

Mix-n-Stain™ Selected References

Key: IF=Immunofluorescence Microscopy; FC=Flow Cytometry; SR=Super-Resolution Microscopy

[Click on the blue link to see journal article](#)

Ahi, Y. S., Vemula, S. V., Hassan, A. O., Costakes, G., Stauffacher, C., and Mittal, S. K. (2015). [Adenoviral L4 33K forms ring-like oligomers and stimulates ATPase activity of IVa2: implications in viral genome packaging](#). *Front Microbiol* 6, 318. 10.3389/fmicb.2015.00318 (*Mix-n-Stain™ CF®350, Mix-n-Stain™ CF®488A; IF*)

Ahi, Y. S., Vemula, S. V., and Mittal, S. K. (2013). [Adenoviral E2 IVa2 Protein Interacts with L4 33K Protein and E2 DNA-Binding Protein](#). *J Gen Virol*. DOI:10.1099/vir.0.049346-0 (*Mix-n-Stain™ CF®350, Mix-n-Stain™ CF®488A; IF*)

Anderson, M. S., Loftus, M. S., and Kedes, D. H. (2014). [Maturation and vesicle-mediated egress of primate gammaherpesvirus rhesus monkey rhadinovirus require inner tegument protein ORF52](#). *J Virol* 88, 9111-9128. 10.1128/JVI.01502-14 (*Mix n Stain™ CF®488A; IF*)

Arul, M., Roslani, A. C., Ng, C. L., and Cheah, S. H. (2013). [Culture of low passage colorectal cancer cells and demonstration of variation in selected tumour marker expression](#). *Cytotechnology*. 10.1007/s10616-013-9600-4 (*Mix n Stain™ CF®488A, Mix n Stain™ CF®594; IF*)

Bolduc, J. F., Hany, L., Barat, C., Ouellet, M., and Tremblay, M. J. (2017). [Epigenetic Metabolite Acetate Inhibits Class I/II Histone Deacetylases, Promotes Histone Acetylation, and Increases HIV-1 Integration in CD4\(+\) T Cells](#). *J Virol* 91. 10.1128/JVI.01943-16 (*Mix n Stain™ CF®405M; FC*)

Breton-Provencher, V., Bakhshetyan, K., Hardy, D., Bammann, R. R., Cavarretta, F., Snapyan, M., Cote, D., Migliore, M., and Saghatelian, A. (2016). [Principal cell activity induces spine relocation of adult-born interneurons in the olfactory bulb](#). *Nat Commun* 7, 12659. 10.1038/ncomms12659 (*Mix-n-Stain™ CF®568; IF*)

Ciapa, B., and Granon, S. (2018). [Expression of Cyclin-D1 in Astrocytes Varies During Aging](#). *Front Aging Neurosci* 10, 104. 10.3389/fnagi.2018.00104 (*Mix-n-Stain™CF®568; IF*)

Darvish-Damavandi, M., Ho, H. K., and Kang, T. S. (2016). [Towards the development of an enzyme replacement therapy for the metabolic disorder propionic acidemia](#). *Mol Genet Metab Rep* 8, 51-60. 10.1016/j.ymgmr.2016.06.009 (*Mix-n-Stain™ CF®488A; IF*)

de la Puente, P., Luderer, M. J., Federico, C., Jin, A., Gilson, R. C., Egbulefu, C., Alhallak, K., Shah, S., Muz, B., Sun, J., et al. (2018). [Enhancing proteasome-inhibitory activity and specificity of bortezomib by CD38 targeted nanoparticles in multiple myeloma](#). *J Control Release* 270, 158-176. 10.1016/j.jconrel.2017.11.045 (*Mix-n-Stain™ biotin*)

Deo, D. I., Gautrot, J. E., Sukhorukov, G. B., and Wang, W. (2014). [Biofunctionalization of PEGylated microcapsules for exclusive binding to protein substrates](#). *Biomacromolecules* 15, 2555-2562. 10.1021/bm500412d (*Mix-n-Stain™ CF®555; IF*)

