
TrueBlack® Lipofuscin Autofluorescence Quencher References

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

Contents:

The use of TrueBlack® Lipofuscin Autofluorescence Quencher has been published in over a hundred journal articles. The full list of publications using TrueBlack® is organized below by the tissue type TrueBlack® was used with.

Brain Tissue

- [Human Brain Tissue](#) p. 2
- [Mouse Brain Tissue](#) p. 4
- [Monkey Brain Tissue](#) p. 6
- [Rat Brain Tissue](#) p. 6
- [Other Brain Tissue](#) p. 7

Retina Tissue

- [Human Retina Tissue](#) p. 7
- [Mouse Retina Tissue](#) p. 7
- [Monkey Retina Tissue](#) p. 8

Lung Tissue

- [Human Lung Tissue](#) p. 8
- [Mouse Lung Tissue](#) p. 8
- [Rat Lung Tissue](#) p. 8

Nasal Tissue

- [Human Nasal Tissue](#) p. 8
- [Mouse Nasal Tissue](#) p. 9

Heart Tissue

- [Human Heart Tissue](#) p. 9
- [Mouse Heart Tissue](#) p. 9

Liver Tissue

- [Human Liver Tissue](#) p. 9
- [Mouse Liver Tissue](#) p. 9

Intestine Tissue

- [Human Intestine Tissue](#) p. 9
- [Mouse Intestine Tissue](#) p. 10

Kidney Tissue

- [Mouse Kidney Tissue](#) p. 10

[Spinal Cord Tissue](#) p. 10

[Red Blood Cells](#) p. 11

[Thick Floating Sections](#) p. 11

[Other Tissue](#) p. 12

Human Brain Tissue

- 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* 235(3), 651-669, (2019), DOI: 10.1111/joa.13068 (human, macaque brain)
- Bernstock, J. D., et al. [Cerebral Ischemia Increases Small Ubiquitin-Like Modifier Conjugation within Human Penumbra Tissue: Radiological-Pathological Correlation](#). *Front Neurol* 8, 738, (2017), DOI: 10.3389/fneur.2017.00738 (human brain)
- Bian, Z., et al. [Accelerated accumulation of fibrinogen peptide chains with Abeta deposition in Alzheimer's disease \(AD\) mice and human AD brains](#). *Brain Res* 1767, 147569, (2021), DOI: 10.1016/j.brainres.2021.147569 (AD mouse brain, AD human brain)
- Bouche, M., et al. [Novel Treatment for Glioblastoma Delivered by a Radiation Responsive and Radiopaque Hydrogel](#). *ACS Biomater Sci Eng* 7(7), 3209-3220, (2021), DOI: 10.1021/acsbomaterials.1c00385 (human GBM brain tissue)
- Chen, W. T., et al. [Spatial Transcriptomics and In Situ Sequencing to Study Alzheimer's Disease](#). *Cell* 182(4), 976-991 e919, (2020), DOI: 10.1016/j.cell.2020.06.038 (human brain in situ sequencing)
- Codron, P., et al. [STochastic Optical Reconstruction Microscopy \(STORM\) reveals the nanoscale organization of pathological aggregates in human brain](#). *Neuropathol Appl Neurobiol* 47(1), 127-142, (2021), DOI: 10.1111/nan.12646 (human brain, STORM)
- De Blasio, D., et al. [Human brain trauma severity is associated with lectin complement pathway activation](#). *J Cereb Blood Flow Metab* 39(5), 794-807, (2019), DOI: 10.1177/0271678X18758881 (human brain)
- Dominy, S. S., et al. [Porphyromonas gingivalis in Alzheimer's disease brains: Evidence for disease causation and treatment with small-molecule inhibitors](#). *Sci Adv* 5(1), eaau3333, (2019), DOI: 10.1126/sciadv.aau3333 (human brain)
- Farmer, K. M., et al. [P53 aggregation, interactions with tau, and impaired DNA damage response in Alzheimer's disease](#). *Acta Neuropathol Commun* 8(1), 132, (2020), DOI: 10.1186/s40478-020-01012-6 (human, mouse brain, pretreatment)
- Fosso, M. Y., et al. [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(2), 171-176, (2016), DOI: 10.1021/acscchemneuro.5b00266 (human brain)
- Fullard, J. F., et al. [Single-nucleus transcriptome analysis of human brain immune response in patients with severe COVID-19](#). *Genome Med* 13(1), 118, (2021), DOI: 10.1186/s13073-021-00933-8 (human brain)
- 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), 21-39, (2018), DOI: 10.1093/jnen/nlx097 (human brain)
- Guttenplan, K. A., et al. [Knockout of reactive astrocyte activating factors slows disease progression in an ALS mouse model](#). *Nat Commun* 11(1), 3753, (2020), DOI: 10.1038/s41467-020-17514-9 (human brain, spinal cord FFPE)

