopoulou, A., et al. [Denial or receipt of expected reward through maternal contact during the neonatal period differentially affect the development of the rat amygdala and program its function in adulthood in a sex-dimorphic way](#). Psychoneuroendocrinology 38(9), 1757, (2013), DOI: DOI:10.1016/j.psyneuen.2013.02.012 (Mix-n-Stain CF488A, Mix-N-Stain CF555; IHC)

Farias, G. G., et al. [Imaging the Polarized Sorting of Proteins from the Golgi Complex in Live Neurons](#). Methods Mol Biol 1496, 13-30, (2016), DOI: 10.1007/978-1-4939-6463-5_2 (CF488A, CF555 & CF640R Mix-n-Stain IF)

Gan, E. S., et al. [Jumonji demethylases moderate precocious flowering at elevated temperature via regulation of FLC in Arabidopsis](#). Nat Commun 5, 5098, (2014), DOI: ncomms6098 [pii] 10.1038/ncomms6098 (CF488A goat anti-rabbit, CF555 goat anti-mouse; ICC)

Giannakopoulou, A., et al. [Acute inflammation alters adult hippocampal neurogenesis in a multiple sclerosis mouse model](#). J Neurosci Res 91(7), 890-900, (2013), DOI: DOI:10.1002/jnr.23226 (CF555 streptavidin; TUNEL, IHC)

Gonzaga, A. C. R., et al. [Profile of cell proliferation and apoptosis activated by the intrinsic and extrinsic pathways in the prostate of aging rats](#). Prostate 77(9), 937-948, (2017), DOI: 10.1002/pros.23349 (CF555 goat anti-mouse)

Gregorieff, A., et al. [Yap-dependent reprogramming of Lgr5\(+\) stem cells drives intestinal regeneration and cancer](#). Nature 526(7575), 715-718, (2015), DOI: 10.1038/nature15382 (CF555 donkey anti-mouse & CF647 donkey anti-rabbit)

He, G. L., et al. [Inhibition of HSP90beta by ganetespib blocks the microglial signalling of evoked pro-inflammatory responses to heat shock](#). Int J Biochem Cell Biol 106, 35-45, (2019), DOI: 10.1016/j.biocel.2018.11.003 (CF633 cytokeratin anti-chicken CF555 donkey anti-mouse)

He, L., et al. [C9orf72 functions in the nucleus to regulate DNA damage repair](#). Cell Death Differ 30, 716-730, (2023), DOI: 10.1038/s41418-022-01074-0 (CF555 alpha-bungarotoxin, fluorescence microscopy (STED))

Held, R. G., et al. [Synapse and Active Zone Assembly in the Absence of Presynaptic Ca\(2+\) Channels and Ca\(2+\) Entry](#). Neuron 107(4), 667-683 e669, (2020), DOI: 10.1016/j.neuron.2020.05.032 (CF555 donkey anti-mouse 3D STORM)

Hu, M., et al. [Parkinson's disease-risk protein TMEM175 is a proton-activated proton channel in lysosomes](#). Cell 185, 2292-2308 e2220, (2022), DOI: 10.1016/j.cell.2022.05.021 (CF555 dextran 10K, fluorescence microscopy (STED))

Ikemoto, H., et al. [Hyaluronan-binding peptide for targeting peritoneal carcinomatosis](#). Tumour Biol 39(5), 1010428317701628, (2017), DOI: 10.1177/1010428317701628 (CF555SE)

Ise, H., et al. [Engulfment and clearance of apoptotic cells based on a GlcNAc-binding lectin-like property of surface vimentin](#). Glycobiology 22(6), 788, (2012), DOI: DOI:10.1093/glycob/cws052 (CF488A goat anti-mouse IgG, CF488A donkey anti-rat IgG, CF555 donkey anti-rabbit IgG; ICC, IHC)

Jiang, L., et al. [A bacterial extracellular vesicle-based intranasal vaccine against SARS-CoV-2 protects against disease and elicits neutralizing antibodies to wild-type and Delta variants](#). bioRxiv, (2022), DOI: 10.1101/2021.06.28.450181 (CF488A SE, CF555 SE, CF647 SE (STED))

Kawashima, Y., et al. [Detection of DNA double-strand breaks by pulsed-field gel electrophoresis](#). Genes Cells 22(1), 84-93, (2017), DOI: 10.1111/gtc.12457 (CF555 donkey anti-rabbit & CF633 donkey anti-mouse)

Kesik, M., et al. [A multi-functional fluorescent scaffold as a multi-colour probe: design and application in targeted cell imaging](#). RSC Advances 5(101), 83361-83367, (2015), DOI: 10.1039/c5ra16600d (CF555 Mix-n-Stain)

Kim, J., et al. [Basolateral to Central Amygdala Neural Circuits for Appetitive Behaviors](#). Neuron 93(6), 1464-1479 e1465, (2017), DOI: 10.1016/j.neuron.2017.02.034 (CF555 streptavidin)

Kiratitanaporn, W., et al. [3D printing a biocompatible elastomer for modeling muscle regeneration after volumetric muscle loss](#). Biomater Adv 142, 213171, (2022), DOI: 10.1016/j.bioadv.2022.213171 (CF555 GAMs, CF555 GARb, CF488A phalloidin, fluorescence microscopy (STED))

Klimas, A., et al. [Magnify is a universal molecular anchoring strategy for expansion microscopy](#). Nat Biotechnol 41, 858-869, (2023) DOI: 10.1038/s41587-022-01546-1 (CF555, super-resolution, ExM, STED)

Kumaran, G. K. and Hanukoglu, I. [Identification and classification of epithelial cells in nephron segments by actin cytoskeleton patterns](#). FEBS J 287(6), 1176-1194, (2019), DOI: <https://doi.org/10.1111/febs.15088> (CF488A phalloidin & CF555 wheat germ agglutinin)

Lehmann, M., et al. [Novel organic dyes for multicolor localization-based super-resolution microscopy](#). J Biophotonics 9(1-2), 161-170, (2016), DOI: 10.1002/jbio.201500119 (CF555, CF568, CF647, CF680, STORM)

Lin, Z. P., et al. [Macrophages Actively Transport Nanoparticles in Tumors After Extravasation](#). ACS Nano 16, 6080-6092, (2022), DOI: 10.1021/acsnano.1c11578 (CF555 GARt, CF647 DARb, 3D microscopy (STED))

Lindberg, O. R., et al. [Epidermal growth factor treatment of the adult brain subventricular zone leads to focal microglia/macrophage accumulation and angiogenesis](#). Stem Cell Reports 2(4), 440-448, (2014), DOI: 10.1016/j.stemcr.2014.02.003 S2213-6711(14)00055-1 [pii] (CF488A donkey anti-mouse CF555 donkey anti-rabbit, donkey anti-goat CF633 donkey anti-goat; IHC)

Lindberg, O. R., et al. [EGF-Induced Expansion of Migratory Cells in the Rostral Migratory Stream](#). PLoS One 7(9), e46380, (2012), DOI: DOI:10.1371/journal.pone.0046380 (CF488A donkey anti-mouse IgG; CF555 donkey anti-goat IgG; CF555 donkey anti-rabbit IgG; CF633 donkey anti-goat IgG; IHC)

Lourbopoulos, A., et al. [Changes in thyroid hormone receptors after permanent cerebral ischemia in male rats](#). J Mol Neurosci 54(1), 78-91, (2014), DOI: 10.1007/s12031-014-0253-3 (CF488A goat anti-rabbit, CF488A donkey anti-rabbit, CF555 goat anti-mouse, CF555 donkey anti-goat, EverBrite; IHC)

Lozada, A. F., et al. [Glutamatergic Synapse Formation is Promoted by alpha7-Containing Nicotinic Acetylcholine Receptors](#). J Neurosci 32(22), 7651-7661, (2012), DOI: DOI:10.1523/JNEUROSCI.6246-11.2012 (CF555 donkey anti-rabbit, CF633 goat anti-guinea pig; IHC)

Martins-Santos, E., et al. [Persistent testicular structural and functional alterations after exposure of adult rats to atrazine](#). Reprod Toxicol 73, 201-213, (2017), DOI: 10.1016/j.reprotox.2017.08.010 (CF555 goat anti-mouse)

Mikosik, A., et al. [Expression of calpain-calpastatin system \(CCS\) member proteins in human lymphocytes of young and elderly individuals: pilot baseline data for the CALPACENT project](#). Immun Ageing 10(1), 27, (2013), DOI: DOI:10.1186/1742-4933-10-27 (Mix-n-Stain CF555; FC)

Miyagishima, S. Y., et al. [DipM is required for peptidoglycan hydrolysis during chloroplast division](#). BMC Plant Biol 14, 57, (2014), DOI: 1471-2229-14-57 [pii] 10.1186/1471-2229-14-57 (CF488A Mix-n-Stain, CF555 Mix-n-Stain; ICC)

Nagornyykh, A. M., et al. [Protocol for chronic hepatitis B virus infection mouse model development by patient-derived orthotopic xenografts](#). PLoS One 17, e0264266, (2022), DOI: 10.1371/journal.pone.0264266 (CF555, Mix-n-Stain)

Nakashima, H., et al. [Canonical TGF-beta Signaling Negatively Regulates Neuronal Morphogenesis through TGIF/Smad Complex-Mediated CRMP2 Suppression](#). J Neurosci 38(20), 4791-4810, (2018), DOI: 10.1523/JNEUROSCI.2423-17.2018 (CF488A, CF555, CF647 secondaries)

Nakashima, H., et al. [MeCP2 controls neural stem cell fate specification through miR-199a-mediated inhibition of BMP-Smad signaling](#). Cell Rep 35(7), 109124, (2021), DOI: 10.1016/j.celrep.2021.109124 (CF488A donkey anti-rabbit CF488A donkey anti-goat CF555 donkey anti-mouse CF555 donkey anti-rabbit CF555 donkey anti-guinea pig CF647 donkey anti-mouse and CF647 donkey anti-rat)

Okada, S., et al. [Localization of heat shock protein 110 in canine mammary gland tumors](#). Vet Immunol Immunopathol 167(3-4), 139-146, (2015), DOI: 10.1016/j.vetimm.2015.07.012 (CF555 goat anti-rabbit)

Pang, H. B., et al. [An endocytosis pathway initiated through neuropilin-1 and regulated by nutrient availability](#). Nat Commun 5, 4904, (2014), DOI: ncomms5904 [pii] 10.1038/ncomms5904 (CF555 succinimidyl ester; nanoparticles)

Persson, A., et al. [Radixin expression in microglia after cortical stroke lesion](#). Glia 61(5), 790-799, (2013), DOI: DOI:10.1002/glia.22473 (CF555 donkey anti-goat IgG; IHC)

Peixoto-Neves, D., et al. [KV7.1 channel blockade inhibits neonatal renal autoregulation triggered by a step decrease in arterial pressure](#). Am J Physiol Renal Physiol 322(2), F197-F207, (2022), DOI: 10.1152/ajprenal.00568.2020 (CF488A donkey anti-rabbit & CF555 donkey anti-goat)

Pillai, A. G., et al. [Delayed effects of corticosterone on slow after-hyperpolarization potentials in mouse hippocampal versus prefrontal cortical pyramidal neurons](#). PLoS One 9(6), e99208, (2014), DOI: 10.1371/journal.pone.0099208 PONE-D-14-01195 [pii] (CF555 hydrazide; fluorescence microscopy)

Platt, J. L., et al. [C3d regulates immune checkpoint blockade and enhances antitumor immunity](#). JCI Insight 2(9), (2017), DOI: 10.1172/jci.insight.90201 (CF488A donkey anti-rat and CF555 donkey anti-goat)

Robajac, D., et al. [Sensitive glycoprofiling of insulin-like growth factor receptors isolated from colon tissue of patients with colorectal carcinoma using lectin-based protein microarray](#). Int J Biol Macromol 144, 932-937, (2020), DOI: 10.1016/j.ijbiomac.2019.09.170 (CF555 and CF647 streptavidin)

Rodkin, S., et al. [The Localization of p53 in the Crayfish Mechanoreceptor Neurons and Its Role in Axotomy-Induced Death of Satellite Glial Cells Remote from the Axon Transection Site](#). J Mol Neurosci 70(4), 532-541, (2020), DOI: 10.1007/s12031-019-01453-2 (CF555 goat anti-mouse IgG1)

Salazar-Cavazos, E., et al. [Multisite EGFR phosphorylation is regulated by adaptor protein abundances and dimer lifetimes](#). Mol Biol Cell 31(7), 695-708, (2020), DOI: 10.1091/mbc.E19-09-0548 (CF555 CF640R SIM)

Schultz, I., et al. [Pulmonary Insulin-like Growth Factor I Delivery from Trehalose and Silk-Fibroin Microparticles](#). ACS Biomaterials Science & Engineering 1(2), 119-129, (2015), DOI: 10.1021/ab500101c (CF555 goat anti-rabbit)

Smith, P. L., et al. [Neonatal peripheral immune challenge activates microglia and inhibits neurogenesis in the developing murine hippocampus](#). Dev Neurosci 36(2), 119-131, (2014), DOI: 000359950 [pii] 10.1159/000359950 (CF555 donkey anti-goat, CF488A donkey anti-chicken & CF647 donkey anti-mouse; IHC)

Song, Y., et al. [Activation of FOXO3a reverses 5-Fluorouracil resistance in human breast cancer cells](#). Exp Mol Pathol 105(1), 57-62, (2018), DOI: 10.1016/j.yexmp.2018.05.013 (CF555 goat anti-rabbit)

Starunov, V. V., et al. [A metameric origin for the annelid pygidium?](#). BMC Evol Biol 15, 25, (2015), DOI: 10.1186/s12862-015-0299-z 10.1186/s12862-015-0299-z [pii] (CF633 goat anti-mouse & CF555 goat anti-rabbit)

Sugahara, K. N., et al. [Tumor-penetrating iRGD peptide inhibits metastasis](#). Mol Cancer Ther 14(1), 120-128, (2015), DOI: 10.1158/1535-7163.MCT-14-0366 (CF555 succinimidyl ester)

Sunderic, M., et al. [Lectin-based protein microarray analysis of differences in serum alpha-2-macroglobulin glycosylation between patients with colorectal cancer and persons without cancer](#). Biotechnol Appl Biochem 63(4), 457-464, (2016), DOI: 10.1002/bab.1407 (CF555 streptavidin)

Tai, L. H., et al. [Transient stimulation of distinct subpopulations of striatal neurons mimics changes in action value](#). Nat Neurosci 15(9), 1281-1289, (2012), DOI: DOI:10.1038/nn.3188 (CF555 donkey anti-mouse IgG, CF647 donkey anti-goat IgG; IHC)

Takakura, M., et al. [Circulating tumour cells detected by a novel adenovirus-mediated system may be a potent therapeutic marker in gynaecological cancers](#). Br J Cancer 107(3), 448, (2012), DOI: DOI:10.1038/bjc.2012.276 (CF555 goat anti-mouse IgG; ICC)

Takeichi, T., et al. [The effects of low-dose methamphetamine pretreatment on endoplasmic reticulum stress and methamphetamine neurotoxicity in the rat midbrain](#). Leg Med (Tokyo) 14(2), 69-77, (2012), DOI: DOI:10.1016/j.legalmed.2011.12.004 (CF555 goat anti-rabbit IgG; IHC)

Theotokis, P., et al. [Connexin43 and connexin47 alterations after neural precursor cells transplantation in experimental autoimmune encephalomyelitis](#). Glia 63(10), 1772-1783, (2015), DOI: 10.1002/glia.22843 (CF555 secondaries)

Uno, T., et al. [Characterization of Rab-interacting lysosomal protein in the brain of Bombyx mori](#). Histochem Cell Biol 141(3), 311-320, (2014), DOI: 10.1007/s00418-013-1160-3 (CF555 donkey anti-mouse; IHC)

Wada, M., et al. [Epitope-Dependent Pathogenicity of Antibodies Targeting a Major Bullous Pemphigoid Autoantigen Collagen XVII/BP180](#). J Invest Dermatol 136(5), 938-946, (2016), DOI: 10.1016/j.jid.2015.11.030 (CF555 Mix-n-Stain time lapse imaging)

Wei, Y., et al. [Synergistic Entry of Individual Nanoparticles into Mammalian Cells Driven by Free Energy Decline and Regulated by Their Sizes](#). ACS Nano 16, 5885-5897, (2022), DOI: 10.1021/acsnano.1c11068 (CF488A, CF555, CF647 SE, nanoparticle labeling, fluorescence microscopy (STED))

West, A. B., et al. [Differential LRRK2 expression in the cortex, striatum, and substantia nigra in transgenic and nontransgenic rodents](#). J Comp Neurol 522(11), 2465-2480, (2014), DOI: 10.1002/cne.23583 (CF555 donkey anti-rabbit & CF488A donkey anti-mouse; IHC)

Willmore, A. M., et al. [Targeted silver nanoparticles for ratiometric cell phenotyping](#). Nanoscale 8(17), 9096-9101, (2016), DOI: 10.1039/c5nr07928d (CF555 succinimidyl ester)

Wu, S., et al. [AUF1 is recruited to the stress granules induced by coxsackievirus B3](#). Virus Res 192, 52-61, (2014), DOI: S0168-1702(14)00331-1 [pii] 10.1016/j.virusres.2014.08.003 (CF488A goat anti-rabbit, CF555 goat anti-mouse & CF555 goat anti-rabbit; ICC)

Wu, S., et al. [Protease 2A induces stress granule formation during coxsackievirus B3 and enterovirus 71 infections](#). Virol J 11, 192, (2014), DOI: 10.1186/s12985-014-0192-1 s12985-014-0192-1 [pii] (CF555 goat anti-mouse & CF555 goat anti-rabbit)

Xu, Y., et al. [Arabidopsis MRG domain proteins bridge two histone modifications to elevate expression of flowering genes](#). Nucleic Acids Res 42(17), 10960-10974, (2014), DOI: 10.1093/nar/gku781 (CF488A goat anti-rabbit & CF555 goat anti-mouse)

Xu, Z., et al. [Mucosal immunoglobulins at respiratory surfaces mark an ancient association that predates the emergence of tetrapods](#). Nat Commun 7, 10728, (2016), DOI: 10.1038/ncomms10728 (CF488A, CF555 & CF640R Mix-n-Stain)

Xu, Z., et al. [Teleost skin, an ancient mucosal surface that elicits gut-like immune responses](#). Proc Natl Acad Sci U S A 110(32), 13097-13102, (2013), DOI: 1304319110 [pii] 10.1073/pnas.1304319110 (CF488A Mix-n-Stain, CF555, CF640R; IHC)

Yang, Y., et al. [Paraoxonase 2 decreases renal reactive oxygen species production, lowers blood pressure, and mediates dopamine D\(2\) receptor-induced inhibition of NADPH oxidase](#). Free Radic Biol Med 53(3), 437, (2012), DOI: DOI:10.1016/j.freeradbiomed.2012.05.015 (Mix-n-Stain CF555; IHC)

Yu, L., et al. [Gemcitabine eliminates double minute chromosomes from human ovarian cancer cells](#). PLoS One 8(8), e71988, (2013), DOI: 10.1371/journal.pone.0071988 PONE-D-13-11229 [pii] (CF555 goat anti-mouse; ICC)

Zhang, Y., et al. [Deep-brain magnetic stimulation promotes adult hippocampal neurogenesis and alleviates stress-related behaviors in mouse models for neuropsychiatric disorders](#). Mol Brain 7, 11, (2014), DOI: 1756-6606-7-11 [pii] 10.1186/1756-6606-7-11 (CF488A goat anti-mouse & CF555 donkey anti-mouse; IHC)

Zhou, H. and Liu, G. [Regulation of density of functional presynaptic terminals by local energy supply](#). Mol Brain 8, 42, (2015), DOI: 10.1186/s13041-015-0132-z (CF488A, CF555 & CF640R secondaries, MitoView 633)

CF®568

Ahi, Y. S., et al. [Adenoviral E2 IVa2 Protein Interacts with L4 33K Protein and E2 DNA-Binding Protein](#). J Gen Virol 94(6), 1325, (2013), DOI: DOI:10.1099/vir.0.049346-0 (Mix-n-Stain CF350, Mix-n-Stain CF488A, CF568 goat anti-mouse IgG; ICC)

Almahayni, K., et al. [Simple, Economic, and Robust Rail-Based Setup for Super-Resolution Localization Microscopy](#). J Phys Chem A 127, 4553-4560, (2023), DOI: 10.1021/acs.jpca.3c01351 (CF568, super-resolution, SMLM, TIRF)

Alshafie, W., et al. [Regulated resurfacing of a somatostatin receptor storage compartment fine-tunes pituitary secretion](#). Journal of Cell Biology 219(1), (2020), DOI: 10.1083/jcb.201904054 (CF568 goat anti-rat STORM)

Bartok, A., et al. [IP3 receptor isoforms differently regulate ER-mitochondrial contacts and local calcium transfer](#). Nat Commun 10(1), 3726, (2019), DOI: 10.1038/s41467-019-11646-3 (CF568 goat anti-rabbit STORM)

Bennett, H. W., et al. [Novel fibrillar structure in the inversin compartment of primary cilia revealed by 3D single-molecule superresolution microscopy](#). Mol Biol Cell 31(7), 619-639, (2020), DOI: 10.1091/mbc.E19-09-0499 (CF568 STORM)

Breton-Provencher, V., et al. [Principal cell activity induces spine relocation of adult-born interneurons in the olfactory bulb](#). Nat Commun 7, 12659, (2016), DOI: 10.1038/ncomms12659 (CF568 Mix-n-Stain IF)

Brock, V. J., et al. [P2X4 and P2X7 are essential players in basal T cell activity and Ca\(2+\) signaling milliseconds after T cell activation](#). Sci Adv 8(5), eabl9770, (2022), DOI: 10.1126/sciadv.abl9770 (CF568)

Catsburg, L. A. E., et al. [Dynamics and nanoscale organization of the postsynaptic endocytic zone at excitatory synapses](#). bioRxiv, (2021), DOI: 10.1101/2021.02.18.431766 (CF568 STED)

Cespedes, P. F., et al. [T-cell trans-synaptic vesicles are distinct and carry greater effector content than constitutive extracellular vesicles](#). Nat Commun 13, 3460, (2022), DOI: 10.1038/s41467-022-31160-3 (CF568, WGA, fluorescence microscopy, TIRF)

Chen, K., et al. [Excitation spectral microscopy for highly multiplexed fluorescence imaging and quantitative biosensing](#). bioRxiv, (2021), DOI: 10.1101/2021.04.13.439576 (CF532 wheat germ agglutinin CF514 succinimidyl ester and CF568 succinimidyl ester)

Ciapa, B. and Granon, S. [Expression of Cyclin-D1 in Astrocytes Varies During Aging](#). Front Aging Neurosci 10, 104, (2018), DOI: 10.3389/fnagi.2018.00104 (CF568 Mix-n-Stain IF)

Cohen, Y., et al. [The yeast p5 type ATPase, spf1, regulates manganese transport into the endoplasmic reticulum](#). PLoS One 8(12), e85519, (2013), DOI: 10.1371/journal.pone.0085519, PONE-D-13-37918 [pii] (CF568 goat anti-rabbit; ICC)

Deng, C., et al. [Impaired dynamic interaction of axonal endoplasmic reticulum and ribosomes contributes to defective stimulus-response in spinal muscular atrophy](#). Transl Neurodegener 11, 31, (2022)., DOI: 10.1186/s40035-022-00304-2 (CF568 DARb, super-resolution, dSTORM)

Davies, B. S., et al. [Assessing mechanisms of GPIHBP1 and lipoprotein lipase movement across endothelial cells](#). J Lipid Res 53(12), 2690, (2012), DOI: DOI:10.1194/jlr.M031559 (CF568 succinimidyl ester; in vivo injection, IHC)

Engdahl, A. K., et al. [Large Field-of-View Super-Resolution Optical Microscopy Based on Planar Polymer Waveguides](#). ACS Photonics 8(7), 1944-1950, (2021), DOI: 10.1021/acsphotonics.1c00914 (CF568 phalloidin TIRF)

Engdahl, A. K., et al. [Switching of Organic Fluorophores by Glycerol-sulfite Interactions for Single-Molecule Super-Resolution Microscopy](#). ChemPhotoChem 7, (2023), DOI: 10.1002/cptc.202200305 (CF647, CF568, Super-Resolution, dSTORM)

Engel, B. J., et al. [Caspase-3 Substrates for Noninvasive Pharmacodynamic Imaging of Apoptosis by PET/CT](#). Bioconjug Chem 29(9), 3180-3195, (2018), DOI: 10.1021/acs.bioconjchem.8b00514 (CF568 Annexin V)

Fernandez, M. A., et al. [High-Dimensional Immunophenotyping with 37-Color Panel Using Full-Spectrum Cytometry](#). Methods Mol Biol 2386, 43-60, (2022), DOI: 10.1007/978-1-0716-1771-7_4 (CF568)

Fu, X., et al. [In Vivo Single-Molecule Detection of Nanoparticles for Multiphoton Fluorescence Correlation Spectroscopy to Quantify Cerebral Blood Flow](#). Nano Lett 20(8), 6135-6141, (2020), DOI: 10.1021/acs.nanolett.0c02280 (CF488A dextran 10K 70K 250K CF568 dextran 10K)

Futatsuki, T., et al. [Involvement of orexin neurons in fasting- and central adenosine-induced hypothermia](#). Sci Rep 8(1), 2717, (2018), DOI: 10.1038/s41598-018-21252-w (CF568 donkey anti-goat and CF647 donkey anti-goat)

Gao, N., et al. [A Role of Lamin A/C in Preventing Neuromuscular Junction Decline in Mice](#). J Neurosci 40(38), 7203-7215, (2020), DOI: 10.1523/JNEUROSCI.0443-20.2020 (CF568 alpha-bungarotoxin)

Garcia, J. D., et al. [Stepwise disassembly of GABAergic synapses during pathogenic excitotoxicity](#). Cell Rep 37, 110142, (2021), DOI: 10.1016/j.celrep.2021.110142 (CF568 GAMs, TIRF)

Gong, Y.-N., et al. [Biological events and molecular signaling following MLKL activation during necroptosis](#). Cell Cycle, 1-13, (2017), DOI: 10.1080/15384101.2017.1371889 (CF568, CF647 STORM)

Gorur, A., et al. [COPII-coated membranes function as transport carriers of intracellular procollagen I](#). The Journal of Cell Biology 216(6), 1745-1759, (2017), DOI: 10.1083/jcb.201702135 (CF568, CF680 STORM)

Guo, D., et al. [Endosomal trafficking of nanoformulated antiretroviral therapy facilitates drug particle carriage and HIV clearance](#). J Virol 88(17), 9504-9513, (2014), DOI: JVI.01557-14 [pii] 10.1128/JVI.01557-14 (CF568 succinimidyl ester labeled atazanavir; fluorescence microscopy)

Guo, X., et al. [Mitochondrial stress is relayed to the cytosol by an OMA1-DELE1-HRI pathway](#). Nature 579(7799), 427-432, (2020), DOI: 10.1038/s41586-020-2078-2 (CF568 donkey anti-rabbit CF680 donkey anti-rabbit STORM)

Haas, K. T., et al. [Single-molecule localization microscopy reveals molecular transactions during RAD51 filament assembly at cellular DNA damage sites](#). Nucleic Acids Res, (2018), DOI: 4788348 [pii] 10.1093/nar/gkx1303 (CF568, STORM)

Haas, K. T., et al. [Multitarget Immunohistochemistry for Confocal and Super-resolution Imaging of Plant Cell Wall Polysaccharides](#). Bio Protoc 10(19), e3783, (2020), DOI: 10.21769/BioProtoc.3783 (CF568)

Hadassa Y. Holzapfel, and Marc R. Birtwistle. [Creating Complex Fluorophore spectra on Antibodies Through Combinatorial Labeling](#). Transl Sci, 1-10, (2016). (CF405M, CF488A, CF568, CF640R, FRET, MnS)

Hamel, Y., et al. [Compromised mitochondrial quality control triggers lipin1-related rhabdomyolysis](#). Cell Rep Med 2, 100370, (2021), DOI: 10.1016/j.xcrm.2021.100370 (CF568 CTX, cholera toxin uptake)

Haselmann, H., et al. [Human Autoantibodies against the AMPA Receptor Subunit GluA2 Induce Receptor Reorganization and Memory Dysfunction](#). Neuron 100(1), 91-105 e109, (2018), DOI: 10.1016/j.neuron.2018.07.048 (CF568 goat anti-guinea pig STORM)

Hauser, M., et al. [The Spectrin-Actin-Based Periodic Cytoskeleton as a Conserved Nanoscale Scaffold and Ruler of the Neural Stem Cell Lineage](#). Cell Rep 24(6), 1512-1522, (2018), DOI: 10.1016/j.celrep.2018.07.005 (CF568 STORM)

He, J., et al. [Cucurbitacin IIa induces caspase-3-dependent apoptosis and enhances autophagy in lipopolysaccharide-stimulated RAW 264.7 macrophages](#). Int Immunopharmacol, (2013), DOI: DOI:10.1016/j.intimp.2013.03.013 (CF488A goat anti-mouse IgG, CF568 goat anti-rabbit IgG; ICC)

Heller, J. [Exploring Nanoscale Organisation of Synapses with Super-Resolution Microscopy](#). OM&P 3(2), 48-58, (2017), DOI: doi:10.20388/omp2017.002.0045 (CF568 STORM)

Henneberger, C., et al. [LTP Induction Boosts Glutamate Spillover by Driving Withdrawal of Perisynaptic Astroglia](#). Neuron 108(5), 919-936 e911, (2020), DOI: 10.1016/j.neuron.2020.08.030 (CF568 DAMs 3-color STORM)

Hirano, K. and Namihira, M. [LSD1 Mediates Neuronal Differentiation of Human Fetal Neural Stem Cells by Controlling the Expression of a Novel Target Gene, HEYL](#). Stem Cells 34(7), 1872-1882, (2016), DOI: 10.1002/stem.2362 (CF488A, CF568, CF647 donkey secondaries)

Hofmann, D., et al. [A small molecule screen identifies novel inhibitors of mechanosensory nematocyst discharge in Hydra](#). Sci Rep 11(1), 20627, (2021), DOI: 10.1038/s41598-021-99974-7 (CF568 ARb)

Home, P., et al. [Altered subcellular localization of transcription factor TEAD4 regulates first mammalian cell lineage commitment](#). Proc Natl Acad Sci U S A 109(19), 7362-7367, (2012), DOI: DOI:10.1073/pnas.1201595109 (Mix-n-Stain CF488A, Mix-n-Stain CF568; ICC)

Hu, B., et al. [Orthogonally engineering matrix topography and rigidity to regulate multicellular morphology](#). Adv Mater 26(33), 5786-5793, (2014), DOI: 10.1002/adma.201402489 (CF568 phalloidin; ICC)

Huang, M. Y., et al. [Chemotherapeutic agent CPT-11 eliminates peritoneal resident macrophages by inducing apoptosis](#). Apoptosis 21(2), 130-142, (2016), DOI: 10.1007/s10495-015-1193-2 (CF568 goat anti-rabbit)

Hui, T. K., et al. [Ablation of Lrp4 in Schwann Cells Promotes Peripheral Nerve Regeneration in Mice](#). Biology (Basel) 10(6), (2021), DOI: 10.3390/biology10060452 (CF568 CF488A alpha-bungarotoxin)

Hui, T., et al. [Neuromuscular junction-specific genes screening by deep RNA-seq analysis](#). Cell Biosci 11(1), 81, (2021), DOI: 10.1186/s13578-021-00590-9 (CF568 bungarotoxin)

Inui, K., et al. [Nasal TRPA1 mediates irritant-induced bradypnea in mice](#). Physiol Rep 4(24), (2016), DOI: 10.14814/phy2.13098 (CF568 donkey anti-guinea pig)

Irie, T. [Essential Role of Somatic Kv2 Channels in High-Frequency Firing in Cartwheel Cells of the Dorsal Cochlear Nucleus](#). eNeuro, (2021), DOI: 10.1523/ENEURO.0515-20.2021 (CF405S IgG2a CF568 IgG2b)

Jaiswal, D. and Brown, J. L. [Nanofiber diameter-dependent MAPK activity in osteoblasts](#). J Biomed Mater Res A 100(11), 2921, (2012), DOI: DOI:10.1002/jbm.a.34234 (CF568 phalloidin; ICC)

J. Kim and M. Wojcik, Y. W., K. Xu, X. Zhang. [Oblique-sectional single-molecule microscopy](#). IEEE Xplore, 1-2, (2018), DOI: 10.1109/IPCon.2018.8527149 (CF568 obSTORM)

Jorgens, D. M., et al. [Deep nuclear invaginations are linked to cytoskeletal filaments - integrated bioimaging of epithelial cells in 3D culture](#). J Cell Sci 130(1), 177-189, (2017), DOI: jcs.190967 [pii] 10.1242/jcs.190967 (CF568 STORM)

Jung, S. J., et al. [Novel cerebello-amygdala connections provide missing link between cerebellum and limbic system](#). (2022), DOI: 10.1101/2022.02.07.479043 (CF568 CTB, confocal microscopy)

Karagounis, I. V., et al. [Repression of the autophagic response sensitises lung cancer cells to radiation and chemotherapy](#). Br J Cancer 115(3), 312-321, (2016), DOI: 10.1038/bjc.2016.202 (CF488A & CF568 secondaries)

Karanasios, E., et al. [Autophagy initiation by ULK complex assembly on ER tubulovesicular regions marked by ATG9 vesicles](#). Nat Commun 7, 12420, (2016), DOI: ncomms12420 [pii] 10.1038/ncomms12420 (CF568 STORM)

Katona, M., et al. [Capture at the ER-mitochondrial contacts licenses IP\(3\) receptors to stimulate local Ca\(2+\) transfer and oxidative metabolism](#). Nat Commun 13, 6779, (2022) DOI: 10.1038/s41467-022-34365-8 (CF568 GARb, super-resolution, STORM)