Diamantopoulou, A., Raftogianni, A., Stamatakis, A., Tzanoulina, S., Oitzl, M. S., and Stylianopoulou, F. (2013). [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. DOI:10.1016/j.psyneuen.2013.02.012 (*Mix-n-Stain™ CF®488A, Mix-n-Stain™ CF®555; IF*)

Donskow-Lysoniewska, K.,Majewski, P.,Brodaczewska, K.,Jozwicka, K., and Doligalska, M. (2012).

[Heligmosmoides polygyrus fourth stages induce protection against DSS- induced colitis and change opioid expression in the intestine.](#) Parasite Immunol. DOI:10.1111/pim.12003 (*Mix-n-Stain™ CF®488A; IF*)

Ersek, B.,Molnar, V.,Balogh, A.,Matko, J.,Cope, A. P.,Buzas, E. I.,Falus, A., and Nagy, G. (2012). [CD3zeta-Chain Expression of Human T Lymphocytes Is Regulated by TNF via Src-like Adaptor Protein-Dependent Proteasomal Degradation.](#) J Immunol. DOI:10.4049/jimmunol.1102365 (*Mix-n-Stain™ CF®488A; IF*)

Farias, G. G.,Britt, D. J., and Bonifacino, J. S. (2016). [Imaging the Polarized Sorting of Proteins from the Golgi Complex in Live Neurons.](#) Methods Mol Biol 1496, 13-30. 10.1007/978-1-4939-6463-5_2 (*Mix-n-Stain™ CF®488A, Mix-n-Stain™ CF®555, Mix-n-Stain™ CF®640R; IF*)

Galletti, G.,Sung, M. S.,Vahdat, L. T.,Shah, M. A.,Santana, S. M.,Altavilla, G.,Kirby, B. J., and Giannakakou, P. (2014). [Isolation of breast cancer and gastric cancer circulating tumor cells by use of an anti HER2-based microfluidic device.](#) Lab Chip 14, 147-156. 10.1039/c3lc51039e (*Mix-n-Stain™ CF®594; IF*)

Geherin, S. A.,Gomez, D.,Glabman, R. A.,Ruthel, G.,Hamann, A., and Debes, G. F. (2016). [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, 2514-2525. 10.4049/jimmunol.1403246 (*Mix-n-Stain™ CF®488A, Mix-n-Stain™ CF®594; IF*)

Gennet, N.,Tamburini, C.,Nan, X., and Li, M. (2016). [FolR1: a novel cell surface marker for isolating midbrain dopamine neural progenitors and nascent dopamine neurons.](#) 10.1038/srep32488 (*Mix-n-Stain™ CF®640R; FC*)

Gonzalez, A.,Gaines, M.,Gallegos, L. Y.,Guevara, R., and Gomez, F. A. (2018). [Enzyme-linked immunosorbent assays \(ELISA\) based on thread, paper, and fabric.](#) Electrophoresis 39, 476-484. 10.1002/elps.201700354 (*Mix-n-Stain™ biotin*)

Gorham, JB, Williams, BA, Gidley, MJ, and Mikkelsen, D.(2016). [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. (*Mix-n-Stain™ CF®660R labeling of enzyme*)

Grzelak, C. A.,Sigglekow, N. D.,Tirnitz-Parker, J. E.,Hamson, E. J.,Warren, A.,Maneck, B.,Chen, J.,Patkunanathan, B.,Boland, J.,Cheng, R., et al. (2017). [Widespread GLI expression but limited canonical hedgehog signaling restricted to the ductular reaction in human chronic liver disease.](#) PLoS One 12, e0171480. 10.1371/journal.pone.0171480 (*Mix-n-Stain™ CF®488A; IF*)

