

TrueBlack® References
[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*)
- 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*)
- 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*)
- 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*)
- 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*)
- 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*)
- 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 (*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*)
- 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*)
- 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*)
- 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*)
- 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*)