Kawase, A., et al. [Decreased radixin function for ATP-binding cassette transporters in liver in adjuvant-induced arthritis rats](#). J Pharm Sci 103(12), 4058-4065, (2014), DOI: 10.1002/jps.24210, S0022-3549(15)30300-2 [pii] (CF488A goat anti-rat & CF568 goat anti-mouse)

Kawase, A., et al. [Effects of ginsenosides on the expression of cytochrome P450s and transporters involved in cholesterol metabolism](#). J Nat Med 68(2), 395, (2013), DOI: DOI:10.1007/s11418-013-0791-y (CF568 goat anti-mouse IgG; IHC)

Ketel, K., et al. [A phosphoinositide conversion mechanism for exit from endosomes](#). Nature 529(7586), 408-412, (2016), DOI: 10.1038/nature16516 (CF568 Cholera toxin)

Kim, J. M., et al. [Regulation of sclerostin by the SIRT1 stabilization pathway in osteocytes](#). Cell Death Differ 29, 1625-1638, (2022), DOI: 10.1038/s41418-022-00952-x (CF568 phalloidin)

Köhler, S., et al. [The interaction of crossover formation and the dynamic architecture of the synaptonemal complex during meiosis](#). bioRxiv, (2020), DOI: 10.1101/2020.02.16.947804 (CF568 succinimidyl ester STORM)

Kraus, F., et al. [Quantitative 3D structured illumination microscopy of nuclear structures](#). Nat Protoc 12(5), 1011-1028, (2017), DOI: nprot.2017.020 [pii] 10.1038/nprot.2017.020 (CF568 azide, CF405M azide, SIM)

Krey, J. F., et al. [ANKRD24 organizes TRIOBP to reinforce stereocilia insertion points](#). J Cell Biol 221, (2022), DOI: 10.1083/jcb.202109134 (CF405M phalloidin, CF568 phalloidin, CF568 WGA, Everbrite, fluorescence microscopy)

Kudou, K., et al. [The requirement of Mettl3-promoted MyoD mRNA maintenance in proliferative myoblasts for skeletal muscle differentiation](#). Open Biol 7(9), (2017), DOI: 10.1098/rsob.170119 (CF568 goat anti-mouse)

Lehmann, M., et al. [Novel organic dyes for multicolor localization-based super-resolution microscopy](#). J Biophotonics 9(1-2), 161-170, (2016), DOI: 10.1002/jbio.201500119 (CF555, CF568, CF647, CF680, STORM)

Li, C. G., et al. [Berberine augments ATP-induced inflammasome activation in macrophages by enhancing AMPK signaling](#). Oncotarget 8(1), 95-109, (2017), DOI: 10.18632/oncotarget.13921 (CF568 goat anti-rabbit)

Liedmann, S., et al. [Localization of a TORC1-eIF4F translation complex during CD8\(+\) T cell activation drives divergent cell fate](#). Mol Cell 82, 2401-2414 e2409, (2022), DOI: 10.1016/j.molcel.2022.04.016 (CF488A DARt, CF568 DAGt, CF568 DARt, super-resolution, STORM)

Lim, A., et al. [Two kinesins drive anterograde neuropeptide transport](#). Mol Biol Cell, (2017), DOI: mbc.E16-12-0820 [pii] 10.1091/mbc.E16-12-0820 (Mix-n-Stain CF568, Direct + indirect IF, CF568 3D SIM)

Liu, K. P., et al. [LC3B-II deacetylation by histone deacetylase 6 is involved in serum-starvation-induced autophagic degradation](#). Biochem Biophys Res Commun 441(4), 970-975, (2013), DOI: S0006-291X(13)01880-9 [pii] 10.1016/j.bbrc.2013.11.007 (CF488A goat anti-mouse CF568 goat anti-rabbit; ICC)

Lukonin, I., et al. [Phenotypic landscape of intestinal organoid regeneration](#). Nature 586(7828), 275-280, (2020), DOI: 10.1038/s41586-020-2776-9 (CF568 Mix-n-Stain)

Luo, F., et al. [Neurexins regulate presynaptic GABAB-receptors at central synapses](#). Nat Commun 12(1), 2380, (2021), DOI: 10.1038/s41467-021-22753-5 (CF568 secondary)

Lycas, M. D., et al. [Nanoscopic dopamine transporter distribution and conformation are inversely regulated by excitatory drive and D2-autoreceptor activity](#). bioRxiv, (2021), DOI: 10.1101/2021.03.09.434538 (CF568 STORM)

Ma, X., et al. [Histone chaperone CAF-1 promotes HIV-1 latency by leading the formation of phase-separated suppressive nuclear bodies](#). EMBO J 40(10), e106632, (2021), DOI: 10.15252/embj.2020106632 (CF568 donkey anti-rabbit CF568 donkey anti-mouse STORM)

Markaki, Y., et al. (2021). [Xist nucleates local protein gradients to propagate silencing across the X chromosome](#). Cell 184, 6174-6192 e6132, (2021), DOI: 10.1016/j.cell.2021.10.022 (CF568 DARb, CF568 azide, TIRF)

Meitinger, F., et al. [TRIM37 prevents formation of condensate-organized ectopic spindle poles to ensure mitotic fidelity](#). J Cell Biol 220(7), (2021), DOI: 10.1083/jcb.202010180 (CF647 CF568)

Meyer, K., et al. [A Predictive 3D Multi-Scale Model of Biliary Fluid Dynamics in the Liver Lobule](#). Cell Syst 4(3), 277-290 e279, (2017), DOI: 10.1016/j.cels.2017.02.008 (CF568 donkey anti-rat)

Mezache, L., et al. [Vascular endothelial growth factor promotes atrial arrhythmias by inducing acute intercalated disk remodeling](#). Sci Rep 10(1), 20463, (2020), DOI: 10.1038/s41598-020-77562-5 (CF568 2-color STORM)

Michaluk, P., et al. [Rapid recycling of glutamate transporters on the astroglial surface](#). Elife 10, (2021), DOI: 10.7554/eLife.64714 (CF568 DAMs)

Mofarrahi, M., et al. [Angiopoietin-1 enhances skeletal muscle regeneration in mice](#). American Journal of Physiology - Regulatory, Integrative and Comparative Physiology 308(7), R576-R589, (2015), DOI: 10.1152/ajpregu.00267.2014 (CF568 Mix-n-Stain)

Neumann, K., et al. [Clec12a is an inhibitory receptor for uric acid crystals that regulates inflammation in response to cell death](#). Immunity 40(3), 389-399, (2014), DOI: S1074-7613(14)00069-7 [pii] 10.1016/j.immuni.2013.12.015 (CF568 succinimidyl ester; fluorescence microscopy)

Niec, R. E., et al. [Lymphatics act as a signaling hub to regulate intestinal stem cell activity](#). Cell Stem Cell 29, 1067-1082 e1018, (2022), DOI: 10.1016/j.stem.2022.05.007 (CF568, Mix-n-Stain)

Noguchi, H., et al. [Expression of DNMT1 in neural stem/precursor cells is critical for survival of newly generated neurons in the adult hippocampus](#). Neurosci Res 95, 1-11, (2015), DOI: S0168-0102(15)00033-4 [pii] 10.1016/j.neures.2015.01.014 (CF488A donkey anti-mouse, CF543 donkey anti-rabbit, CF568 donkey anti-rat, CF647 donkey anti-goat, CF647 donkey anti-rabbit)

Olasz, F., et al. [Immunological and biochemical characterisation of 7ap, a short protein translated from an alternative frame of ORF7 of PRRSV](#). Acta Vet Hung 64(2), 273-287, (2016), DOI: 10.1556/004.2016.027 (CF568 goat anti-swine & CF594 goat anti-mouse)

Ouyang, D. Y., et al. [Autophagy is differentially induced in prostate cancer LNCaP, DU145 and PC-3 cells via distinct splicing profiles of ATG5](#). Autophagy 9(1), 20-32, (2013), DOI: DOI:10.4161/auto.22397 (CF568 goat anti-rabbit IgG; ICC)

Ouyang, D. Y., et al. [Piperine inhibits the proliferation of human prostate cancer cells via induction of cell cycle arrest and autophagy](#). Food Chem Toxicol 60, 424-430, (2013), DOI: S0278-6915(13)00549-8 [pii] 10.1016/j.fct.2013.08.007 (CF568 goat anti-rabbit; ICC)

Pang, J., et al. [Step-by-step preparation of mouse eye sections for routine histology, immunofluorescence, and RNA in situ hybridization multiplexing](#). STAR Protoc 2, 100879, (2021), DOI: 10.1016/j.xpro.2021.100879 (CF488A DAMs, CF568 DARb, CF488A DAGt)

Park, K. S., et al. [A crucial requirement for Hedgehog signaling in small cell lung cancer](#). Nat Med 17(11), 1504-1508, (2011), DOI: DOI:10.1038/nm.2473 (Mix-n-Stain CF488A, Mix-n-Stain CF568, and Mix-n-Stain CF647; ICC)

Petrelli, A., et al. [Efficient Conjugation of Oligosaccharides to Polymer Particles through Furan/Maleimide Diels-Alder Reaction: Application to the Capture of Carbohydrate-Binding Proteins](#). Chembiochem 18(2), 206-212, (2017), DOI: 10.1002/cbic.201600509 (CF488A, CF568 Mix-n-Stain)

Poddar, A., et al. [Effects of individual base-pairs on in vivo target search and destruction kinetics of small RNA](#). bioRxiv, (2020), DOI: 10.1101/2020.07.24.215095 (CF568 succinimidyl ester)

Pons, M., et al. [Phosphorylation of filamin A regulates chemokine receptor CCR2 recycling](#). J Cell Sci 130(2), 490-501, (2017), DOI: 10.1242/jcs.193821 (CF488A, CF568 Mix-n-Stain)

Rahman, A., et al. [GMAP is an Atg8a-interacting protein that regulates Golgi turnover in Drosophila](#). Cell Rep 39, 110903, (2022), DOI: 10.1016/j.celrep.2022.110903 (CF488A, CF568 secondaries, fluorescence microscopy)

Rantataro, S., et al. [Nanoscale geometry determines mechanical biocompatibility of vertically aligned nanofibers](#). Acta Biomater 146, 235-247, (2022), DOI: 10.1016/j.actbio.2022.04.032 (CF568 phalloidin)

Ren, S., et al. [Anti-proliferative effect of 23,24-dihydrocucurbitacin F on human prostate cancer cells through induction of actin aggregation and cofilin-actin rod formation](#). Cancer Chemother Pharmacol 70(3), 415, (2012), DOI: DOI:10.1007/s00280-012-1921-z (CF488A goat anti-mouse, CF568 goat anti-rabbit; ICC)

Robertson, T. F., et al. [Lymphocyte egress signal sphingosine-1-phosphate promotes ERM-guided, bleb-based migration](#). J Cell Biol 220, (2021), DOI: 10.1083/jcb.202007182 (CF568, Mix-n-Stain)

Rojas-Sánchez, L., et al. [Synthesis and intracellular tracing surface-functionalized calcium phosphate nanoparticles by super-resolution microscopy \(STORM\)](#). Materialia 12, (2020), DOI: 10.1016/j.mtla.2020.100773 (CF568 wheat germ agglutinin STORM)

Sabbah, S., et al. [A retinal code for motion along the gravitational and body axes](#). Nature 546(7659), 492-497, (2017), DOI: 10.1038/nature22818 (CF568 Cholera Toxin retrograde tracing)

Saliba, D. G., et al. [Composition and structure of synaptic ectosomes exporting antigen receptor linked to functional CD40 ligand from helper T-cells](#). (2019), DOI: 10.1101/600551 (CF568 wheat germ agglutinin STORM)

Samarasinghe, A. E., et al. [Eosinophils Promote Antiviral Immunity in Mice Infected with Influenza A Virus](#). J Immunol 198(8), 3214-3226, (2017), DOI: 10.4049/jimmunol.1600787 (CF568 donkey anti-mouse)

Sando, R., et al. [Engineered synaptic tools reveal localized cAMP signaling in synapse assembly](#). J Cell Biol 221, (2022), DOI: 10.1083/jcb.202109111 (CF568, AMs)

Sato-Yamada, Y., et al. [A SARM1-mitochondrial feedback loop drives neuropathogenesis in a Charcot-Marie-Tooth disease type 2A rat model](#). J Clin Invest 132, (2022), DOI: 10.1172/JCI161566 (CF568 bungarotoxin, microscopy)

Schmerl, B., et al. [The synaptic scaffold protein MPP2 interacts with GABAA receptors at the periphery of the postsynaptic density of glutamatergic synapses](#). PLoS Biol 20, e3001503, (2022), DOI: 10.1371/journal.pbio.3001503 (CF568 SE, CF680 SE, super-resolution, dSTORM,)

Schmidt, C., et al. [The Dual Role of a Polyvalent IgM/IgA-Enriched Immunoglobulin Preparation in Activating and Inhibiting the Complement System](#). Biomedicines 9(7), (2021), DOI: 10.3390/biomedicines9070817 (CF488A DAGt CF568 DAM WB)

Schmitner, N., et al. [Differential Responses of Neural Retina Progenitor Populations to Chronic Hyperglycemia](#). Cells 10, (2021), DOI: 10.3390/cells10113265 (CF568 PNA)

Seo, J., et al. [PICASSO allows ultra-multiplexed fluorescence imaging of spatially overlapping proteins without reference spectra measurements](#). Nat Commun 13, 2475, (2022), DOI: 10.1038/s41467-022-30168-z (CF405S, CF405M, CF405L, CF488A, CF568, CF633, CF660R, CF680R)

Shi, L., et al. [Non-thermal plasma induces a stress response in mesothelioma cells resulting in increased endocytosis, lysosome biogenesis and autophagy](#). Free Radic Biol Med 108, 904-917, (2017), DOI: 10.1016/j.freeradbiomed.2017.04.368 (CF488A & CF568 secondaries)

Shibata, T., et al. [Crosslinking of a Peritrophic Matrix Protein Protects Gut Epithelia from Bacterial Exotoxins](#). PLoS Pathog 11(10), e1005244, (2015), DOI: 10.1371/journal.ppat.1005244 (CF488A goat anti-mouse, CF568 goat anti-mouse and CF568 goat anti-rabbit)

Shiheido, H., et al. [A phthalimide derivative that inhibits centrosomal clustering is effective on multiple myeloma](#). PLoS One 7(6), e38878, (2012), DOI: DOI:10.1371/journal.pone.0038878 (CF568 anti-rabbit IgG; ICC)

Shimizu, F., et al. [Glucose-regulated protein 78 autoantibody associates with blood-brain barrier disruption in neuromyelitis optica](#). Sci Transl Med 9(397), (2017), DOI: 10.1126/scitranslmed.aai9111 (CF568 Mix-n-Stain)

Sonoda, Y., et al. [Applications of cell resealing to reconstitute microRNA loading to extracellular vesicles](#). Sci Rep 11(1), 2900, (2021), DOI: 10.1038/s41598-021-82452-5 (CF568 Dextran 10K exosomes STORM)

Sørensen, G., et al. [Disrupted PDZ-domain binding motif of the dopamine transporter uniquely alters nanoscale distribution, dopamine homeostasis and reward motivation](#). BioRxiv, (2021), DOI: 10.1101/2021.01.06.425643 (CF568 succinimidyl ester 2-color STORM)

Sullivan, P. M., et al. [The ALS/FTLD associated protein C9orf72 associates with SMCR8 and WDR41 to regulate the autophagy-lysosome pathway](#). Acta Neuropathol Commun 4(1), 51, (2016), DOI: 10.1186/s40478-016-0324-5 (CF568 donkey anti-mouse)

Tashiro, S., et al. [Odour-induced analgesia mediated by hypothalamic orexin neurons in mice](#). Sci Rep 6, 37129, (2016), DOI: 10.1038/srep37129 (CF568 donkey anti-goat)

Turkowyd, B., et al. [From single molecules to life: microscopy at the nanoscale](#). Anal Bioanal Chem 408(25), 6885-6911, (2016), DOI: 10.1007/s00216-016-9781-8, 10.1007/s00216-016-9781-8 [pii] (CF568, CF647, CF660C, CF680 super res review)

Uchida, T., et al. [CUL2-mediated clearance of misfolded TDP-43 is paradoxically affected by VHL in oligodendrocytes in ALS](#). Sci Rep 6, 19118, (2016), DOI: 10.1038/srep19118 (CF568 goat anti-rabbit)

Uesaka, T., et al. [Neuronal Differentiation in Schwann Cell Lineage Underlies Postnatal Neurogenesis in the Enteric Nervous System](#). J Neurosci 35(27), 9879-9888, (2015), DOI: 10.1523/JNEUROSCI.1239-15.2015 (CF568 & CF640R secondaries)

Varriale, A., et al. [On the possibility of ephedrine detection: time-resolved fluorescence resonance energy transfer \(FRET\)-based approach](#). Anal Bioanal Chem 408, 6329-6336, (2016), DOI: 10.1007/s00216-016-9738-y (CF568, CF647, FRET)

Wang, H., et al. [Functional Hydrogels with Chondroitin Sulfate Release Properties Regulate the Angiogenesis Behaviors of Endothelial Cells](#). Gels 8, (2022), DOI: 10.3390/gels8050261 (CF568 phalloidin)

Wang, Y., et al. [Cucurbitacin IIb exhibits anti-inflammatory activity through modulating multiple cellular behaviors of mouse lymphocytes](#). PLoS One 9(2), e89751, (2014), DOI: 10.1371/journal.pone.0089751, PONE-D-13-38697 [pii] (CF488A goat anti-mouse, CF568 goat anti-rabbit; ICC)

Watson, J. F., et al. [AMPA receptor anchoring at CA1 synapses is determined by N-terminal domain and TARP gamma8 interactions](#). Nat Commun 12(1), 5083, (2021), DOI: 10.1038/s41467-021-25281-4 (CF568 succinimidyl ester)

Willems, J., and MacGillavry, H. D. [A coordinate-based co-localization index to quantify and visualize spatial associations in single-molecule localization microscopy](#). Sci Rep 12, 4676, (2022), DOI: 10.1038/s41598-022-08746-4 (CF568 secondaries, TIRF)

Wisnovsky, S., et al. [Genome-wide CRISPR screens reveal a specific ligand for the glycan-binding immune checkpoint receptor Siglec-7](#). Proc Natl Acad Sci U S A 118(5), (2021), DOI: 10.1073/pnas.2015024118 (CF568 Mix-n-Stain STORM)

Wolf, K. J., et al. [A mode of cell adhesion and migration facilitated by CD44-dependent microtentacles](#). Proc Natl Acad Sci U S A 117(21), 11432-11443, (2020), DOI: 10.1073/pnas.1914294117 (CF568 STORM)

Wu, L. and Lambert, J. D. [A serpin is required for ectomesoderm, a hallmark of spiralian development](#). Dev Biol 469, 172-181, (2021), DOI: 10.1016/j.ydbio.2020.10.011 (CF568 10KDa Dextran)

Xu, H., et al. [Molecular organization of mammalian meiotic chromosome axis revealed by expansion STORM microscopy](#). Proc Natl Acad Sci U S A 116(37), 18423-18428, (2019), DOI: 10.1073/pnas.1902440116 (CF568 ExSTORM)

Yamamoto, S., et al. [Contribution of calumin to embryogenesis through participation in the endoplasmic reticulum-associated degradation activity](#). Dev Biol 393(1), 33-43, (2014), DOI: S0012-1606(14)00320-0 [pii] 10.1016/j.ydbio.2014.06.024 (CF488A Mix-n-Stain, CF568 Mix-n-Stain; IHC)

Yang, H., et al. [Role of the sigma-1 receptor chaperone in rod and cone photoreceptor degenerations in a mouse model of retinitis pigmentosa](#). Mol Neurodegener 12(1), 68, (2017), DOI: 10.1186/s13024-017-0202-z (CF568 tyramide)

Yan, R., et al. [SURF4-induced tubular ERGIC selectively expedites ER-to-Golgi transport](#). Dev Cell, (2022), DOI: 10.1016/j.devcel.2021.12.018 (CF568 succinimidyl ester)

Yang, W., et al. [Regulation of Meristem Morphogenesis by Cell Wall Synthases in Arabidopsis](#). Curr Biol 26(11), 1404-1415, (2016), DOI: 10.1016/j.cub.2016.04.026 (CF488A & CF568 secondaries)

Yang, Z., et al. [USP18 negatively regulates NF-kappaB signaling by targeting TAK1 and NEMO for deubiquitination through distinct mechanisms](#). Sci Rep 5, 12738, (2015), DOI: 10.1038/srep12738 (CF568 goat anti-mouse)

Yin, C. and Li, P. [Growth Suppression of Glioma Cells Using HDAC6 Inhibitor, Tubacin](#). Open Med (Wars) 13, 221-226, (2018), DOI: 10.1515/med-2018-0034 (CF488A goat anti-rabbit CF568 goat anti-mouse)

Zanetti-Domingues, L. C., et al. [Hydrophobic fluorescent probes introduce artifacts into single molecule tracking experiments due to non-specific binding](#). PLoS One 8(9), e74200, (2013), DOI: 10.1371/journal.pone.0074200, PONE-D-13-17594 [pii] (CF488A, CF568, CF633 & CF640R maleimide, TIRF)

Zha, Q. B., et al. [ATP-Induced Inflammasome Activation and Pyroptosis Is Regulated by AMP-Activated Protein Kinase in Macrophages](#). Front Immunol 7, 597, (2016), DOI: 10.3389/fimmu.2016.00597 (CF568 goat anti-rabbit)

Zha, Q. B., et al. [Cucurbitacin E Induces Autophagy via Downregulating mTORC1 Signaling and Upregulating AMPK Activity](#). PLoS One 10(5), e0124355, (2015), DOI: 10.1371/journal.pone.0124355 (CF488A goat anti-mouse & CF568 goat anti-rabbit)

Zhang, M., et al. [Translocation of interleukin-1beta into a vesicle intermediate in autophagy-mediated secretion](#). eLife 4, (2015), DOI: 10.7554/eLife.11205 (CF568 STORM)

Zhang, X., et al. [Teneurins assemble into presynaptic nanoclusters that promote synapse formation via postsynaptic non-teneurin ligands](#). Nat Commun 13, 2297, (2022), DOI: 10.1038/s41467-022-29751-1 (CF568 secondary, super-resolution, dSTORM)

Zhang, Y., et al. [Endothelial Cell Migration Regulated by Surface Topography of Poly\(epsilon-caprolactone\) Nanofibers](#). ACS Biomater Sci Eng 7, 4959-4970, (2021), DOI: 10.1021/acsbiomaterials.1c00951 (CF568, phalloidin)

Zhang, Y. T., et al. [Formation of cofilin-actin rods following cucurbitacin-B-induced actin aggregation depends on slingshot homolog 1-mediated cofilin hyperactivation](#). J Cell Biochem 114(10), 2415, (2013), DOI: DOI:10.1002/jcb.24587 (CF488A goat anti-mouse, CF568 goat anti-rabbit; ICC)

Zhang, Y. T., et al. [VASP activation via the Galpha13/RhoA/PKA pathway mediates cucurbitacin-B-induced actin aggregation and cofilin-actin rod formation](#). PLoS One 9(4), e93547, (2014), DOI: 10.1371/journal.pone.0093547, PONE-D-13-34310 [pii] (CF488A goat anti-mouse, CF568 goat anti-rabbit; ICC)

Zhao, K., et al. [Muscle Yap Is a Regulator of Neuromuscular Junction Formation and Regeneration](#). J Neurosci 37(13), 3465-3477, (2017), DOI: 10.1523/JNEUROSCI.2934-16.2017 (CF568 bungarotoxin)

Zhu, J. S., et al. [Cucurbitacin B induces cell cycle arrest, apoptosis and autophagy associated with G actin reduction and persistent activation of cofilin in Jurkat cells](#). Pharmacology 89(5-6), 348-346, (2012), DOI: DOI:10.1159/000338757 (CF488A goat anti-mouse, CF568 goat anti-rabbit; ICC)

CF®583R

Chen, K., et al. [Super-multiplexing excitation spectral microscopy with multiple fluorescence bands](#). Biomed Opt Express 13, 6048-6060, (2022), DOI: 10.1364/BOE.473241 (CF583R, CF620R, CF514, LipidSpot488, super-resolution, Excitation spectral microscopy)

Wang, B., et al. [Transforming Rhodamine Dyes for \(d\)STORM Super-Resolution Microscopy via 1,3-Disubstituted Imidazolium Substitution](#). Angew Chem Int Ed Engl, e202113612, (2021), DOI: 10.1002/anie.202113612 (CF597R, CF583R, dSTORM)

CF®594

Ahi, Y. S., et al. [Functional Interplay Between Murine Leukemia Virus Glycag, Serinc5, and Surface Glycoprotein Governs Virus Entry, with Opposite Effects on Gammaretroviral and Ebolavirus Glycoproteins](#). mBio 7(6), (2016), DOI: 10.1128/mBio.01985-16 (CF488A goat anti-rabbit & CF594 donkey anti-mouse)

Akhmetshina, A., et al. [Activation of canonical Wnt signalling is required for TGF-beta-mediated fibrosis](#). Nat Commun 3, 735, (2012), DOI: DOI:10.1038/ncomms1734 (CF488A and CF594 streptavidin; IHC, ICC)

Arul, M., et al. [Culture of low passage colorectal cancer cells and demonstration of variation in selected tumour marker expression](#). Cytotechnology 66(3), 481, (2013), DOI: 10.1007/s10616-013-9600-4 (CF488A Mix-n-Stain, CF594 Mix-n-Stain, CF488A goat anti-mouse; ICC)

Agustinus, A., et al. [Epigenetic dysregulation from chromosomal transit in micronuclei](#). (2022), DOI: 10.1101/2022.01.12.475944 (CF594 tyramide)

Azzedine, H., et al. [PLEKHG5 deficiency leads to an intermediate form of autosomal-recessive Charcot-Marie-Tooth disease](#). Hum Mol Genet 22(20), 4224, (2013), DOI: DOI:10.1093/hmg/ddt274 (CF594 a-bungarotoxin; IHC)

Babu, R., et al. [Interphase Chromosome Profiling: A Method for Conventional Banded Chromosome Analysis Using Interphase Nuclei](#). Arch Pathol Lab Med 142(2), 213-228, (2018), DOI: 10.5858/arpa.2016-0621-OA (CF594 dUTP)

Balez, R., et al. [Neuroprotective effects of apigenin against inflammation, neuronal excitability and apoptosis in an induced pluripotent stem cell model of Alzheimer's disease](#). Sci Rep 6, 31450, (2016), DOI: 10.1038/srep31450 (CF594 Annexin V, NucView primary neurons)

Bauer, S., et al. [Functional coupling of human pancreatic islets and liver spheroids on-a-chip: Towards a novel human ex vivo type 2 diabetes model](#). Sci Rep 7(1), 14620, (2017), DOI: 10.1038/s41598-017-14815-w (CF594 goat anti-guinea pig, CF594 goat anti-mouse and CF594 goat anti-rabbit)

Bernard-Marissal, N., et al. [Dysfunction in endoplasmic reticulum-mitochondria crosstalk underlies SIGMAR1 loss of function mediated motor neuron degeneration](#). Brain 138(Pt 4), 875-890, (2015), DOI: 10.1093/brain/awv008 (CF594 bungarotoxin)

Bielska, E., et al. [Pathogen-derived extracellular vesicles mediate virulence in the fatal human pathogen Cryptococcus gattii](#). Nat Commun 9(1), 1556, (2018), DOI: 10.1038/s41467-018-03991-6 (CF594 goat anti-mouse)

Boratko, A. and Csortos, C. [NHERF2 is crucial in ERM phosphorylation in pulmonary endothelial cells](#). Cell Commun Signal 11, 99, (2013), DOI: 1478-811X-11-99 [pii] 10.1186/1478-811X-11-99 (CF594 phalloidin; ICC)

Chew, W. L., et al. [A multifunctional AAV-CRISPR-Cas9 and its host response](#). Nat Methods 13(10), 868-874, (2016), DOI: nmeth.3993 [pii] 10.1038/nmeth.3993 (CF633 goat anti-mouse & CF594 goat anti-mouse)

Daly, D. T. and Ariel, M. [A novel cerebellar commissure and other myelinated axons in the Purkinje cell layer of a pond turtle \(*Trachemys scripta elegans*\)](#). J Comp Neurol 526(17), 2802-2823, (2018), DOI: 10.1002/cne.24528 (CF488A and CF594 10K dextran)

Elting, M. W., et al. [Single-molecule fluorescence imaging of processive myosin with enhanced background suppression using linear zero-mode waveguides \(ZMWs\) and convex lens induced confinement \(CLIC\)](#). Opt Express 21(1), 1189-1202, (2013), DOI: 248311 [pii] (CF594 anti-GFP; single molecule imaging)

Felix-Ortiz, A. C., et al. [BLA to vHPC inputs modulate anxiety-related behaviors](#). Neuron 79(4), 658-664, (2013), DOI: S0896-6273(13)00532-1 [pii] 10.1016/j.neuron.2013.06.016 (CF594 Streptavidin; IHC)

Feng, X., et al. [Receptor-interacting protein 140 attenuates endoplasmic reticulum stress in neurons and protects against cell death](#). Nat Commun 5, 4487, (2014), DOI: ncomms5487 [pii] 10.1038/ncomms5487 (CF594-dUTP; tunel assay)

Feng, X., et al. [beta-arrestin protects neurons by mediating endogenous opioid arrest of inflammatory microglia](#). Cell Death Differ 21(3), 397-406, (2014), DOI: cdd2013152 [pii] 10.1038/cdd.2013.152 (CF594-dUTP; tunel assay)

Frontera, J. L., et al. [Bidirectional control of fear memories by the cerebellum through the ventrolateral periaqueductal grey](#). (2020), DOI: 10.1101/2020.02.19.956375 (CF594 cholera toxin B)

Galletti, G., et al. [Isolation of breast cancer and gastric cancer circulating tumor cells by use of an anti HER2-based microfluidic device](#). Lab Chip 14(1), 147-156, (2014), DOI: 10.1039/c3lc51039e (CF594 Mix-n-Stain; ICC)

Geherin, S. A., et al. [IL-10+ Innate-like B Cells Are Part of the Skin Immune System and Require alpha4beta1 Integrin To Migrate between the Peritoneum and Inflamed Skin](#). J Immunol 196(6), 2514-2525, (2016), DOI: 10.4049/jimmunol.1403246 (CF488A CF594 Mix-n-Stain)

Grix, T., et al. [Bioprinting Perfusion-Enabled Liver Equivalents for Advanced Organ-on-a-Chip Applications](#). Genes (Basel) 9(4), (2018), DOI: 10.3390/genes9040176 (CF594 goat anti-mouse & CF594 goat anti-rabbit)

Gonzalez Pisfil, M., et al. [Stimulated emission depletion microscopy with a single depletion laser using five fluorochromes and fluorescence lifetime phasor separation](#). Sci Rep 12, 14027, (2022), DOI: 10.1038/s41598-022-17825-5 (CF594 WGA, CF680R DAMs, super-resolution, confocal, FLIM, fluorescence microscopy, 2-Photon)

Harris, M. J., et al. [Quantifying persistence in the T-cell signaling network using an optically controllable antigen receptor](#). Mol Syst Biol 17(5), e10091, (2021), DOI: 10.15252/msb.202010091 (CF594)

Hashimoto, Y. and Tanaka, H. [Ongoing replication forks delay nuclear envelope breakdown upon mitotic entry](#). J Biol Chem, (2020), DOI: 10.1074/jbc.RA120.015142 (CF594 dUTP)

Higa, M., et al. [TRF2 recruits ORC through TRFH domain dimerization](#). *Biochim Biophys Acta Mol Cell Res* 1864(1), 191-201, (2017), DOI: 10.1016/j.bbamcr.2016.11.004 (CF488A goat anti-rabbit, CF594 donkey anti-rat, CF594 goat anti-mouse)

Hindul, N. L., et al. [Construction of a human hTERT RPE-1 cell line with inducible Cre for editing of endogenous genes](#). *Biol Open* 11(2), (2022), DOI: 10.1242/bio.059056 (CF594 goat anti-mouse IgG1)

Holton, A. B., et al. [Microfluidic Biopsy Trapping Device for the Real-Time Monitoring of Tumor Microenvironment](#). *PLoS One* 12(1), e0169797, (2017), DOI: 10.1371/journal.pone.0169797 (CF594 Annexin V)

Hwang, J., et al. [Aptamer-conjugated live human immune cell based biosensors for the accurate detection of C-reactive protein](#). *Sci Rep* 6, 34778, (2016), DOI: 10.1038/srep34778 (CF594 Mix-n-Stain)

Hwang, R. D., et al. [The neuroprotective effect of human uncoupling protein 2 \(hUCP2\) requires cAMP-dependent protein kinase in a toxin model of Parkinson's disease](#). *Neurobiol Dis* 69, 180-191, (2014), DOI: S0969-9961(14)00150-8 [pii] 10.1016/j.nbd.2014.05.032 (CF594 Annexin V; IHC)

Inutsuka, A., et al. [Concurrent and robust regulation of feeding behaviors and metabolism by orexin neurons](#). *Neuropharmacology* 85, 451-460, (2014), DOI: S0028-3908(14)00237-8 [pii] 10.1016/j.neuropharm.2014.06.015 (CF488 CF594 CF647; IHC)

Inutsuka, A., et al. [The integrative role of orexin/hypocretin neurons in nociceptive perception and analgesic regulation](#). *Sci Rep* 6, 29480, (2016), DOI: 10.1038/srep29480 (CF488A, CF594 & CF647 secondaries)