Hirtz, M, Feng, W, Fuchs, H, and Levkin, PA. (2016). [Click-Chemistry Immobilized 3D-Infused Microarrays in Nanoporous Polymer Substrates.](#) Advanced Materials Interfaces. DOI: 10.1002/admi.201500469 (*Mix-n-Stain™ CF®488A; microarray*)

Homann, S., Hofmann, C., Gorin, A. M., Nguyen, H. C. X., Huynh, D., Hamid, P., Maithel, N., Yacoubian, V., Mu, W., Kossyvakis, A., et al. (2017). [A novel rapid and reproducible flow cytometric method for optimization of transfection efficiency in cells](#). *PLoS One* 12, e0182941. 10.1371/journal.pone.0182941 (*Mix-n-Stain™ CF®647; FC*)

Home, P., Saha, B., Ray, S., Dutta, D., Gunewardena, S., Yoo, B., Pal, A., Vivian, J. L., Larson, M., Petroff, M., et al. (2012). [Altered subcellular localization of transcription factor TEAD4 regulates first mammalian cell lineage commitment](#). *Proc Natl Acad Sci U S A* 109, 7362-7367. DOI:10.1073/pnas.1201595109 (*Mix-n-Stain™ CF®488A, Mix-n-Stain™ CF®568; IF*)

Hwang, J., Seo, Y., Jo, Y., Son, J., and Choi, J. (2016). [Aptamer-conjugated live human immune cell based biosensors for the accurate detection of C-reactive protein](#). *Sci Rep* 6, 34778. 10.1038/srep34778 (*Mix-n-Stain™ CF®594; IF*)

Iwamoto, T., Okamoto, A., Ishinaga, H., Shimizu, K., Gayle, A. A., Arai, N., Takeuchi, K., and Okuda, M. (2016). [A novel approach to predict cetuximab-induced hypersensitivity reaction: detection of drug-specific IgE on basophils](#). *Cancer Med* 5, 1004-1012. 10.1002/cam4.658 (*Mix-n-Stain™ CF®405S; FC*)

Jankowska, K. I., Williamson, E. K., Roy, N. H., Blumenthal, D., Chandra, V., Baumgart, T., and Burkhardt, J. K. (2018). [Integrins Modulate T Cell Receptor Signaling by Constraining Actin Flow at the Immunological Synapse](#). *Front Immunol* 9, 25. 10.3389/fimmu.2018.00025 (*Mix-n-Stain™ CF®488A; IF*)

Kesik, M., Demir, B., Barlas, FB, Geyik, C, Cevher, SC, Odaci Demirkol, D, Timur, S, Cirpan, A, and Toppare, L. (2015). [A multi-functional fluorescent scaffold as a multi-colour probe: design and application in targeted cell imaging](#). *RSC Advances* 5: 83361-83367. (*Mix-n-Stain™ CF®555; FC*)

Lim, A., Rechtsteiner, A., and Saxton, W. M. (2017). [Two kinesins drive anterograde neuropeptide transport](#). *Mol Biol Cell* 28, 3542-3553. 10.1091/mbc.E16-12-0820 (*Mix-n-Stain™ CF®568; IF; SR*)

Lim, A. K., Lorthongpanich, C., Chew, T. G., Tan, C. W., Shue, Y. T., Balu, S., Gounko, N., Kuramochi-Miyagawa, S., Matzuk, M. M., Chuma, S., et al. (2013). [The nuage mediates retrotransposon silencing in mouse primordial ovarian follicles](#). *Development* 140, 3819-3825. 10.1242/dev.099184 (*Mix-n-Stain™ CF®488A; IF*)

Lin, T-B, Hsieh, M-C, Lai, C-Y, Cheng, J-K, Wang, H-H, Chau, Y-P, Chen, G-D, and Peng, H-Y. (2016). [Melatonin relieves neuropathic allodynia through spinal MT2-enhanced PP2Ac and downstream HDAC4 shuttling-dependent epigenetic modification of hmgb1 transcription](#). *Journal of Pineal Research* DOI: 10.1111/jpi.12307 (*Mix-n-Stain™; IF*)