- Hirai, S., et al. [Brain Angiopathy and Impaired Glucose Logistics in Mice with Psychosis-related Higher Brain Dysfunction](#). bioRxiv, 2020.2002.2014.939546, (2020), DOI: 10.1101/2020.02.14.939546 (human brain FFPE)
- Huang, M., et al. [Network analysis of the progranulin-deficient mouse brain proteome reveals pathogenic mechanisms shared in human frontotemporal dementia caused by GRN mutations](#). Acta Neuropathol Commun 8(1), 163, (2020), DOI: 10.1186/s40478-020-01037-x (human, mouse brain FFPE)
- Ising, C., et al. [NLRP3 inflammasome activation drives tau pathology](#). Nature 575(7784), 669-673, (2019), DOI: 10.1038/s41586-019-1769-z 10.1038/s41586-019-1769-z [pii] (human frontal temporal dementia brain)
- Jarvela, T. S., et al. [The neural chaperone proSAAS blocks alpha-synuclein fibrillation and neurotoxicity](#). Proc Natl Acad Sci U S A 113(32), E4708-4715, (2016), DOI: 1601091113 [pii] 10.1073/pnas.1601091113 (human brain)
- Jolly, S., et al. [Single-Cell Quantification of mRNA Expression in The Human Brain](#). Sci Rep 9(1), 12353, (2019), DOI: 10.1038/s41598-019-48787-w (human brain, RNAscope)
- 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 79(1), 3-21, (2020), DOI: 10.1093/jnen/nlz116 (human Alzheimer disease brain)
- Krienen, F. M., et al. [Innovations present in the primate interneuron repertoire](#). Nature 586(7828), 262-269, (2020), DOI: 10.1038/s41586-020-2781-z (primate brain 120 um sections)
- 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 Relat Disord 64, 66-72, (2019), DOI: 10.1016/j.parkreldis.2019.03.003 (human brain)
- Lackie, R. E., et al. [Increased levels of Stress-inducible phosphoprotein-1 accelerates amyloid-beta deposition in a mouse model of Alzheimer's disease](#). Acta Neuropathol Commun 8(1), 143, (2020), DOI: 10.1186/s40478-020-01013-5 (human brain FFPE)
- Lee, J. H., et al. [Astrocytes phagocytose adult hippocampal synapses for circuit homeostasis](#). Nature 590(7847), 612-617, (2021), DOI: 10.1038/s41586-020-03060-3 (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), 1337-1353 e1335, (2018), DOI: 10.1016/j.neuron.2018.10.031 (human brain)
- Miskic, T., et al. [Adult Upper Cortical Layer Specific Transcription Factor CUX2 Is Expressed in Transient Subplate and Marginal Zone Neurons of the Developing Human Brain](#). Cells 10(2), (2021), DOI: 10.3390/cells10020415 (human brain FFPE)
- Palasz, A., et al. [The first identification of nesfatin-1-expressing neurons in the human bed nucleus of the stria terminalis](#). J Neural Transm (Vienna) 126(3), 349-355, (2019), DOI: 10.1007/s00702-019-01984-3 (human brain FFPE)
- Palasz, A., et al. [Spexin-expressing neurons in the magnocellular nuclei of the human hypothalamus](#). J Chem Neuroanat 111, 101883, (2021), DOI: 10.1016/j.jchemneu.2020.101883 (human brain IF and IHC)

Pfisterer, U., et al. [Identification of epilepsy-associated neuronal subtypes and gene expression underlying epileptogenesis](#). Nat Commun 11(1), 5038, (2020), DOI: 10.1038/s41467-020-18752-7 (human brain RNAScope)

Reinert, L. S., et al. [Brain immune cells undergo cGAS/STING-dependent apoptosis during herpes simplex virus type 1 infection to limit type I IFN production](#). J Clin Invest 131(1), (2021), DOI: 10.1172/JCI136824 (human brain FFPE)

Riggle, B. A., et al. [CD8+ T cells target cerebrovasculature in children with cerebral malaria](#). J Clin Invest 130(3), 1128-1138, (2020), DOI: 10.1172/JCI133474 (human brain)

Rocco, B. R., et al. [Fluorescence-based cell-specific detection for laser-capture microdissection in human brain](#). Sci Rep 7(1), 14213, (2017), DOI: 10.1038/s41598-017-14484-9 (human brain)

Senatorov, V. V., et al. [Blood-brain barrier dysfunction in aging induces hyper-activation of TGF-beta signaling and chronic yet reversible neural dysfunction](#). bioRxiv, 537431, (2019), DOI: 10.1101/537431 (human, aged mouse brain)

Shahidehpour, R. K., et al. [Dystrophic microglia are associated with neurodegenerative disease and not healthy aging in the human brain](#). Neurobiol Aging 99, 19-27, (2021), DOI: 10.1016/j.neurobiolaging.2020.12.003 (human brain FFPE)

Tanaka, M., et al. [Aggregation of scaffolding protein DISC1 dysregulates phosphodiesterase 4 in Huntington's disease](#). J Clin Invest 127(4), 1438-1450, (2017), DOI: 10.1172/JCI85594 (human brain)

Watanabe, R., et al. [Intracellular dynamics of Ataxin-2 in the human brains with normal and frontotemporal lobar degeneration with TDP-43 inclusions](#). Acta Neuropathol Commun 8(1), 176, (2020), DOI: 10.1186/s40478-020-01055-9 (human brain)