Jalewa, J., et al. [Neural circuit interactions between the dorsal raphe nucleus and the lateral hypothalamus: an experimental and computational study](#). *PLoS One* 9(2), e88003, (2014), DOI: 10.1371/journal.pone.0088003 PONE-D-13-34080 [pii] (CF594 donkey anti-chicken CF633 donkey anti-guinea pig; IHC)

Jeon, E. J., et al. [rgs-CaM Detects and Counteracts Viral RNA Silencing Suppressors in Plant Immune Priming](#). *J Virol* 91(19), (2017), DOI: 10.1128/JVI.00761-17 (CF594 goat anti-rabbit)

Kaneda, M., et al. [Synaptic localisation of SRF coactivators, MKL1 and MKL2, and their role in dendritic spine morphology](#). *Sci Rep* 8(1), 727, (2018), DOI: 10.1038/s41598-017-18905-7 (CF488A goat anti-rabbit, CF488A goat anti-mouse and CF594 goat anti-guinea pig)

Kapoor, N., et al. [Human Granuloma In Vitro Model, for TB Dormancy and Resuscitation](#). *PLoS One* 8(1), e53657, (2013), DOI: DOI:10.1371/journal.pone.0053657 (CF594 TUNEL Assay Apoptosis Detection Kit; TUNEL assay)

Kijani, S., et al. [Filter-Dense Multicolor Microscopy](#). *PLoS One* 10(3), e0119499, (2015), DOI: 10.1371/journal.pone.0119499 PONE-D-14-20661 [pii] (CF488A, CF543, CF594)

Koban, R., et al. [A novel three-dimensional cell culture method enhances antiviral drug screening in primary human cells](#). *Antiviral Res* 150, 20-29, (2018), DOI: 10.1016/j.antiviral.2017.12.005 (CF594 wheat germ agglutinin)

Krolewski, D. M., et al. [Quantitative validation of immunofluorescence and lectin staining using reduced CLARITY acrylamide formulations](#). *Brain Struct Funct* 223(2), 987-999, (2018), DOI: 10.1007/s00429-017-1583-z (CF594 CF647)

Kume, K. and Nishizuka, S. S. [Colony Lysate Arrays for Proteomic Profiling of Drug-Tolerant Persisters of Cancer Cell](#). Anal Chem 89(17), 8626-8631, (2017), DOI: 10.1021/acs.analchem.7b01215 (CF488A goat anti-mouse & CF594 goat anti-rabbit)

Lannin, T. B., et al. [Comparison and optimization of machine learning methods for automated classification of circulating tumor cells](#). Cytometry A 89(10), 922-931, (2016), DOI: 10.1002/cyto.a.22993 (CF543 pan anti-cytokeratin & CF594 Mix-n-Stain)

Lavinsky, D., et al. [Nondamaging Retinal Laser Therapy: Rationale and Applications to the Macula](#). Invest Ophthalmol Vis Sci 57(6), 2488-2500, (2016), DOI: 10.1167/iovs.15-18981 (CF594 goat anti-mouse IHF)

Maenohara, S., et al. [Role of UHRF1 in de novo DNA methylation in oocytes and maintenance methylation in preimplantation embryos](#). PLoS Genet 13(10), e1007042, (2017), DOI: 10.1371/journal.pgen.1007042 (CF488A donkey anti-rabbit, CF488A donkey anti-rat and CF594 donkey anti-rabbit)

Moller, R., et al. [IgG trafficking in the adult pig small intestine: one- or bidirectional transfer across the enterocyte brush border?](#). Histochem Cell Biol 147(3), 399-411, (2017), DOI: 10.1007/s00418-016-1492-x (CF594 hydrazide)

Nabeshima, Y., et al. [Calpain 1 inhibitor BDA-410 ameliorates alpha-klotho-deficiency phenotypes resembling human aging-related syndromes](#). Sci Rep 4, 5847, (2014), DOI: srep05847 [pii] 10.1038/srep05847 (CF594 goat anti-rat & CF488 goat anti-chicken; IHC)

Ohira, M., et al. [A novel anti-microtubule agent with carbazole and benzohydrazide structures suppresses tumor cell growth in vivo](#). Biochim Biophys Acta 1850(9), 1676-1684, (2015), DOI: 10.1016/j.bbagen.2015.04.013 (CF488A goat anti-rabbit, CF594 goat anti-mouse)

Olasz, F., et al. [Immunological and biochemical characterisation of 7ap, a short protein translated from an alternative frame of ORF7 of PRRSV](#). Acta Vet Hung 64(2), 273-287, (2016), DOI: 10.1556/004.2016.027 (CF568 goat anti-swine CF594 goat anti-mouse)

Orr, J. S., et al. [CC-chemokine receptor 7 \(CCR7\) deficiency alters adipose tissue leukocyte populations in mice](#). Physiol Rep 4(18), (2016), DOI: 10.14814/phy2.12971 (CF594 succinimidyl ester)

Perez Bay, A. E., et al. [Galectin-4-mediated transcytosis of transferrin receptor](#). J Cell Sci 127(20), 4457, (2014), DOI: jcs.153437 [pii] 10.1242/jcs.153437 (CF594 succinimidyl ester labeled transferrin; fluorescence microscopy)

Perez Bay, A. E., et al. [The kinesin KIF16B mediates apical transcytosis of transferrin receptor in AP-1B-deficient epithelia](#). EMBO J 32(15), 2125-2139, (2013), DOI: emboj2013130 [pii] 10.1038/emboj.2013.130 (CF594 succinimidyl ester labeled transferrin; fluorescence microscopy)

Pignoloni, B., et al. [Distinct Roles for Human Cytomegalovirus Immediate Early Proteins IE1 and IE2 in the Transcriptional Regulation of MICA and PVR/CD155 Expression](#). J Immunol 197(10), 4066-4078, (2016), DOI: 10.4049/jimmunol.1502527 (CF594 rabbit anti-mouse)

Pontes-Quero, S., et al. [Dual ifgMosaic: A Versatile Method for Multispectral and Combinatorial Mosaic Gene-Function Analysis](#). Cell 170(4), 800-814 e818, (2017), DOI: 10.1016/j.cell.2017.07.031 (CF594 Rabbit anti-RFP)

Prapa, E., et al. [Effect of Anti-Mullerian hormone \(AMH\) and bone morphogenetic protein 15 \(BMP-15\) on steroidogenesis in primary-cultured human luteinizing granulosa cells through Smad5 signalling](#). J Assist

Reprod Genet 32(7), 1079-1088, (2015), DOI: 10.1007/s10815-015-0494-2 (CF488A donkey anti-rabbit & CF594 donkey anti-mouse)

Qu, X., et al. [Relative impact of uniaxial alignment vs. form-induced stress on differentiation of human adipose derived stem cells](#). Biomaterials 34(38), 9812-9818, (2013), DOI: S0142-9612(13)01092-2 [pii] 10.1016/j.biomaterials.2013.09.009 (CF594 goat anti-rabbit; fluorescence microscopy)

Reichhart, N., et al. [Rab27a GTPase modulates L-type Ca₂₊ channel function via interaction with the II-III linker of CaV1.3 subunit](#). Cell Signal 27(11), 2231-2240, (2015), DOI: 10.1016/j.cellsig.2015.07.023 (CF594 donkey anti-goat)

Sawano, S., et al. [A One-Step Immunostaining Method to Visualize Rodent Muscle Fiber Type within a Single Specimen](#). PLoS One 11(11), e0166080, (2016), DOI: 10.1371/journal.pone.0166080 PONE-D-16-33181 [pii] (CF594 streptavidin)

Sheu, J. R., et al. [A novel indication of platonin, a therapeutic immunomodulating medicine, on neuroprotection against ischemic stroke in mice](#). Sci Rep 7, 42277, (2017), DOI: 10.1038/srep42277 (CF488A donkey anti-mouse, CF488A donkey anti-rabbit, CF594 donkey anti-mouse and CF594 donkey anti-rabbit)

Stoltzfus, C. R., et al. [CytoMAP: A Spatial Analysis Toolbox Reveals Features of Myeloid Cell Organization in Lymphoid Tissues](#). Cell Rep 31(3), 107523, (2020), DOI: 10.1016/j.celrep.2020.107523 (CF750, CF594, CF514, CF633 Mix-n-Stain)

Trofimenko, E., et al. [The endocytic pathway taken by cationic substances requires Rab14 but not Rab5 and Rab7](#). Cell Rep 37, 109945, (2021), DOI: 10.1016/j.celrep.2021.109945 (CF405M, CF488A, CF594, CF647, Mix-n-Stain, fluorescence microscopy)

Trofimenko, E., et al. [Identification of a new, Rab14-dependent, endo-lysosomal pathway](#). bioRxiv, (2021), DOI: 10.1101/2021.03.25.436964 (CF405M, CF594, Mix-n-Stain)

Wagner, M. C., et al. [Mechanism of increased clearance of glycated albumin by proximal tubule cells](#). Am J Physiol Renal Physiol 310(10), F1089-1102, (2016), DOI: 10.1152/ajprenal.00605.2015 (CF594 2-photon)

Wang, Q., et al. [The hepatokine TSK maintains myofiber integrity and exercise endurance and contributes to muscle regeneration](#). JCI Insight, (2022), DOI: 10.1172/jci.insight.154746 (CF594 bungarotoxin)

Wang, X., et al. [The somite-secreted factor Maeg promotes zebrafish embryonic angiogenesis](#). Oncotarget 7, 77749-77763, (2016), DOI: <https://doi.org/10.18632/oncotarget.12793> (CF594 donkey anti-rat)

Watari, K., et al. [Tumor-derived interleukin-1 promotes lymphangiogenesis and lymph node metastasis through M2-type macrophages](#). PLoS One 9(6), e99568, (2014), DOI: 10.1371/journal.pone.0099568 PONE-D-13-45479 [pii] (CF488A DARat & CF594 donkey anti-rabbit; IHC)

Yamashita, A., et al. [Increased metabolite levels of glycolysis and pentose phosphate pathway in rabbit atherosclerotic arteries and hypoxic macrophage](#). PLoS One 9(1), e86426, (2014), DOI: 10.1371/journal.pone.0086426 PONE-D-13-24660 [pii] (CF488A Mix-n-Stain, CF594 Mix-n-Stain; IHC)

Yates, J. L., et al. [T Cell-Dependent IgM Memory B Cells Generated during Bacterial Infection Are Required for IgG Responses to Antigen Challenge](#). *J Immunol* 191(3), 1240-1249, (2013), DOI: DOI:10.4049/jimmunol.1300062 (PE-CF594 conjugated anti-CD19 (BD Biosciences); IHC)

Yu, C. H., et al. [Integrin-matrix clusters form podosome-like adhesions in the absence of traction forces](#). *Cell Rep* 5(5), 1456-1468, (2013), DOI: S2211-1247(13)00639-6 [pii] 10.1016/j.celrep.2013.10.040 (CF594 phalloidin; ICC)

Zaichik, S., et al. [Development and In Vitro Characterization of Transferrin-Decorated Nanoemulsion Utilizing Hydrophobic Ion Pairing for Targeted Cellular Uptake](#). *Journal of Pharmaceutical Innovation*, (2021), DOI: 10.1007/s12247-021-09549-2 (CF594 Transferrin)

CF®597R

Wang, B., et al. [Transforming Rhodamine Dyes for \(d\)STORM Super-Resolution Microscopy via 1,3-Disubstituted Imidazolium Substitution](#). *Angew Chem Int Ed Engl*, e202113612, (2021), DOI: 10.1002/anie.202113612 (CF597R, CF583R dSTORM)

CF®620R

Chen, K., et al. [Super-multiplexing excitation spectral microscopy with multiple fluorescence bands](#). *Biomed Opt Express* 13, 6048-6060, (2022), DOI: 10.1364/BOE.473241 (CF583R, CF620R, CF514, LipidSpot488, super-resolution, Excitation spectral microscopy)

CF®633

Ache, J. M., et al. [A direct descending pathway informing locomotor networks about tactile sensor movement](#). *J Neurosci* 35(9), 4081-4091, (2015), DOI: 10.1523/JNEUROSCI.3350-14.2015 (CF633 Streptavidin)

Balogh, A., et al. [Sex hormone-binding globulin provides a novel entry pathway for estradiol and influences subsequent signaling in lymphocytes via membrane receptor](#). *Sci Rep* 9(1), 4, (2019), DOI: 10.1038/s41598-018-36882-3 (CF633 Mix-n-Stain)

Blomeke, L., et al. [Quantitative detection of alpha-Synuclein and Tau oligomers and other aggregates by digital single particle counting](#). *NPJ Parkinsons Dis* 8, 68, (2022), DOI: 10.1038/s41531-022-00330-x (CF488A SE, CF633 SE, TIRF)

Chew, W. L., et al. [A multifunctional AAV-CRISPR-Cas9 and its host response](#). *Nat Methods* 13(10), 868-874, (2016), DOI: nmeth.3993 [pii] 10.1038/nmeth.3993 (CF633 goat anti-mouse & CF594 goat anti-mouse)

Cutler, C. P., et al. [Aquaporin \(AQP\) channels in the spiny dogfish, *Squalus acanthias* II: Localization of AQP3, AQP4 and AQP15 in the kidney](#). *Comp Biochem Physiol B Biochem Mol Biol* 258, 110701, (2022), DOI: 10.1016/j.cbpb.2021.110701 (CF633)

Egashira, Y., et al. [Spontaneously Recycling Synaptic Vesicles Constitute Readily Releasable Vesicles in Intact Neuromuscular Synapses](#). *J Neurosci* 42, 3523-3536, (2022), DOI: 10.1523/JNEUROSCI.2005-21.2022 (CF633 alpha-bungarotoxin, fluorescence microscopy)

Gaudreau-Lapierre, A., et al. [Expansion microscopy-based imaging of nuclear structures in cultured cells](#). STAR Protocols 2(3), (2021), DOI: 10.1016/j.xpro.2021.100630 (CF633 goat anti-mouse & CF633 goat anti-rabbit)

Grob, R., et al. [Rotation of skylight polarization during learning walks is necessary to trigger neuronal plasticity in Cataglyphis ants](#). Proc Biol Sci 289(1967), 20212499, (2022), DOI: 10.1098/rspb.2021.2499 (CF633 phalloidin)

Huang, T., et al. [Simultaneous Multicolor Single-Molecule Tracking with Single-Laser Excitation via Spectral Imaging](#). Biophys J 114(2), 301-310, (2018), DOI: S0006-3495(17)31242-0 [pii] 10.1016/j.bpj.2017.11.013 (CF633, CF660C, CF680, CF680R STORM)

Jalewa, J., et al. [Neural circuit interactions between the dorsal raphe nucleus and the lateral hypothalamus: an experimental and computational study](#). PLoS One 9(2), e88003, (2014), DOI: 10.1371/journal.pone.0088003 PONE-D-13-34080 [pii] (CF594 donkey anti-chicken CF633 donkey anti-guinea-pig; IHC)

Jung, Y., et al. [CD45 pre-exclusion from the tips of T cell microvilli prior to antigen recognition](#). Nat Commun 12(1), 3872, (2021), DOI: 10.1038/s41467-021-23792-8 (CF633)

Kaljas, Y., et al. [Human adenosine deaminases ADA1 and ADA2 bind to different subsets of immune cells](#). Cell Mol Life Sci 74(3), 555-570, (2017), DOI: 10.1007/s00018-016-2357-0 (CF633 succinimidyl ester)

Kawaguchi, S. Y. and Sakaba, T. [Fast Ca\(2+\) Buffer-Dependent Reliable but Plastic Transmission at Small CNS Synapses Revealed by Direct Bouton Recording](#). Cell Rep 21(12), 3338-3345, (2017), DOI: 10.1016/j.celrep.2017.11.072 (CF633 succinimidyl ester neuronal tracing)

Kawashima, Y., et al. [Detection of DNA double-strand breaks by pulsed-field gel electrophoresis](#). Genes Cells 22(1), 84-93, (2017), DOI: 10.1111/gtc.12457 (CF555 donkey anti-rabbit & CF633 donkey anti-mouse)

Kim, H. J. and Selvin, P. R. [Fluorescence Imaging with One Nanometer Accuracy](#). in: SpringerReference, Springer, (2013), (CF633; Fluorescence Imaging with One Nanometer Accuracy (FIONA))

Klimas, A., et al. [Nanoscopic Imaging of Human Tissue Sections via Physical and Isotropic Expansion](#). J Vis Exp (151), (2019), DOI: 10.3791/60195 (CF488A CF633 goat anti-chicken)

Kohara, K., et al. [Cell type-specific genetic and optogenetic tools reveal hippocampal CA2 circuits](#). Nat Neurosci 17(2), 269-279, (2014), DOI: nn.3614 [pii] 10.1038/nn.3614 (CF633 goat anti-mouse, CF633 goat anti-rabbit, CF633 SA, CF405S goat anti-mouse; IHC)

Kunasegaran, K., et al. [Transcriptional repressor Tbx3 is required for the hormone-sensing cell lineage in mammary epithelium](#). PLoS One 9(10), e110191, (2014), DOI: 10.1371/journal.pone.0110191 (CF633 donkey anti-rat)

Li, T., et al. [Magnetic resonance imaging of folic acid-coated magnetite nanoparticles reflects tissue biodistribution of long-acting antiretroviral therapy](#). Int J Nanomedicine 10, 3779-3790, (2015), DOI: 10.2147/IJN.S83279 (CF633 succinimidyl ester)

Liang, Y., et al. [Biomimetic Tubular Matrix Induces Periodontal Ligament Principal Fiber Formation and Inhibits Osteogenic Differentiation of Periodontal Ligament Stem Cells](#). ACS Appl Mater Interfaces 14, 36451-36461, (2022), DOI: 10.1021/acsami.2c09420 (CF633 phalloidin, fluorescence microscopy)

Lindberg, O. R., et al. [Epidermal growth factor treatment of the adult brain subventricular zone leads to focal microglia/macrophage accumulation and angiogenesis](#). Stem Cell Reports 2(4), 440-448, (2014), DOI: 10.1016/j.stemcr.2014.02.003 S2213-6711(14)00055-1 [pii] (CF488A donkey anti-mouse CF555 donkey anti-rabbit, donkey anti-goat CF633 donkey anti-goat; IHC)

Lindberg, O. R., et al. [EGF-Induced Expansion of Migratory Cells in the Rostral Migratory Stream](#). PLoS One 7(9), e46380, (2012), DOI: DOI:10.1371/journal.pone.0046380 (CF488A donkey anti-mouse IgG; CF555 donkey anti-goat IgG; CF555 donkey anti-rabbit IgG; CF633 donkey anti-goat IgG; IHC)

Lozada, A. F., et al. [Glutamatergic Synapse Formation is Promoted by alpha7-Containing Nicotinic Acetylcholine Receptors](#). J Neurosci 32(22), 7651-7661, (2012), DOI: DOI:10.1523/JNEUROSCI.6246-11.2012 (CF555 donkey anti-rabbit, CF633 goat anti-guinea pig; IHC)

Lu, K., et al. [Progressive stretch enhances growth and maturation of 3D stem-cell-derived myocardium](#). Theranostics 11(13), 6138-6153, (2021), DOI: 10.7150/thno.54999 (CF633 wheat germ agglutinin)

Ma, L., et al. [Cryopreserved dental pulp tissues of exfoliated deciduous teeth is a feasible stem cell resource for regenerative medicine](#). PLoS One 7(12), e51777, (2012), DOI: DOI:10.1371/journal.pone.0051777 (CF633 secondary antibodies; ICC, IHC)

Martinez-Skinner, A. L., et al. [Cellular Responses and Tissue Depots for Nanoformulated Antiretroviral Therapy](#). PLoS One 10, (2015), DOI: 10.1371/journal.pone.0145966 (CF633 SE)

Miller, S. J. [Targeted detection of murine colonic dysplasia in vivo with flexible multispectral scanning fiber endoscopy](#). Journal of Biomedical Optics 17(2), (2012), DOI: DOI:10.1117/1.JBO.17.2.021103 (CF633 succinimidyl ester-labeled peptide, endoscopy)

Muoio, F., et al. [Development of a Biodegradable Microcarrier for the Cultivation of Human Adipose Stem Cells \(hASCs\) with a Defined Xeno- and Serum-Free Medium](#). Applied Sciences 11(3), (2021), DOI: 10.3390/app11030925 (CF633)

Okumura, R., et al. [Lypd8 promotes the segregation of flagellated microbiota and colonic epithelia](#). Nature 532(7597), 117-121, (2016), DOI: 10.1038/nature17406 (CF633 succinimidyl ester)

Oliveira, V. E. M., et al. [Oxytocin and vasopressin within the ventral and dorsal lateral septum modulate aggression in female rats](#). Nat Commun 12(1), 2900, (2021), DOI: 10.1038/s41467-021-23064-5 (CF633 streptavidin)

Orme, M. H., et al. [The unconventional myosin CRINKLED and its mammalian orthologue MYO7A regulate caspases in their signalling roles](#). Nat Commun 7, 10972, (2016), DOI: 10.1038/ncomms10972 (CF488A donkey anti-mouse, CF633 donkey anti-rabbit & CF633 donkey anti-goat)

Park, C. E., et al. [Super-Resolution Three-Dimensional Imaging of Actin Filaments in Cultured Cells and the Brain via Expansion Microscopy](#). ACS Nano 14(11), 14999-15010, (2020), DOI: 10.1021/acsnano.0c04915 (CF633 secondary expansion microscopy)

Pena-Hernandez, R., et al. [Genome-wide targeting of the epigenetic regulatory protein CTCF to gene promoters by the transcription factor TFII-I](#). Proc Natl Acad Sci U S A 112(7), E677-686, (2015), DOI: 10.1073/pnas.1416674112 (CF633 Annexin V flow)

Polackwich, R. J., et al. [A Novel 3D Fibril Force Assay Implicates Src in Tumor Cell Force Generation in Collagen Networks](#). PLoS One 8(3), e58138, (2013), DOI: DOI:10.1371/journal.pone.0058138 (CF633 succinimidyl ester (used to label collagen); 3D Fibril Force Assay, ICC)

Puligujja, P., et al. [Macrophage Folate Receptor-Targeted Antiretroviral Therapy Facilitates Drug Entry, Retention, Antiretroviral Activities and Biodistribution for Reduction of Human Immunodeficiency Virus Infections](#). Nanomedicine 9(8), 1263, (2013), DOI: DOI:10.1016/j.nano.2013.05.003 (CF633 succinimidyl ester nanoparticles; ICC)

Quinn, K. E., et al. [Activation of the CXCL12/CXCR4 signaling axis may drive vascularization of the ovine placenta](#). Domest Anim Endocrinol 47, 11-21, (2014), DOI: S0739-7240(13)00159-8 [pii] 10.1016/j.domaniend.2013.12.004 (CF633 goat anti-rabbit; IHC)

Ren, J., et al. [Sirt1 Protects Subventricular Zone-Derived Neural Stem Cells from DNA Double-Strand Breaks and Contributes to Olfactory Function Maintenance in Aging Mice](#). Stem Cells 40, 493-507, (2022), DOI: 10.1093/stmcls/sxac008 (CF488A, CF543, CF633 secondaries. fluorescence microscopy)

Reed, A. P., et al. [Dominant-Negative Effect of a Missense Variant in the TASK-2 \(KCNK5\) K⁺ Channel Associated with Balkan Endemic Nephropathy](#). PLoS One 11(5), e0156456, (2016), DOI: 10.1371/journal.pone.0156456 (CF633 wheat germ agglutinin, xenopus oocytes IF)

Roche, S. L., et al. [Loss of glial neurofascin155 delays developmental synapse elimination at the neuromuscular junction](#). J Neurosci 34(38), 12904-12918, (2014), DOI: 10.1523/JNEUROSCI.1725-14.2014 (CF633 bungarotoxin)

Romero, G., et al. [Localized Excitation of Neural Activity via Rapid Magnetothermal Drug Release](#). Advanced Functional Materials 26(35), 6471-6478, (2016), DOI: 10.1002/adfm.201602189 (CF633 amine)

Roy, D. S., et al. [Memory retrieval by activating engram cells in mouse models of early Alzheimer's disease](#). Nature 531(7595), 508-512, (2016), DOI: 10.1038/nature17172 (CF633 streptavidin)

Schmitt, F., et al. [Neuropeptides in the desert ant *Cataglyphis fortis*: Mass spectrometric analysis, localization, and age-related changes](#). J Comp Neurol 525(4), 901-918, (2017), DOI: 10.1002/cne.24109 (CF633 goat anti-mouse)

Senatore, A., et al. [Neuropeptidergic integration of behavior in Trichoplax adhaerens, an animal without synapses](#). J Exp Biol 220(Pt 18), 3381-3390, (2017), DOI: 10.1242/jeb.162396 (CF633 succinimidyl ester)

Seo, J., et al. [PICASSO: Ultra-multiplexed fluorescence imaging of biomolecules through single-round imaging and blind source unmixing](#). bioRxiv, (2021), DOI: 10.1101/2021.01.27.428247 (CF633, CF660R, CF680R)

Seo, J., et al. [PICASSO allows ultra-multiplexed fluorescence imaging of spatially overlapping proteins without reference spectra measurements](#). Nat Commun 13, 2475, (2022), DOI: 10.1038/s41467-022-30168-z (CF405S, CF405M, CF405L, CF488A, CF568, CF633, CF660R, CF680R)

Sevenich, M., et al. [Development of an alpha-synuclein fibril and oligomer specific tracer for diagnosis of Parkinson's disease, dementia with Lewy bodies and multiple system atrophy](#). Neurochem Int 161, 105422, (2022), DOI: 10.1016/j.neuint.2022.105422 (CF488A SE, CF633 maleimide)

Shiozaki, H. M. and Kazama, H. [Parallel encoding of recent visual experience and self-motion during navigation in Drosophila](#). Nat Neurosci 20(10), 1395-1403, (2017), DOI: 10.1038/nn.4628 (CF488A anti-rat, CF633 anti-mouse and CF633 anti-rabbit)

Simonson, P. D., et al. [Single-molecule-based super-resolution images in the presence of multiple fluorophores](#). Nano Lett 11(11), 5090-5096, (2011), DOI: DOI:10.1021/nl203560r (CF633 goat anti-mouse IgG; ICC, gSHRImP)

Starunov, V. V., et al. [A metameric origin for the annelid pygidium?](#). BMC Evol Biol 15, 25, (2015), DOI: 10.1186/s12862-015-0299-z 10.1186/s12862-015-0299-z [pii] (CF633 goat anti-mouse & CF555 goat anti-rabbit)

Stempel, H., et al. [Strain-specific Loss of Formyl Peptide Receptor 3 in the Murine Vomeronasal and Immune Systems](#). J Biol Chem 291(18), 9762-9775, (2016), DOI: 10.1074/jbc.M116.714493 (CF633 donkey anti-rat)

Stoltzfus, C. R., et al. [CytoMAP: A Spatial Analysis Toolbox Reveals Features of Myeloid Cell Organization in Lymphoid Tissues](#). Cell Rep 31(3), 107523, (2020), DOI: 10.1016/j.celrep.2020.107523 (CF750, CF594, CF514, CF633 Mix-n-Stain)

Sawanwongse, K. and Shabarek, N. [Clinical features and outcome of HIV/SARS-CoV-2 coinfected patients in The Bronx, New York city](#). J Med Virol 92(11), 2387-2389, (2020), DOI: 10.1002/jmv.26077 (CF633 goat anti-rabbit)

Tian, T., et al. [Dynamics of exosome internalization and trafficking](#). J Cell Physiol 228(7), 1487-1495, (2013), DOI: 10.1002/jcp.24304 (CF633 goat anti-rabbit; ICC)

Truckenbrodt, S., et al. [A practical guide to optimization in X10 expansion microscopy](#). Nat Protoc 14(3), 832-863, (2019), DOI: 10.1038/s41596-018-0117-3 (CF633 donkey anti-mouse expansion microscopy)

Vlahava, V. M., et al. [CD40 ligand exhibits a direct antiviral effect on Herpes Simplex Virus type-1 infection via a PI3K-dependent, autophagy-independent mechanism](#). Cell Signal 27(6), 1253-1263, (2015), DOI: S0898-6568(15)00078-9 [pii] 10.1016/j.cellsig.2015.03.002 (CF633 FC)

von Walden, F., et al. [Mechanical loading induces the expression of a Pol I regulon at the onset of skeletal muscle hypertrophy](#). Am J Physiol Cell Physiol 302(10), C1523-1530, (2012), DOI: DOI:10.1152/ajpcell.00460.2011 (CF488A goat anti-mouse IgG, CF633 goat anti-rabbit IgG; ICC, IHC)

Wakisaka, N., et al. [An Adenosine Receptor for Olfaction in Fish](#). Curr Biol 27(10), 1437-1447 e1434, (2017), DOI: 10.1016/j.cub.2017.04.014 (CF633 goat anti-mouse)

Wang, H., et al. [Cellular nanovesicles with bioorthogonal targeting enhance photodynamic/photothermal therapy in psoriasis](#). Acta Biomater 134, 674-685, (2021), DOI: 10.1016/j.actbio.2021.07.068 (CF633 succinimidyl ester)

Wang, W., et al. [A light- and calcium-gated transcription factor for imaging and manipulating activated neurons](#). Nat Biotechnol 35(9), 864-871, (2017), DOI: 10.1038/nbt.3909 (CF633 streptavidin)

Witteveldt, J., et al. [The influence of viral RNA secondary structure on interactions with innate host cell defences](#). Nucleic Acids Res 42(5), 3314-3329, (2014), DOI: gkt1291 [pii] 10.1093/nar/gkt1291 (CF633 anti-biotin; ICC)

Xia, S. H., et al. [Cortical and Thalamic Interaction with Amygdala-to-Accumbens Synapses](#). J Neurosci 40(37), 7119-7132, (2020), DOI: 10.1523/JNEUROSCI.1121-20.2020 (CF633 hydrazide)

Xie, H., et al. [TLR4 is involved in the pathogenic effects observed in a murine model of antiphospholipid syndrome](#). Clin Immunol 160(2), 198-210, (2015), DOI: 10.1016/j.clim.2015.05.017 (CF633 succinimidyl ester)

Yang, X., et al. [Dihomo-gamma-linolenic acid inhibits growth of xenograft tumors in mice bearing human pancreatic cancer cells \(BxPC-3\) transfected with delta-5-desaturase shRNA](#). Redox Biol 20, 236-246, (2019), DOI: 10.1016/j.redox.2018.10.001 (CF633 goat anti-rabbit CF633 goat anti-mouse)

Yu, Y., et al. [Expression, Localization and Functional Activity of the Major Na\(+\)/H\(+\) Exchange Isoforms Expressed in the Intestinal Cell Line Caco-2BBe](#). Cell Physiol Biochem 52(5), 1017-1038, (2019), DOI: 10.33594/000000070 (CF633 wheat germ agglutinin)

Zambon, P., Palani, S., Jadhav, S. S., Gayathri, P., Balasubramanian, M. K. . [Genetic Suppression of Defective Profilin by Attenuated Myosin II Reveals a Potential Role for Myosin II in Actin Dynamics in vivo in fission yeast](#). Mol Biol Cell 31(19), 2107-2114, (2020), DOI: 10.1091/mbc.E20-04-0224 (CF633 phalloidin fission yeast)

Zanetti-Domingues, L. C., et al. [Hydrophobic fluorescent probes introduce artifacts into single molecule tracking experiments due to non-specific binding](#). PLoS One 8(9), e74200, (2013), DOI: 10.1371/journal.pone.0074200 PONE-D-13-17594 [pii] (CF488A, CF568, CF633 & CF640R maleimide, TIRF)

Zhao, Y., et al. [Nanoscale imaging of clinical specimens using pathology-optimized expansion microscopy](#). Nat Biotechnol 35(8), 757-764, (2017), DOI: 10.1038/nbt.3892 (CF633 goat anti-mouse)

Zolnerciks, J. K., et al. [The Q loops of the human multidrug resistance transporter ABCB1 are necessary to couple drug binding to the ATP catalytic cycle](#). FASEB J 28(10), 4335-4346, (2014), DOI: fj.13-245639 [pii] 10.1096/fj.13-245639 (CF633 Mix-n-Stain; ICC)

CF®640R

Adolfsson, K., et al. [Ingestion of gallium phosphide nanowires has no adverse effect on Drosophila tissue function](#). Nanotechnology 24(28), 285101, (2013), DOI: DOI:10.1088/0957-4484/24/28/285101 (CF640R wheat germ agglutinin; ICC)

Aggarwal, N., et al. [Secreted osteopontin from CD4\(+\) T cells limits acute graft-versus-host disease](#). Cell Rep 37, 110170, (2021), DOI: 10.1016/j.celrep.2021.110170 (CF640R)

Agrawal, A., et al. [Manipulation of coacervate droplets with an electric field](#). Proc Natl Acad Sci U S A 119, e2203483119, (2022), DOI: 10.1073/pnas.2203483119 (CF488A BSA, CF640R BSA, fluorescence microscopy)

Akin, E. J., et al. [Single-Molecule Imaging of Nav1.6 on the Surface of Hippocampal Neurons Reveals Somatic Nanoclusters](#). Biophys J 111(6), 1235-1247, (2016), DOI: 10.1016/j.bpj.2016.08.016 (CF640R in TIRF)

Ara, A., et al. [Mechanism of Enhanced HIV Restriction by Virion Coencapsidated Cytidine Deaminases APOBEC3F and APOBEC3G](#). J Virol 91(3), (2017), DOI: 10.1128/JVI.02230-16 (CF543 Rabbit anti-V5 & CF640R Rabbit anti-HA)

Bosch, P. J., et al. [Evaluation of fluorophores to label SNAP-tag fused proteins for multicolor single-molecule tracking microscopy in live cells](#). Biophys J 107(4), 803-814, (2014), DOI: S0006-3495(14)00686-9 [pii] 10.1016/j.bpj.2014.06.040 (CF633 succinimidyl ester, CF640R succinimidyl ester; TIRF)