Loftus, M. S., Verville, N., and Kedes, D. H. (2017). [A Conserved Leucine Zipper Motif in Gammaherpesvirus ORF52 Is Critical for Distinct Microtubule Rearrangements](#). *J Virol* 91. 10.1128/JVI.00304-17 (*Mix-n-Stain™ CF®488A; IF*)

Mikosik, A., Foerster, J., Jasiulewicz, A., Frackowiak, J., Colonna-Romano, G., Bulati, M., Buffa, S., Martorana, A., Caruso, C., Bryl, E., et al. (2013). [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, 27. DOI:10.1186/1742-4933-10-27 (*Mix-n-Stain™ CF®555; FC*)

Miyagishima, S. Y., Kabeya, Y., Sugita, C., Sugita, M., and Fujiwara, T. (2014). [DipM is required for peptidoglycan hydrolysis during chloroplast division](#). *BMC Plant Biol* 14, 57. 10.1186/1471-2229-14-57 (Mix-n-Stain™ CF®488A, Mix-n-Stain™ CF®555; IF)

Mofarrahi, M., McClung, JM, Kontos, CD, Davis, EC, Tappuni, B, Moroz, N, Pickett, AE, Huck, L, Harel, S, Danialou, G, et al. (2015). [Angiopoietin-1 enhances skeletal muscle regeneration in mice](#). *American Journal of Physiology - Regulatory, Integrative and Comparative Physiology* 308: R576-R589 (Mix-n-Stain™ CF®568; IF)

Montesinos-Rongen, M, Purschke, FG, Brunn, A, May, C, Nordhoff, E, Marcus, K, and Deckert, M. (2015). [Primary Central Nervous System \(CNS\) Lymphoma B Cell Receptors Recognize CNS Proteins](#). *J Immunol* 195: 1312-1319. (Mix-n-Stain™ CF®488A; IF)

Okumura, R., Kurakawa, T., Nakano, T., Kayama, H., Kinoshita, M., Motooka, D., Gotoh, K., Kimura, T., Kamiyama, N., Kusu, T., et al. (2016). [Lypd8 promotes the segregation of flagellated microbiota and colonic epithelia](#). *Nature* 532, 117-121. 10.1038/nature17406 (Mix-n-Stain™ CF®633; IF)

Osteikoetxea, X, Balogh, A, Szabó-Taylor, K, Németh, A, Szabó, TG, Pálóczi, K, Sódar, B, Kittel, Á, György, B, Pállinger, É, et al. (2015). [Improved Characterization of EV Preparations Based on Protein to Lipid Ratio and Lipid Properties](#). *PLoS One* 10: e0121184 (Mix-n-Stain™ CF®488A; FC)

Park, K. S., Martelotto, L. G., Peifer, M., Sos, M. L., Karnezis, A. N., Mahjoub, M. R., Bernard, K., Conklin, J. F., Szczepny, A., Yuan, J., et al. (2011). [A crucial requirement for Hedgehog signaling in small cell lung cancer](#). *Nat Med* 17, 1504-1508. DOI:10.1038/nm.2473 (Mix-n-Stain™ CF®488A, Mix-n-Stain™ CF®568, Mix-n-Stain™ CF®647; IF)

Paz, H., Joo, E. J., Chou, C. H., Fei, F., Mayo, K. H., Abdel-Azim, H., Ghazarian, H., Groffen, J., and Heisterkamp, N. (2018). [Treatment of B-cell precursor acute lymphoblastic leukemia with the Galectin-1 inhibitor PTX008](#). *J Exp Clin Cancer Res* 37, 67. 10.1186/s13046-018-0721-7 (Mix-n-Stain™ CF®647; FC)

