

Selected References for TrueBlack® Lipofuscin Autofluorescence Quencher

Specimen/sample type is shown in italics

Click on a link below to see journal article

- Aldrich, A., et al. [Efficacy of phosphodiesterase-4 inhibitors in juvenile Batten disease \(CLN3\)](#). *Annals of Neurology* 80, 909-923 (2016). doi:10.1002/ana.24815 (*mouse brain 30 um floating sections*)
- Aparicio, J.G. et al. [Temporal expression of CD184\(CXCR4\) and CD171\(L1CAM\) identifies distinct early developmental stages of human retinal ganglion cells in embryonic stem cell derived retina](#). *Exp Eye Res.* S0014-4835(16)30468-7 (2016). (*human retina*)
- Arhatta, M. [TMEM33 regulates intracellular calcium homeostasis in renal tubular epithelial cells](#). *Nature Communications* 10:2024 (2019). (*mouse kidney*)
- Axelrod, H.D., et al. [Optimization of Immunofluorescent Detection of Bone Marrow Disseminated Tumor Cells](#). *Biological Procedures Online* 20, 13 (2018). (*human bone marrow*)
- Badders, N.M., et al. [Selective modulation of the androgen receptor AF2 domain rescues degeneration in spinal bulbar muscular atrophy](#). *Nature Medicine* 24, 427 (2018). (*mouse spinal cord, skeletal muscle*)
- Banovac, I., et al. [Somato-dendritic morphology and axon origin site specify von Economo neurons as a subclass of modified pyramidal neurons in the human anterior cingulate cortex](#). *J. Anat* (2019) <https://doi.org/10.1111/joa.13068> (*human, macaque brain*)
- Barrenas, F., et al. [Macrophage-associated wound healing contributes to African green monkey SIV pathogenesis control](#). *Nat Commun* 10, 5101 (2019). doi:10.1038/s41467-019-12987-9 (*green monkey rectum*)
- Bosch, M.E., et al. [Self-Complementary AAV9 Gene Delivery Partially Corrects Pathology Associated with Juvenile Neuronal Ceroid Lipofuscinosis \(CLN3\)](#). *J Neurosci* 36, 9669-9682 (2016). (*mouse brain 30 um floating sections*)
- Chan, T., et al. [Human organic anion transporting polypeptide 1A2 \(OATP1A2\) mediates cellular uptake of all-trans-retinol in human retinal pigmented epithelial cells](#). *Br J Pharmacol* 172, 2343-2353 (2015). (*human retina*)
- De Blasio, D., et al. [Human brain trauma severity is associated with lectin complement pathway activation](#). *J Cereb Blood Flow Metab.* 39(5): 794 (2019). (*human brain*)
- Ellsworth J.L., et al. [Clade F AAVHSCs cross the blood brain barrier and transduce the central nervous system in addition to peripheral tissues following intravenous administration in nonhuman primates](#). *PLoS ONE* 14(11) e0225582 (2019). <https://doi.org/10.1371/journal.pone.0225582> (*cynomolgous brain, spinal chord, dorsal root ganglia 40 um floating sections*)
- Fosso, M.Y., [Differential Effects of Structural Modifications on the Competition of Chalcones for the PIB Amyloid Imaging Ligand-Binding Site in Alzheimer's Disease Brain and Synthetic Abeta Fibrils](#). *ACS Chem Neurosci* 7, 171-176 (2016). 10.1021/acscchemneuro.5b00266 (*human brain*)
- Futia, G. L., et al. [Quantitative image cytometry measurements of lipids, DNA, CD45 and cytokeratin for circulating tumor cell identification in a model system](#). *SPIE BiOS.* 15 (2016). (*polycarbonate filters*)
- Gal, J., et al. [Detergent Insoluble Proteins and Inclusion Body-Like Structures Immunoreactive for PRKDC/DNA-PK/DNA-PKcs, FTL, NNT, and AIFM1 in the Amygdala of Cognitively Impaired Elderly Persons](#). *J. Neuropathol Exp Neurol* 77(1), 21 (2018). (*human brain*)