Woolf, Z., et al. [Single-cell image analysis reveals a protective role for microglia in glioblastoma](#). Neurooncol Adv 3(1), vdab031, (2021), DOI: 10.1093/noajnl/vdab031 (human glioblastoma tissue)

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 1655, 168-175, (2017), DOI: 10.1016/j.brainres.2016.10.030 (human brain)

Zhao, J., et al. [Apolipoprotein E regulates lipid metabolism and alpha-synuclein pathology in human iPSC-derived cerebral organoids](#). Acta Neuropathol, (2021), DOI: 10.1007/s00401-021-02361-9 (human cerebral organoids)

Zhou, X., et al. [Mature myelin maintenance requires Qki to coactivate PPARbeta-RXRalpha-mediated lipid metabolism](#). J Clin Invest 130(5), 2220-2236, (2020), DOI: 10.1172/JCI131800 (human brain)

Mouse Brain Tissue

Aldrich, A., et al. [Efficacy of phosphodiesterase-4 inhibitors in juvenile Batten disease \(CLN3\)](#). Ann Neurol 80(6), 909-923, (2016), DOI: 10.1002/ana.24815 (mouse brain 30 um floating sections)

Bian, Z., et al. [Accelerated accumulation of fibrinogen peptide chains with Abeta deposition in Alzheimer's disease \(AD\) mice and human AD brains](#). Brain Res 1767, 147569, (2021), DOI: 10.1016/j.brainres.2021.147569 (AD mouse brain, AD human brain)

- Bosch, M. E., et al. [Self-Complementary AAV9 Gene Delivery Partially Corrects Pathology Associated with Juvenile Neuronal Ceroid Lipofuscinosis \(CLN3\)](#). J Neurosci 36(37), 9669-9682, (2016), DOI: 10.1523/JNEUROSCI.1635-16.2016 (mouse brain 30 um floating sections)
- Demarest, T. G., et al. [Biological sex and DNA repair deficiency drive Alzheimer's disease via systemic metabolic remodeling and brain mitochondrial dysfunction](#). Acta Neuropathol 140(1), 25-47, (2020), DOI: 10.1007/s00401-020-02152-8 (AD mouse brain)
- Everard, A., et al. [Intestinal epithelial N-acylphosphatidylethanolamine phospholipase D links dietary fat to metabolic adaptations in obesity and steatosis](#). Nat Commun 10(1), 457, (2019), DOI: 10.1038/s41467-018-08051-7 (mouse brain)
- Farmer, K. M., et al. [P53 aggregation, interactions with tau, and impaired DNA damage response in Alzheimer's disease](#). Acta Neuropathol Commun 8(1), 132, (2020), DOI: 10.1186/s40478-020-01012-6 (human, mouse brain, pretreatment)
- Feng, T., et al. [Loss of TMEM106B and PGRN leads to severe lysosomal abnormalities and neurodegeneration in mice](#). EMBO Rep 21(10), e50219, (2020), DOI: 10.15252/embr.202050219 (mouse brain)
- Huang, M., et al. [Network analysis of the progranulin-deficient mouse brain proteome reveals pathogenic mechanisms shared in human frontotemporal dementia caused by GRN mutations](#). Acta Neuropathol Commun 8(1), 163, (2020), DOI: 10.1186/s40478-020-01037-x (human, mouse brain FFPE)
- Kaczmarczyk, L., et al. [Slc1a3-2A-CreERT2 mice reveal unique features of Bergmann glia and augment a growing collection of Cre drivers and effectors in the 129S4 genetic background](#). Sci Rep 11(1), 5412, (2021), DOI: 10.1038/s41598-021-84887-2 (mouse brain FFPE)
- Mayagoitia, K., et al. [Loss of APP in mice increases thigmotaxis and is associated with elevated brain expression of IL-13 and IP-10/CXCL10](#). Physiol Behav 240, 113533, (2021), DOI: 10.1016/j.physbeh.2021.113533 (AD mouse brain tissue)
- Mikulka, C. R., et al. [Cell-autonomous expression of the acid hydrolase galactocerebrosidase](#). Proc Natl Acad Sci U S A 117(16), 9032-9041, (2020), DOI: 10.1073/pnas.1917675117 (mouse brain)
- Plummer, N. W., et al. [Two Subpopulations of Noradrenergic Neurons in the Locus Coeruleus Complex Distinguished by Expression of the Dorsal Neural Tube Marker Pax7](#). Front Neuroanat 11, 60, (2017), DOI: 10.3389/fnana.2017.00060 (mouse brain 40 um floating sections)
- Polyzos, A. A., et al. [Metabolic Reprogramming in Astrocytes Distinguishes Region-Specific Neuronal Susceptibility in Huntington Mice](#). Cell Metab 29(6), 1258-1273 e1211, (2019), DOI: 10.1016/j.cmet.2019.03.004 (mouse brain)
- Rey, N. L., et al. [Widespread transneuronal propagation of alpha-synucleinopathy triggered in olfactory bulb mimics prodromal Parkinson's disease](#). J Exp Med 213(9), 1759-1778, (2016), DOI: 10.1084/jem.20160368 (mouse brain 30 um sections)
- Sadhukhan, T., et al. [In a mouse model of INCL reduced S-palmitoylation of cytosolic thioesterase APT1 contributes to microglia proliferation and neuroinflammation](#). J Inherit Metab Dis, (2021), DOI: 10.1002/jimd.12379 (mouse brain FFPE)