Bazzurro, V., et al. [Involvement of GABA\(A\) receptors containing alpha\(6\) subtypes in antsecretory factor activity on rat cerebellar granule cells studied by two-photon uncaging](#). Eur J Neurosci 56, 4505-4513, (2022), DOI: 10.1111/ejn.15775 (CF640R biocytin, fluorescence microscopy, 2-photon)

Bzowska, M., et al. [Involvement of cell surface 90 kDa heat shock protein \(HSP90\) in pattern recognition by human monocyte-derived macrophages](#). J Leukoc Biol 102(3), 763-774, (2017), DOI: 10.1189/jlb.2MA0117-019R (CF640R anti-biotin)

Dasgupta, P., et al. [Rapid assessment of G protein signaling of four opioid receptors using a real-time fluorescence-based membrane potential assay](#). Eur J Pharmacol 890, 173640, (2021), DOI: 10.1016/j.ejphar.2020.173640 (CF488A Anti-6X His CF640R Anti-FLAG)

De Pace, R., et al. [Synaptic Vesicle Precursors and Lysosomes Are Transported by Different Mechanisms in the Axon of Mammalian Neurons](#). Cell Rep 31(11), 107775, (2020), DOI: 10.1016/j.celrep.2020.107775 (CF640R Mix-n-Stain)

Engel, U. [Four-Channel Super-Resolution Imaging by 3-D Structured Illumination](#). Methods Mol Biol 1663, 79-94, (2017), DOI: 10.1007/978-1-4939-7265-4_7 (CF640R SIM)

Farias, G. G., et al. [Feedback-Driven Mechanisms between Microtubules and the Endoplasmic Reticulum Instruct Neuronal Polarity](#). Neuron 102(1), 184-201 e188, (2019), DOI: 10.1016/j.neuron.2019.01.030 (CF640R Mix-n-Stain)

Farias, G. G., et al. [Imaging the Polarized Sorting of Proteins from the Golgi Complex in Live Neurons](#). Methods Mol Biol 1496, 13-30, (2016), DOI: 10.1007/978-1-4939-6463-5_2 (CF488A, CF555 & CF640R Mix-n-Stain IF)

Forero-Quintero, L. S., et al. [Live-cell imaging reveals the spatiotemporal organization of endogenous RNA polymerase II phosphorylation at a single gene](#). Nat Commun 12(1), 3158, (2021), DOI: 10.1038/s41467-021-23417-0 (CF640R succinimidyl ester)

Gennet, N., et al. [FolR1: a novel cell surface marker for isolating midbrain dopamine neural progenitors and nascent dopamine neurons](#). Sci Rep 6, 32488, (2016), DOI: 10.1038/srep32488 (CF640R Mix-n-Stain FACS)

Gilboa, T., et al. [Single-Molecule DNA Methylation Quantification Using Electro-optical Sensing in Solid-State Nanopores](#). ACS Nano 10(9), 8861-8870, (2016), DOI: 10.1021/acsnano.6b04748 (CF640R succinimidyl ester)

Giraldo-Vela, J. P., et al. [Single-cell detection of mRNA expression using nanofountain-probe electroporated molecular beacons](#). Small 11(20), 2386-2391, (2015), DOI: 10.1002/smll.201401137 (CF640R succinimidyl ester)

Goldsmith, J. R., et al. [TNFAIP8 controls murine intestinal stem cell homeostasis and regeneration by regulating microbiome-induced Akt signaling](#). Nat Commun 11, 2591, (2020), DOI: 10.1038/s41467-020-16379-2 (CF640R WGA, fluorescence microscopy)

Hadassa Y. Holzapfel, and Marc R. Birtwistle. [Creating Complex Fluorophore spectra on Antibodies Through Combinatorial Labeling](#). Transl Sci, 1-10, (2016). (CF405M, CF488A, CF568, CF640R, FRET, MnS)

Herrera Estrada, L., et al. [Bioengineering Bacterially Derived Immunomodulants: A Therapeutic Approach to Inflammatory Bowel Disease](#). ACS Nano 11(10), 9650-9662, (2017), DOI: 10.1021/acsnano.7b03239 (CF640R transferrin, CF640R BSA and CF640R dextran 10,000 Mt)

Jeffet, J., et al. [Multi-modal Single-molecule Imaging with continuously Controlled Spectral-resolution \(CoCoS\) Microscopy](#). bioRxiv, (2020), DOI: 10.1101/2020.10.13.330910 (CF640R BCN)

Karwelat, D., et al. [A rodent thyroid-liver chip to capture thyroid toxicity on organ function level](#). ALTEX 40, 83-102, (2023), DOI: 10.14573/altex.2108262 (CF640R)

Keren-Kaplan, T., et al. [RUFY3 and RUFY4 are ARL8 effectors that promote coupling of endolysosomes to dynein-dynactin](#). Nat Commun 13, 1506, (2022), DOI: 10.1038/s41467-022-28952-y (CF640R, Mix-n-Stain)

Kolahi, K. S., et al. [Cytotrophoblast, Not Syncytiotrophoblast, Dominates Glycolysis and Oxidative Phosphorylation in Human Term Placenta](#). Sci Rep 7, 42941, (2017), DOI: 10.1038/srep42941 (CF640R wheat germ agglutinin)

Lee, H. J., et al. [ATP synthase inhibitory factor 1 \(IF1\), a novel myokine, regulates glucose metabolism by AMPK and Akt dual pathways](#). FASEB J 33(12), 14825-14840, (2019), DOI: 10.1096/fj.201901440RR (CF640R wheat germ agglutinin)

Loh, L. N., et al. [Dissecting Bacterial Cell Wall Entry and Signaling in Eukaryotic Cells: an Actin-Dependent Pathway Parallels Platelet-Activating Factor Receptor-Mediated Endocytosis](#). MBio 8(1), (2017), DOI: mBio.02030-16 [pii] 10.1128/mBio.02030-16 (CF405M wheat germ agglutinin, CF488A succinimidyl ester, CF640R succinimidyl ester; CF640R SIM)

Martin-Fernandez, M. L., et al. [A 'pocket guide' to total internal reflection fluorescence](#). J Microsc 252(1), 16-22, (2013), DOI: 10.1111/jmi.12070 (CF640R; TIRF)

Matsuzaki, T., et al. [Generation and characterization of a mouse line carrying Reck-CreERT2 knock-in allele](#). Genesis 56(4), e23099, (2018), DOI: 10.1002/dvg.23099 (CF640R goat anti-mouse)

Mattiazzi Usaj, M., et al. [Systematic genetics and single-cell imaging reveal widespread morphological pleiotropy and cell-to-cell variability](#). Mol Syst Biol 16(2), e9243, (2020), DOI: 10.15252/msb.20199243 (CF640R wheat germ agglutinin)

Motozono, C., et al. [An emerging SARS-CoV-2 mutant evading cellular immunity and increasing viral infectivity](#). bioRxiv, (2021), DOI: 10.1101/2021.04.02.438288 (CF640R succinimidyl ester)

Needham, S. R., et al. [EGFR oligomerization organizes kinase-active dimers into competent signalling platforms](#). Nat Commun 7, 13307, (2016), DOI: ncomms13307 [pii] 10.1038/ncomms13307 (CF640R FLImP)

Needham, S. R., et al. [Determining the geometry of oligomers of the human epidermal growth factor family on cells with <10 nm resolution](#). Biochem Soc Trans 43(3), 309-314, (2015), DOI: BST20140318 [pii] 10.1042/BST20140318 (CF640R FLImP)

Orosz, L., et al. [IL-17A and IL-17F induce autophagy in RAW 264.7 macrophages](#). Biomed Pharmacother 77, 129-134, (2016), DOI: 10.1016/j.biopha.2015.12.020 (CF640R goat anti-rabbit)

Parreno, J., et al. [Microstructural determinants of lens stiffness in rat versus guinea pig lenses](#). (2021), DOI: 10.1101/2021.02.15.431302 (CF640R wheat germ agglutinin, Hoechst)

Sainski, A. M., et al. [Casp8p41 generated by HIV protease kills CD4 T cells through direct Bak activation](#). J Cell Biol 207(1), 159, (2014), DOI: jcb.20140505109252014c [pii] 10.1083/jcb.20140505109252014c (CF640R Mix-n-Stain; FC)

Salazar-Cavazos, E., et al. [Multisite EGFR phosphorylation is regulated by adaptor protein abundances and dimer lifetimes](#). Mol Biol Cell 31(7), 695-708, (2020), DOI: 10.1091/mbc.E19-09-0548 (CF555 CF640R SIM)

Samec, N., Jovcevska, I., Stojan, J., Zottel, A., Liovic, M., Myers, M. P., Muyldermans, S., Šribar, J., Kržaj, I., Komel, R. . [Glioblastoma-specific anti-TUFM nanobody for in-vitro immunoimaging and cancer stem cell targeting](#). Oncotarget 9, 17282-17299, (2018), DOI: https://doi.org/10.18632/oncotarget.24629 (CF640R goat anti-mouse Fab)

Sanmarco, L. M., et al. [Gut-licensed IFNgamma\(+\) NK cells drive LAMP1\(+\)TRAIL\(+\) anti-inflammatory astrocytes](#). Nature 590(7846), 473-479, (2021), DOI: 10.1038/s41586-020-03116-4 (CF640R wheat germ agglutinin)

Sawaguchi, S., et al. [O-GlcNAc on NOTCH1 EGF repeats regulates ligand-induced Notch signaling and vascular development in mammals](#). eLife 6, (2017), DOI: 10.7554/eLife.24419 (CF488A donkey anti-goat, CF640R donkey anti-rat and CF488A streptavidin)

Shapira, S., et al. [Selective eradication of cancer cells by delivery of adenovirus-based toxins](#). Oncotarget 8(24), 28581-38591, (2017), DOI: 10.18632/oncotarget.16934 (CF640R Annexin RedDot2)

Singh, A. K., et al. [Structure and Sialyllactose Binding of the Carboxy-Terminal Head Domain of the Fibre from a Siadenovirus, Turkey Adenovirus 3](#). PLoS One 10(9), e0139339, (2015), DOI: 10.1371/journal.pone.0139339 (CF640R anti-6XHis)

Song, Y., et al. [Real-time imaging of single engineered RNA transcripts in living cells using ratiometric bimolecular beacons](#). J Vis Exp 6(90), e51544, (2014), DOI: 10.3791/51544 (CF640R succinimidyl ester molecular beacons RNA; fluorescence microscopy)

Stasevich, T. J., et al. [Quantifying histone and RNA polymerase II post-translational modification dynamics in mother and daughter cells](#). Methods 70(2-3), 77, (2014), DOI: S1046-2023(14)00260-6 [pii] 10.1016/jymeth.2014.08.002 (CF640R succinimidyl ester; live cell imaging)

Tokunaga, M., et al. [Hereditary spastic paraparesis protein spartin is an FK506-binding protein identified by mRNA display](#). Chem Biol 20(7), 935-942, (2013), DOI: S1074-5521(13)00236-6 [pii] 10.1016/j.chembiol.2013.05.011 (CF640R goat anti-mouse; ICC)

Uesaka, T., et al. [Neuronal Differentiation in Schwann Cell Lineage Underlies Postnatal Neurogenesis in the Enteric Nervous System](#). J Neurosci 35(27), 9879-9888, (2015), DOI: 10.1523/JNEUROSCI.1239-15.2015 (CF568 & CF640R secondaries)

Utratna, M., et al. [Rapid screening for specific glycosylation and pathogen interactions on a 78 species avian egg white glycoprotein microarray](#). Sci Rep 7(1), 6477, (2017), DOI: 10.1038/s41598-017-06797-6 (CF640R anti-6X His)

Valent, A. M., et al. [Hyperglycemia and gestational diabetes suppress placental glycolysis and mitochondrial function and alter lipid processing](#). FASEB J 35(3), e21423, (2021), DOI: 10.1096/fj.202000326RR (CF640R wheat germ agglutinin)

Wang, X., et al. [Disrupted Coordination of Hypoglossal Motor Control in a Mouse Model of Pediatric Dysphagia in DiGeorge/22q11.2 Deletion Syndrome](#). eNeuro 7(5), (2020), DOI: 10.1523/ENEURO.0520-19.2020 (CF640R biocytin)

Webb, S. E., et al. [Nanometric molecular separation measurements by single molecule photobleaching](#). Methods 88, 76-80, (2015), DOI: 10.1016/j.ymeth.2015.05.009 (CF640R succinimidyl ester)

Wu, M., et al. [Unbalanced Regulation of alpha7 nAChRs by Ly6h and NACHO Contributes to Neurotoxicity in Alzheimer's Disease](#). J Neurosci 41(41), 8461-8474, (2021), DOI: 10.1523/JNEUROSCI.0494-21.2021 (CF640R alpha-bungarotoxin)

Xu, Z., et al. [Teleost skin, an ancient mucosal surface that elicits gut-like immune responses](#). Proc Natl Acad Sci U S A 110(32), 13097-13102, (2013), DOI: 10.1073/pnas.1304319110 [pii] 10.1073/pnas.1304319110 (CF488A Mix-n-Stain, CF555, CF640R; IHC)

Yamada, S., et al. [Swelling-activated CIC-3 activity regulates prostaglandin E2 release in human OUMS-27 chondrocytes](#). Biochem Biophys Res Commun 537, 29-35, (2021), DOI: 10.1016/j.bbrc.2020.12.068 (CF640R wheat germ agglutinin)

Yu, Y., et al. [Single-phagosome imaging reveals that homotypic fusion impairs phagosome degradative function](#). Biophys J 121, 459-469, (2022), DOI: 10.1016/j.bpj.2021.12.032 (CF488A CTB, CF640R amine, TIRF)

Zahradnik, J., et al. [SARS-CoV-2 variant prediction and antiviral drug design are enabled by RBD in vitro evolution](#). Nat Microbiol 6(9), 1188-1198, (2021), DOI: 10.1038/s41564-021-00954-4 (CF640R succinimidyl ester)

Zanetti-Domingues, L. C., et al. [Determining the geometry of oligomers of the human epidermal growth factor family on cells with 7 nm resolution](#). Prog Biophys Mol Biol 118(3), 139-152, (2015), DOI: S0079-6107(15)00047-4 [pii] 10.1016/j.pbiomolbio.2015.04.002 (CF640R FLImP)

Zanetti-Domingues, L. C., et al. [Hydrophobic fluorescent probes introduce artifacts into single molecule tracking experiments due to non-specific binding](#). PLoS One 8(9), e74200, (2013), DOI: 10.1371/journal.pone.0074200 PONE-D-13-17594 [pii] (CF488A, CF568, CF633 & CF640R maleimide)

Zhang, R., et al. [Structural insight into TPX2-stimulated microtubule assembly](#). eLife 6, (2017), DOI: 10.7554/eLife.30959

Zhang, X., et al. [Quantitative assessment of ratiometric bimolecular beacons as a tool for imaging single engineered RNA transcripts and measuring gene expression in living cells](#). Nucleic Acids Res 41(15), e152, (2013), DOI: DOI:10.1093/nar/gkt561 (CF640R succinimidyl ester; RNA probes for use in living cells (Ratiometric Bimolecular Beacon (RBMB); ICC)

Zhou, H. and Liu, G. [Regulation of density of functional presynaptic terminals by local energy supply](#). Mol Brain 8, 42, (2015), DOI: 10.1186/s13041-015-0132-z (CF488A, CF555 & CF640R secondaries, MitoView 633)

Zurla, C., et al. [Characterizing mRNA interactions with RNA granules during translation initiation inhibition](#). PLoS One 6(5), e19727, (2011), DOI: DOI:10.1371/journal.pone.0019727 (CF640R tetravalent RNA imaging probe; FISH)

CF®647

Ahmed, S. M., et al. [KIF14 negatively regulates Rap1a-Radil signaling during breast cancer progression](#). J Cell Biol 199(6), 951-967, (2012), DOI: DOI:10.1083/jcb.201206051 (CF647 phalloidin; ICC)

Amann, R., et al. [A new rabies vaccine based on a recombinant Orf virus \(Parapoxvirus\) expressing the rabies virus glycoprotein](#). J Virol 87(3), 1618, (2012), DOI: DOI:10.1128/JVI.02470-12 (CF647 phalloidin; ICC)

Barbirou, M., et al. [Single Circulating-Tumor-Cell-Targeted Sequencing to Identify Somatic Variants in Liquid Biopsies in Non-Small-Cell Lung Cancer Patients](#). Current Issues in Molecular Biology 44(2), 750-763, (2022), DOI: 10.3390/cimb44020052 (CF405 succinimidyl ester & CF647 succinimidyl ester)

Belicky, S., et al. [Label-free chronopotentiometric glycoprofiling of prostate specific antigen using sialic acid recognizing lectins](#). Bioelectrochemistry 117, 89-94, (2017), DOI: 10.1016/j.bioelechem.2017.06.005 (CF647 streptavidin)

Bertok, T., et al. [Analysis of serum glycome by lectin microarrays for prostate cancer patients - a search for aberrant glycoforms](#). Glycoconj J 37(6), 703-711, (2020), DOI: 10.1007/s10719-020-09958-4 (CF647 streptavidin)

Cai, L., et al. [Optimal Poly\(L-lysine\) Grafting Density in Hydrogels for Promoting Neural Progenitor Cell Functions](#). Biomacromolecules 13(5), 1663, (2012), DOI: DOI:10.1021/bm300381d (CF647 goat anti-rabbit IgG; ICC)

Cantuti-Castelvetri, L., et al. [Neuropilin-1 facilitates SARS-CoV-2 cell entry and infectivity](#). Science 370(6518), 856-860, (2020), DOI: 10.1126/science.abd2985 (CF488A, CF555, CF647)

Castellani, S., et al. [Nanoparticle delivery of grape seed-derived proanthocyanidins to airway epithelial cells dampens oxidative stress and inflammation](#). J Transl Med 16(1), 140, (2018), DOI: 10.1186/s12967-018-1509-4 (CF647 Annexin V)

Coustets, M., et al. [Development of a near infrared protein nanoprobe targeting Thomsen-Friedenreich antigen for intraoperative detection of submillimeter nodules in an ovarian peritoneal carcinomatosis mouse model](#). Biomaterials 241, 119908, (2020), DOI: 10.1016/j.biomaterials.2020.119908 (CF647 protein labeling kit)

El Kojok, H., et al. [Fluorescence polarization assay to detect the presence of traces of ciprofloxacin](#). Sci Rep 10(1), 4550, (2020), DOI: 10.1038/s41598-020-61395-3 (CF647 succinimidyl ester toxin-receptor complex labeling)

Engdahl, A. K., et al. (2023). [Switching of Organic Fluorophores by Glycerol-sulfite Interactions for Single-Molecule Super-Resolution Microscopy](#). ChemPhotoChem 7, (2023), DOI: 10.1002/cptc.202200305 (CF647, CF568, Super-Resolution, dSTORM)

Estrada, A. L., et al. [Extracellular vesicle secretion is tissue-dependent ex vivo and skeletal muscle myofiber extracellular vesicles reach the circulation in vivo](#). Am J Physiol Cell Physiol 322, C246-C259, (2022), DOI: 10.1152/ajpcell.00580.2020 (CF647 CD63, extracellular vesicles, imaging immunocapture)

Futatsuki, T., et al. [Involvement of orexin neurons in fasting- and central adenosine-induced hypothermia](#). Sci Rep 8(1), 2717, (2018), DOI: 10.1038/s41598-018-21252-w (CF568 donkey anti-goat and CF647 donkey anti-goat)

Gabr, M. T. and Gambhir, S. S. [Discovery and Optimization of Small-Molecule Ligands for V-Domain Ig Suppressor of T-Cell Activation \(VISTA\)](#). J Am Chem Soc 142(38), 16194-16198, (2020), DOI: 10.1021/jacs.0c07276 (CF647 unknown)

Gomes de Castro, M. A., et al. [Differential organization of tonic and chronic B cell antigen receptors in the plasma membrane](#). Nat Commun 10(1), 820, (2019), DOI: 10.1038/s41467-019-08677-1 (CF647 maleimide)

Gong, Y.-N., et al. [Biological events and molecular signaling following MLKL activation during necroptosis](#). Cell Cycle, 1-13, (2017), DOI: 10.1080/15384101.2017.1371889 (CF568, CF647 STORM)

Gregorieff, A., et al. [Yap-dependent reprogramming of Lgr5\(+\) stem cells drives intestinal regeneration and cancer](#). Nature 526(7575), 715-718, (2015), DOI: 10.1038/nature15382 (CF555 donkey anti-mouse & CF647 donkey anti-rabbit)

Hess, D., et al. [High-throughput, quantitative enzyme kinetic analysis in microdroplets using stroboscopic epifluorescence imaging](#). Anal Chem 87(9), 4965-4972, (2015), DOI: 10.1021/acs.analchem.5b00766 (CF647 hydrazide-fluidics tracer)

Hirano, K. and Namihira, M. [LSD1 Mediates Neuronal Differentiation of Human Fetal Neural Stem Cells by Controlling the Expression of a Novel Target Gene, HEYL](#). Stem Cells 34(7), 1872-1882, (2016), DOI: 10.1002/stem.2362 (CF488A, CF568, CF647 donkey secondaries)

Homann, S., et al. [A novel rapid and reproducible flow cytometric method for optimization of transfection efficiency in cells](#). PLoS One 12(9), e0182941, (2017), DOI: 10.1371/journal.pone.0182941 (CF647 Mix-n-Stain)

Huh, D., et al. [Microfabrication of human organs-on-chips](#). Nat Protoc 8(11), 2135-2157, (2013), DOI: nprot.2013.137 [pii] 10.1038/nprot.2013.137 (CF647 phalloidin; ICC)

Inutsuka, A., et al. [Concurrent and robust regulation of feeding behaviors and metabolism by orexin neurons](#). Neuropharmacology 85, 451-460, (2014), DOI: S0028-3908(14)00237-8 [pii] 10.1016/j.neuropharm.2014.06.015 (CF488 CF594 CF647; IHC)

Inutsuka, A., et al. [The integrative role of orexin/hypocretin neurons in nociceptive perception and analgesic regulation](#). Sci Rep 6, 29480, (2016), DOI: 10.1038/srep29480 (CF488A, CF594 & CF647 secondaries)

Jain, D., et al. [Prognostic significance of NT5E/CD73 in neuroblastoma and its function in CSC stemness maintenance](#). Cell Biol Toxicol, (2021), DOI: 10.1007/s10565-021-09658-1 (CF647)

Jiang, L., et al. [A bacterial extracellular vesicle-based intranasal vaccine against SARS-CoV-2 protects against disease and elicits neutralizing antibodies to wild-type and Delta variants](#). bioRxiv, (2022), DOI: 10.1101/2021.06.28.450181 (CF488A SE, CF555 SE, CF647 SE, Single-particle interferometric reflectance imaging sensing)

Kanzaki, H., et al. [A-Disintegrin and Metalloproteinase \(ADAM\) 17 Enzymatically Degrades Interferon-gamma](#). Sci Rep 6, 32259, (2016), DOI: 10.1038/srep32259 (CF647 succinimidyl ester)

Kim, H. J. and Ingber, D. E. [Gut-on-a-Chip microenvironment induces human intestinal cells to undergo villus differentiation](#). Integr Biol (Camb) 5(9), 1130, (2013), DOI: DOI:10.1039/c3ib40126j (CF647 phalloidin; ICC)

Kong, D., et al. [Prolonged mitosis results in structurally aberrant and over-elongated centrioles](#). J Cell Biol 219(6), (2020), DOI: 10.1083/jcb.201910019 (CF647 goat anti-rabbit F(ab)2 STORM)

Koo, S. J., et al. [Vesicular Synaptobrevin/VAMP2 Levels Guarded by AP180 Control Efficient Neurotransmission](#). Neuron 88(2), 330-344, (2015), DOI: 10.1016/j.neuron.2015.08.034 (CF647 succinimidyl ester)

Krolewski, D. M., et al. [Quantitative validation of immunofluorescence and lectin staining using reduced CLARITY acrylamide formulations](#). Brain Struct Funct 223(2), 987-999, (2018), DOI: 10.1007/s00429-017-1583-z (CF594 CF647)

Kroll, A. V., et al. [Nanoparticulate Delivery of Cancer Cell Membrane Elicits Multiantigenic Antitumor Immunity](#). Adv Mater 29(47), (2017), DOI: 10.1002/adma.201703969 (CF647 succinimidyl ester)

Lee, K. S., et al. [Functional Synaptic Architecture of Callosal Inputs in Mouse Primary Visual Cortex](#). Neuron 101(3), 421-428 e425, (2019), DOI: 10.1016/j.neuron.2018.12.005 (CF405M CF647 expansion microscopy)

Lehmann, M., et al. [Nanoscale coupling of endocytic pit growth and stability](#). Sci Adv 5(11), eaax5775, (2019), DOI: 10.1126/sciadv.aax5775 (CF647 goat anti-mouse CF680 goat anti-rabbit STORM)

Lehmann, M., et al. [Novel organic dyes for multicolor localization-based super-resolution microscopy](#). J Biophotonics 9(1-2), 161-170, (2016), DOI: 10.1002/jbio.201500119 (CF555, CF568, CF647, CF680, STORM)

Liebel, M., et al. [3D tracking of extracellular vesicles by holographic fluorescence imaging](#). Sci Adv 6(45), (2020), DOI: 10.1126/sciadv.abc2508 (CF647 amine)

Lin, Z. P., et al. [Macrophages Actively Transport Nanoparticles in Tumors After Extravasation](#). ACS Nano 16, 6080-6092, (2022), DOI: 10.1021/acsnano.1c11578 (CF555 GARt, CF647 DARb, 3D microscopy)

Linhoff, M. W., et al. [A high-resolution imaging approach to investigate chromatin architecture in complex tissues](#). Cell 163(1), 246-255, (2015), DOI: S0092-8674(15)01122-8 [pii] 10.1016/j.cell.2015.09.002 (CF488A GAMs2a CF488A goat anti-chicken CF647 goat anti-mouse)

Liu, T., et al. [Quantification of antibody avidities and accurate detection of SARS-CoV-2 antibodies in serum and saliva on plasmonic substrates](#). Nat Biomed Eng 4, 1188-1196, (2020), DOI: 10.1038/s41551-020-00642-4 (CF647, antibody avidity assay, streptavidin)

Marenda, M., et al. [Parameter-free molecular super-structures quantification in single-molecule localization microscopy](#). J Cell Biol 220(5), (2021), DOI: 10.1083/jcb.202010003 (CF647)

Mattsson, L., et al. [Modified carbon black as label in a colorimetric on-chip immunoassay for histamine](#). Sensors and Actuators B: Chemical 246, 1092-1099, (2017), DOI: 10.1016/j.snb.2016.11.141 (CF647 goat anti-rabbit)

Meitinger, F., et al. [TRIM37 prevents formation of condensate-organized ectopic spindle poles to ensure mitotic fidelity](#). J Cell Biol 220(7), (2021), DOI: 10.1083/jcb.202010180 (CF647 CF568)

Merched, A. J., et al. [Specific autoantigens in experimental autoimmunity-associated atherosclerosis](#). FASEB J 30(6), 2123-2134, (2016), DOI: 10.1096/fj.201500131 (CF647 goat anti-mouse)

Merklinger, E., et al. [No Evidence for Spontaneous Lipid Transfer at ER-PM Membrane Contact Sites](#). J Membr Biol 249(1-2), 41-56, (2016), DOI: 10.1007/s00232-015-9845-2 (CF647 rabbit anti-goat)

Monaghan, C. E., et al. [REST corepressors RCOR1 and RCOR2 and the repressor INSM1 regulate the proliferation-differentiation balance in the developing brain](#). Proc Natl Acad Sci U S A 114(3), E406-E415, (2017), DOI: 10.1073/pnas.1620230114 (CF647 goat anti-mouse & CF647 goat anti-mouse IgG1)

Murao, N., et al. [Np95/Uhrf1 regulates tumor suppressor gene expression of neural stem/precursor cells, contributing to neurogenesis in the adult mouse brain](#). Neurosci Res 143, 31-43, (2019), DOI: 10.1016/j.neures.2018.05.007 (CF Donkey secondaries CF647 goat anti-chicken)

Nakamura, S., et al. [Dual role for DOCK7 in tangential migration of interneuron precursors in the postnatal forebrain](#). J Cell Biol 216(12), 4313-4330, (2017), DOI: 10.1083/jcb.201704157 (CF647 donkey anti-goat)

Nakashima, H., et al. [Canonical TGF-beta Signaling Negatively Regulates Neuronal Morphogenesis through TGIF/Smad Complex-Mediated CRMP2 Suppression](#). J Neurosci 38(20), 4791-4810, (2018), DOI: 10.1523/JNEUROSCI.2423-17.2018 (CF488A, CF555, CF647 secondaries)

Nakashima, H., et al. [MeCP2 controls neural stem cell fate specification through miR-199a-mediated inhibition of BMP-Smad signaling](#). Cell Rep 35(7), 109124, (2021), DOI: 10.1016/j.celrep.2021.109124 (CF488A donkey anti-rabbit CF488A donkey anti-goat CF555 donkey anti-mouse CF555 donkey anti-rabbit CF555 donkey anti-guinea pig CF647 donkey anti-mouse and CF647 donkey anti-rat)

Noguchi, H., et al. [Expression of DNMT1 in neural stem/precursor cells is critical for survival of newly generated neurons in the adult hippocampus](#). Neurosci Res 95, 1-11, (2015), DOI: S0168-0102(15)00033-4 [pii] 10.1016/j.neures.2015.01.014 (CF488A donkey anti-mouse, CF543 donkey anti-rabbit, CF568 donkey anti-rat, CF647 donkey anti-goat, CF647 donkey anti-rabbit)

Olivier, N., et al. [Simple buffers for 3D STORM microscopy](#). Biomed Opt Express 4(6), 885-899, (2013), DOI: 10.1364/BOE.4.000885 187153 [pii] (CF647; 3D STORM)

Park, K. S., et al. [A crucial requirement for Hedgehog signaling in small cell lung cancer](#). Nat Med 17(11), 1504-1508, (2011), DOI: DOI:10.1038/nm.2473 (Mix-n-Stain CF488A, Mix-n-Stain CF568, and Mix-n-Stain CF647; ICC)

Paz, H., et al. [Treatment of B-cell precursor acute lymphoblastic leukemia with the Galectin-1 inhibitor PTX008](#). J Exp Clin Cancer Res 37(1), 67, (2018), DOI: 10.1186/s13046-018-0721-7 (CF647 Mix-n-Stain)

Pennacchio, A., et al. [A novel fluorescence polarization assay for determination of penicillin G in milk](#). Food Chem 190, 381-385, (2016), DOI: S0308-8146(15)00871-7 [pii] 10.1016/j.foodchem.2015.05.127 (CF647 succinimidyl ester)

Raynes, J. K., et al. [Coaggregation of kappa-Casein and beta-Lactoglobulin Produces Morphologically Distinct Amyloid Fibrils](#). Small 13(14), (2017), DOI: 10.1002/smll.201603591 (CF488A maleimide CF647 maleimide)

Robajac, D., et al. [Sensitive glycoprofiling of insulin-like growth factor receptors isolated from colon tissue of patients with colorectal carcinoma using lectin-based protein microarray](#). Int J Biol Macromol 144, 932-937, (2020), DOI: 10.1016/j.ijbiomac.2019.09.170 (CF555 and CF647 streptavidin)

Robajac, D., et al. [Screening for the best detergent for the isolation of placental membrane proteins](#). Int J Biol Macromol 102, 431-437, (2017), DOI: 10.1016/j.ijbiomac.2017.04.046 (CF647 streptavidin)

Rohde, J., et al. [New Orf virus \(Parapoxvirus\) recombinant expressing H5 hemagglutinin protects mice against H5N1 and H1N1 influenza A virus](#). PLoS One 8(12), e83802, (2013), DOI: 10.1371/journal.pone.0083802 PONE-D-13-40047 [pii] (CF647 phalloidin; ICC)

Rohm, M., et al. [Data of rational process optimization for the production of a full IgG and its Fab fragment from hybridoma cells](#). Data Brief 8, 426-435, (2016), DOI: 10.1016/j.dib.2016.05.067 (CF647 Mix-n-Stain)

Sabath, D. E., et al. [Clinical Validation of a Circulating Tumor Cell Assay Using Density Centrifugation and Automated Immunofluorescence Microscopy](#). Am J Clin Pathol 158, 270-276, (2022), DOI: 10.1093/ajcp/aqac040 (CF647 Fluorescence Microscopy)

Samson, A. L., et al. [Physicochemical properties that control protein aggregation also determine whether a protein is retained or released from necrotic cells](#). Open Biol 6(11), (2016), DOI: 10.1098/rsob.160098 (CF647 donkey anti-rabbit)

Santacatterina, F., et al. [Pyruvate kinase M2 and the mitochondrial ATPase Inhibitory Factor 1 provide novel biomarkers of dermatomyositis: a metabolic link to oncogenesis](#). J Transl Med 15(1), 29, (2017), DOI: 10.1186/s12967-017-1136-5 (CF647 goat anti-mouse)

Schuhenn, J., et al. [Differential interferon-alpha subtype induced immune signatures are associated with suppression of SARS-CoV-2 infection](#). Proc Natl Acad Sci U S A 119(8), (2022), DOI: 10.1073/pnas.2111600119 (CF647 phalloidin)

Skindersoe, M. E., et al. [A novel and rapid apoptosis assay based on thiol redox status](#). Cytometry A 81(5), 430, (2012), DOI: DOI:10.1002/cyto.a.22032 (CF647 Annexin V; ICC, FC)

Smith, P. L., et al. [Neonatal peripheral immune challenge activates microglia and inhibits neurogenesis in the developing murine hippocampus](#). Dev Neurosci 36(2), 119-131, (2014), DOI: 000359950 [pii] 10.1159/000359950 (CF555 donkey anti-goat , CF488A donkey anti-chicken & CF647 donkey anti-mouse; IHC)