- Hanael, E., et al. [Blood-brain barrier dysfunction in canine epileptic seizures detected by dynamic contrast-enhanced magnetic resonance imaging](#). *Epilepsia* 60 (5), 1005 (2019). (*canine brain*)
- Harun-Or-Rashid, M., et al. [Structural and functional rescue of chronic metabolically stressed optic nerves through respiration](#). *J. Neurosci* 38 (22), 5122 (2018). (*human retina*)
- Inoue, Y., et al. [RS9, a novel Nrf2 activator, attenuates light-induced death of cells of photoreceptor cells and Muller glia cells](#). *J Neurochem* 141, 750-765 (2017). doi:10.1111/jnc.14029 (*mouse retina*)
- Ising, C., et al. [NLRP3 inflammasome activation drives tau pathology](#). *Nature* 575, 669–673 (2019). doi:10.1038/s41586-019-1769-z (*human frontal temporal dementia brain*)
- Isonaka, R., et al. [Alpha-Synuclein Deposition Within Sympathetic Noradrenergic Neurons Is Associated With Myocardial Noradrenergic Deficiency in Neurogenic Orthostatic Hypotension](#). *Hypertension* 73:910–918 (2019). (*human skin biopsy*)
- Ito, N., et al. [Kososan, a Kampo medicine, prevents a social avoidance behavior and attenuates neuroinflammation in socially defeated mice](#). *J Neuroinflammation* 14, 98 (2017). doi:10.1186/s12974-017-0876-8 (*50 um mouse brain floating sections*)
- Jarvela, T.S. et al. [The neural chaperone proSAAS blocks alpha-synuclein fibrillation and neurotoxicity](#). *Proc Natl Acad Sci U S A* 113, E4708-4715 (2016). (*human brain*)
- Jolly, S., et al. [Single-Cell Quantification of mRNA Expression in The Human Brain](#). *Scientific Reports* 9, 12353 (2019). (*human brain, RNAscope*)
- Kageyama, Y., et al. [Amyloid \$\beta\$ toxic conformer has dynamic localization in the human inferior parietal cortex in absence of amyloid plaques](#). *Scientific Reports* 8, 16895 (2018). (*human brain*)
- Katsumata, Y., et al. [Alzheimer Disease Pathology-Associated Polymorphism in a Complex Variable Number of Tandem Repeat Region Within the MUC6 Gene, Near the AP2A2 Gene](#). *J. Neuropathol Exp Neurol*, nlz116 (2019). <https://doi.org/10.1093/jnen/nlz116> (*human Alzheimer disease brain*)
- Kuzkina, A., et al. [The aggregation state of \$\alpha\$ -synuclein deposits in dermal nerve fibers of patients with Parkinson's disease resembles that in the brain](#). *Parkinsonism & Related Disorders* 64, 66 (2019). (*human brain*)
- Lenz N., et al. [Evaluation of antivirals against tick borne encephalitis virus in organotypic brain slices of rat cerebellum](#). *PLoS ONE* 13(10), e0205294 (2018). (*rat cerebellum*)
- LeVine, H., III et al. [Down syndrome: age-dependence of PiB binding in postmortem frontal cortex across the lifespan](#). *Neurobiology of Aging* 54, 163-169 (2017). doi:10.1016/j.neurobiolaging.2017.03.005 (*human brain*)
- Litvinchuk, A., et al. [Complement C3aR Inactivation Attenuates Tau Pathology and Reverses an Immune Network Deregulated in Tauopathy Models and Alzheimer's Disease](#). *Neuron* 100 (6), 1337 (2018). (*human brain*)
- Maejima, Y., et al. [The hypothalamus to brainstem circuit suppresses late-onset body weight gain](#). *Sci Rep* 9, 18360 (2019). doi:10.1038/s41598-019-54870-z (aged rat brains 40 um floating sections, 8 min. pre-treatment)
- Nakagawa, H., et al. [Treatment With the Neutralizing Antibody Against Repulsive Guidance Molecule-a Promotes Recovery From Impaired Manual Dexterity in a Primate Model of Spinal Cord Injury](#). *Cerebral Cortex* 29(2), 561 (2019). (*rhesus monkey brain 50 um sections, spinal cord 40 um sections*)
- O'Hara, J.M., et al. [Generation of protective pneumococcal-specific nasal resident memory CD4+ T cells via parenteral immunization](#). *Mucosal Immunol* (2019). doi:10.1038/s41385-019-0218-5 (*mouse nasal mucosa*)