Sala Frigerio, C., et al. [The Major Risk Factors for Alzheimer's Disease: Age, Sex, and Genes Modulate the Microglia Response to Abeta Plaques](#). Cell Rep 27(4), 1293-1306 e1296, (2019), DOI: 10.1016/j.celrep.2019.03.099 (mouse brain)

Sarkar, C., et al. [Cln1-mutations suppress Rab7-RILP interaction and impair autophagy contributing to neuropathology in a mouse model of infantile neuronal ceroid lipofuscinosis](#). J Inherit Metab Dis 43(5), 1082-1101, (2020), DOI: 10.1002/jimd.12242 (mouse brain FFPE)

Senatorov, V. V., et al. [Blood-brain barrier dysfunction in aging induces hyper-activation of TGF-beta signaling and chronic yet reversible neural dysfunction](#). bioRxiv, 537431, (2019), DOI: 10.1101/537431 (human, aged mouse brain)

Tanriover, G., et al. [Prominent microglial inclusions in transgenic mouse models of alpha-synucleinopathy that are distinct from neuronal lesions](#). Acta Neuropathol Commun 8(1), 133, (2020), DOI: 10.1186/s40478-020-00993-8 (mouse brain)

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, 241-250, (2019), DOI: 10.1016/j.intimp.2019.03.039 (mouse brain)

Yamagata, T., et al. [CRISPR/dCas9-based Scn1a gene activation in inhibitory neurons ameliorates epileptic and behavioral phenotypes of Dravet syndrome model mice](#). Neurobiol Dis 141, 104954, (2020), DOI: 10.1016/j.nbd.2020.104954 (mouse brain FFPE)

Monkey Brain Tissue

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 235(3), 651-669, (2019), DOI: 10.1111/joa.13068 (human, macaque brain)

Beckman, D., et al. [Oligomeric Abeta in the monkey brain impacts synaptic integrity and induces accelerated cortical aging](#). Proc Natl Acad Sci U S A, (2019), DOI: 10.1073/pnas.1902301116 (monkey brain 50 um floating sections)

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), DOI: 10.1371/journal.pone.0225582
PONE-D-19-21158 [pii] (cynomolgous brain, spinal cord, DRG 40 um floating sections)

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](#). Cereb Cortex 29(2), 561-572, (2019), DOI: 10.1093/cercor/bhx338 (rhesus monkey brain 50 um sections, spinal cord 40 um sections)

Rat Brain Tissue

De Miranda, B. R., et al. [The industrial solvent trichloroethylene induces LRRK2 kinase activity and dopaminergic neurodegeneration in a rat model of Parkinson's disease](#). Neurobiol Dis 153, 105312, (2021), DOI: 10.1016/j.nbd.2021.105312 (rat brain 35 um floating sections)

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), DOI: 10.1371/journal.pone.0205294 (rat cerebellum)

Maejima, Y., et al. [The hypothalamus to brainstem circuit suppresses late-onset body weight gain](#). Sci Rep 9(1), 18360, (2019), DOI: 10.1038/s41598-019-54870-z
10.1038/s41598-019-54870-z [pii] (aged rat brains 40 um floating sections, 8 min. pre-treatment)

Other Brain Tissue

Hanael, E., et al. [Blood-brain barrier dysfunction in canine epileptic seizures detected by dynamic contrast-enhanced magnetic resonance imaging](#). Epilepsia 60(5), 1005-1016, (2019), DOI: 10.1111/epi.14739 (canine brain)

Human Retina Tissue

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 154, 177-189, (2017), DOI: 10.1016/j.exer.2016.11.013 (human retina)

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(9), 2343-2353, (2015), DOI: 10.1111/bph.13060 (human retina)

Dopper, H., et al. [Differentiation Protocol for 3D Retinal Organoids, Immunostaining and Signal Quantitation](#). Curr Protoc Stem Cell Biol 55(1), e120, (2020), DOI: 10.1002/cpsc.120 (3D retinal organoids)

Harun-Or-Rashid, M., et al. [Structural and Functional Rescue of Chronic Metabolically Stressed Optic Nerves through Respiration](#). J Neurosci 38(22), 5122-5139, (2018), DOI: 10.1523/JNEUROSCI.3652-17.2018 (human retina)

Liu, Z., et al. [Submacular integration of hESC-RPE monolayer xenografts in a surgical non-human primate model](#). Stem Cell Res Ther 12(1), 423, (2021), DOI: 10.1186/s13287-021-02395-6 (human retinal pigment epithelial cells)

Voigt, A. P., et al. [Bulk and single-cell gene expression analyses reveal aging human choriocapillaris has pro-inflammatory phenotype](#). Microvasc Res 131, 104031, (2020), DOI: 10.1016/j.mvr.2020.104031 (human retina)