Petrelli, A., Samain, E., Pradeau, S., Halila, S., and Fort, S. (2016). [Efficient Conjugation of Oligosaccharides to Polymer Particles by Furan/Maleimide Diels-Alder Reaction: Application to the Capture of Carbohydrate-Binding Proteins](#). *Chembiochem*. 10.1002/cbic.201600509 (Mix-n-Stain™ CF®488A, Mix-n-Stain™ CF®568; IF)

Pons, M., Izquierdo, I., Andreu-Carbo, M., Garrido, G., Planaguma, J., Muriel, O., Del Pozo, M. A., Geli, M. I., and Aragay, A. M. (2017). [Phosphorylation of filamin A regulates chemokine receptor CCR2 recycling](#). *J Cell Sci* 130, 490-501. 10.1242/jcs.193821 (Mix-n-Stain™ CF®488A, Mix-n-Stain™ CF®568; FC)

Rohm, M., Handl, A., Konig, M., Mavoungou, C., Handrick, R., and Schindowski, K. (2016). [Data of rational process optimization for the production of a full IgG and its Fab fragment from hybridoma cells](#). *Data Brief* 8, 426-435. 10.1016/j.dib.2016.05.067 (Mix-n-Stain™ CF®647; FC)

Ruger, M., Ackermann, M., and Reichl, U. (2014). [Species-specific viability analysis of Pseudomonas aeruginosa, Burkholderia cepacia and Staphylococcus aureus in mixed culture by flow cytometry](#). *BMC Microbiol* 14, 56. 10.1186/1471-2180-14-56 (CF®405S Mix-n-Stain™; FC)