Tai, L. H., et al. [Transient stimulation of distinct subpopulations of striatal neurons mimics changes in action value](#). Nat Neurosci 15(9), 1281-1289, (2012), DOI: DOI:10.1038/nn.3188 (CF555 donkey anti-mouse IgG, CF647 donkey anti-goat IgG; IHC)

Takenaka, N., et al. [A critical role of the small GTPase Rac1 in Akt2-mediated GLUT4 translocation in mouse skeletal muscle](#). FEBS J 281(5), 1493-1504, (2014), DOI: 10.1111/febs.12719 (CF488A, CF543, CF647 goat anti-mouse, goat anti-rabbit, goat anti-rat; ICC)

Takenaka, N., et al. [Rac1 Activation Caused by Membrane Translocation of a Guanine Nucleotide Exchange Factor in Akt2-Mediated Insulin Signaling in Mouse Skeletal Muscle](#). PLoS One 11(5), e0155292, (2016), DOI: 10.1371/journal.pone.0155292 (CF405, CF488A, CF543 & CF647 IHF)

Takenaka, N., et al. [Role of the guanine nucleotide exchange factor in Akt2-mediated plasma membrane translocation of GLUT4 in insulin-stimulated skeletal muscle](#). Cell Signal 26(11), 2460-2469, (2014), DOI: S0898-6568(14)00221-6 [pii] 10.1016/j.cellsig.2014.07.002 (CF350, CF543 & CF647; ICC)

Trofimenko, E., et al. [The endocytic pathway taken by cationic substances requires Rab14 but not Rab5 and Rab7](#). Cell Rep 37, 109945, (2021), DOI: 10.1016/j.celrep.2021.109945 (CF405M, CF488A, CF594, CF647, Mix-n-Stain, fluorescence microscopy)

Turkowyd, B., et al. [From single molecules to life: microscopy at the nanoscale](#). Anal Bioanal Chem 408(25), 6885-6911, (2016), DOI: 10.1007/s00216-016-9781-8, 10.1007/s00216-016-9781-8 [pii] (CF568, CF647, CF660C, CF680 super res review)

Varriale, A., et al. [A Fluorescence Polarization Assay To Detect Steroid Hormone Traces in Milk](#). J Agric Food Chem 63(41), 9159-9164, (2015), DOI: 10.1021/acs.jafc.5b03689 (CF647 succinimidyl ester)

Varriale, A., et al. [On the possibility of ephedrine detection: time-resolved fluorescence resonance energy transfer \(FRET\)-based approach](#). Anal Bioanal Chem 408, 6329-6336, (2016), DOI: 10.1007/s00216-016-9738-y (CF568, CF647, FRET)

Virant, D., et al. [Unraveling the kinetochore nanostructure in *Schizosaccharomyces pombe* using multi-color single-molecule localization microscopy](#). (2022), DOI: 10.1101/2021.12.01.469981 (CF647, SMLM)

Wang, H. Y., et al. [Synthesis and single-molecule imaging reveal stereospecific enhancement of binding kinetics by the antitumour eEF1A antagonist SR-A3](#). Nat Chem 14, 1443-1450, (2022), DOI: 10.1038/s41557-022-01039-3 (CF647 azide, flow cytometry)

Wei, Y., et al. [Synergistic Entry of Individual Nanoparticles into Mammalian Cells Driven by Free Energy Decline and Regulated by Their Sizes](#). ACS Nano 16, 5885-5897, (2022), DOI: 10.1021/acsnano.1c11068 (CF488A, CF555, CF647 SE, nanoparticle labeling, fluorescence microscopy)

Xu, N., et al. [Neonatal Pharmacokinetics and Biodistribution of Polymeric Nanoparticles and Effect of Surfactant](#). Pharmaceutics 15, (2023), DOI: 10.3390/pharmaceutics15041176 (CF647 SE, in vivo)

Yuan, B., et al. [Plasmon-enhanced fluorescence imaging with silicon-based silver chips for protein and nucleic acid assay](#). Anal Chim Acta 955, 98-107, (2017), DOI: 10.1016/j.aca.2016.11.059 (CF647 streptavidin)

CF®660C

Diekmann, R., et al. [Optimizing imaging speed and excitation intensity for single-molecule localization microscopy](#). Nat Methods 17(9), 909-912, (2020), DOI: 10.1038/s41592-020-0918-5 (CF660C, CF680)

Gu, L., et al. [Molecular-scale axial localization by repetitive optical selective exposure](#). Nat Methods 18(4), 369-373, (2021), DOI: 10.1038/s41592-021-01099-2 (STORM CF660C donkey anti-mouse)

Gwosch, K. C., et al. [MINFLUX nanoscopy delivers 3D multicolor nanometer resolution in cells](#). Nat Methods 17(2), 217-224, (2020), DOI: 10.1038/s41592-019-0688-0 (CF660C succinimidyl ester CF680 wheat germ agglutinin)

Huang, T., et al. [Simultaneous Multicolor Single-Molecule Tracking with Single-Laser Excitation via Spectral Imaging](#). Biophys J 114(2), 301-310, (2018), DOI: S0006-3495(17)31242-0 [pii] 10.1016/j.bpj.2017.11.013 (CF633, CF660C, CF680, CF680R STORM)

Li, Y., et al. [Global fitting for high-accuracy multi-channel single-molecule localization](#). Nat Commun 13, 3133, (2022), DOI: 10.1038/s41467-022-30719-4 (CF660C GARb, CF680 WGA, SMLM, fluorescence microscopy)

Nahidazar, L. and Harkes, R. [Multicolor Localization-Based Super Resolution Microscopy](#). Methods Mol Biol 2350, 69-76, (2021), DOI: 10.1007/978-1-0716-1593-5_5 (CF488A, CF680, CF660C, CF514 SMLM)

Pape, J. K., et al. [Multicolor 3D MINFLUX nanoscopy of mitochondrial MICOS proteins](#). Proc Natl Acad Sci U S A 117(34), 20607-20614, (2020), DOI: 10.1073/pnas.2009364117 (CF660C succinimidyl ester CF680 succinimidyl ester)

Turkowyd, B., et al. [From single molecules to life: microscopy at the nanoscale](#). Anal Bioanal Chem 408(25), 6885-6911, (2016), DOI: 10.1007/s00216-016-9781-8 10.1007/s00216-016-9781-8 [pii] (CF568, CF647, CF660C, CF680 super res review)

Wu, W., et al. [Tetra-color superresolution microscopy based on excitation spectral demixing](#). Light Sci Appl 12, 9, (2023), DOI: 10.1038/s41377-022-01054-6 (CF660C, CF680, super-resolution, ExR STORM)

Wu, Y. L., et al. [Maximum-likelihood model fitting for quantitative analysis of SMLM data](#). Nat Methods 20, 139-148, (2023), DOI: 10.1038/s41592-022-01676-z (CF660C, CF680 secondaries, super-resolution, SMLM)

Zhang, Y., et al. [Multicolor super-resolution imaging using spectroscopic single-molecule localization microscopy with optimal spectral dispersion](#). Appl Opt 58(9), 2248-2255, (2019), DOI: 10.1364/AO.58.002248 (CF660C CF680 sSMLM)

Zhang, Y., et al. [Nanoscale subcellular architecture revealed by multicolor three-dimensional salvaged fluorescence imaging](#). Nat Methods 17(2), 225-231, (2020), DOI: 10.1038/s41592-019-0676-4 (CF660C & CF680 single labeled, CF680-wheat germ agglutinin)

Zhang, Y., et al. [Minimizing Molecular Misidentification in Imaging Low-Abundance Protein Interactions Using Spectroscopic Single-Molecule Localization Microscopy](#). Anal Chem 94, 13834-13841, (2022), DOI: 10.1021/acs.analchem.2c02417 (CF660C, CF680 secondaries, super-resolution, SMLM)

Zhang, Z., et al. [Ultrahigh-throughput single-molecule spectroscopy and spectrally resolved super-resolution microscopy](#). Nat Methods 12(10), 935-938, (2015), DOI: nmeth.3528 [pii] 10.1038/nmeth.3528 (CF660C, CF680 STORM)

CF®660R

Civitci, F., et al. [Fast and multiplexed superresolution imaging with DNA-PAINT-ERS](#). Nat Commun 11(1), 4339, (2020), DOI: 10.1038/s41467-020-18181-6 (CF660R SMLM)

Gorham, J. B., et al. [Visualization of microbe-dietary remnant interactions in digesta from pigs, by fluorescence in situ hybridization and staining methods; effects of a dietary arabinoxylan-rich wheat fraction](#). Food Hydrocolloids 52, 952-962, (2016), DOI: http://dx.doi.org/10.1016/j.foodhyd.2015.09.011 (660R, Mix-n-Stain)

Heidarsson, P. O., et al. [Disordered Proteins Enable Histone Chaperoning on the Nucleosome](#). (2020), DOI: 10.1101/2020.04.17.046243 (CF660R FRET)

MacDonald, A. B. [Lyme ELISA seronegative patients with borrelia antibody coated borrelia microbes in their blood](#). Med Clin Res 7, (2022), ISSN: 2577 - 8005 (CF405S, CF660R, Mix-n-Stain)

Seo, J., et al. [PICASSO: Ultra-multiplexed fluorescence imaging of biomolecules through single-round imaging and blind source unmixing](#). bioRxiv, (2021), DOI: 10.1101/2021.01.27.428247 (CF633, CF660R, CF680R)

Seo, J., et al. [PICASSO allows ultra-multiplexed fluorescence imaging of spatially overlapping proteins without reference spectra measurements](#). Nat Commun 13, 2475, (2022), DOI: 10.1038/s41467-022-30168-z (CF405S, CF405M, CF405L, CF488A, CF568, CF633, CF660R, CF680R)

Shirasaki, Y., et al. [Real-time single-cell imaging of protein secretion](#). Sci Rep 4, 4736, (2014), DOI: srep04736 [pii] 10.1038/srep04736 (CF660R; TIRF)

Schubert, J., et al. [Two-colour single-molecule photoinduced electron transfer fluorescence imaging microscopy of chaperone dynamics](#). Nat Commun 12, 6964, (2021), DOI: 10.1038/s41467-021-27286-5 (CF660R maleimide, PET)

CF®680

Avelar, G. M., et al. [A rhodopsin-guanylyl cyclase gene fusion functions in visual perception in a fungus](#). Curr Biol 24(11), 1234-1240, (2014), DOI: S0960-9822(14)00409-6 [pii] 10.1016/j.cub.2014.04.009 (CF680; WB; LI-COR Odyssey)

Bonnaud, E. M., et al. [Borna disease virus phosphoprotein modulates epigenetic signaling in neurons to control viral replication](#). J Virol 89(11), 5996-6008, (2015), DOI: 10.1128/JVI.00454-15 (CF680 & CF770 WB)

Cho, H., et al. [Study of osteoarthritis treatment with anti-inflammatory drugs: cyclooxygenase-2 inhibitor and steroids](#). Biomed Res Int 2015, 595273, (2015), DOI: 10.1155/2015/595273 (CF680SE in vivo)

Diekmann, R., et al. [Optimizing imaging speed and excitation intensity for single-molecule localization microscopy](#). Nat Methods 17(9), 909-912, (2020), DOI: 10.1038/s41592-020-0918-5 (CF660C, CF680)

Ducheix, S., et al. [Dietary oleic acid regulates hepatic lipogenesis through a liver X receptor-dependent signaling](#). PLoS One 12(7), e0181393, (2017), DOI: 10.1371/journal.pone.0181393 (CF680 & CF770 secondaries WB)

Frank, C. L., et al. [Regulation of chromatin accessibility and Zic binding at enhancers in the developing cerebellum](#). Nat Neurosci 18(5), 647-656, (2015), DOI: 10.1038/nn.3995 (CF680 goat anti-mouse & CF770 goat anti-rabbit WB)

Fruh, S. M., et al. [Molecular architecture of native fibronectin fibrils](#). Nat Commun 6, 7275, (2015), DOI: ncomms8275 [pii] 10.1038/ncomms8275 (CF680 STORM)

Gauthier, T., et al. [Deoxynivalenol impairs the immune functions of neutrophils](#). Mol Nutr Food Res 57(6), 1026, (2013), DOI: DOI:10.1002/mnfr.201200755 (CF680 goat anti-rabbit IgG, CF770 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Glebov, O. O., et al. [Nanoscale Structural Plasticity of the Active Zone Matrix Modulates Presynaptic Function](#). Cell Rep 18(11), 2715-2728, (2017), DOI: S2211-1247(17)30279-6 [pii] 10.1016/j.celrep.2017.02.064 (CF680 STORM)

Gon, S., et al. [Fit alphabeta T-cell receptor suppresses leukemogenesis of Pten-deficient thymocytes](#). Haematologica 103(6), 999-1007, (2018), DOI: 10.3324/haematol.2018.188359 (CF680 goat anti-mouse and CF770 goat anti-rabbit WB)

Gorur, A., et al. [COPII-coated membranes function as transport carriers of intracellular procollagen I](#). The Journal of Cell Biology 216(6), 1745-1759, (2017), DOI: 10.1083/jcb.201702135 (CF568, CF680 STORM)

Graber, T. E., et al. [UPF1 Governs Synaptic Plasticity through Association with a STAU2 RNA Granule](#). J Neurosci 37(38), 9116-9131, (2017), DOI: 10.1523/JNEUROSCI.0088-17.2017 (CF680 & CF770 secondaries WB)

Guo, X., et al. [Mitochondrial stress is relayed to the cytosol by an OMA1-DELE1-HRI pathway](#). Nature 579(7799), 427-432, (2020), DOI: 10.1038/s41586-020-2078-2 (CF568 donkey anti-rabbit CF680 donkey anti-rabbit STORM)

Gwosch, K. C., et al. [MINFLUX nanoscopy delivers 3D multicolor nanometer resolution in cells](#). Nat Methods 17(2), 217-224, (2020), DOI: 10.1038/s41592-019-0688-0 (CF660C succinimidyl ester CF680 wheat germ agglutinin)

Hapuarachchige, S., et al. [Bioorthogonal two-component drug delivery in HER2\(+\) breast cancer mouse models](#). Sci Rep 6, 24298, (2016), DOI: 10.1038/srep24298 (CF680 succinimidyl ester & CF750 succinimidyl ester, in vivo)

Hilbers, F. S., et al. [Functional Analysis of Missense Variants in the Putative Breast Cancer Susceptibility Gene XRCC2](#). Hum Mutat 37(9), 914-925, (2016), DOI: 10.1002/humu.23019 (CF680 goat anti-rabbit & CF770 goat anti-mouse WB)

Hoess, P., et al. [Dual-Color and 3D Super-Resolution Microscopy of Multi-protein Assemblies](#). Methods Mol Biol 1764, 237-251, (2018), DOI: 10.1007/978-1-4939-7759-8_14 (CF680 goat anti-rabbit 2-color STORM)

Huang, T., et al. [Simultaneous Multicolor Single-Molecule Tracking with Single-Laser Excitation via Spectral Imaging](#). Biophys J 114(2), 301-310, (2018), DOI: S0006-3495(17)31242-0 [pii] 10.1016/j.bpj.2017.11.013 (CF633, CF660C, CF680, CF680R STORM)

Jing, H., et al. [Combination treatment with hypofractionated radiotherapy plus IL-2/anti-IL-2 complexes and its theranostic evaluation](#). J Immunother Cancer 7(1), 55, (2019), DOI: 10.1186/s40425-019-0537-9 (CF680 VivoBrite)

Klevanski, M., et al. [Automated highly multiplexed super-resolution imaging of protein nano-architecture in cells and tissues](#). Nat Commun 11(1), 1552, (2020), DOI: 10.1038/s41467-020-15362-1 (CF680 multi-round STORM)

Kohler, C., et al. [Extracytoplasmic function \(ECF\) sigma factor sigmaF is involved in Caulobacter crescentus response to heavy metal stress](#). BMC Microbiol 12, 210, (2012), DOI: DOI:10.1186/1471-2180-12-210 (CF680 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Lee, S., et al. [Localization of near-infrared labeled antibodies to the central nervous system in experimental autoimmune encephalomyelitis](#). PLoS One 14(2), e0212357, (2019), DOI: 10.1371/journal.pone.0212357 (CF680 in vivo)

Lehmann, M., et al. [Nanoscale coupling of endocytic pit growth and stability](#). Sci Adv 5(11), eaax5775, (2019), DOI: 10.1126/sciadv.aax5775 (CF647 goat anti-mouse CF680 goat anti-rabbit STORM)

Lehmann, M., et al. [Novel organic dyes for multicolor localization-based super-resolution microscopy](#). J Biophotonics 9(1-2), 161-170, (2016), DOI: 10.1002/jbio.201500119 (CF555, CF568, CF647, CF680, STORM)

Li, C. G., et al. [Berberine augments ATP-induced inflammasome activation in macrophages by enhancing AMPK signaling](#). Oncotarget 8(1), 95-109, (2017), DOI: 10.18632/oncotarget.13921 (CF568 goat anti-rabbit)

Li, Y., et al. [Global fitting for high-accuracy multi-channel single-molecule localization](#). Nat Commun 13, 3133, (2022), DOI: 10.1038/s41467-022-30719-4 (CF660C GARb, CF680 WGA, SMLM, fluorescence microscopy)

Nahidazar, L. and Harkes, R. [Multicolor Localization-Based Super Resolution Microscopy](#). Methods Mol Biol 2350, 69-76, (2021), DOI: 10.1007/978-1-0716-1593-5_5 (CF488A, CF680, CF660C, CF514 SMLM)

Najimi, M., et al. [Human liver mesenchymal stem/progenitor cells inhibit hepatic stellate cell activation: in vitro and in vivo evaluation](#). Stem Cell Res Ther 8(1), 131, (2017), DOI: 10.1186/s13287-017-0575-5 (CF680 donkey anti-rabbit & CF790 donkey anti-mouse WB)

Newman, Z. L., et al. [Determinants of synapse diversity revealed by super-resolution quantal transmission and active zone imaging](#). Nat Commun 13, 229, (2022), DOI: 10.1038/s41467-021-27815-2 (CF680 SE STORM)

Oliveira, A. F. and Yasuda, R. [Neurofibromin is the major ras inactivator in dendritic spines](#). J Neurosci 34(3), 776-783, (2014), DOI: 34/3/776 [pii] 10.1523/JNEUROSCI.3096-13.2014 (CF680 goat anti-mouse, CF770 goat anti-rabbit; WB; LI-COR Odyssey)

Ottenheijm, C. A., et al. [Deleting exon 55 from the nebulin gene induces severe muscle weakness in a mouse model for nemaline myopathy](#). Brain 136(Pt 6), 1718-1731, (2013), DOI: awt113 [pii] 10.1093/brain/awt113 (CF680 goat anti-rabbit; WB; LI-COR Odyssey)

Pape, J. K., et al. [Multicolor 3D MINFLUX nanoscopy of mitochondrial MICOS proteins](#). Proc Natl Acad Sci U S A 117(34), 20607-20614, (2020), DOI: 10.1073/pnas.2009364117 (CF660C succinimidyl ester CF680 succinimidyl ester)

Platonova, E., et al. [A simple method for GFP- and RFP-based dual color single-molecule localization microscopy](#). ACS Chem Biol 10(6), 1411-1416, (2015), DOI: 10.1021/acscchembio.5b00046 (CF680 STORM)

Platonova, E., et al. [Single-molecule microscopy of molecules tagged with GFP or RFP derivatives in mammalian cells using nanobody binders](#). Methods 88, 89-97, (2015), DOI: http://dx.doi.org/10.1016/j.ymeth.2015.06.018 (CF680 STORM)

Rashid, F. M., et al. [Hi-NESS: a family of genetically encoded DNA labels based on a bacterial nucleoid-associated protein](#). Nucleic Acids Res 50, e10, (2022), DOI: 10.1093/nar/gkab993 (CF680 GARb, CF770 GAMs)

Rojas-Sánchez, L., et al. [Synthesis and intracellular tracing surface-functionalized calcium phosphate nanoparticles by super-resolution microscopy \(STORM\)](#). Materialia 12, (2020), DOI: 10.1016/j.mtla.2020.100773 (CF568 wheat germ agglutinin STORM)

Rother, M. B., et al. [The Human Thymus Is Enriched for Autoreactive B Cells](#). J Immunol 197(2), 441-448, (2016), DOI: 10.4049/jimmunol.1501992 (CF680 GAHu & CF770 goat anti-mouse WB)

Salvador-Gallego, R., et al. [Bax assembly into rings and arcs in apoptotic mitochondria is linked to membrane pores](#). EMBO J 35(4), 389-401, (2016), DOI: 10.15252/embj.201593384 (CF680 goat anti-mouse STORM)

Sarveswaran, K., et al. [Synthetic Capillaries to Control Microscopic Blood Flow](#). Sci Rep 6, 21885, (2016), DOI: 10.1038/srep21885 (CF680 succinimidyl ester IF)

Schimmel, J., et al. [Mutational signatures of non-homologous and polymerase theta-mediated end-joining in embryonic stem cells](#). EMBO J 36(24), 3634-3649, (2017), DOI: 10.15252/embj.201796948 (CF680 goat anti-rabbit & CF770 GARMs WB)

Shrestha, R. L., et al. [Aurora-B kinase pathway controls the lateral to end-on conversion of kinetochore-microtubule attachments in human cells](#). Nat Commun 8(1), 150, (2017), DOI: 10.1038/s41467-017-00209-z 10.1038/s41467-017-00209-z [pii] (CF680 DaHu STORM)

Schmerl, B., et al. [The synaptic scaffold protein MPP2 interacts with GABA_A receptors at the periphery of the postsynaptic density of glutamatergic synapses](#). PLoS Biol 20, e3001503, (2022), DOI: 10.1371/journal.pbio.3001503 (CF568 SE, CF680 SE, super-resolution, dSTORM,)

Sun, Y., et al. [Single tumor imaging with multiple antibodies targeting different antigens](#). Biotechniques 52(4), 1-3, (2012), DOI: DOI:10.2144/000113855 (XenoLight CF680, XenoLight CF770; in vivo imaging; PerkinElmer IVIS Lumina)

Titova, E. and Obenaus, A. [Immunological Response to Experimental Intracerebral Hemorrhage: Morphological Assessments](#). In: Chen, J., et al. (Eds.), Animal Models of Acute Neurological Injuries II. Humana Press, pp. 625-652, (2012), DOI: DOI:10.1007/978-1-61779-576-3_48 (CF680 and CF770 secondary antibody conjugates; IHC in floating brain sections; LI-COR Odyssey)

Titova, E. M., et al. [The late phase of post-stroke neurorepair in aged rats is reflected by MRI-based measures](#). Neuroscience, (2014), DOI: S0306-4522(14)00770-2 [pii] 10.1016/j.neuroscience.2014.09.028 (CF680 goat anti-mouse & CF770 goat anti-rat; IHC; LI-COR Odyssey)

Thiele, J. C., et al. [Isotropic three-dimensional dual-color super-resolution microscopy with metal-induced energy transfer](#). Science Advances 8, eabo2506, (2022), DOI: 10.1126/sciadv.abo2506 (CF680 SE, MIET-dSTORM, super-resolution microscopy)

Turkowyd, B., et al. [From single molecules to life: microscopy at the nanoscale](#). Anal Bioanal Chem 408(25), 6885-6911, (2016), DOI: 10.1007/s00216-016-9781-8, 10.1007/s00216-016-9781-8 [pii] (CF568, CF647, CF660C, CF680 super res review)

van den Heuvel, D., et al. [A CSB-PAF1C axis restores processive transcription elongation after DNA damage repair](#). Nat Commun 12(1), 1342, (2021), DOI: 10.1038/s41467-021-21520-w (CF680 goat anti-rabbit CF770 goat anti-mouse WB Odyssey)

van der Weegen, Y., et al. [The sequential and cooperative action of CSB, CSA, and UVSSA targets the TFIH complex to DNA damage-stalled RNA polymerase II](#). bioRxiv, (2019), DOI: 10.1101/707216 (CF680 donkey anti-goat CF680 goat anti-rabbit CF770 goat anti-mouse WB)

Venkataramani, V., et al. [SuReSim: simulating localization microscopy experiments from ground truth models](#). Nat Methods 13(4), 319-321, (2016), DOI: nmeth.3775 [pii] 10.1038/nmeth.3775

Winterflood, C. M., et al. [Dual-color 3D superresolution microscopy by combined spectral-demixing and biplane imaging](#). Biophys J 109(1), 3-6, (2015), DOI: S0006-3495(15)00538-X [pii] 10.1016/j.bpj.2015.05.026 (CF680 3D SMLM)

Wu, C. L., et al. [Interplay between Cell Migration and Neurite Outgrowth Determines SH2B1beta-Enhanced Neurite Regeneration of Differentiated PC12 Cells](#). PLoS One 7(4), e34999, (2012), DOI: DOI:10.1371/journal.pone.0034999 (CF680 goat anti-mouse IgG; WB; LI-COR Odyssey)

Wu, W., et al. [Tetra-color superresolution microscopy based on excitation spectral demixing](#). Light Sci Appl 12, 9, (2023), DOI: 10.1038/s41377-022-01054-6 (CF660C, CF680, super-resolution, ExR STORM)

Wu, Y. L., et al. [Maximum-likelihood model fitting for quantitative analysis of SMLM data](#). Nat Methods 20, 139-148, (2023), DOI: 10.1038/s41592-022-01676-z (CF660C, CF680 secondaries, super-resolution, SMLM)

Zhang, Y., et al. [Multicolor super-resolution imaging using spectroscopic single-molecule localization microscopy with optimal spectral dispersion](#). Appl Opt 58(9), 2248-2255, (2019), DOI: 10.1364/AO.58.002248 (CF660C CF680 sSMLM)

Zhang, Y., et al. [Nanoscale subcellular architecture revealed by multicolor three-dimensional salvaged fluorescence imaging](#). Nat Methods 17(2), 225-231, (2020), DOI: 10.1038/s41592-019-0676-4 (CF660C & CF680 single labeled, CF680-wheat germ agglutinin)

Zhang, Y., et al. [Minimizing Molecular Misidentification in Imaging Low-Abundance Protein Interactions Using Spectroscopic Single-Molecule Localization Microscopy](#). Anal Chem 94, 13834-13841, (2022), DOI: 10.1021/acs.analchem.2c02417 (CF660C, CF680 secondaries, super-resolution, SMLM)

Zhang, Z., et al. [Ultrahigh-throughput single-molecule spectroscopy and spectrally resolved super-resolution microscopy](#). Nat Methods 12(10), 935-938, (2015), DOI: nmeth.3528 [pii] 10.1038/nmeth.3528 (CF660C, CF680 STORM)

Zijlstra, N., et al. [Rapid Microfluidic Dilution for Single-Molecule Spectroscopy of Low-Affinity Biomolecular Complexes](#). Angew Chem Int Ed Engl 56(25), 7126-7129, (2017), DOI: 10.1002/anie.201702439 (CF680 3 color FRET)

CF®680R

Conley, G. M., et al. [Jamming and overpacking fuzzy microgels: Deformation, interpenetration, and compression](#). Sci Adv 3(10), e1700969, (2017), DOI: 10.1126/sciadv.1700969 1700969 [pii] (CF680R, STORM)

Gao, F., et al. [Ultrasmall \[\(64\)Cu\]Cu nanoclusters for targeting orthotopic lung tumors using accurate positron emission tomography imaging](#). ACS Nano 9(5), 4976-4986, (2015), DOI: 10.1021/nn507130k (CF680R)

Gonzalez Pisfil, M., et al. [Stimulated emission depletion microscopy with a single depletion laser using five fluorochromes and fluorescence lifetime phasor separation](#). Sci Rep 12, 14027, (2022), DOI: 10.1038/s41598-022-17825-5 (CF594 WGA, CF680R DAMs, super-resolution, confocal, FLIM, fluorescence microscopy, 2-Photon)

Gorlitz, F., et al. [A STED Microscope Designed for Routine Biomedical Applications](#). Progress In Electromagnetics Research 147, 57-68, (2014), DOI: http://dx.doi.org/10.2528/PIER14042708 (CF680R STED)

Gourbeyre, P., et al. [Pattern recognition receptors in the gut: analysis of their expression along the intestinal tract and the crypt/villus axis](#). Physiol Rep 3(2), (2015), DOI: 10.14814/phy2.12225 (CF680R goat anti-rabbit)

Harmsen, S., et al. [A dual-modal PET/near infrared fluorescent nanotag for long-term immune cell tracking](#). Biomaterials 269, 120630, (2021), DOI: 10.1016/j.biomaterials.2020.120630 (CF680R maleimide)

Huang, T., et al. [Simultaneous Multicolor Single-Molecule Tracking with Single-Laser Excitation via Spectral Imaging](#). Biophys J 114(2), 301-310, (2018), DOI: S0006-3495(17)31242-0 [pii] 10.1016/j.bpj.2017.11.013 (CF680R STORM)

Konig, I., et al. [Impact of incell and invitro crowding on the conformations and dynamics of an intrinsically disordered protein](#). Angew Chem Int Ed Engl 60(19), 10724-10729, (2021), DOI: 10.1002/anie.202016804 (CF680R maleimide) (FCF680R)

Konig, I., et al. [Single-molecule spectroscopy of protein conformational dynamics in live eukaryotic cells](#). Nat Methods 12(8), 773-779, (2015), DOI: nmeth.3475 [pii] 10.1038/nmeth.3475 (CF680R single molecule spectroscopy)

Moran, I., et al. [Memory B cells are reactivated in subcapsular proliferative foci of lymph nodes](#). Nat Commun 9(1), 3372, (2018), DOI: 10.1038/s41467-018-05772-7 (CF680R 2-photon)

Seo, J., et al. [PICASSO: Ultra-multiplexed fluorescence imaging of biomolecules through single-round imaging and blind source unmixing](#). bioRxiv, (2021), DOI: 10.1101/2021.01.27.428247 (CF633, CF660R, CF680R)

Seo, J., et al. [PICASSO allows ultra-multiplexed fluorescence imaging of spatially overlapping proteins without reference spectra measurements](#). Nat Commun 13, 2475, (2022), DOI: 10.1038/s41467-022-30168-z (CF405S, CF405M, CF405L, CF488A, CF568, CF633, CF660R, CF680R)

Sun, X., et al. [Longitudinal monitoring of tumor antiangiogenic therapy with near-infrared fluorophore-labeled agents targeted to integrin alphavbeta3 and vascular endothelial growth factor](#). Eur J Nucl Med Mol Imaging 41(7), 1428-1439, (2014), DOI: 10.1007/s00259-014-2702-1 (Xenolight CF680 CF770; in vivo imaging)

Valdoz, J. C., et al. [Soluble ECM promotes organotypic formation in lung alveolar model](#). Biomaterials 283, 121464, (2022), DOI: 10.1016/j.biomaterials.2022.121464 (CF680R streptavidin, fluorescence microscopy, 3D culture matrigel)

Winter, F. R., et al. [Multicolour nanoscopy of fixed and living cells with a single STED beam and hyperspectral detection](#). Sci Rep 7, 46492, (2017), DOI: srep46492 [pii] 10.1038/srep46492 (CF680R STED)

Yamashiro, S. and Watanabe, N. [An Infrared Actin Probe for Deep-Cell Electroporation-Based Single-Molecule Speckle \(eSiMS\) Microscopy](#). Sensors (Basel) 17(7), (2017), DOI: 10.3390/s17071545 (CF680R succinimidyl ester, CF750 succinimidyl ester & CF770 succinimidyl ester)

Yu, C. H., et al. [Integrin-beta3 clusters recruit clathrin-mediated endocytic machinery in the absence of traction force](#). Nat Commun 6, 8672, (2015), DOI: 10.1038/ncomms9672 (CF680R maleimide and CF405 phalloidin)

CF®750

Alawieh, A., et al. [Modulation of post-stroke degenerative and regenerative processes and subacute protection by site-targeted inhibition of the alternative pathway of complement](#). Journal of

Neuroinflammation 12(1), 1-15, (2015), DOI: 10.1186/s12974-015-0464-8 (CF750 succinimidyl ester in vivo)

Aredo, B., et al. [Differences in the distribution, phenotype and gene expression of subretinal microglia/macrophages in C57BL/6N \(Crb1 rd8/rd8\) versus C57BL6/J \(Crb1 wt/wt\) mice](#). J Neuroinflammation 12, 6, (2015), DOI: 10.1186/s12974-014-0221-4, s12974-014-0221-4 [pii] (CF750 donkey anti-rabbit)

Cai, H., et al. [Positron Emission Tomography Imaging of Platelet-Derived Growth Factor Receptor beta in Colorectal Tumor Xenograft Using Zirconium-89 Labeled Dimeric Affibody Molecule](#). Mol Pharm 16(5), 1950-1957, (2019), DOI: 10.1021/acs.molpharmaceut.8b01317 (CF750SE in vivo)

Chen, X. and Zhang, W. [Diamond nanostructures for drug delivery, bioimaging, and biosensing](#). Chem Soc Rev 46(3), 734-760, (2017), DOI: 10.1039/c6cs00109b (CF750SE in vivo)

Darniot, M., et al. [RNA interference in vitro and in vivo using DsiRNA targeting the nucleocapsid N mRNA of human metapneumovirus](#). Antiviral Res 93(3), 364-373, (2012), DOI: 10.1016/j.antiviral.2012.01.004 (CF750 succinimidyl ester (to label siRNA); in vivo imaging; Visen Medical FMT 2500 imaging system)

Fan, Q., et al. [Biological evaluation of 131I- and CF750-labeled Dmab\(scFv\)-Fc antibodies for xenograft imaging of CD25-positive tumors](#). Biomed Res Int 2014, 459676, (2014), DOI: 10.1155/2014/459676 (CF750; in vivo imaging; Caliper IVIS)

Guo, K., et al. [Engineering the first chimeric antibody in targeting intracellular PRL-3 oncoprotein for cancer therapy in mice](#). Oncotarget 3(2), 158-171, (2012), (XenoLight CF750 NIR Fluorescent Labeling Kit; in vivo imaging; PerkinElmer IVIS Spectrum imaging system)