Mouse Retina Tissue

Harun-Or-Rashid, M., et al. [MCT2 overexpression rescues metabolic vulnerability and protects retinal ganglion cells in two models of glaucoma](#). Neurobiol Dis 141, 104944, (2020), DOI: 10.1016/j.nbd.2020.104944 (mouse 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(5), 750-765, (2017), DOI: 10.1111/jnc.14029 (mouse retina)

Massengill, M. T., et al. [Sectoral activation of glia in an inducible mouse model of autosomal dominant retinitis pigmentosa](#). Sci Rep 10(1), 16967, (2020), DOI: 10.1038/s41598-020-73749-y (mouse retina)

Takahashi, K., et al. [Platelet-Derived Growth Factor-BB Lessens Light-Induced Rod Photoreceptor Damage in Mice](#). Invest Ophthalmol Vis Sci 58(14), 6299-6307, (2017), DOI: 10.1167/iovs.17-22556 (mouse retina)

Monkey Retina Tissue

Gnanaguru, G., et al. [Targeting of miR-33 ameliorates phenotypes linked to age-related macular degeneration](#). Mol Ther, (2021), DOI: 10.1016/j.ymthe.2021.03.014 (primate retina)

Hughes, S., et al. [Prolonged intraocular residence and retinal tissue distribution of a fourth-generation compstatin-based C3 inhibitor in non-human primates](#). Clin Immunol 214, 108391, (2020), DOI: 10.1016/j.clim.2020.108391 (primate retina)

Human Lung Tissue

Cassandras, M., et al. [Gli1\(+\) mesenchymal stromal cells form a pathological niche to promote airway progenitor metaplasia in the fibrotic lung](#). Nat Cell Biol 22(11), 1295-1306, (2020), DOI: 10.1038/s41556-020-00591-9 (human, mouse lung)

Dora, D., et al. [Characterization of Tumor-Associated Macrophages and the Immune Microenvironment in Limited-Stage Neuroendocrine-High and -Low Small Cell Lung Cancer](#). Biology (Basel) 10(6), (2021), DOI: 10.3390/biology10060502 (human lung FFPE tissue)

Evren, E., et al. [Distinct developmental pathways from blood monocytes generate human lung macrophage diversity](#). Immunity 54(2), 259-275 e257, (2021), DOI: 10.1016/j.immuni.2020.12.003 (human lung)

Masso-Silva, J. A., et al. [Increased peripheral blood neutrophil activation phenotypes and NETosis in critically ill COVID-19 patients](#). medRxiv, 2021.2001.2014.21249831, (2021), DOI: 10.1101/2021.01.14.21249831 (COVID lung)

Tsitoura, E., et al. [Accumulation of damaged mitochondria in alveolar macrophages with reduced OXPPOS related gene expression in IPF](#). Respir Res 20(1), 264, (2019), DOI: 10.1186/s12931-019-1196-6 10.1186/s12931-019-1196-6 [pii] (human broncho-alveolar lavage cells)

Mouse Lung Tissue

Cassandras, M., et al. [Gli1\(+\) mesenchymal stromal cells form a pathological niche to promote airway progenitor metaplasia in the fibrotic lung](#). Nat Cell Biol 22(11), 1295-1306, (2020), DOI: 10.1038/s41556-020-00591-9 (human, mouse lung)

Young, R. E., et al. [Smooth Muscle Differentiation Is Essential for Airway Size, Tracheal Cartilage Segmentation, but Dispensable for Epithelial Branching](#). Dev Cell 53(1), 73-85 e75, (2020), DOI: 10.1016/j.devcel.2020.02.001 (mouse embryonic lung sections and whole mount)

Rat Lung Tissue

Andreu, I., et al. [The force loading rate drives cell mechanosensing through both reinforcement and fluidization](#). bioRxiv, 2021.2003.2008.434428, (2021), DOI: 10.1101/2021.03.08.434428 (rat lung 70 um frozen)

Human Nasal Tissue

Patel, N. N., et al. Fungal extracts stimulate solitary chemosensory cell expansion in noninvasive fungal rhinosinusitis. Int Forum Allergy Rhinol 9(7), 730-737, (2019), DOI: 10.1002/alr.22334 (human sinonasal tissue)

Mouse Nasal Tissue

Bloom, M. L., et al. [Renewal and Differentiation of GCD Necklace Olfactory Sensory Neurons](#). Chem Senses 45(5), 333-346, (2020), DOI: 10.1093/chemse/bjaa027 (mouse olfactory bulb)

O'Hara, J. M., et al. [Generation of protective pneumococcal-specific nasal resident memory CD4\(+\) T cells via parenteral immunization](#). Mucosal Immunol 13(1), 172-182, (2020), DOI: 10.1038/s41385-019-0218-5 (mouse nasal mucosa)

Human Heart Tissue

Miyosawa, K., et al. [Enhanced monocyte migratory activity in the pathogenesis of structural remodeling in atrial fibrillation](#). PLoS One 15(10), e0240540, (2020), DOI: 10.1371/journal.pone.0240540 (human heart)