- Patel, N.N., et al. [Fungal extracts stimulate solitary chemosensory cell expansion in noninvasive fungal rhinosinusitis](#). International Forum of Allergy & Rhinology (2019). doi: 10.1002/alr.22334 (*human sinonasal tissue*)
- Petrasek, J., et al. [Protective Effect of SMAD-Specific E3 Ubiquitin Protein Ligase 1 in Alcoholic Steatohepatitis in Mice](#). Hepatology Communications doi:10.1002/hep4.1427 (*human liver*)
- Pieknik, J.R., et al. [Herpes Simplex Virus 2 in Autonomic Ganglia: Evidence for Spontaneous Reactivation](#). J. Virol. (2019). doi: 10.1128/JVI.00227-19 (*guinea pig dorsal root ganglia*)
- Plummer, N.W. [Two Subpopulations of Noradrenergic Neurons in the Locus Coeruleus Complex Distinguished by Expression of the Dorsal Neural Tube Marker Pax7](#). Front Neuroanat. (2017) 11, 60. doi: 10.3389/fnana.2017.00060 (*40 um mouse brain floating sections*)
- Polyzos, A.A., et al. [Metabolic Reprogramming in Astrocytes Distinguishes Region-Specific Neuronal Susceptibility in Huntington Mice](#). Cell Metab 29 (6), 1258 (2019). (*mouse brain*)
- Rey, N., et al. [Widespread transneuronal propagation of alpha-synucleinopathy triggered in olfactory bulb mimics prodromal Parkinson's disease](#). J Exp Med 213, 1759-1778 (2016). (*mouse brain 30 um sections*)
- Sala Frigerio, C., et al. [The Major Risk Factors for Alzheimer's Disease: Age, Sex, and Genes Modulate the Microglia Response to A \$\beta\$ Plaques](#). Cell Reports 27 (4), 1293 (2019). (*mouse brain*)
- Sorrelle, N. [Improved Multiplex Immunohistochemistry for Immune Microenvironment Evaluation of Mouse Formalin-Fixed, Paraffin-Embedded Tissues](#). J. Immunol (2018). doi:10.4049/jimmunol.1800878 (*mouse spleen, tumor xenograft*)
- Sun, Q., et al. [SEMA6D regulates perinatal cardiomyocyte proliferation and maturation in mice](#). Developmental Biology 452 (1), 1-7 (2019). (*mouse embryo*)
- Takahashi, K., et al. [Platelet-Derived Growth Factor-BB Lessens Light-Induced Rod Photoreceptor Damage in Mice](#). IOVS 58, 6299 (2017). (*mouse retina*)
- Tanaka, M., et al. [Aggregation of scaffolding protein DISC1 dysregulates phosphodiesterase 4 in Huntington's disease](#). J Clin Invest 127, 1438-1450 (2017). doi:85594 [pii]10.1172/JCI85594 (*human brain*)
- Tsitoura, E., et al. [Accumulation of damaged mitochondria in alveolar macrophages with reduced OXPHOS related gene expression in IPF](#). Respiratory Research (2019). 20:264 <https://doi.org/10.1186/s12931-019-1196-6> (*human bronchoalveolar lavage cells*)
- vonderEmbse, A.N., et al. [Dysfunctional microglia:neuron interactions with significant female bias in a developmental gene x environment rodent model of Alzheimer's disease](#). Int. Immunopharmacol. 71, 241 (2019). (*mouse brain*)
- Whittington, N.C. and Wray, S. [Suppression of Red Blood Cell Autofluorescence for Immunocytochemistry on Fixed Embryonic Mouse Tissue](#). Curr Protoc Neurosci 81, 2.28.1 (2017). (*mouse embryo*)
- Yamamoto, T., et al. [Expression of secreted phosphoprotein 1 \(osteopontin\) in human sensorimotor cortex and spinal cord: Changes in patients with amyotrophic lateral sclerosis](#). Brain Res. S0006-8993(16)30727-2 (2016). (*human brain*)
- Yang, J., et al. [Quenching autofluorescence in tissue immunofluorescence](#). Wellcome Open Res, 2:79 (2017). doi: 10.12688/wellcomeopenres.12251.1 (*mouse kidney, mouse placenta, human teratoma*)
- You, A.Y.F., et al. [Raman spectroscopy imaging reveals interplay between atherosclerosis and medial calcification in the human aorta](#). Sci. Adv. 3, e1701156 (2017). (*human aorta*)