Hapuarachchige, S., et al. [Bioorthogonal two-component drug delivery in HER2\(+\) breast cancer mouse models](#). Sci Rep 6, 24298, (2016), DOI: 10.1038/srep24298 (CF680 succinimidyl ester & CF750 succinimidyl ester, in vivo)

Hashimoto, Y., et al. [Claudin-targeted drug development using anti-claudin monoclonal antibodies to treat hepatitis and cancer](#). Ann N Y Acad Sci 1397(1), 5-16, (2017), DOI: 10.1111/nyas.13337 (CF750SE)

Herz, C., et al. [The isothiocyanate erucin abrogates telomerase in hepatocellular carcinoma cells in vitro and in an orthotopic xenograft tumour model of HCC](#). J Cell Mol Med, (2014), DOI: 10.1111/jcmm.12412 (Xenolight CF750; in vivo imaging; Bruker Image Station)

Ito, A., et al. [New whole-body multimodality imaging of gastric cancer peritoneal metastasis combining fluorescence imaging with ICG-labeled antibody and MRI in mice](#). Gastric Cancer 17(3), 497-507, (2014), DOI: 10.1007/s10120-013-0316-0 (Xenolight CF750; in vivo imaging; IVIS Lumina II)

Kimbrough, C. W., et al. [Orthotopic pancreatic tumors detected by optoacoustic tomography using Syndecan-1](#). J Surg Res 193(1), 246-254, (2015), DOI: 10.1016/j.jss.2014.06.045 (CF750SE)

Li, X., et al. [Development of an anti-claudin-3 and -4 bispecific monoclonal antibody for cancer diagnosis and therapy](#). J Pharmacol Exp Ther 351(1), 206-213, (2014), DOI: jpet.114.216911 [pii] 10.1124/jpet.114.216911 (Xenolight CF750; in vivo imaging)

Li, X., et al. [Tissue distribution and safety evaluation of a claudin-targeting molecule, the C-terminal fragment of Clostridium perfringens enterotoxin](#). Eur J Pharm Sci 52, 132-137, (2014), DOI: S0928-0987(13)00425-9 [pii] 10.1016/j.ejps.2013.10.018 (Xenolight CF750; FC, in vivo imaging)

Matsuo, H., et al. [Engineered hepatitis B virus surface antigen L protein particles for in vivo active targeting of splenic dendritic cells](#). Int J Nanomedicine 7, 3341-3350, (2012), DOI: DOI:10.2147/IJN.S32813 (CF750 succinimidyl ester; bionanocapsules (BNCs); in vivo imaging; Olympus OV-100 imaging system)

Moore, L., et al. [Diamond-Lipid Hybrids Enhance Chemotherapeutic Tolerance and Mediate Tumor Regression](#). Adv Mater 25(26), 3532, (2013), DOI: DOI:10.1002/adma.201300343 (XenoLight CF750 in Nanodiamond-lipid hybrid particles (NDLPs); in vivo imaging; PerkinElmer IVIS Spectrum imaging system)

Stoltzfus, C. R., et al. [CytoMAP: A Spatial Analysis Toolbox Reveals Features of Myeloid Cell Organization in Lymphoid Tissues](#). Cell Rep 31(3), 107523, (2020), DOI: 10.1016/j.celrep.2020.107523 (CF750, CF594, CF514, CF633 Mix-n-Stain)

Sun, X., et al. [Longitudinal monitoring of tumor antiangiogenic therapy with near-infrared fluorophore-labeled agents targeted to integrin alphavbeta3 and vascular endothelial growth factor](#). Eur J Nucl Med Mol Imaging 41(7), 1428-1439, (2014), DOI: 10.1007/s00259-014-2702-1 (Xenolight CF680 CF770; in vivo imaging)

Wei, D., et al. [CF750-A33scFv-Fc-Based Optical Imaging of Subcutaneous and Orthotopic Xenografts of GPA33-Positive Colorectal Cancer in Mice](#). BioMed Research International 2015, 11, (2015), DOI: 10.1155/2015/505183 (CF750 succinimidyl ester in vivo)

Wu, J., et al. [Optical time-stretch microscopy at visible wavelengths](#). Frontiers in Optics, (2016), DOI: <https://doi.org/10.1364/FIO.2016.FF2A.3> (CF750 succinimidyl ester)

Yamashiro, S. and Watanabe, N. [An Infrared Actin Probe for Deep-Cell Electroporation-Based Single-Molecule Speckle \(eSiMS\) Microscopy](#). Sensors (Basel) 17(7), (2017), DOI: 10.3390/s17071545 (CF680R succinimidyl ester, CF750 succinimidyl ester & CF770 succinimidyl ester)

Zeiderman, M. R., et al. [Targeting of BRAF resistant melanoma via extracellular matrix metalloproteinase inducer receptor](#). J Surg Res 190(1), 111-118, (2014), DOI: S0022-4804(14)00165-6 [pii] 10.1016/j.jss.2014.02.021 (CF750 succinimidyl ester; in vivo imaging, FC)

Zhang, Y., et al. [One-Shot Immunomodulatory Nanodiamond Agents for Cancer Immunotherapy](#). Adv Mater 28(14), 2699-2708, (2016), DOI: 10.1002/adma.201506232 (CF750SE in vivo)

CF®770

Abe, S., et al. [A novel targeting therapy of malignant mesothelioma using anti-podoplanin antibody](#). J Immunol 190(12), 6239-6249, (2013), DOI: DOI:10.4049/jimmunol.1300448 (XenoLight CF770; in vivo imaging; PerkinElmer IVIS Spectrum imaging system)

Ait-Belgnaoui, A., et al. [Probiotic gut effect prevents the chronic psychological stress-induced brain activity abnormality in mice](#). Neurogastroenterol Motil 26(4), 510-520, (2014), DOI: 10.1111/nmo.12295 (CF770; WB; LI-COR Odyssey)

Alfonso-Loeches, S., et al. [Toll-like receptor 4 participates in the myelin disruptions associated with chronic alcohol abuse](#). Glia 60(6), 948-964, (2012), DOI: DOI:10.1002/glia.22327 (XenoLight CF770; in vivo imaging (mouse brain); PerkinElmer IVIS 200 imaging system)

Alfonso-Loeches, S., et al. [Gender differences in alcohol-induced neurotoxicity and brain damage](#). Toxicology 311(1-2), 27, (2013), DOI: DOI:10.1016/j.tox.2013.03.001 (XenoLight CF770; in vivo imaging (mouse brain); PerkinElmer IVIS 200 imaging system)

Alfonso-Loeches, S., et al. [Ethanol-Induced TLR4/NLRP3 Neuroinflammatory Response in Microglial Cells Promotes Leukocyte Infiltration Across the BBB](#). Neurochem Res 41(1-2), 193-209, (2016), DOI: 10.1007/s11064-015-1760-5 (CF770 neg con in vivo imaging)

Andriamihaja, M., et al. [The deleterious metabolic and genotoxic effects of the bacterial metabolite p-cresol on colonic epithelial cells](#). Free Radical Biology and Medicine 85, 219-227, (2015), DOI: http://dx.doi.org/10.1016/j.freeradbiomed.2015.04.004 (CF770 RedDot1 in cell western)

Annahazi, A., et al. [Luminal Cysteine-Proteases Degrade Colonic Tight Junction Structure and Are Responsible for Abdominal Pain in Constipation-Predominant IBS](#). Am J Gastroenterol 108(8), 1392, (2013), DOI: DOI:10.1038/ajg.2013.152 (CF770 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Audebert, M., et al. [Comparative potency approach based on H2AX assay for estimating the genotoxicity of polycyclic aromatic hydrocarbons](#). Toxicol Appl Pharmacol 260(1), 58, (2012), DOI: DOI:10.1016/j.taap.2012.01.022 (CF770 goat anti-rabbit IgG; ICW; LI-COR Odyssey)

Azhdarinia, A., et al. [Dual-labeling strategies for nuclear and fluorescence molecular imaging: a review and analysis](#). Mol Imaging Biol 14(3), 261-276, (2012), DOI: 10.1007/s11307-011-0528-9 (CF770, CF790; in vivo imaging)

Babelova, A., et al. [Activation of Rac-1 and RhoA contributes to podocyte injury in chronic kidney disease](#). PLoS One 8(11), e80328, (2013), DOI: 10.1371/journal.pone.0080328, PONE-D-13-15419 [pii] (CF770 goat anti-rabbit; WB; LI-COR Odyssey)

Beaumont, M., et al. [Detimental effects for colonocytes of an increased exposure to luminal hydrogen sulfide: The adaptive response](#). Free Radic Biol Med 93, 155-164, (2016), DOI: S0891-5849(16)00042-3 [pii] 10.1016/j.freeradbiomed.2016.01.028 (CF770 goat anti-rabbit, RedDot2 ICW)

Bonnaud, E. M., et al. [Borna disease virus phosphoprotein modulates epigenetic signaling in neurons to control viral replication](#). J Virol 89(11), 5996-6008, (2015), DOI: 10.1128/JVI.00454-15 (CF680 & CF770 WB)

Budin, C., et al. [Versicolorin A enhances the genotoxicity of aflatoxin B1 in human liver cells by inducing the transactivation of the Ah-receptor](#). Food Chem Toxicol 153, 112258, (2021), DOI: 10.1016/j.fct.2021.112258 (CF770 goat anti-rat RedDot2)

Chevreau, M., et al. [Role of human sulfotransferase 1A1 and N-acetyltransferase 2 in the metabolic activation of 16 heterocyclic amines and related heterocyclics to genotoxins in recombinant V79 cells](#). Arch Toxicol 91(9), 3175-3184, (2017), DOI: 10.1007/s00204-017-1935-8 (CF770 secondary RedDot ICW)

Comor, L., et al. [Joining the in vitro immunization of alpaca lymphocytes and phage display: rapid and cost effective pipeline for sdAb synthesis](#). Microb Cell Fact 16(1), 13, (2017), DOI: 10.1186/s12934-017-0630-z (CF770 anti-GFP)

Dabek, M., et al. [Intracolonic infusion of fecal supernatants from ulcerative colitis patients triggers altered permeability and inflammation in mice: role of cathepsin G and protease-activated receptor-4](#). Inflamm Bowel Dis 17(6), 1409-1414, (2011), DOI: DOI:10.1002/ibd.21454 (CF770 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Degrelle, S. A., et al. [Annexin-A5 organized in 2D-network at the plasmalemma eases human trophoblast fusion](#). Sci Rep 7, 42173, (2017), DOI: 10.1038/srep42173 (CF770 AnnexinV blot overlay)

Ducheix, S., et al. [Dietary oleic acid regulates hepatic lipogenesis through a liver X receptor-dependent signaling](#). PLoS One 12(7), e0181393, (2017), DOI: 10.1371/journal.pone.0181393 (CF680 & CF770 secondaries WB)

Frank, C. L., et al. [Regulation of chromatin accessibility and Zic binding at enhancers in the developing cerebellum](#). Nat Neurosci 18(5), 647-656, (2015), DOI: 10.1038/nn.3995 (CF680 goat anti-mouse & CF770 goat anti-rabbit WB)

Gauthier, T., et al. [Deoxynivalenol impairs the immune functions of neutrophils](#). Mol Nutr Food Res 57(6), 1026, (2013), DOI: DOI:10.1002/mnfr.201200755 (CF680 goat anti-rabbit IgG, CF770 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Gon, S., et al. [Fit alphabeta T-cell receptor suppresses leukemogenesis of Pten-deficient thymocytes](#). Haematologica 103(6), 999-1007, (2018), DOI: 10.3324/haematol.2018.188359 (CF680 goat anti-mouse and CF770 goat anti-rabbit WB)

Graber, T. E., et al. [UPF1 Governs Synaptic Plasticity through Association with a STAU2 RNA Granule](#). J Neurosci 37(38), 9116-9131, (2017), DOI: 10.1523/JNEUROSCI.0088-17.2017 (CF680 & CF770 secondaries WB)

Hilbers, F. S., et al. [Functional Analysis of Missense Variants in the Putative Breast Cancer Susceptibility Gene XRCC2](#). Hum Mutat 37(9), 914-925, (2016), DOI: 10.1002/humu.23019 (CF680 goat anti-rabbit & CF770 goat anti-mouse WB)

Huc, L., et al. [Low concentrations of bisphenol A induce lipid accumulation mediated by the production of reactive oxygen species in the mitochondria of HepG2 cells](#). Toxicol In Vitro 26(5), 709-717, (2012), DOI: DOI:10.1016/j.tiv.2012.03.017 (CF770 goat anti-mouse IgG; WB; LI-COR Odyssey)

Jamin, E. L., et al. [Combined Genotoxic Effects of a Polycyclic Aromatic Hydrocarbon \(B\(a\)P\) and an Heterocyclic Amine \(PhIP\) in Relation to Colorectal Carcinogenesis](#). PLoS One 8(3), e58591, (2013), DOI: DOI:10.1371/journal.pone.0058591 (CF770 goat anti-rabbit IgG; ICW; LI-COR Odyssey)

Khoury, L., et al. [Validation of high-throughput genotoxicity assay screening using gammaH2AX in-cell western assay on HepG2 cells](#). Environ Mol Mutagen 54(9), 737-746, (2013), DOI: 10.1002/em.21817 (CF770; ICW)

Khoury, L., et al. [Evaluation of four human cell lines with distinct biotransformation properties for genotoxic screening](#). Mutagenesis 31(1), 83-96, (2016), DOI: 10.1093/mutage/gev058 (CF770 goat anti-rabbit, RedDot2, ICW)

Kim, J. Y., et al. [Efficacy of gastro-retentive forms of ecabet sodium in the treatment of gastric ulcer in rats](#). Arch Pharm Res 37(8), 1053-1062, (2014), DOI: 10.1007/s12272-013-0278-0 (CF770; in vivo imaging)

Lai, S. R., et al. [In vitro anti-tubulin effects of mebendazole and fenbendazole on canine glioma cells](#). Vet Comp Oncol 15(4), 1445-1454, (2017), DOI: 10.1111/vco.12288 (CF770 goat anti-mouse WB)

Leskoske, K. [Modulation of TOR Complex 2 signaling and maintenance of plasma membrane homeostasis in *Saccharomyces cerevisiae*](#). in: Graduate Division, PhD in Molecular and Cell Biology, UC Berkeley, 106, (2017), (CF770 goat anti-mouse WB)

Leskoske, K. L., Roelants, F. M., Martinez Marshall, M. N., Hill, J. M., Thorner, J. . [The stress-sensing TORC2 complex activates yeast AGC-family protein kinase Ypk1 at multiple novel sites](#). Genetics 207(1), 179-195, (2017), DOI: <https://doi.org/10.1534/genetics.117.1124> (CF770 goat anti-mouse)

Liang, Y., et al. [The lymphoid lineage-specific actin-uncapping protein Rltp is essential for costimulation via CD28 and the development of regulatory T cells](#). Nat Immunol 14(8), 858-866, (2013), DOI: [ni.2634 \[pii\]](https://doi.org/10.1038/ni.2634) 10.1038/ni.2634 (CF770; WB)

Lucioli, J., et al. [The food contaminant deoxynivalenol activates the mitogen activated protein kinases in the intestine: Interest of ex vivo models as an alternative to in vivo experiments](#). Toxicon 66, 31, (2013), DOI: [DOI:10.1016/j.toxicon.2013.01.024](https://doi.org/10.1016/j.toxicon.2013.01.024) (CF770 goat anti-mouse IgG, CF770 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Martinez Marshall, M. N., et al. [Analysis of the roles of phosphatidylinositol-4,5-bisphosphate and individual subunits in assembly, localization, and function of *Saccharomyces cerevisiae* target of rapamycin complex 2](#). Mol Biol Cell 30(12), 1555-1574, (2019), DOI: [10.1091/mbc.E18-10-0682](https://doi.org/10.1091/mbc.E18-10-0682) (CF770 goat anti-mouse WB)

Martinsen, A., et al. [Rho kinase regulation of vasopressin-induced calcium entry in vascular smooth muscle cell: Comparison between rat isolated aorta and cultured aortic cells](#). Cell Calcium 52(6), 413, (2012), DOI: [DOI:10.1016/j.ceca.2012.07.002](https://doi.org/10.1016/j.ceca.2012.07.002) (CF770 goat anti-mouse IgG, CF770 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Mazzocco, C., et al. [In vivo imaging of prostate cancer using an anti-PSMA scFv fragment as a probe](#). Sci Rep 6, 23314, (2016), DOI: [10.1038/srep23314](https://doi.org/10.1038/srep23314) (XenoLite CF770)

Moussa, L., et al. [A new soy germ fermented ingredient displays estrogenic and protease inhibitor activities able to prevent irritable bowel syndrome-like symptoms in stressed female rats](#). Clin Nutr 32(1), 51, (2012), DOI: [DOI:10.1016/j.clnu.2012.05.021](https://doi.org/10.1016/j.clnu.2012.05.021) (CF770 anti-rabbit IgG; WB; LI-COR Odyssey)

Nebot-Vivinus, M., et al. [Multispecies probiotic protects gut barrier function in experimental models](#). World J Gastroenterol 20(22), 6832-6843, (2014), DOI: [10.3748/wjg.v20.i22.6832](https://doi.org/10.3748/wjg.v20.i22.6832) (CF770 goat anti-rabbit WB Licor)

Nguyen, T. M. D., et al. [Inhibition by fluoxetine of LH-stimulated cyclic AMP synthesis in tumor Leydig cells partly involves AMPK activation](#). PLoS One 14(6), e0217519, (2019), DOI: [10.1371/journal.pone.0217519](https://doi.org/10.1371/journal.pone.0217519) (CF770 anti-rabbit WB)

Norris, J. D., et al. [Androgen receptor antagonism drives cytochrome P450 17A1 inhibitor efficacy in prostate cancer](#). J Clin Invest 127(6), 2326-2338, (2017), DOI: [10.1172/JCI87328](https://doi.org/10.1172/JCI87328) (CF770 goat anti-rabbit In Cell Western)

Oliveira, A. F. and Yasuda, R. [Neurofibromin is the major ras inactivator in dendritic spines](#). J Neurosci 34(3), 776-783, (2014), DOI: [34/3/776 \[pii\] 10.1523/JNEUROSCI.3096-13.2014](https://doi.org/10.1523/JNEUROSCI.3096-13.2014) (CF680 goat anti-mouse, CF770 goat anti-rabbit; WB; LI-COR Odyssey)

Olivier, I., et al. [Is Crohn's creeping fat an adipose tissue?](#). Inflamm Bowel Dis 17(3), 747-757, (2011), DOI: [DOI:10.1002/ibd.21413](https://doi.org/10.1002/ibd.21413) (CF770 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Pollock, J. A., et al. [Inhibiting androgen receptor nuclear entry in castration-resistant prostate cancer](#). Nat Chem Biol 12(10), 795-801, (2016), DOI: [10.1038/nchembio.2131](https://doi.org/10.1038/nchembio.2131) (CF770 goat anti-rabbit In Cell Western)

Pons, B. J., et al. [Cell transfection of purified cytolethal distending toxin B subunits allows comparing their nuclease activity while plasmid degradation assay does not](#). PLoS One 14(3), e0214313, (2019), DOI: 10.1371/journal.pone.0214313 (CF770 goat anti-mouse WB)

Pulzova, L., et al. [Identification of B-cell epitopes of Borrelia burgdorferi outer surface protein C by screening a phage-displayed gene fragment library](#). Microbiol Immunol 60(10), 669-677, (2016), DOI: 10.1111/1348-0421.12438 (CF770SE)

Rashid, F. M., et al. [HI-NESS: a family of genetically encoded DNA labels based on a bacterial nucleoid-associated protein](#). Nucleic Acids Res 50, e10, (2022), DOI: 10.1093/nar/gkab993 (CF680 GARb, CF770 GAMs)

Rother, M. B., et al. [The Human Thymus Is Enriched for Autoreactive B Cells](#). J Immunol 197(2), 441-448, (2016), DOI: 10.4049/jimmunol.1501992 (CF680 GAHu & CF770 goat anti-mouse WB)

Schimmel, J., et al. [Mutational signatures of non-homologous and polymerase theta-mediated end-joining in embryonic stem cells](#). EMBO J 36(24), 3634-3649, (2017), DOI: 10.15252/embj.201796948 (CF680 goat anti-rabbit & CF770 goat anti-mouse WB)

Suemizu, H., et al. [A versatile technique for the in vivo imaging of human tumor xenografts using near-infrared fluorochrome-conjugated macromolecule probes](#). PLoS One 8(12), e82708, (2013), DOI: 10.1371/journal.pone.0082708 PONE-D-13-13248 [pii] (Xenolight CF770; in vivo imaging)

Sun, Y., et al. [Single tumor imaging with multiple antibodies targeting different antigens](#). Biotechniques 52(4), 1-3, (2012), DOI: DOI:10.2144/000113855 (XenoLight CF680, XenoLight CF770; in vivo imaging; PerkinElmer IVIS Lumina)

Theumer, M. G., et al. [Genotoxicity of aflatoxins and their precursors in human cells](#). Toxicol Lett 287, 100-107, (2018), DOI: 10.1016/j.toxlet.2018.02.007 (CF770 & RedDot ICW)

Titova, E. and Obenaus, A. [Immunological Response to Experimental Intracerebral Hemorrhage: Morphological Assessments](#). In: Chen, J., et al. (Eds.), Animal Models of Acute Neurological Injuries II. Humana Press, pp. 625-652, (2012), DOI: DOI:10.1007/978-1-61779-576-3_48 (CF680 and CF770 secondary antibody conjugates; IHC in floating brain sections; LI-COR Odyssey)

Titova, E. M., et al. [The late phase of post-stroke neurorepair in aged rats is reflected by MRI-based measures](#). Neuroscience, (2014), DOI: S0306-4522(14)00770-2 [pii] 10.1016/j.neuroscience.2014.09.028 (CF680 goat anti-mouse, CF770 goat anti-rat; IHC; LI-COR Odyssey)

van den Heuvel, D., et al. [A CSB-PAF1C axis restores processive transcription elongation after DNA damage repair](#). Nat Commun 12(1), 1342, (2021), DOI: 10.1038/s41467-021-21520-w (CF680 goat anti-rabbit CF770 goat anti-mouse WB Odyssey)

van der Weegen, Y., et al. [The sequential and cooperative action of CSB, CSA, and UVSSA targets the TFIIB complex to DNA damage-stalled RNA polymerase II](#). bioRxiv, (2019), DOI: 10.1101/707216 (CF680 donkey anti-goat CF680 goat anti-rabbit CF770 goat anti-mouse WB)

Yamashiro, S. and Watanabe, N. [An Infrared Actin Probe for Deep-Cell Electroporation-Based Single-Molecule Speckle \(eSiMS\) Microscopy](#). Sensors (Basel) 17(7), (2017), DOI: 10.3390/s17071545 (CF680R succinimidyl ester, CF750 succinimidyl ester, CF770 succinimidyl ester)

CF®790

Azhdarinia, A., et al. [Dual-labeling strategies for nuclear and fluorescence molecular imaging: a review and analysis](#). Mol Imaging Biol 14(3), 261-276, (2012), DOI: 10.1007/s11307-011-0528-9 (CF770, CF790; in vivo imaging)

Ghosh, S. C., et al. [Multimodal chelation platform for near-infrared fluorescence/nuclear imaging](#). J Med Chem 56(2), 406-416, (2013), DOI: 10.1021/jm300906g (CF790 in vivo imaging)

Najimi, M., et al. [Human liver mesenchymal stem/progenitor cells inhibit hepatic stellate cell activation: in vitro and in vivo evaluation](#). Stem Cell Res Ther 8(1), 131, (2017), DOI: 10.1186/s13287-017-0575-5 (CF680 donkey anti-rabbit & CF790 donkey anti-mouse WB)

Takeyama, H., et al. [A novel endoscopic fluorescent clip visible with near-infrared imaging during laparoscopic surgery in a porcine model](#). Surg Endosc 28(6), 1984-1990, (2014), DOI: 10.1007/s00464-014-3423-z (CF790; in vivo imaging)

CF® Dye References by Application

CF® Dye Mix-n-Stain™ Antibody Labeling Kits

Ahi, Y. S., et al. [Adenoviral E2 IVa2 Protein Interacts with L4 33K Protein and E2 DNA-Binding Protein](#). J Gen Virol 94(6), 1325, (2013), DOI: DOI:10.1099/vir.0.049346-0 (Mix-n-Stain CF350, Mix-n-Stain CF488A, CF568 goat anti-mouse IgG; ICC)

Anderson, M. S., et al. [Maturation and vesicle-mediated egress of primate gammaherpesvirus rhesus monkey rhadinovirus require inner tegument protein ORF52](#). J Virol 88(16), 9111-9128, (2014), DOI: JVI.01502-14 [pii] 10.1128/JVI.01502-14 (CF488A Mix n Stain; ICC)

Aravindan, S., et al. [Polyphenols from marine brown algae target radiotherapy-coordinated EMT and stemness-maintenance in residual pancreatic cancer](#). Stem Cell Res Ther 6, 182, (2015), DOI: 10.1186/s13287-015-0173-3 (Mix-n-Stain CF488A)

Arul, M., et al. [Culture of low passage colorectal cancer cells and demonstration of variation in selected tumour marker expression](#). Cytotechnology 66(3), 481, (2013), DOI: 10.1007/s10616-013-9600-4 (CF488A Mix-n-Stain, CF594 Mix-n-Stain, CF488A goat anti-mouse; ICC)

Balogh, A., et al. [Sex hormone-binding globulin provides a novel entry pathway for estradiol and influences subsequent signaling in lymphocytes via membrane receptor](#). Sci Rep 9(1), 4, (2019), DOI: 10.1038/s41598-018-36882-3 (CF633 Mix-n-Stain)

Bolduc, J. F., et al. [Epigenetic Metabolite Acetate Inhibits Class I/II Histone Deacetylases, Promotes Histone Acetylation, and Increases HIV-1 Integration in CD4\(+\) T Cells](#). J Virol 91(16), (2017), DOI: 10.1128/JVI.01943-16 (CF4505M Mix-n-Stain flow)

Breton-Provencher, V., et al. [Principal cell activity induces spine relocation of adult-born interneurons in the olfactory bulb](#). Nat Commun 7, 12659, (2016), DOI: 10.1038/ncomms12659 (CF568 Mix-n-Stain IF)

Ciapa, B. and Granon, S. [Expression of Cyclin-D1 in Astrocytes Varies During Aging](#). Front Aging Neurosci 10, 104, (2018), DOI: 10.3389/fnagi.2018.00104 (CF568 Mix-n-Stain IF)

Darvish-Damavandi, M., et al. [Towards the development of an enzyme replacement therapy for the metabolic disorder propionic acidemia](#). Mol Genet Metab Rep 8, 51-60, (2016), DOI: 10.1016/j.ymgmr.2016.06.009 S2214-4269(16)30048-9 [pii] (CF488A Mix-n-Stain)

De Pace, R., et al. [Synaptic Vesicle Precursors and Lysosomes Are Transported by Different Mechanisms in the Axon of Mammalian Neurons](#). Cell Rep 31(11), 107775, (2020), DOI: 10.1016/j.celrep.2020.107775 (CF640R Mix-n-Stain)

Deo, D. I., et al. [Biofunctionalization of PEGylated microcapsules for exclusive binding to protein substrates](#). Biomacromolecules 15(7), 2555-2562, (2014), DOI: 10.1021/bm500412d (CF555 Mix-n-Stain microcapsules; fluorescence microscopy)

Diamantopoulou, A., et al. [Denial or receipt of expected reward through maternal contact during the neonatal period differentially affect the development of the rat amygdala and program its function in adulthood in a sex-dimorphic way](#). Psychoneuroendocrinology 38(9), 1757, (2013), DOI: DOI:10.1016/j.psyneuen.2013.02.012 (Mix-n-Stain CF488A, Mix-N-Stain CF555; IHC)

Donskow-Lysoniewska, K., et al. [Heligmosmoides polygyrus fourth stages induce protection against DSS-induced colitis and change opioid expression in the intestine](#). Parasite Immunol 34(11), 536, (2012), DOI: DOI:10.1111/pim.12003 (Mix-n-Stain CF488A; IHC)

Ersek, B., et al. [CD3zeta-Chain Expression of Human T Lymphocytes Is Regulated by TNF via Src-like Adaptor Protein-Dependent Proteasomal Degradation](#). J Immunol 189(4), 1602, (2012), DOI: DOI:10.4049/jimmunol.1102365 (Mix-n-Stain CF488A; ICC)

Farias, G. G., et al. [Feedback-Driven Mechanisms between Microtubules and the Endoplasmic Reticulum Instruct Neuronal Polarity](#). Neuron 102(1), 184-201 e188, (2019), DOI: 10.1016/j.neuron.2019.01.030 (CF640R Mix-n-Stain)

Farias, G. G., et al. [Imaging the Polarized Sorting of Proteins from the Golgi Complex in Live Neurons](#). Methods Mol Biol 1496, 13-30, (2016), DOI: 10.1007/978-1-4939-6463-5_2 (CF488A, CF555 & CF640R Mix-n-Stain IF)

Galletti, G., et al. [Isolation of breast cancer and gastric cancer circulating tumor cells by use of an anti HER2-based microfluidic device](#). Lab Chip 14(1), 147-156, (2014), DOI: 10.1039/c3lc51039e (CF594 Mix-n-Stain; ICC)

Geherin, S. A., et al. [IL-10+ Innate-like B Cells Are Part of the Skin Immune System and Require alpha4beta1 Integrin To Migrate between the Peritoneum and Inflamed Skin](#). J Immunol 196(6), 2514-2525, (2016), DOI: 10.4049/jimmunol.1403246 (CF488A CF594 Mix-n-Stain)

Gennet, N., et al. [FolR1: a novel cell surface marker for isolating midbrain dopamine neural progenitors and nascent dopamine neurons](#). Sci Rep 6, 32488, (2016), DOI: 10.1038/srep32488 (CF640R Mix-n-Stain FACS)

Gorham, J. B., et al. [Visualization of microbe-dietary remnant interactions in digesta from pigs, by fluorescence in situ hybridization and staining methods; effects of a dietary arabinoxylan-rich wheat fraction](#). Food Hydrocolloids 52, 952-962, (2016), DOI: http://dx.doi.org/10.1016/j.foodhyd.2015.09.011 (660R, Mix-n-Stain)

Hadassa Y. Holzapfel, and Marc R. Birtwistle. [Creating Complex Fluorophore spectra on Antibodies Through Combinatorial Labeling](#). Transl Sci, 1-10, (2016). (CF405M, CF488A, CF568, CF640R, FRET, MnS)

Hirtz, M., et al. [Click-Chemistry Immobilized 3D-Infused Microarrays in Nanoporous Polymer Substrates](#). Advanced Materials Interfaces, n/a-n/a, (2016), DOI: 10.1002/admi.201500469 (CF488A Mix-n-Stain)

Homann, S., et al. [A novel rapid and reproducible flow cytometric method for optimization of transfection efficiency in cells](#). PLoS One 12(9), e0182941, (2017), DOI: 10.1371/journal.pone.0182941 (CF647 Mix-n-Stain)

Home, P., et al. [Altered subcellular localization of transcription factor TEAD4 regulates first mammalian cell lineage commitment](#). Proc Natl Acad Sci U S A 109(19), 7362-7367, (2012), DOI: DOI:10.1073/pnas.1201595109 (Mix-n-Stain CF488A, Mix-n-Stain CF568; ICC)

Hwang, J., et al. [Aptamer-conjugated live human immune cell based biosensors for the accurate detection of C-reactive protein](#). Sci Rep 6, 34778, (2016), DOI: 10.1038/srep34778 (CF594 Mix-n-Stain)

Hwang, J., et al. [Aptamer-conjugated live human immune cell based biosensors for the accurate detection of C-reactive protein](#). Sci Rep 6, 34778, (2016), DOI: 10.1038/srep34778 (CF594 Mix-n-Stain)

Jankowska, K. I., et al. [Integrins Modulate T Cell Receptor Signaling by Constraining Actin Flow at the Immunological Synapse](#). Front Immunol 9, 25, (2018), DOI: 10.3389/fimmu.2018.00025 (CF488A Mix-n-Stain, CF405 phalloidin)

Keren-Kaplan, T., et al. [RUFY3 and RUFY4 are ARL8 effectors that promote coupling of endolysosomes to dynein-dynactin](#). Nat Commun 13, 1506, (2022), DOI: 10.1038/s41467-022-28952-y (CF640R, Mix-n-Stain)

Kesik, M., et al. [A multi-functional fluorescent scaffold as a multi-colour probe: design and application in targeted cell imaging](#). RSC Advances 5(101), 83361-83367, (2015), DOI: 10.1039/c5ra16600d (CF555 Mix-n-Stain)

Lim, A., et al. [Two kinesins drive anterograde neuropeptide transport](#). Mol Biol Cell, (2017), DOI: mbc.E16-12-0820 [pii] 10.1091/mbc.E16-12-0820 (Mix-n-Stain CF568, Direct + indirect IF, CF568 3D SIM)

Lim, A. K., et al. [The nuage mediates retrotransposon silencing in mouse primordial ovarian follicles](#). Development 140(18), 3819-3825, (2013), DOI: dev.099184 [pii] 10.1242/dev.099184 (CF488A Mix-n-Stain; IHC)

Lin, T.-B., et al. [Melatonin relieves neuropathic allodynia through spinal MT2-enhanced PP2Ac and downstream HDAC4 shuttling-dependent epigenetic modification of hmgb1 transcription](#). Journal of Pineal Research, n/a-n/a, (2016), DOI: 10.1111/jpi.12307 (CF Dyes Mix-n-Stain)