You, A. Y. F., et al. [Raman spectroscopy imaging reveals interplay between atherosclerosis and medial calcification in the human aorta](#). Sci Adv 3(12), e1701156, (2017), DOI: 10.1126/sciadv.1701156 (human aorta)

Mouse Heart Tissue

O'Brien, M., et al. [Pressure overload generates a cardiac-specific profile of inflammatory mediators](#). Am J Physiol Heart Circ Physiol 319(2), H331-H340, (2020), DOI: 10.1152/ajpheart.00274.2020 (mouse myocardial sections)

Weckbach, L. T., et al. [Midkine drives cardiac inflammation by promoting neutrophil trafficking and NETosis in myocarditis](#). J Exp Med 216(2), 350-368, (2019), DOI: 10.1084/jem.20181102 (mouse heart red blood cells)

Human Liver Tissue

Ortega-Ribera, M., et al. [The Hepatic Sinusoid in Aging and Disease: Update and Advances From the 20th Liver Sinusoid Meeting](#). Hepatol Commun 4(7), 1087-1098, (2020), DOI: 10.1002/hep4.1517 (human liver sinusoids)

Sabin Bhandari, R. L., et al. [Effect of glucocorticoids on liver sinusoidal endothelial cells in vitro](#). 20th International Symposium On Cells Of The Hepatic Sinusoid, (2019), (human liver sinusoid cells)

Mouse Liver Tissue

Hannemann, C., et al. [Deficiency of inactive rhomboid protein 2 \(iRhom2\) attenuates diet-induced hyperlipidemia and early atherogenesis](#). Cardiovasc Res, (2021), DOI: 10.1093/cvr/cvab041 (mouse liver)

Lindquist, M. E., et al. [Exploring Crimean-Congo Hemorrhagic Fever Virus-Induced Hepatic Injury Using Antibody-Mediated Type I Interferon Blockade in Mice](#). J Virol 92(21), (2018), DOI: 10.1128/JVI.01083-18 (mouse liver)

Human Intestine Tissue

Rao, A., et al. [Cytokines regulate the antigen-presenting characteristics of human circulating and tissue-resident intestinal ILCs](#). Nat Commun 11(1), 2049, (2020), DOI: 10.1038/s41467-020-15695-x (human intestine)

Mouse Intestine Tissue

Guzman, M., et al. [An integrin alphaEbeta7-dependent mechanism of IgA transcytosis requires direct plasma cell contact with intestinal epithelium](#). Mucosal Immunol, (2021), DOI: 10.1038/s41385-021-00439-x (mouse intestinal epithelium)

Janowitz, C., et al. [Disruption of Intestinal Homeostasis and Intestinal Microbiota During Experimental Autoimmune Uveitis](#). Invest Ophthalmol Vis Sci 60(1), 420-429, (2019), DOI: 10.1167/iovs.18-24813 (mouse intestine)

Jarret, A., et al. [Enteric Nervous System-Derived IL-18 Orchestrates Mucosal Barrier Immunity](#). Cell 180(1), 50-63 e12, (2020), DOI: 10.1016/j.cell.2019.12.016 (mouse colon smFISH)

Mouse Kidney Tissue

Nespoux, J., et al. [Gene knockout of the Na\(+\)-glucose cotransporter SGLT2 in a murine model of acute kidney injury induced by ischemia-reperfusion](#). Am J Physiol Renal Physiol 318(5), F1100-F1112, (2020), DOI: 10.1152/ajprenal.00607.2019 (mouse kidney FFPE)

Walter, D. L., et al. [Coxsackievirus B4 Exposure Results in Variable Pattern Recognition Response in the Kidneys of Female Non-Obese Diabetic Mice Before Establishment of Diabetes](#). Viral Immunol 33(7), 494-506, (2020), DOI: 10.1089/vim.2019.0188 (mouse kidney FFPE)

Yang, J., et al. [Quenching autofluorescence in tissue immunofluorescence \[version 1; referees: 1 approved with reservations\]](#). Wellcome Open Res 2(79), (2017), DOI: doi: 10.12688/wellcomeopenres.12251.1 (mouse kidney, mouse placenta, human teratoma)

Spinal Cord Tissue

Badders, N. M., et al. [Selective modulation of the androgen receptor AF2 domain rescues degeneration in spinal bulbar muscular atrophy](#). Nat Med 24(4), 427-437, (2018), DOI: 10.1038/nm.4500 (mouse spinal cord, skeletal muscle)

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), DOI: 10.1371/journal.pone.0225582
PONE-D-19-21158 [pii] (cynomolgous brain, spinal cord, DRG 40 um floating sections)

Guttenplan, K. A., et al. [Knockout of reactive astrocyte activating factors slows disease progression in an ALS mouse model](#). Nat Commun 11(1), 3753, (2020), DOI: 10.1038/s41467-020-17514-9 (human brain, spinal cord FFPE)

MacLean, M., et al. [Microglia RAGE exacerbates the progression of neurodegeneration within the SOD1\(G93A\) murine model of amyotrophic lateral sclerosis in a sex-dependent manner](#). J Neuroinflammation 18(1), 139, (2021), DOI: 10.1186/s12974-021-02191-2 (mouse gastrocnemius tissue, human spinal cord)