- Ruger, M., Bensch, G., Tungler, R., and Reichl, U. (2012). [A flow cytometric method for viability assessment of Staphylococcus aureus and Burkholderia cepacia in mixed culture](#). *Cytometry A*. DOI:10.1002/cyto.a.22219 (CF®405S Mix-n-Stain™ labeling of wheat germ agglutinin; FC)
- Sainski, A. M., Dai, H., Natesampillai, S., Pang, Y. P., Bren, G. D., Cummins, N. W., Correia, C., Meng, X. W., Tarara, J. E., Ramirez-Alvarado, M., et al. (2014). [Casp8p41 generated by HIV protease kills CD4 T cells through direct Bak activation](#). *J Cell Biol* 207, 159. 10.1083/jcb.20140505109252014c (Mix-n-Stain™ CF®640R; FC)
- Shi, Y. Z., Xiong, S., Zhang, Y., Chin, L. K., Chen, Y., Zhang, J. B., Zhang, T. H., Ser, W., Larson, A., Hoi, L. S., et al. (2018). [Sculpting nanoparticle dynamics for single-bacteria-level screening and direct binding-efficiency measurement](#). *Nat Commun* 9, 815. 10.1038/s41467-018-03156-5 (Mix-n-Stain™ biotin; antibody affinity measurement)
- Shimizu, F., Schaller, K. L., Owens, G. P., Coteleur, A. C., Kellner, D., Takeshita, Y., Obermeier, B., Kryzer, T. J., Sano, Y., Kanda, T., et al. (2017). [Glucose-regulated protein 78 autoantibody associates with blood-brain barrier disruption in neuromyelitis optica](#). 10.1126/scitranslmed.aai9111 (Mix-n-Stain™ CF®568; IF)
- Tschumper, R. C., Dispenzieri, A., Abraham, R. S., Henderson, K. J., and Jelinek, D. F. (2013). [Molecular analysis of immunoglobulin genes reveals frequent clonal relatedness in double monoclonal gammopathies](#). *Blood Cancer J* 3, e112. DOI:10.1038/bcj.2013.12 (Mix-n-Stain™ CF®350; IF)
- Wada, M., Nishie, W., Ujiie, H., Izumi, K., Iwata, H., Natsuga, K., Nakamura, H., Kitagawa, Y., and Shimizu, H. (2016). [Epitope-Dependent Pathogenicity of Antibodies Targeting a Major Bullous Pemphigoid Autoantigen Collagen XVII/BP180](#). *J Invest Dermatol* 136, 938-946. 10.1016/j.jid.2015.11.030 (Mix-n-Stain™ CF®555; IF)
- Wienand, K., and Shires, K. (2015). [The Use of MAGE C1 and Flow Cytometry to Determine the Malignant Cell Type in Multiple Myeloma](#). *PLoS ONE* 10(3): e0120734. doi:10.1371/journal.pone.0120734 (Mix-n-Stain™ CF®488A; FC)
- Xu, Z., Parra, D., Gomez, D., Salinas, I., Zhang, Y. A., von Gersdorff Jorgensen, L., Heinecke, R. D., Buchmann, K., LaPatra, S., and Sunyer, J. O. (2013). [Teleost skin, an ancient mucosal surface that elicits gut-like immune responses](#). *Proc Natl Acad Sci U S A* 110, 13097-13102. 10.1073/pnas.1304319110 (Mix-n-Stain™ CF®488A, Mix-n-Stain™ CF®555, Mix-n-Stain™ CF®640R; IF)
- Xu, Z., Takizawa, F., Parra, D., Gomez, D., von Gersdorff Jorgensen, L., LaPatra, S. E., and Sunyer, J. O. (2016). [Mucosal immunoglobulins at respiratory surfaces mark an ancient association that predates the emergence of tetrapods](#). *Nat Commun* 7, 10728. 10.1038/ncomms10728 (Mix-n-Stain™ CF®488A, Mix-n-Stain™ CF®555, Mix-n-Stain™ CF®640R; IF)
- Yamamoto, S., Yamazaki, T., Komazaki, S., Yamashita, T., Osaki, M., Matsubayashi, M., Kidoya, H., Takakura, N., Yamazaki, D., and Kakizawa, S. (2014). [Contribution of calumin to embryogenesis through participation in the endoplasmic reticulum-associated degradation activity](#). *Dev Biol* 393, 33-43. 10.1016/j.ydbio.2014.06.024 (Mix-n-Stain™ CF®488A, Mix-n-Stain™ CF®568; IF)
- Yamashita, A., Zhao, Y., Matsuura, Y., Yamasaki, K., Moriguchi-Goto, S., Sugita, C., Iwakiri, T., Okuyama, N., Koshimoto, C., Kawai, K., et al. (2014). [Increased metabolite levels of glycolysis and pentose phosphate](#)

[pathway in rabbit atherosclerotic arteries and hypoxic macrophage](#). PLoS One 9, e86426.
10.1371/journal.pone.0086426 (*Mix-n-Stain™ CF®488A, Mix-n-Stain™ CF®594; IF*)

Yang, S., Yang, Y., Yu, P., Yang, J., Jiang, X., Villar, V. A., Sibley, D. R., Jose, P. A., and Zeng, C. (2015). [Dopamine D1 and D5 receptors differentially regulate oxidative stress through paraoxonase 2 in kidney cells](#). Free Radic Res 49, 397-410. 10.3109/10715762.2015.1006215 (*Mix-n-Stain™*)

Yang, Y., Zhang, Y., Cuevas-Gonzalez, S., Villar, V. A., Escano, C., Asico, L., Yu, P., Grandy, D. K., Felder, R. A., Armando, I., et al. (2012). [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. DOI:10.1016/j.freeradbiomed.2012.05.015 (*Mix-n-Stain™ CF®555; IF*)

Zolnerciks, J. K., Akkaya, B. G., Snippe, M., Chiba, P., Seelig, A., and Linton, K. J. (2014). [The Q loops of the human multidrug resistance transporter ABCB1 are necessary to couple drug binding to the ATP catalytic cycle](#). FASEB J 28, 4335-4346. 10.1096/fj.13-245639 (*Mix-n-Stain™ CF®633; IF*)