Loftus, M. S., et al. [A Conserved Leucine Zipper Motif in Gammaherpesvirus ORF52 Is Critical for Distinct Microtubule Rearrangements](#). J Virol 91(17), (2017), DOI: 10.1128/JVI.00304-17 (CF488A Mix-n-Stain)

Lukonin, I., et al. [Phenotypic landscape of intestinal organoid regeneration](#). Nature 586(7828), 275-280, (2020), DOI: 10.1038/s41586-020-2776-9 (CF568 Mix-n-Stain)

MacDonald, A. B. [Lyme ELISA seronegative patients with borrelia antibody coated borrelia microbes in their blood](#). Med Clin Res 7, (2022), ISSN: 2577 - 8005 (CF405S, CF660R, Mix-n-Stain)

Mikosik, A., et al. [Expression of calpain-calpastatin system \(CCS\) member proteins in human lymphocytes of young and elderly individuals; pilot baseline data for the CALPACENT project](#). Immun Ageing 10(1), 27, (2013), DOI: DOI:10.1186/1742-4933-10-27 (Mix-n-Stain CF555; FC)

Miyagishima, S. Y., et al. [DipM is required for peptidoglycan hydrolysis during chloroplast division](#). BMC Plant Biol 14, 57, (2014), DOI: 1471-2229-14-57 [pii] 10.1186/1471-2229-14-57 (CF488A Mix-n-Stain, CF555 Mix-n-Stain; ICC)

Mofarrahi, M., et al. [Angiopoietin-1 enhances skeletal muscle regeneration in mice](#). American Journal of Physiology - Regulatory, Integrative and Comparative Physiology 308(7), R576-R589, (2015), DOI: 10.1152/ajpregu.00267.2014 (CF568 Mix-n-Stain)

Montesinos-Rongen, M., et al. [Primary Central Nervous System \(CNS\) Lymphoma B Cell Receptors Recognize CNS Proteins](#). J Immunol 195(3), 1312-1319, (2015), DOI: jimmunol.1402341 [pii] 10.4049/jimmunol.1402341 (Mix-n-Stain CF488A)

Nagornikh, A. M., et al. [Protocol for chronic hepatitis B virus infection mouse model development by patient-derived orthotopic xenografts](#). PLoS One 17, e0264266, (2022), DOI: 10.1371/journal.pone.0264266 (CF555, Mix-n-Stain)

Niec, R. E., et al. [Lymphatics act as a signaling hub to regulate intestinal stem cell activity](#). Cell Stem Cell 29, 1067-1082 e1018, (2022), DOI: 10.1016/j.stem.2022.05.007 (CF568, Mix-n-Stain)

Park, K. S., et al. [A crucial requirement for Hedgehog signaling in small cell lung cancer](#). Nat Med 17(11), 1504-1508, (2011), DOI: DOI:10.1038/nm.2473 (Mix-n-Stain CF488A, Mix-n-Stain CF568, and Mix-n-Stain CF647; ICC)

Paz, H., et al. [Treatment of B-cell precursor acute lymphoblastic leukemia with the Galectin-1 inhibitor PTX008](#). J Exp Clin Cancer Res 37(1), 67, (2018), DOI: 10.1186/s13046-018-0721-7 (CF647 Mix-n-Stain)

Petrelli, A., et al. [Efficient Conjugation of Oligosaccharides to Polymer Particles through Furan/Maleimide Diels-Alder Reaction: Application to the Capture of Carbohydrate-Binding Proteins](#). Chembiochem 18(2), 206-212, (2017), DOI: 10.1002/cbic.201600509 (CF488A, CF568 Mix-n-Stain)

Pons, M., et al. [Phosphorylation of filamin A regulates chemokine receptor CCR2 recycling](#). J Cell Sci 130(2), 490-501, (2017), DOI: 10.1242/jcs.193821 (CF488A, CF568 Mix-n-Stain)

Robertson, T. F., et al. [Lymphocyte egress signal sphingosine-1-phosphate promotes ERM-guided, bleb-based migration](#). J Cell Biol 220, (2021), DOI: 10.1083/jcb.202007182 (CF568, Mix-n-Stain)

Rohm, M., et al. [Data of rational process optimization for the production of a full IgG and its Fab fragment from hybridoma cells](#). Data Brief 8, 426-435, (2016), DOI: 10.1016/j.dib.2016.05.067 (CF647 Mix-n-Stain)

Ruger, M., et al. [Species-specific viability analysis of Pseudomonas aeruginosa, Burkholderia cepacia and Staphylococcus aureus in mixed culture by flow cytometry](#). BMC Microbiol 14, 56, (2014), DOI: 1471-2180-14-56 [pii] 10.1186/1471-2180-14-56 (CF405S Mix-n-Stain; FC)

Ruger, M., et al. [A flow cytometric method for viability assessment of Staphylococcus aureus and Burkholderia cepacia in mixed culture](#). Cytometry A 81(12), 1055, (2012), DOI: DOI:10.1002/cyto.a.22219 (CF405S Mix-n-Stain labeling of wheat germ agglutinin; FC)

Sainski, A. M., et al. [Casp8p41 generated by HIV protease kills CD4 T cells through direct Bak activation](#). J Cell Biol 207(1), 159, (2014), DOI: jcb.20140505109252014c [pii] 10.1083/jcb.20140505109252014c (CF640R Mix-n-Stain; FC)

Shimizu, F., et al. [Glucose-regulated protein 78 autoantibody associates with blood-brain barrier disruption in neuromyelitis optica](#). Sci Transl Med 9(397), (2017), DOI: 10.1126/scitranslmed.aai9111 (CF568 Mix-n-Stain)

Stoltzfus, C. R., et al. [CytoMAP: A Spatial Analysis Toolbox Reveals Features of Myeloid Cell Organization in Lymphoid Tissues](#). Cell Rep 31(3), 107523, (2020), DOI: 10.1016/j.celrep.2020.107523 (CF750, CF594, CF514, CF633 Mix-n-Stain)

Trofimenko, E., et al. [The endocytic pathway taken by cationic substances requires Rab14 but not Rab5 and Rab7](#). Cell Rep 37, 109945, (2021), DOI: 10.1016/j.celrep.2021.109945 (CF405M, CF488A, CF594, CF647, Mix-n-Stain)

Trofimenko, E., et al. [Identification of a new, Rab14-dependent, endo-lysosomal pathway](#). bioRxiv, (2021), DOI: 10.1101/2021.03.25.436964 (CF405M, CF594, Mix-n-Stain)

Tschumper, R. C., et al. [Molecular analysis of immunoglobulin genes reveals frequent clonal relatedness in double monoclonal gammopathies](#). Blood Cancer J 3, e112, (2013), DOI: DOI:10.1038/bcj.2013.12 (Mix-n-Stain CF350; ICC)

Wada, M., et al. [Epitope-Dependent Pathogenicity of Antibodies Targeting a Major Bullous Pemphigoid Autoantigen Collagen XVII/BP180](#). J Invest Dermatol 136(5), 938-946, (2016), DOI: 10.1016/j.jid.2015.11.030 (CF555 Mix-n-Stain time lapse imaging)

Wisnovsky, S., et al. [Genome-wide CRISPR screens reveal a specific ligand for the glycan-binding immune checkpoint receptor Siglec-7](#). Proc Natl Acad Sci U S A 118(5), (2021), DOI: 10.1073/pnas.2015024118 (CF568 Mix-n-Stain STORM)

Xu, Z., et al. [Mucosal immunoglobulins at respiratory surfaces mark an ancient association that predates the emergence of tetrapods](#). Nat Commun 7, 10728, (2016), DOI: 10.1038/ncomms10728 (CF488A, CF555 & CF640R Mix-n-Stain)

Xu, Z., et al. [Teleost skin, an ancient mucosal surface that elicits gut-like immune responses](#). Proc Natl Acad Sci U S A 110(32), 13097-13102, (2013), DOI: 1304319110 [pii] 10.1073/pnas.1304319110 (CF488A Mix-n-Stain, CF555, CF640R; IHC)

Yamamoto, S., et al. [Contribution of calumin to embryogenesis through participation in the endoplasmic reticulum-associated degradation activity](#). Dev Biol 393(1), 33-43, (2014), DOI: S0012-1606(14)00320-0 [pii] 10.1016/j.ydbio.2014.06.024 (CF488A Mix-n-Stain, CF568 Mix-n-Stain; IHC)

Yamashita, A., et al. [Increased metabolite levels of glycolysis and pentose phosphate pathway in rabbit atherosclerotic arteries and hypoxic macrophage](#). PLoS One 9(1), e86426, (2014), DOI: 10.1371/journal.pone.0086426 PONE-D-13-24660 [pii] (CF488A Mix-n-Stain, CF594 Mix-n-Stain; IHC)

Yang, Y., et al. [Paraoxonase 2 decreases renal reactive oxygen species production, lowers blood pressure, and mediates dopamine D\(2\) receptor-induced inhibition of NADPH oxidase](#). Free Radic Biol Med 53(3), 437, (2012), DOI: DOI:10.1016/j.freeradbiomed.2012.05.015 (Mix-n-Stain CF555; IHC)

Zolnerciks, J. K., et al. [The Q loops of the human multidrug resistance transporter ABCB1 are necessary to couple drug binding to the ATP catalytic cycle](#). FASEB J 28(10), 4335-4346, (2014), DOI: fj.13-245639 [pii] 10.1096/fj.13-245639 (CF633 Mix-n-Stain; ICC)

Flow Cytometry

Bolduc, J. F., et al. [Epigenetic Metabolite Acetate Inhibits Class I/II Histone Deacetylases, Promotes Histone Acetylation, and Increases HIV-1 Integration in CD4\(+\) T Cells](#). J Virol 91(16), (2017), DOI: 10.1128/JVI.01943-16 (CF4505M Mix-n-Stain flow)

Endo, H., et al. [Ibuprofen enhances the anticancer activity of cisplatin in lung cancer cells by inhibiting the heat shock protein 70](#). Cell Death Dis 5, e1027, (2014), DOI: cddis2013550 [pii] 10.1038/cddis.2013.550 (CF488A Annexin V; FC)

Gautam, N., et al. [Preclinical Pharmacokinetics and Tissue Distribution of Long-Acting Nanoformulated Antiretroviral Therapy](#). Antimicrob Agents Chemother 57(7), 3110, (2013), DOI: DOI:10.1128/AAC.00267-13 (CF633 succinimidyl ester; nanoparticles; ICC; FC)

Gielen, P. R., et al. [Connexin43 confers Temozolomide resistance in human glioma cells by modulating the mitochondrial apoptosis pathway](#). Neuropharmacology, (2013), DOI: DOI:10.1016/j.neuropharm.2013.05.002 (CF405M Annexin V; FC)

Giuntini, S., et al. [Combined roles of human IgG subclass, alternative complement pathway activation, and epitope density in the bactericidal activity of antibodies to meningococcal factor h binding protein](#). Infect Immun 80(1), 187-194, (2012), DOI: DOI:10.1128/IAI.05956-11 (CF488A goat anti-human IgG; FC)

Hwang, J., et al. [Aptamer-conjugated live human immune cell based biosensors for the accurate detection of C-reactive protein](#). Sci Rep 6, 34778, (2016), DOI: 10.1038/srep34778 (CF594 Mix-n-Stain)

Kesik, M., et al. [A multi-functional fluorescent scaffold as a multi-colour probe: design and application in targeted cell imaging](#). RSC Advances 5(101), 83361-83367, (2015), DOI: 10.1039/c5ra16600d (CF555 Mix-n-Stain)

Li, X., et al. [Tissue distribution and safety evaluation of a claudin-targeting molecule, the C-terminal fragment of Clostridium perfringens enterotoxin](#). Eur J Pharm Sci 52, 132-137, (2014), DOI: S0928-0987(13)00425-9 [pii] 10.1016/j.ejps.2013.10.018 (Xenolight CF750; FC, in vivo imaging)

Mahassni, S. H. and Al-Reemi, R. M. [Apoptosis and necrosis of human breast cancer cells by an aqueous extract of garden cress \(*Lepidium sativum*\) seeds](#). Saudi Journal of Biological Sciences 20(2), 131-139, (2013), DOI: DOI: http://dx.doi.org/10.1016/j.sjbs.2012.12.002 (CF488A Annexin V; ICC, FC)

McCausland, M., et al. [With great power comes great responsibility: high-dimensional spectral flow cytometry to support clinical trials](#). Bioanalysis 13, 1597-1616, (2021), DOI: 10.4155/bio-2021-0201 (Spectral flow)

Mikosik, A., et al. [Expression of calpain-calpastatin system \(CCS\) member proteins in human lymphocytes of young and elderly individuals: pilot baseline data for the CALPACENT project](#). Immun Ageing 10(1), 27, (2013), DOI: DOI:10.1186/1742-4933-10-27 (Mix-n-Stain CF555; FC)

Pena-Hernandez, R., et al. [Genome-wide targeting of the epigenetic regulatory protein CTCF to gene promoters by the transcription factor TFIID](#). Proc Natl Acad Sci U S A 112(7), E677-686, (2015), DOI: 10.1073/pnas.1416674112 (CF633 Annexin V flow)

Ruger, M., et al. [Species-specific viability analysis of *Pseudomonas aeruginosa*, *Burkholderia cepacia* and *Staphylococcus aureus* in mixed culture by flow cytometry](#). BMC Microbiol 14, 56, (2014), DOI: 1471-2180-14-56 [pii] 10.1186/1471-2180-14-56 (CF405S Mix-n-Stain; FC)

Ruger, M., et al. [A flow cytometric method for viability assessment of *Staphylococcus aureus* and *Burkholderia cepacia* in mixed culture](#). Cytometry A 81(12), 1055, (2012), DOI: DOI:10.1002/cyto.a.22219 (CF405S Mix-n-Stain labeling of wheat germ agglutinin; FC)

Sainski, A. M., et al. [Casp8p41 generated by HIV protease kills CD4 T cells through direct Bak activation](#). J Cell Biol 207(1), 159, (2014), DOI: jcb.20140505109252014c [pii] 10.1083/jcb.20140505109252014c (CF640R Mix-n-Stain; FC)

Skindersoe, M. E., et al. [A novel and rapid apoptosis assay based on thiol redox status](#). Cytometry A 81(5), 430, (2012), DOI: DOI:10.1002/cyto.a.22032 (CF647 Annexin V; ICC, FC)

Sutherland, D. R., et al. [High-sensitivity 5-, 6-, and 7-color PNH WBC assays for both Canto II and Navios platforms](#). Cytometry B Clin Cytom 94(4), 637-651, (2018), DOI: 10.1002/cyto.b.21626 (CF405M succinimidyl ester flow)

Vlahava, V. M., et al. [CD40 ligand exhibits a direct antiviral effect on Herpes Simplex Virus type-1 infection via a PI3K-dependent, autophagy-independent mechanism](#). Cell Signal 27(6), 1253-1263, (2015), DOI: S0898-6568(15)00078-9 [pii] 10.1016/j.cellsig.2015.03.002 (CF633 FC)

Wang, H. Y., et al. [Synthesis and single-molecule imaging reveal stereospecific enhancement of binding kinetics by the antitumour eEF1A antagonist SR-A3](#). Nat Chem 14, 1443-1450, (2022), DOI: 10.1038/s41557-022-01039-3 (CF647 azide, flow cytometry)

Zeiderman, M. R., et al. [Targeting of BRAF resistant melanoma via extracellular matrix metalloproteinase inducer receptor](#). J Surg Res 190(1), 111-118, (2014), DOI: S0022-4804(14)00165-6 [pii] 10.1016/j.jss.2014.02.021 (CF750 succinimidyl ester; in vivo imaging, FC)

Western Blotting

Ait-Belgnaoui, A., et al. [Probiotic gut effect prevents the chronic psychological stress-induced brain activity abnormality in mice](#). Neurogastroenterol Motil 26(4), 510-520, (2014), DOI: 10.1111/nmo.12295 (CF770; WB; LI-COR Odyssey)

Andriamihaja, M., et al. [The deleterious metabolic and genotoxic effects of the bacterial metabolite p-cresol on colonic epithelial cells](#). Free Radical Biology and Medicine 85, 219-227, (2015), DOI: http://dx.doi.org/10.1016/j.freeradbiomed.2015.04.004 (CF770 RedDot1 in cell western)

Annahazi, A., et al. [Luminal Cysteine-Proteases Degrade Colonic Tight Junction Structure and Are Responsible for Abdominal Pain in Constipation-Predominant IBS](#). Am J Gastroenterol 108(8), 1392, (2013), DOI: DOI:10.1038/ajg.2013.152 (CF770 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Audebert, M., et al. [Comparative potency approach based on H2AX assay for estimating the genotoxicity of polycyclic aromatic hydrocarbons](#). Toxicol Appl Pharmacol 260(1), 58, (2012), DOI: DOI:10.1016/j.taap.2012.01.022 (CF770 goat anti-rabbit IgG; ICW; LI-COR Odyssey)

Avelar, G. M., et al. [A rhodopsin-guanylyl cyclase gene fusion functions in visual perception in a fungus](#). Curr Biol 24(11), 1234-1240, (2014), DOI: S0960-9822(14)00409-6 [pii] 10.1016/j.cub.2014.04.009 (CF680; WB; LI-COR Odyssey)

Babelova, A., et al. [Activation of Rac-1 and RhoA contributes to podocyte injury in chronic kidney disease](#). PLoS One 8(11), e80328, (2013), DOI: 10.1371/journal.pone.0080328 PONE-D-13-15419 [pii] (CF770 goat anti-rabbit; WB; LI-COR Odyssey)

Dabek, M., et al. [Intracolonic infusion of fecal supernatants from ulcerative colitis patients triggers altered permeability and inflammation in mice: role of cathepsin G and protease-activated receptor-4](#). Inflamm Bowel Dis 17(6), 1409-1414, (2011), DOI: DOI:10.1002/ibd.21454 (CF770 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Ducheix, S., et al. [Dietary oleic acid regulates hepatic lipogenesis through a liver X receptor-dependent signaling](#). PLoS One 12(7), e0181393, (2017), DOI: 10.1371/journal.pone.0181393 (CF680 & CF770 secondaries WB)

Gauthier, T., et al. [Deoxynivalenol impairs the immune functions of neutrophils](#). Mol Nutr Food Res 57(6), 1026, (2013), DOI: DOI:10.1002/mnfr.201200755 (CF680 goat anti-rabbit IgG, CF770 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Gon, S., et al. [Fit alphabeta T-cell receptor suppresses leukemogenesis of Pten-deficient thymocytes](#). Haematologica 103(6), 999-1007, (2018), DOI: 10.3324/haematol.2018.188359 (CF680 goat anti-mouse and CF770 goat anti-rabbit WB)

Graber, T. E., et al. [UPF1 Governs Synaptic Plasticity through Association with a STAU2 RNA Granule](#). J Neurosci 37(38), 9116-9131, (2017), DOI: 10.1523/JNEUROSCI.0088-17.2017 (CF680 & CF770 secondaries WB)

Hilbers, F. S., et al. [Functional Analysis of Missense Variants in the Putative Breast Cancer Susceptibility Gene XRCC2](#). Hum Mutat 37(9), 914-925, (2016), DOI: 10.1002/humu.23019 (CF680 goat anti-rabbit & CF770 goat anti-mouse WB)

Huc, L., et al. [Low concentrations of bisphenol A induce lipid accumulation mediated by the production of reactive oxygen species in the mitochondria of HepG2 cells](#). Toxicol In Vitro 26(5), 709-717, (2012), DOI: DOI:10.1016/j.tiv.2012.03.017 (CF770 goat anti-mouse IgG; WB; LI-COR Odyssey)

Jamin, E. L., et al. [Combined Genotoxic Effects of a Polycyclic Aromatic Hydrocarbon \(B\(a\)P\) and an Heterocyclic Amine \(PhIP\) in Relation to Colorectal Carcinogenesis](#). PLoS One 8(3), e58591, (2013), DOI: DOI:10.1371/journal.pone.0058591 (CF770 goat anti-rabbit IgG; ICW; LI-COR Odyssey)

Kohler, C., et al. [Extracytoplasmic function \(ECF\) sigma factor sigmaF is involved in Caulobacter crescentus response to heavy metal stress](#). BMC Microbiol 12, 210, (2012), DOI: DOI:10.1186/1471-2180-12-210 (CF680 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Lai, S. R., et al. [In vitro anti-tubulin effects of mebendazole and fenbendazole on canine glioma cells](#). Vet Comp Oncol 15(4), 1445-1454, (2017), DOI: 10.1111/vco.12288 (CF770 goat anti-mouse WB)

Leskoske, K. [Modulation of TOR Complex 2 signaling and maintenance of plasma membrane homeostasis in Saccharomyces cerevisiae](#). in: Graduate Division, PhD in Molecular and Cell Biology, UC Berkeley, 106, (2017), (CF770 goat anti-mouse WB)

Liang, Y., et al. [The lymphoid lineage-specific actin-uncapping protein Rltp is essential for costimulation via CD28 and the development of regulatory T cells](#). Nat Immunol 14(8), 858-866, (2013), DOI: ni.2634 [pii] 10.1038/ni.2634 (CF770; WB)

Lucioli, J., et al. [The food contaminant deoxynivalenol activates the mitogen activated protein kinases in the intestine: Interest of ex vivo models as an alternative to in vivo experiments](#). Toxicon 66, 31, (2013), DOI: DOI:10.1016/j.toxicon.2013.01.024 (CF770 goat anti-mouse IgG, CF770 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Martinez Marshall, M. N., et al. [Analysis of the roles of phosphatidylinositol-4,5-bisphosphate and individual subunits in assembly, localization, and function of *Saccharomyces cerevisiae* target of rapamycin complex 2](#). Mol Biol Cell 30(12), 1555-1574, (2019), DOI: 10.1091/mbc.E18-10-0682 (CF770 goat anti-mouse WB)

Martinsen, A., et al. [Rho kinase regulation of vasopressin-induced calcium entry in vascular smooth muscle cell: Comparison between rat isolated aorta and cultured aortic cells](#). Cell Calcium 52(6), 413, (2012), DOI: DOI:10.1016/j.ceca.2012.07.002 (CF770 goat anti-mouse IgG, CF770 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Moussa, L., et al. [A new soy germ fermented ingredient displays estrogenic and protease inhibitor activities able to prevent irritable bowel syndrome-like symptoms in stressed female rats](#). Clin Nutr 32(1), 51, (2012), DOI: DOI:10.1016/j.clnu.2012.05.021 (CF770 anti-rabbit IgG; WB; LI-COR Odyssey)

Najimi, M., et al. [Human liver mesenchymal stem/progenitor cells inhibit hepatic stellate cell activation: in vitro and in vivo evaluation](#). Stem Cell Res Ther 8(1), 131, (2017), DOI: 10.1186/s13287-017-0575-5 (CF680 donkey anti-rabbit & CF790 donkey anti-mouse WB)

Nguyen, T. M. D., et al. [Inhibition by fluoxetine of LH-stimulated cyclic AMP synthesis in tumor Leydig cells partly involves AMPK activation](#). PLoS One 14(6), e0217519, (2019), DOI: 10.1371/journal.pone.0217519 (CF770 anti-rabbit WB)

Oliveira, A. F. and Yasuda, R. [Neurofibromin is the major ras inactivator in dendritic spines](#). J Neurosci 34(3), 776-783, (2014), DOI: 34/3/776 [pii] 10.1523/JNEUROSCI.3096-13.2014 (CF680 goat anti-mouse, CF770 goat anti-rabbit; WB; LI-COR Odyssey)

Olivier, I., et al. [Is Crohn's creeping fat an adipose tissue?](#). Inflamm Bowel Dis 17(3), 747-757, (2011), DOI: DOI:10.1002/ibd.21413 (CF770 goat anti-rabbit IgG; WB; LI-COR Odyssey)

Ottenheijm, C. A., et al. [Deleting exon 55 from the nebulin gene induces severe muscle weakness in a mouse model for nemaline myopathy](#). Brain 136(Pt 6), 1718-1731, (2013), DOI: awt113 [pii] 10.1093/brain/awt113 (CF680 goat anti-rabbit; WB; LI-COR Odyssey)

Pons, B. J., et al. [Cell transfection of purified cytolethal distending toxin B subunits allows comparing their nuclease activity while plasmid degradation assay does not](#). PLoS One 14(3), e0214313, (2019), DOI: 10.1371/journal.pone.0214313 (CF770 goat anti-mouse WB)

Rother, M. B., et al. [The Human Thymus Is Enriched for Autoreactive B Cells](#). J Immunol 197(2), 441-448, (2016), DOI: 10.4049/jimmunol.1501992 (CF680 goat anti-human & CF770 goat anti-mouse WB)

Schimmel, J., et al. [Mutational signatures of non-homologous and polymerase theta-mediated end-joining in embryonic stem cells](#). EMBO J 36(24), 3634-3649, (2017), DOI: 10.15252/embj.201796948 (CF680 goat anti-rabbit & CF770 goat anti-mouse WB)

Schmidt, C., et al. [The Dual Role of a Polyvalent IgM/IgA-Enriched Immunoglobulin Preparation in Activating and Inhibiting the Complement System](#). Biomedicines 9(7), (2021), DOI: 10.3390/biomedicines9070817 (CF488A DAGt CF568 donkey anti-mouse, western blotting)

Titova, E. and Obenaus, A. [Immunological Response to Experimental Intracerebral Hemorrhage: Morphological Assessments](#). In: Chen, J., et al. (Eds.), Animal Models of Acute Neurological Injuries II. Humana Press, pp. 625-652, (2012), DOI: DOI:10.1007/978-1-61779-576-3_48 (CF680- and CF770- secondary antibody conjugates; IHC in floating brain sections; LI-COR Odyssey)

Titova, E. M., et al. [The late phase of post-stroke neurorepair in aged rats is reflected by MRI-based measures](#). Neuroscience, (2014), DOI: S0306-4522(14)00770-2 [pii] 10.1016/j.neuroscience.2014.09.028 (CF680 goat anti-mouse & CF770 goat anti-rat; IHC; LI-COR Odyssey)

van den Heuvel, D., et al. [A CSB-PAF1C axis restores processive transcription elongation after DNA damage repair](#). Nat Commun 12(1), 1342, (2021), DOI: 10.1038/s41467-021-21520-w (CF680 goat anti-rabbit CF770 goat anti-mouse WB Odyssey)

van der Weegen, Y., et al. [The sequential and cooperative action of CSB, CSA, and UVSSA targets the TFIIH complex to DNA damage-stalled RNA polymerase II](#). bioRxiv, (2019), DOI: 10.1101/707216 (CF680 donkey anti-goat CF680 goat anti-rabbit CF770 goat anti-mouse WB)

Wu, C. L., et al. [Interplay between Cell Migration and Neurite Outgrowth Determines SH2B1beta-Enhanced Neurite Regeneration of Differentiated PC12 Cells](#). PLoS One 7(4), e34999, (2012), DOI: DOI:10.1371/journal.pone.0034999 (CF680 goat anti-mouse IgG; WB; LI-COR Odyssey)

In-Cell Western®

Andriamihaja, M., et al. [The deleterious metabolic and genotoxic effects of the bacterial metabolite p-cresol on colonic epithelial cells](#). Free Radical Biology and Medicine 85, 219-227, (2015), DOI: http://dx.doi.org/10.1016/j.freeradbiomed.2015.04.004 (CF770 RedDot1 in cell western)

Audebert, M., et al. [Comparative potency approach based on H2AX assay for estimating the genotoxicity of polycyclic aromatic hydrocarbons](#). Toxicol Appl Pharmacol 260(1), 58, (2012), DOI: DOI:10.1016/j.taap.2012.01.022 (CF770 goat anti-rabbit IgG; ICW; LI-COR Odyssey)

Chevereau, M., et al. [Role of human sulfotransferase 1A1 and N-acetyltransferase 2 in the metabolic activation of 16 heterocyclic amines and related heterocyclics to genotoxins in recombinant V79 cells](#). Arch Toxicol 91(9), 3175-3184, (2017), DOI: 10.1007/s00204-017-1935-8 (CF770 secondary RedDot ICW)

Jamin, E. L., et al. [Combined Genotoxic Effects of a Polycyclic Aromatic Hydrocarbon \(B\(a\)P\) and an Heterocyclic Amine \(PhIP\) in Relation to Colorectal Carcinogenesis](#). PLoS One 8(3), e58591, (2013), DOI: DOI:10.1371/journal.pone.0058591 (CF770 goat anti-rabbit IgG; ICW; LI-COR Odyssey)

Khoury, L., et al. [Validation of high-throughput genotoxicity assay screening using gammaH2AX in-cell western assay on HepG2 cells](#). Environ Mol Mutagen 54(9), 737-746, (2013), DOI: 10.1002/em.21817 (CF770; ICW)

Khoury, L., et al. [Evaluation of four human cell lines with distinct biotransformation properties for genotoxic screening](#). Mutagenesis 31(1), 83-96, (2016), DOI: 10.1093/mutage/gev058 (CF770 goat anti-rabbit, RedDot2, ICW)

Miyagishima, S. Y., et al. [DipM is required for peptidoglycan hydrolysis during chloroplast division](#). BMC Plant Biol 14, 57, (2014), DOI: 1471-2229-14-57 [pii] 10.1186/1471-2229-14-57 (CF488A Mix-n-Stain, CF555 Mix-n-Stain; ICC)

Norris, J. D., et al. [Androgen receptor antagonism drives cytochrome P450 17A1 inhibitor efficacy in prostate cancer](#). J Clin Invest 127(6), 2326-2338, (2017), DOI: 10.1172/JCI87328 (CF770 goat anti-rabbit In Cell Western)

Pollock, J. A., et al. [Inhibiting androgen receptor nuclear entry in castration-resistant prostate cancer](#). Nat Chem Biol 12(10), 795-801, (2016), DOI: 10.1038/nchembio.2131 (CF770 goat anti-rabbit In Cell Western)

Theumer, M. G., et al. [Genotoxicity of aflatoxins and their precursors in human cells](#). Toxicol Lett 287, 100-107, (2018), DOI: 10.1016/j.toxlet.2018.02.007 (CF770 & RedDot ICW)

In Vivo Imaging

Abe, S., et al. [A novel targeting therapy of malignant mesothelioma using anti-podoplanin antibody](#). J Immunol 190(12), 6239-6249, (2013), DOI: DOI:10.4049/jimmunol.1300448 (XenoLight CF770; in vivo imaging; PerkinElmer IVIS Spectrum imaging system)

Alawieh, A., et al. [Modulation of post-stroke degenerative and regenerative processes and subacute protection by site-targeted inhibition of the alternative pathway of complement](#). Journal of Neuroinflammation 12(1), 1-15, (2015), DOI: 10.1186/s12974-015-0464-8 (CF750 succinimidyl ester in vivo)

Alfonso-Loeches, S., et al. [Toll-like receptor 4 participates in the myelin disruptions associated with chronic alcohol abuse](#). Glia 60(6), 948-964, (2012), DOI: DOI:10.1002/glia.22327 (XenoLight CF770; in vivo imaging (mouse brain); PerkinElmer IVIS 200 imaging system)

Alfonso-Loeches, S., et al. [Gender differences in alcohol-induced neurotoxicity and brain damage](#). Toxicology 311(1-2), 27, (2013), DOI: DOI:10.1016/j.tox.2013.03.001 (XenoLight CF770; in vivo imaging (mouse brain); PerkinElmer IVIS 200 imaging system)

Alfonso-Loeches, S., et al. [Ethanol-Induced TLR4/NLRP3 Neuroinflammatory Response in Microglial Cells Promotes Leukocyte Infiltration Across the BBB](#). Neurochem Res 41(1-2), 193-209, (2016), DOI: 10.1007/s11064-015-1760-5 (CF770 neg con in vivo imaging)

Azhdarinia, A., et al. [Dual-labeling strategies for nuclear and fluorescence molecular imaging: a review and analysis](#). Mol Imaging Biol 14(3), 261-276, (2012), DOI: 10.1007/s11307-011-0528-9 (CF770, CF790; in vivo imaging)

Cai, H., et al. [Positron Emission Tomography Imaging of Platelet-Derived Growth Factor Receptor beta in Colorectal Tumor Xenograft Using Zirconium-89 Labeled Dimeric Affibody Molecule](#). Mol Pharm 16(5), 1950-1957, (2019), DOI: 10.1021/acs.molpharmaceut.8b01317 (CF750 succinimidyl ester in vivo)

Chen, X. and Zhang, W. [Diamond nanostructures for drug delivery, bioimaging, and biosensing](#). Chem Soc Rev 46(3), 734-760, (2017), DOI: 10.1039/c6cs00109b (CF750 succinimidyl ester in vivo)

Cho, H., et al. [Study of osteoarthritis treatment with anti-inflammatory drugs: cyclooxygenase-2 inhibitor and steroids](#). Biomed Res Int 2015, 595273, (2015), DOI: 10.1155/2015/595273 (CF680 succinimidyl ester in vivo)

Darniot, M., et al. [RNA interference in vitro and in vivo using DsiRNA targeting the nucleocapsid N mRNA of human metapneumovirus](#). Antiviral Res 93(3), 364-373, (2012), DOI: DOI:10.1016/j.antiviral.2012.01.004 (CF750 succinimidyl ester (to label siRNA); in vivo imaging; Visen Medical FMT 2500 imaging system)

Davies, B. S., et al. [Assessing mechanisms of GPIHBP1 and lipoprotein lipase movement across endothelial cells](#). J Lipid Res 53(12), 2690, (2012), DOI: DOI:10.1194/jlr.M031559 (CF568 succinimidyl ester; in vivo injection, IHC)

Fan, Q., et al. [Biological evaluation of 131I- and CF750-labeled Dmab\(scFv\)-Fc antibodies for xenograft imaging of CD25-positive tumors](#). Biomed Res Int 2014, 459676, (2014), DOI: 10.1155/2014/459676 (CF750; in vivo imaging; Caliper IVIS)