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](#). Cereb Cortex 29(2), 561-572, (2019), DOI: 10.1093/cercor/bhx338 (rhesus monkey brain 50 um sections, spinal cord 40 um sections)

Red Blood Cells

Weckbach, L. T., et al. [Midkine drives cardiac inflammation by promoting neutrophil trafficking and NETosis in myocarditis](#). J Exp Med 216(2), 350-368, (2019), DOI: 10.1084/jem.20181102 (mouse heart red blood cells)

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 21-22 28 12, (2017), DOI: 10.1002/cpns.35 (mouse embryo red blood cells)

Thick Floating Sections

Aldrich, A., et al. [Efficacy of phosphodiesterase-4 inhibitors in juvenile Batten disease \(CLN3\)](#). Ann Neurol 80(6), 909-923, (2016), DOI: 10.1002/ana.24815 (mouse brain 30 um floating sections)

Andreu, I., et al. [The force loading rate drives cell mechanosensing through both reinforcement and fluidization](#). bioRxiv, 2021.2003.2008.434428, (2021), DOI: 10.1101/2021.03.08.434428 (rat lung 70 um frozen)

Beckman, D., et al. [Oligomeric Abeta in the monkey brain impacts synaptic integrity and induces accelerated cortical aging](#). Proc Natl Acad Sci U S A, (2019), DOI: 10.1073/pnas.1902301116 (monkey brain 50 um floating sections)

Bosch, M. E., et al. [Self-Complementary AAV9 Gene Delivery Partially Corrects Pathology Associated with Juvenile Neuronal Ceroid Lipofuscinosis \(CLN3\)](#). J Neurosci 36(37), 9669-9682, (2016), DOI: 10.1523/JNEUROSCI.1635-16.2016 (mouse brain 30 um floating sections)

De Miranda, B. R., et al. [The industrial solvent trichloroethylene induces LRRK2 kinase activity and dopaminergic neurodegeneration in a rat model of Parkinson's disease](#). Neurobiol Dis 153, 105312, (2021), DOI: 10.1016/j.nbd.2021.105312 (rat brain 35 um floating sections)

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), DOI: 10.1371/journal.pone.0225582 PONE-D-19-21158 [pii] (cynomolgous brain, spinal cord, DRG 40 um floating sections)

Ito, N., et al. [Kososan, a Kampo medicine, prevents a social avoidance behavior and attenuates neuroinflammation in socially defeated mice](#). J Neuroinflammation 14(1), 98, (2017), DOI: 10.1186/s12974-017-0876-8 10.1186/s12974-017-0876-8 [pii] (mouse brain 50 um floating sections)

Jimenez, J. A., et al. [Chd8 haploinsufficiency impairs early brain development and protein homeostasis later in life](#). Mol Autism 11(1), 74, (2020), DOI: 10.1186/s13229-020-00369-8 (mouse brain 60 um floating sections)

Krienen, F. M., et al. [Innovations present in the primate interneuron repertoire](#). Nature 586(7828), 262-269, (2020), DOI: 10.1038/s41586-020-2781-z (primate brain 120 um sections)

Maejima, Y., et al. [The hypothalamus to brainstem circuit suppresses late-onset body weight gain](#). Sci Rep 9(1), 18360, (2019), DOI: 10.1038/s41598-019-54870-z 10.1038/s41598-019-54870-z [pii] (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](#). Cereb

Cortex 29(2), 561-572, (2019), DOI: 10.1093/cercor/bhx338 (rhesus monkey brain 50 um sections, spinal cord 40 um sections)

Plummer, N. W., et al. [Two Subpopulations of Noradrenergic Neurons in the Locus Coeruleus Complex Distinguished by Expression of the Dorsal Neural Tube Marker Pax7](#). Front Neuroanat 11, 60, (2017), DOI: 10.3389/fnana.2017.00060 (mouse brain 40 um floating sections)

Rey, N. L., et al. [Widespread transneuronal propagation of alpha-synucleinopathy triggered in olfactory bulb mimics prodromal Parkinson's disease](#). J Exp Med 213(9), 1759-1778, (2016), DOI: 10.1084/jem.20160368 (mouse brain 30 um sections)

Other Tissue

Axelrod, H. D., et al. [Optimization of Immunofluorescent Detection of Bone Marrow Disseminated Tumor Cells](#). Biol Proced Online 20, 13, (2018), DOI: 10.1186/s12575-018-0078-5 (human bone marrow)

Barrenas, F., et al. [Macrophage-associated wound healing contributes to African green monkey SIV pathogenesis control](#). Nat Commun 10(1), 5101, (2019), DOI: 10.1038/s41467-019-12987-9 10.1038/s41467-019-12987-9 [pii] (African green monkey rectum)

Chu, A., et al. [Epithelial membrane protein 2 \(Emp2\) modulates innate immune cell population recruitment at the maternal-fetal interface](#). J Reprod Immunol 145, 103309, (2021), DOI: 10.1016/j.jri.2021.103309 (mouse FFPE tissues)

Cugurra, A., et al. [Skull and vertebral bone marrow are myeloid cell reservoirs for the meninges and CNS parenchyma](#). Science 373(6553), (2021), DOI: 10.1126/science.abf7844 (mouse bone marrow)

Dhouchak, S., et al. [Heparan sulfate proteoglycans in beta cells provide a critical link between endoplasmic reticulum stress, oxidative stress and type 2 diabetes](#). PLoS One 16(6), e0252607, (2021), DOI: 10.1371/journal.pone.0252607 (human and mouse pancreatic islet beta cells, T2D)

Freichel, T., et al. [Sequence-Defined Heteromultivalent Precision Glycomacromolecules Bearing Sulfonated/Sulfated Nonglycosidic Moieties Preferentially Bind Galectin-3 and Delay Wound Healing of a Galectin-3 Positive Tumor Cell Line in an In Vitro Wound Scratch Assay](#). Macromol Biosci 20(9), e2000163, (2020), DOI: 10.1002/mabi.202000163 (HEK-293, MCF-7 cell lines)

Futia, G. L., et al. [Quantitative image cytometry measurements of lipids, DNA, CD45 and cytokeratin for circulating tumor cell identification in a model system](#). Paper presented at: SPIE BiOS, (SPIE), (2016)(polycarbonate filters)

Guccini, I., et al. [Senescence Reprogramming by TIMP1 Deficiency Promotes Prostate Cancer Metastasis](#). Cancer Cell 39(1), 68-82 e69, (2021), DOI: 10.1016/j.ccell.2020.10.012 (prostate cancer xenografts, single molecule FISH)

Isonaka, R., et al. [Alpha-Synuclein Deposition Within Sympathetic Noradrenergic Neurons Is Associated With Myocardial Noradrenergic Deficiency in Neurogenic Orthostatic Hypotension](#). Hypertension 73(4), 910-918, (2019), DOI: 10.1161/HYPERTENSIONAHA.118.12642 (human skin biopsy)

Jiang, C., et al. [The Role of IgM Antibodies in T Cell Lymphoma Protection in a Novel Model Resembling Anaplastic Large Cell Lymphoma](#). J Immunol 206(10), 2468-2477, (2021), DOI: 10.4049/jimmunol.2001279 (mouse T cell lymphocytes)

Jin, H., et al. [Local Delivery of miR-21 Stabilizes Fibrous Caps in Vulnerable Atherosclerotic Lesions](#). Mol Ther 26(4), 1040-1055, (2018), DOI: 10.1016/j.ymthe.2018.01.011 (human atherosclerotic plaques)

- Kanie, K., et al. [Mechanistic insights into immune checkpoint inhibitor-related hypophysitis: a form of paraneoplastic syndrome](#). *Cancer Immunol Immunother*, (2021), DOI: 10.1007/s00262-021-02955-y (human pituitary tissue)
- Li, Y., et al. [Matrix metalloproteinase \(MMP\)-degradable tissue engineered periosteum coordinates allograft healing via early stage recruitment and support of host neurovasculature](#). *Biomaterials* 268, 120535, (2021), DOI: 10.1016/j.biomaterials.2020.120535 (mouse bone FFPE)
- Luo, P., et al. [An autofluorescence-based isolation of Leydig cells for testosterone deficiency treatment](#). *Mol Cell Endocrinol* 535, 111389, (2021), DOI: 10.1016/j.mce.2021.111389 (mouse testicular tissue)
- Pieknik, J. R., et al. [Herpes Simplex Virus 2 in Autonomic Ganglia: Evidence for Spontaneous Reactivation](#). *J Virol* 93(11), (2019), DOI: 10.1128/JVI.00227-19 (guinea pig dorsal root ganglia)
- Rotgers, E., et al. [Constitutive expression of Steroidogenic factor-1 \(NR5A1\) disrupts ovarian functions, fertility, and metabolic homeostasis in female mice](#). *FASEB J* 35(8), e21770, (2021), DOI: 10.1096/fj.202100304R (mouse steroidogenic tissue)
- Sorrelle, N., et al. [Improved Multiplex Immunohistochemistry for Immune Microenvironment Evaluation of Mouse Formalin-Fixed, Paraffin-Embedded Tissues](#). *J Immunol* 202(1), 292-299, (2019), DOI: 10.4049/jimmunol.1800878 (mouse spleen, tumor xenograft)
- Sun, Q., et al. [SEMA6D regulates perinatal cardiomyocyte proliferation and maturation in mice](#). *Dev Biol* 452(1), 1-7, (2019), DOI: 10.1016/j.ydbio.2019.04.013 (mouse embryo)
- Yang, J., et al. [Quenching autofluorescence in tissue immunofluorescence \[version 1; referees: 1 approved with reservations\]](#). *Wellcome Open Res* 2(79), (2017), DOI: doi: 10.12688/wellcomeopenres.12251.1 (mouse kidney, mouse placenta, human teratoma)
- Zhuang, X., et al. [Time- and Stimulus-Dependent Characteristics of Innate Immune Cells in Organ-Cultured Human Corneal Tissue](#). *J Innate Immun*, 1-14, (2021), DOI: 10.1159/000516669 (human corneal tissue)