Guo, K., et al. [Engineering the first chimeric antibody in targeting intracellular PRL-3 oncprotein for cancer therapy in mice](#). Oncotarget 3(2), 158-171, (2012), (XenoLight CF750 NIR Fluorescent Labeling Kit; in vivo imaging; PerkinElmer IVIS Spectrum imaging system)

Hapuarachchige, S., et al. [Bioorthogonal two-component drug delivery in HER2\(+\) breast cancer mouse models](#). Sci Rep 6, 24298, (2016), DOI: 10.1038/srep24298 (CF680 succinimidyl ester & CF750 succinimidyl ester, in vivo)

Herz, C., et al. [The isothiocyanate erucin abrogates telomerase in hepatocellular carcinoma cells in vitro and in an orthotopic xenograft tumour model of HCC](#). J Cell Mol Med, (2014), DOI: 10.1111/jcmm.12412 (Xenolight CF750; in vivo imaging; Bruker Image Station)

Ito, A., et al. [New whole-body multimodality imaging of gastric cancer peritoneal metastasis combining fluorescence imaging with ICG-labeled antibody and MRI in mice](#). Gastric Cancer 17(3), 497-507, (2014), DOI: 10.1007/s10120-013-0316-0 (Xenolight CF750; in vivo imaging; IVIS Lumina II)

Kim, J. Y., et al. [Efficacy of gastro-retentive forms of ecabet sodium in the treatment of gastric ulcer in rats](#). Arch Pharm Res 37(8), 1053-1062, (2014), DOI: 10.1007/s12272-013-0278-0 (CF770; in vivo imaging)

Lee, S., et al. [Localization of near-infrared labeled antibodies to the central nervous system in experimental autoimmune encephalomyelitis](#). PLoS One 14(2), e0212357, (2019), DOI: 10.1371/journal.pone.0212357 (CF680 in vivo)

Li, X., et al. [Development of an anti-claudin-3 and -4 bispecific monoclonal antibody for cancer diagnosis and therapy](#). J Pharmacol Exp Ther 351(1), 206-213, (2014), DOI: jpet.114.216911 [pii] 10.1124/jpet.114.216911 (Xenolight CF750; in vivo imaging)

Li, X., et al. [Tissue distribution and safety evaluation of a claudin-targeting molecule, the C-terminal fragment of Clostridium perfringens enterotoxin](#). Eur J Pharm Sci 52, 132-137, (2014), DOI: S0928-0987(13)00425-9 [pii] 10.1016/j.ejps.2013.10.018 (Xenolight CF750; FC, in vivo imaging)

Ma, X., et al. [Histone chaperone CAF-1 promotes HIV-1 latency by leading the formation of phase-separated suppressive nuclear bodies](#). EMBO J 40(10), e106632, (2021), DOI: 10.15252/embj.2020106632 (CF568 donkey anti-rabbit CF568 donkey anti-mouse STORM)

Matsuo, H., et al. [Engineered hepatitis B virus surface antigen L protein particles for in vivo active targeting of splenic dendritic cells](#). Int J Nanomedicine 7, 3341-3350, (2012), DOI: DOI:10.2147/IJN.S32813 (CF750 succinimidyl ester; bionanocapsules (BNCs); in vivo imaging; Olympus OV-100 imaging system)

Moore, L., et al. [Diamond-Lipid Hybrids Enhance Chemotherapeutic Tolerance and Mediate Tumor Regression](#). Adv Mater 25(26), 3532, (2013), DOI: DOI:10.1002/adma.201300343 (Xenolight CF750 in Nanodiamond-lipid hybrid particles (NDLPs); in vivo imaging; PerkinElmer IVIS Spectrum imaging system)

Stasevich, T. J., et al. [Quantifying histone and RNA polymerase II post-translational modification dynamics in mother and daughter cells](#). Methods 70(2-3), 77, (2014), DOI: S1046-2023(14)00260-6 [pii] 10.1016/jymeth.2014.08.002 (CF640R succinimidyl ester; live cell imaging)

Suemizu, H., et al. [A versatile technique for the in vivo imaging of human tumor xenografts using near-infrared fluorochrome-conjugated macromolecule probes](#). PLoS One 8(12), e82708, (2013), DOI: 10.1371/journal.pone.0082708 PONE-D-13-13248 [pii] (Xenolight CF770; in vivo imaging)

Sun, X., et al. [Longitudinal monitoring of tumor antiangiogenic therapy with near-infrared fluorophore-labeled agents targeted to integrin alphavbeta3 and vascular endothelial growth factor](#). Eur J Nucl Med Mol Imaging 41(7), 1428-1439, (2014), DOI: 10.1007/s00259-014-2702-1 (Xenolight CF680 CF770; in vivo imaging)

Sun, Y., et al. [Single tumor imaging with multiple antibodies targeting different antigens](#). Biotechniques 52(4), 1-3, (2012), DOI: DOI:10.2144/000113855 (XenoLight CF680, XenoLight CF770; in vivo imaging; PerkinElmer IVIS Lumina)

Takeyama, H., et al. [A novel endoscopic fluorescent clip visible with near-infrared imaging during laparoscopic surgery in a porcine model](#). Surg Endosc 28(6), 1984-1990, (2014), DOI: 10.1007/s00464-014-3423-z (CF790; in vivo imaging)

Wei, D., et al. [CF750-A33scFv-Fc-Based Optical Imaging of Subcutaneous and Orthotopic Xenografts of GPA33-Positive Colorectal Cancer in Mice](#). BioMed Research International 2015, 11, (2015), DOI: 10.1155/2015/505183 (CF750 succinimidyl ester in vivo)

Xu, N., et al. [Neonatal Pharmacokinetics and Biodistribution of Polymeric Nanoparticles and Effect of Surfactant](#). Pharmaceutics 15, (2023), DOI: 10.3390/pharmaceutics15041176 (CF647 SE, in vivo)

Zeiderman, M. R., et al. [Targeting of BRAF resistant melanoma via extracellular matrix metalloproteinase inducer receptor](#). J Surg Res 190(1), 111-118, (2014), DOI: S0022-4804(14)00165-6 [pii] 10.1016/j.jss.2014.02.021 (CF750 succinimidyl ester; in vivo imaging, FC)

Zhang, Y., et al. [One-Shot Immunomodulatory Nanodiamond Agents for Cancer Immunotherapy](#). Adv Mater 28(14), 2699-2708, (2016), DOI: 10.1002/adma.201506232 (CF750 succinimidyl ester in vivo)

Super-Resolution Microscopy

Alshafie, W., et al. [Regulated resurfacing of a somatostatin receptor storage compartment fine-tunes pituitary secretion](#). Journal of Cell Biology 219(1), (2020), DOI: 10.1083/jcb.201904054 (CF568 goat anti-rat STORM)

Angelov, B. and Angelova, A. [Nanoscale clustering of the neurotrophin receptor TrkB revealed by super-resolution STED microscopy](#). Nanoscale 9(28), 9797-9804, (2017), DOI: 10.1039/c7nr03454g (CF488A STED)

Bartok, A., et al. [IP3 receptor isoforms differently regulate ER-mitochondrial contacts and local calcium transfer](#). Nat Commun 10(1), 3726, (2019), DOI: 10.1038/s41467-019-11646-3 (CF568 goat anti-rabbit STORM)

Bennett, H. W., et al. [Novel fibrillar structure in the inversin compartment of primary cilia revealed by 3D single-molecule superresolution microscopy](#). Mol Biol Cell 31(7), 619-639, (2020), DOI: 10.1091/mbc.E19-09-0499 (CF568 STORM)

Catsburg, L. A. E., et al. [Dynamics and nanoscale organization of the postsynaptic endocytic zone at excitatory synapses](#). bioRxiv, (2021), DOI: 10.1101/2021.02.18.431766 (CF568 STED)

Civitci, F., et al. [Fast and multiplexed superresolution imaging with DNA-PAINT-ERS](#). Nat Commun 11(1), 4339, (2020), DOI: 10.1038/s41467-020-18181-6 (CF660R SMLM)

Conley, G. M., et al. [Jamming and overpacking fuzzy microgels: Deformation, interpenetration, and compression](#). Sci Adv 3(10), e1700969, (2017), DOI: 10.1126/sciadv.1700969 [pii] (CF680R, STORM)

Demmerle, J., et al. [Strategic and practical guidelines for successful structured illumination microscopy](#). Nat Protoc 12(5), 988-1010, (2017), DOI: nprot.2017.019 [pii] 10.1038/nprot.2017.019 (CF405S and CF405M SIM)

Engel, U. [Four-Channel Super-Resolution Imaging by 3-D Structured Illumination](#). Methods Mol Biol 1663, 79-94, (2017), DOI: 10.1007/978-1-4939-7265-4_7 (CF640R SIM)

Essig, K., et al. [Roquin Suppresses the PI3K-mTOR Signaling Pathway to Inhibit T Helper Cell Differentiation and Conversion of Treg to Tfr Cells](#). Immunity, (2017), DOI: 10.1016/j.immuni.2017.11.008 (CF405S, SIM)

Fruh, S. M., et al. [Molecular architecture of native fibronectin fibrils](#). Nat Commun 6, 7275, (2015), DOI: ncomms8275 [pii] 10.1038/ncomms8275 (CF680 STORM)

Gao, R., et al. [Cortical column and whole-brain imaging with molecular contrast and nanoscale resolution](#). Science 363(6424), (2019), DOI: 10.1126/science.aau8302 (CF633 donkey anti-mouse CF633 goat anti-mouse expansion microscopy)

Gaudreau-Lapierre, A., et al. [Expansion microscopy-based imaging of nuclear structures in cultured cells](#). STAR Protocols 2(3), (2021), DOI: 10.1016/j.xpro.2021.100630 (CF633 goat anti-mouse & CF633 goat anti-rabbit)

Glebov, O. O., et al. [Nanoscale Structural Plasticity of the Active Zone Matrix Modulates Presynaptic Function](#). Cell Rep 18(11), 2715-2728, (2017), DOI: S2211-1247(17)30279-6 [pii] 10.1016/j.celrep.2017.02.064 (CF680 STORM)

Gomes de Castro, M. A., et al. [Differential organization of tonic and chronic B cell antigen receptors in the plasma membrane](#). Nat Commun 10(1), 820, (2019), DOI: 10.1038/s41467-019-08677-1 (CF647 maleimide STORM)

Gong, Y.-N., et al. [Biological events and molecular signaling following MLKL activation during necroptosis](#). Cell Cycle, 1-13, (2017), DOI: 10.1080/15384101.2017.1371889 (CF568, CF647 STORM)

Gorlitz, F., et al. [A STED Microscope Designed for Routine Biomedical Applications](#). Progress In Electromagnetics Research 147, 57-68, (2014), DOI: http://dx.doi.org/10.2528/PIER14042708 (CF680R STED)

Gorur, A., et al. [COPII-coated membranes function as transport carriers of intracellular procollagen I](#). The Journal of Cell Biology 216(6), 1745-1759, (2017), DOI: 10.1083/jcb.201702135 (CF568, CF680 STORM)

Guo, X., et al. [Mitochondrial stress is relayed to the cytosol by an OMA1-DELE1-HRI pathway](#). Nature 579(7799), 427-432, (2020), DOI: 10.1038/s41586-020-2078-2 (CF568 donkey anti-rabbit CF680 donkey anti-rabbit STORM)

Gu, L., et al. [Molecular-scale axial localization by repetitive optical selective exposure](#). Nat Methods 18(4), 369-373, (2021), DOI: 10.1038/s41592-021-01099-2 (STORM CF660C donkey anti-mouse)

Haas, K. T., et al. [Multitarget Immunohistochemistry for Confocal and Super-resolution Imaging of Plant Cell Wall Polysaccharides](#). Bio Protoc 10(19), e3783, (2020), DOI: 10.21769/BioProtoc.3783 (CF568)

Haas, K. T., et al. [Single-molecule localization microscopy reveals molecular transactions during RAD51 filament assembly at cellular DNA damage sites](#). Nucleic Acids Res, (2018), DOI: 4788348 [pii] 10.1093/nar/gkx1303 (CF568, STORM)

Haselmann, H., et al. [Human Autoantibodies against the AMPA Receptor Subunit GluA2 Induce Receptor Reorganization and Memory Dysfunction](#). Neuron 100(1), 91-105 e109, (2018), DOI: 10.1016/j.neuron.2018.07.048 (CF568 goat anti-guinea pig STORM)

Hauser, M., et al. [The Spectrin-Actin-Based Periodic Cytoskeleton as a Conserved Nanoscale Scaffold and Ruler of the Neural Stem Cell Lineage](#). Cell Rep 24(6), 1512-1522, (2018), DOI: 10.1016/j.celrep.2018.07.005 (CF568 STORM)

Held, R. G., et al. [Synapse and Active Zone Assembly in the Absence of Presynaptic Ca\(2+\) Channels and Ca\(2+\) Entry](#). Neuron 107(4), 667-683 e669, (2020), DOI: 10.1016/j.neuron.2020.05.032 (CF555 donkey anti-mouse 3D STORM)

Heller, J. [Exploring Nanoscale Organisation of Synapses with Super-Resolution Microscopy](#). OM&P 3(2), 48-58, (2017), DOI: doi:10.20388/omp2017.002.0045 (CF568 STORM)

Henneberger, C., et al. [LTP Induction Boosts Glutamate Spillover by Driving Withdrawal of Perisynaptic Astroglia](#). Neuron 108(5), 919-936 e911, (2020), DOI: 10.1016/j.neuron.2020.08.030 (CF568 DAMs 3-color STORM)

Hoess, P., et al. [Dual-Color and 3D Super-Resolution Microscopy of Multi-protein Assemblies](#). Methods Mol Biol 1764, 237-251, (2018), DOI: 10.1007/978-1-4939-7759-8_14 (CF680 goat anti-rabbit 2-color STORM)

Huang, T., et al. [Simultaneous Multicolor Single-Molecule Tracking with Single-Laser Excitation via Spectral Imaging](#). Biophys J 114(2), 301-310, (2018), DOI: S0006-3495(17)31242-0 [pii] 10.1016/j.bpj.2017.11.013 (CF633, CF660C, CF680, CF680R STORM)

J. Kim and M. Wojcik, Y. W., K. Xu, X. Zhang. [Oblique-sectional single-molecule microscopy](#). IEEE Xplore, 1-2, (2018), DOI: 10.1109/IPCon.2018.8527149 (CF568 obSTORM)

Jorgens, D. M., et al. [Deep nuclear invaginations are linked to cytoskeletal filaments - integrated bioimaging of epithelial cells in 3D culture](#). J Cell Sci 130(1), 177-189, (2017), DOI: jcs.190967 [pii] 10.1242/jcs.190967 (CF568 STORM)

Karanasios, E., et al. [Autophagy initiation by ULK complex assembly on ER tubulovesicular regions marked by ATG9 vesicles](#). Nat Commun 7, 12420, (2016), DOI: ncomms12420 [pii] 10.1038/ncomms12420 (CF568 STORM)

Kim, H. J. and Selvin, P. R. [Fluorescence Imaging with One Nanometer Accuracy](#). in: SpringerReference, Springer, (2013), (CF633; Fluorescence Imaging with One Nanometer Accuracy (FIONA))

Klevanski, M., et al. [Automated highly multiplexed super-resolution imaging of protein nano-architecture in cells and tissues](#). Nat Commun 11(1), 1552, (2020), DOI: 10.1038/s41467-020-15362-1 (CF680 multi-round STORM)

Köhler, S., et al. [The interaction of crossover formation and the dynamic architecture of the synaptonemal complex during meiosis](#). bioRxiv, (2020), DOI: 10.1101/2020.02.16.947804 (CF568 succinimidyl ester STORM)

Kong, D., et al. [Prolonged mitosis results in structurally aberrant and over-elongated centrioles](#). J Cell Biol 219(6), (2020), DOI: 10.1083/jcb.201910019 (CF647 goat anti-rabbit F(ab)2 STORM)

Konig, I., et al. [Single-molecule spectroscopy of protein conformational dynamics in live eukaryotic cells](#). Nat Methods 12(8), 773-779, (2015), DOI: nmeth.3475 [pii] 10.1038/nmeth.3475 (CF680R single molecule spectroscopy)

Kraus, F., et al. [Quantitative 3D structured illumination microscopy of nuclear structures](#). Nat Protoc 12(5), 1011-1028, (2017), DOI: nprot.2017.020 [pii] 10.1038/nprot.2017.020 (CF568 azide, CF405M azide, SIM)

Lee, K. S., et al. [Functional Synaptic Architecture of Callosal Inputs in Mouse Primary Visual Cortex](#). Neuron 101(3), 421-428 e425, (2019), DOI: 10.1016/j.neuron.2018.12.005 (CF405M CF647 expansion microscopy)

Lehmann, M., et al. [Nanoscale coupling of endocytic pit growth and stability](#). Sci Adv 5(11), eaax5775, (2019), DOI: 10.1126/sciadv.aax5775 (CF647 goat anti-mouse CF680 goat anti-rabbit STORM)

Lehmann, M., et al. [Novel organic dyes for multicolor localization-based super-resolution microscopy](#). J Biophotonics 9(1-2), 161-170, (2016), DOI: 10.1002/jbio.201500119 (CF555, CF568, CF647, CF680, STORM)

Lim, A., et al. [Two kinesins drive anterograde neuropeptide transport](#). Mol Biol Cell, (2017), DOI: mbc.E16-12-0820 [pii] 10.1091/mbc.E16-12-0820 (Mix-n-Stain CF568, Direct + indirect IF, CF568 3D SIM)

Lycas, M. D., et al. [Nanoscopic dopamine transporter distribution and conformation are inversely regulated by excitatory drive and D2-autoreceptor activity](#). bioRxiv, (2021), DOI: 10.1101/2021.03.09.434538 (CF568 STORM)

Marenda, M., et al. [Parameter-free molecular super-structures quantification in single-molecule localization microscopy](#). J Cell Biol 220(5), (2021), DOI: 10.1083/jcb.202010003 (CF647 STORM)

Markaki, Y., et al. [Fluorescence In Situ Hybridization Applications for Super-Resolution 3D Structured Illumination Microscopy](#). Methods Mol Biol 950, 43-64, (2013), DOI: DOI:10.1007/978-1-62703-137-0_4 (CF405M secondary antibodies; super resolution microscopy (3D-SIM); 3D-FISH)

Ma, X., et al. [Histone chaperone CAF-1 promotes HIV-1 latency by leading the formation of phase-separated suppressive nuclear bodies](#). EMBO J 40(10), e106632, (2021), DOI: 10.15252/embj.2020106632 (CF568 donkey anti-rabbit CF568 donkey anti-mouse STORM)

Meitinger, F., et al. [TRIM37 prevents formation of condensate-organized ectopic spindle poles to ensure mitotic fidelity](#). J Cell Biol 220(7), (2021), DOI: 10.1083/jcb.202010180 (CF647 CF568)

Mezache, L., et al. [Vascular endothelial growth factor promotes atrial arrhythmias by inducing acute intercalated disk remodeling](#). Sci Rep 10(1), 20463, (2020), DOI: 10.1038/s41598-020-77562-5 (CF568 2-color STORM)

Miron, E., et al. [In Vivo and In Situ Replication Labeling Methods for Super-resolution Structured Illumination Microscopy of Chromosome Territories and Chromatin Domains](#). Methods Mol Biol 1431, 127-140, (2016), DOI: 10.1007/978-1-4939-3631-1_10 (CF405M SIM)

Nahidazar, L. and Harkes, R. [Multicolor Localization-Based Super Resolution Microscopy](#). Methods Mol Biol 2350, 69-76, (2021), DOI: 10.1007/978-1-0716-1593-5_5 (CF488A, CF680, CF660C, CF514 SMLM)

Needham, S. R., et al. [EGFR oligomerization organizes kinase-active dimers into competent signalling platforms](#). Nat Commun 7, 13307, (2016), DOI: ncomms13307 [pii] 10.1038/ncomms13307 (CF640R FLImP)

Needham, S. R., et al. [Determining the geometry of oligomers of the human epidermal growth factor family on cells with <10 nm resolution](#). Biochem Soc Trans 43(3), 309-314, (2015), DOI: BST20140318 [pii] 10.1042/BST20140318 (CF640R FLImP)

Ohgomori, T., et al. [Differential activation of neuronal and glial STAT3 in the spinal cord of the SOD1G93A mouse model of amyotrophic lateral sclerosis](#). Eur J Neurosci 46(4), 2001-2014, (2017), DOI: 10.1111/ejn.13650 (CF405 streptavidin SIM)

Olivier, N., et al. [Simple buffers for 3D STORM microscopy](#). Biomed Opt Express 4(6), 885-899, (2013), DOI: 10.1364/BOE.4.000885187153 [pii] (CF647; Super Resolution Microscopy)

Park, C. E., et al. [Super-Resolution Three-Dimensional Imaging of Actin Filaments in Cultured Cells and the Brain via Expansion Microscopy](#). ACS Nano 14(11), 14999-15010, (2020), DOI: 10.1021/acsnano.0c04915 (CF633 secondary expansion microscopy)

Platonova, E., et al. [A simple method for GFP- and RFP-based dual color single-molecule localization microscopy](#). ACS Chem Biol 10(6), 1411-1416, (2015), DOI: 10.1021/acscchembio.5b00046 (CF680 STORM)

Platonova, E., et al. [Single-molecule microscopy of molecules tagged with GFP or RFP derivatives in mammalian cells using nanobody binders](#). Methods 88, 89-97, (2015), DOI: http://dx.doi.org/10.1016/j.ymeth.2015.06.018 (CF680 STORM)

Rojas-Sánchez, L., et al. [Synthesis and intracellular tracing surface-functionalized calcium phosphate nanoparticles by super-resolution microscopy \(STORM\)](#). Materialia 12, (2020), DOI: 10.1016/j.mtla.2020.100773 (CF568 wheat germ agglutinin STORM)

Salazar-Cavazos, E., et al. [Multisite EGFR phosphorylation is regulated by adaptor protein abundances and dimer lifetimes](#). Mol Biol Cell 31(7), 695-708, (2020), DOI: 10.1091/mbc.E19-09-0548 (CF555 CF640R SIM)

Saliba, D. G., et al. [Composition and structure of synaptic ectosomes exporting antigen receptor linked to functional CD40 ligand from helper T-cells..](#) (2019), DOI: 10.1101/600551 (CF568 wheat germ agglutinin STORM)

Salvador-Gallego, R., et al. [Bax assembly into rings and arcs in apoptotic mitochondria is linked to membrane pores](#). EMBO J 35(4), 389-401, (2016), DOI: 10.15252/embj.201593384 (CF680 goat anti-mouse STORM)

Schneider, K., et al. [Dissection of cell cycle-dependent dynamics of Dnmt1 by FRAP and diffusion-coupled modeling](#). Nucleic Acids Res 41(9), 4860-4876, (2013), DOI: gkt191 [pii] 10.1093/nar/gkt191 (CF405S goat anti-mouse; FRAP)

Shrestha, R. L., et al. [Aurora-B kinase pathway controls the lateral to end-on conversion of kinetochore-microtubule attachments in human cells](#). Nat Commun 8(1), 150, (2017), DOI: 10.1038/s41467-017-00209-z [pii] (CF680 donkey anti-human STORM)

Simonson, P. D., et al. [Single-molecule-based super-resolution images in the presence of multiple fluorophores](#). Nano Lett 11(11), 5090-5096, (2011), DOI: DOI:10.1021/nl203560r (CF633 goat anti-mouse IgG; ICC, gSHRImP)

Sonoda, Y., et al. [Applications of cell resealing to reconstitute microRNA loading to extracellular vesicles](#). Sci Rep 11(1), 2900, (2021), DOI: 10.1038/s41598-021-82452-5 (CF568 Dextran 10K exosomes STORM)

Sørensen, G., et al. [Disrupted PDZ-domain binding motif of the dopamine transporter uniquely alters nanoscale distribution, dopamine homeostasis and reward motivation](#). BioRxiv, (2021), DOI: 10.1101/2021.01.06.425643 (CF568 succinimidyl ester 2-color STORM)

Turkowyd, B., et al. [From single molecules to life: microscopy at the nanoscale](#). Anal Bioanal Chem 408(25), 6885-6911, (2016), DOI: 10.1007/s00216-016-9781-8 [pii] (CF568, CF647, CF660C, CF680 super res review)

Wang, B., et al. [Transforming Rhodamine Dyes for \(d\)STORM Super-Resolution Microscopy via 1,3-Disubstituted Imidazolium Substitution](#). Angew Chem Int Ed Engl, e202113612, (2021), DOI: 10.1002/anie.202113612 (CF597R, CF583R dSTORM)

Watson, J. F., et al. [AMPA receptor anchoring at CA1 synapses is determined by N-terminal domain and TARP gamma8 interactions](#). Nat Commun 12(1), 5083, (2021), DOI: 10.1038/s41467-021-25281-4 (CF568 succinimidyl ester STORM)

Winter, F. R., et al. [Multicolour nanoscopy of fixed and living cells with a single STED beam and hyperspectral detection](#). Sci Rep 7, 46492, (2017), DOI: srep46492 [pii] 10.1038/srep46492 (CF680R STED)

Wisnovsky, S., et al. [Genome-wide CRISPR screens reveal a specific ligand for the glycan-binding immune checkpoint receptor Siglec-7](#). Proc Natl Acad Sci U S A 118(5), (2021), DOI: 10.1073/pnas.2015024118 (CF568 Mix-n-Stain STORM)

Xu, H., et al. [Molecular organization of mammalian meiotic chromosome axis revealed by expansion STORM microscopy](#). Proc Natl Acad Sci U S A 116(37), 18423-18428, (2019), DOI: 10.1073/pnas.1902440116 (CF568 ExSTORM)

Zanetti-Domingues, L. C., et al. [Determining the geometry of oligomers of the human epidermal growth factor family on cells with 7 nm resolution](#). Prog Biophys Mol Biol 118(3), 139-152, (2015), DOI: S0079-6107(15)00047-4 [pii] 10.1016/j.pbiomolbio.2015.04.002 (CF640R FLImP)

Zhang, M., et al. [Translocation of interleukin-1beta into a vesicle intermediate in autophagy-mediated secretion](#). eLife 4, (2015), DOI: 10.7554/eLife.11205 (CF568 STORM)

Zhang, R., et al. [The mechanisms of dynamin-actin interaction](#). bioRxiv, (2019), DOI: 10.1101/586461 (CF405M goat anti-rabbit, STED)

Zhang, Y., et al. [Multicolor super-resolution imaging using spectroscopic single-molecule localization microscopy with optimal spectral dispersion](#). Appl Opt 58(9), 2248-2255, (2019), DOI: 10.1364/AO.58.002248 (CF660C CF680 sSMLM)

Zhang, Z., et al. [Ultrahigh-throughput single-molecule spectroscopy and spectrally resolved super-resolution microscopy](#). Nat Methods 12(10), 935-938, (2015), DOI: nmeth.3528 [pii] 10.1038/nmeth.3528 (CF660C, CF680 STORM)

TIRF Microscopy

Akin, E. J., et al. [Single-Molecule Imaging of Nav1.6 on the Surface of Hippocampal Neurons Reveals Somatic Nanoclusters](#). Biophys J 111(6), 1235-1247, (2016), DOI: 10.1016/j.bpj.2016.08.016 (CF640R in TIRF)

Almahayni, K., et al. [Simple, Economic, and Robust Rail-Based Setup for Super-Resolution Localization Microscopy](#). J Phys Chem A 127, 4553-4560, (2023), DOI: 10.1021/acs.jpca.3c01351 (CF568, super-resolution, SMLM, TIRF)

Blomeke, L., et al. [Quantitative detection of alpha-Synuclein and Tau oligomers and other aggregates by digital single particle counting](#). NPJ Parkinsons Dis 8, 68, (2022), DOI: 10.1038/s41531-022-00330-x (CF488A SE, CF633 SE, TIRF)

Bosch, P. et al. (2014). [Evaluation of fluorophores to label SNAP-tag fused proteins for multicolor single-molecule tracking microscopy in live cells](#). Biophys J 107, 803-814. 10.1016/j.bpj.2014.06.040 (CF633 succinimidyl ester, CF640R succinimidyl ester; TIRF)

Cespedes, P. F., et al. [T-cell trans-synaptic vesicles are distinct and carry greater effector content than constitutive extracellular vesicles](#). Nat Commun 13, 3460, (2022), DOI: 10.1038/s41467-022-31160-3 (CF568, WGA, fluorescence microscopy, TIRF)

Engdahl, A. K., et al. [Large Field-of-View Super-Resolution Optical Microscopy Based on Planar Polymer Waveguides](#). ACS Photonics 8(7), 1944-1950, (2021), DOI: 10.1021/acsphotonics.1c00914 (CF568 phalloidin TIRF)

Garcia, J. D., et al. [Stepwise disassembly of GABAergic synapses during pathogenic excitotoxicity](#). Cell Rep 37, 110142, (2021), DOI: 10.1016/j.celrep.2021.110142 (CF568 GAMs, TIRF)

Markaki, Y., et al. [Xist nucleates local protein gradients to propagate silencing across the X chromosome](#). Cell 184, 6174-6192 e6132, (2021), DOI: 10.1016/j.cell.2021.10.022 (CF568 DARb, CF568 azide, TIRF)

Martin-Fernandez, M. et al. (2013). [A 'pocket guide' to total internal reflection fluorescence](#). J Microsc 252, 16-22. 10.1111/jmi.12070 (CF640R; TIRF)

Shirasaki, Y., et al. (2014). [Real-time single-cell imaging of protein secretion](#). Sci Rep 4, 4736. 10.1038/srep04736 (CF660R; TIRF)

Sorensen, G., et al. [Disruption of the PDZ domain-binding motif of the dopamine transporter uniquely alters nanoscale distribution, dopamine homeostasis, and reward motivation](#). J Biol Chem 297, 101361, (2021), DOI: 10.1016/j.jbc.2021.101361 (CF568SE, TIRF)

Willems, J., and MacGillavry, H. D. [A coordinate-based co-localization index to quantify and visualize spatial associations in single-molecule localization microscopy](#). Sci Rep 12, 4676, (2022), DOI: 10.1038/s41598-022-08746-4 (CF568 secondaries, TIRF)

Yu, Y., et al. [Single-phagosome imaging reveals that homotypic fusion impairs phagosome degradative function](#). Biophys J 121, 459-469, (2022), DOI: 10.1016/j.bpj.2021.12.032 (CF488A CTB, CF640R amine, TIRF)

Zanetti-Domingues, L. C., et al. [Hydrophobic fluorescent probes introduce artifacts into single molecule tracking experiments due to non-specific binding](#). PLoS One 8(9), e74200, (2013), DOI: 10.1371/journal.pone.0074200, PONE-D-13-17594 [pii] (CF488A, CF568, CF633 & CF640R maleimide, TIRF)

Zhang, R., et al. [Structural insight into TPX2-stimulated microtubule assembly](#). eLife 6, (2017), DOI: 10.7554/eLife.30959 (CF640R TIRFM)

Fluorescence Polarization Assay

Antolín-Urbaneja, J. et al. (2013). [Compact and cost effective instrument for detecting drug precursors in different environments based on fluorescence polarization](#). Paper presented at: Optical Sensors 2013 (Prague, Czech Republic, Proc SPIE). DOI:10.1117/12.2017257 (CF488A succinimidyl ester; fluorescence polarization assay)

Di Giovanni, S., Varriale, et al. (2012). [Determination of benzyl methyl ketone - a commonly used precursor in amphetamine manufacture](#). Analytical Methods 4, 3558-3564. DOI:10.1039/c2ay25772f (CF488A succinimidyl ester; fluorescence polarization assay)

Pennacchio, A, et al. (2016). [A novel fluorescence polarization assay for determination of penicillin G in milk](#). Food Chemistry 190: 381-385. (CF647 succinimidyl ester; fluorescence polarization assay)

Varriale, A, et al. (2015). [A Fluorescence Polarization Assay To Detect Steroid Hormone Traces in Milk](#). Journal of Agricultural and Food Chemistry 63: 9159-9164. (CF647 succinimidyl ester; fluorescence polarization assay)

2-Photon Microscopy

Bazzurro, V., et al. [Involvement of GABA\(A\) receptors containing alpha\(6\) subtypes in antisecretory factor activity on rat cerebellar granule cells studied by two-photon uncaging](#). Eur J Neurosci 56, 4505-4513, (2022), DOI: 10.1111/ejn.15775 (CF640R biocytin, fluorescence microscopy, 2-photon)

Gonzalez Pisfil, M., et al. (2022). [Stimulated emission depletion microscopy with a single depletion laser using five fluorochromes and fluorescence lifetime phasor separation](#). Sci Rep 12, 14027, (2022), DOI: 10.1038/s41598-022-17825-5 (CF594 WGA, CF680R DAMs, super-resolution, confocal, FLIM, fluorescence microscopy, 2-Photon)

Kim, Y. R., et al. [Neutrophils Return to Bloodstream Through the Brain Blood Vessel After Crosstalk With Microglia During LPS-Induced Neuroinflammation](#). Front Cell Dev Biol 8, 613733, (2020), DOI: 10.3389/fcell.2020.613733 (CF405M wheat-germ agglutinin, 2-photon)

Mercier, L., et al. [In vivo imaging of skeletal muscle in mice highlights muscle defects in a model of myotubular myopathy](#). Intravital 5(1), e1168553, (2016), DOI: 10.1080/21659087.2016.1168553 (CF488A alpha-bungarotoxin, 2-photon)

Moran, I., et al. [Memory B cells are reactivated in subcapsular proliferative foci of lymph nodes](#). Nat Commun 9(1), 3372, (2018), DOI: 10.1038/s41467-018-05772-7 (CF680R 2-photon)

Wagner, M. C., et al. [Mechanism of increased clearance of glycated albumin by proximal tubule cells](#). Am J Physiol Renal Physiol 310(10), F1089-1102, (2016), DOI: 10.1152/ajprenal.00605.2015 (CF594 2-photon)