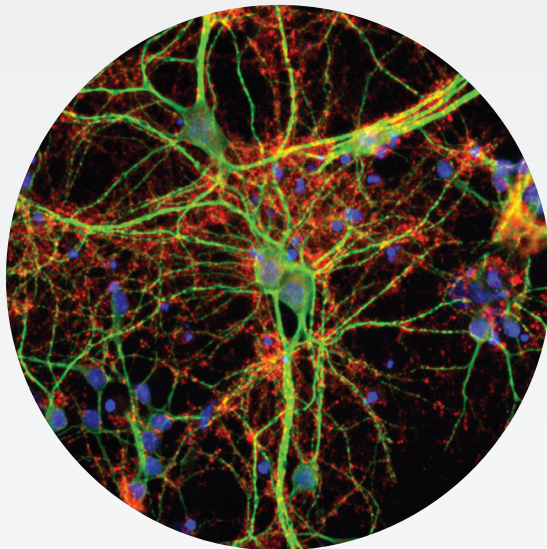
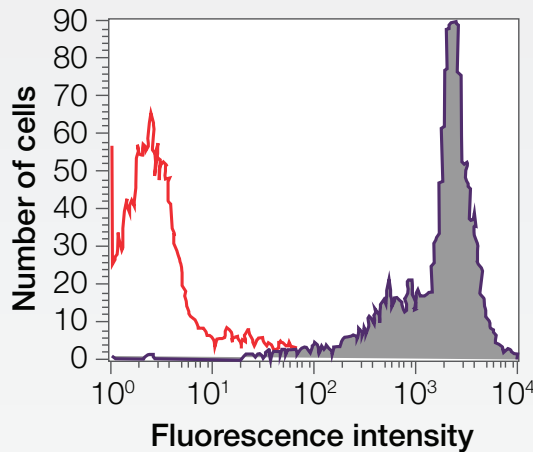


Antibodies for neuroscience research



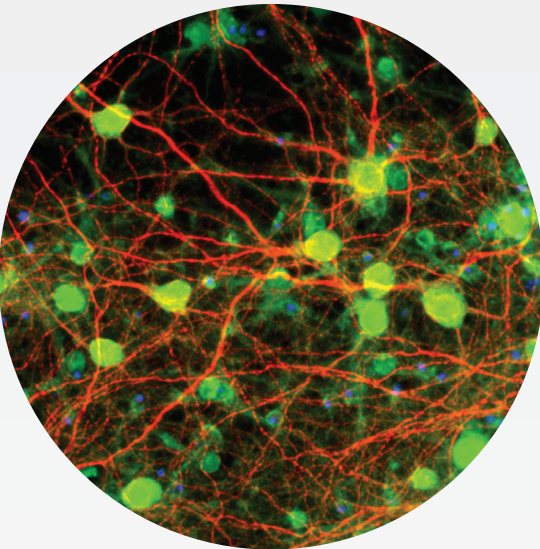
Neuron markers

Cell type	Target	Cat. No. (clone number)
Cholinergic neurons	ChAT	PA1-4710, PA1-4738, PA1-9027, PA1-18313, PA5-29653
Dopaminergic (DA) progenitor cells, DA neurons	LMX1A	710980 (20HCLC)
	Nurr1	MA1-195 (N1404)
	OTX2	701948 (14H14L5), MA5-15854 (1H12C4B5), MA5-15855 (1H12G8B2), PA5-23406, PA5-29914
	PITX3	701181 (5H10L5), 710212 (7M5HCLC), 38-2850
GABA neurons	Tyrosine hydroxylase	P21962, 701949, 710982, MA1-24654 (185)
	GABA	PA5-32241
	GAD65	398200 (ZG002), PA5-32430, PA5-22260, PA5-18369, MA5-18053 (144)
	GAD65/GAD67	PA1-84572
Glutamatergic neurons	Somatostatin	710968 (9HCLC), MA5-16987 (YGT), MA5-17182 (7G5)
	NMDAR1	700685 (1H13L3), 32-0500 (54.1), PA3-102
	NMDAR2B	71-8600, PA3-105, A-6474, MA1-2014 (NR2B), 32-0700 (B3-13B11)
	VGLUT1	48-2400
Motor neuron progenitor cells, motor neurons	VGLUT2	42-7800
	HLXB9	PA5-23407
	HOXB4	PA5-23366
	ISL1	MA5-15515 (1H9), MA5-15516 (1B1), PA5-27789
Neurons	NeuroG2	PA5-34472, PA5-11892
	NKX6.1/6.2	PA5-23070
	Synapsin 1	A6442
	Beta-3 tubulin	MA1-118 (2G10), 480011 (38F4)
	Calretinin	710559 (4HCLC), MA5-11479 (5A5), MA5-12539 (DAK Calret 1), MA514540 (SP13), PA51-6681
	DCX	MA5-17066 (2G5), 48-1200
	HuC/HuD	A21271, A21272
	MAP2	MA5-12823 (AP20), 13-1500 (M13), MA5-12826 (AP18)
	Neurofilament H, M, L, or pan	711025 (4HCLC), MA1-2012 (3G3), MA1-2010 (DA2), MA1-2011 (3H11), 13-0500 (RMO-44), 13-0700 (RMO-270), 34-1000 (RMO 14.9), 13-1300 (RMdO-20)
	PSD93	711031 (7HCLC), PA1-043, 34-4700
	PSD95	700902 (23H23L19), MA1-045 (6G6-1C9), MA1-046 (7E3-1B8), 51-6900



Glial progenitor cell markers

Target	Cat. No. (clone number)
A2B5	43-3110 (105)
CD44	701406 (19H8L4), 710413 (19HCLC), MA5-12394 (5F12), MA5-13887 (156-3C11), MA5-13890 (156-3C11), MHCD4401 (MEM-85), MHCD4404, HM5704 (IM7.8.1)

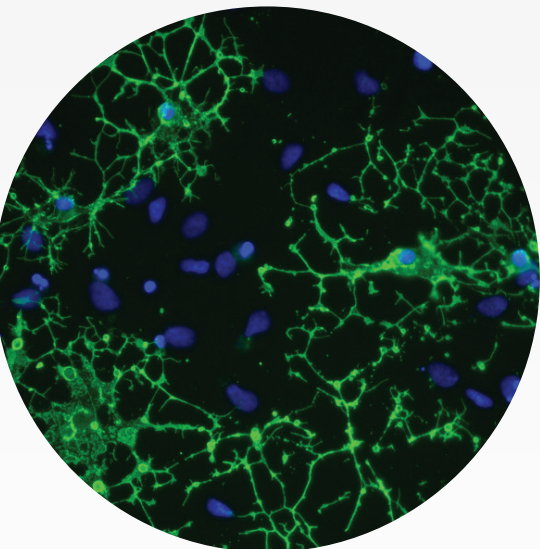


Presynaptic and postsynaptic markers

Cell type	Target	Cat. No. (clone number)
Postsynaptic	HOMER1	PA5-23023, PA5-14188
	PSD95	MA1-045 (6G6-1C9), MA1-046 (7E3-1B8), 700902 (23H23L19), 51-6900
Presynaptic	Synaptophysin	701503 (8H2L12), 710532 (8HCLC), 08-0130, 08-1130, MA5-11575 (SYP02), MA5-14532 (SP11), PA1-1043, PA5-16417
	Synaptotagmin	OSS00022W
	VGLUT1	48-2400, OSV00006W

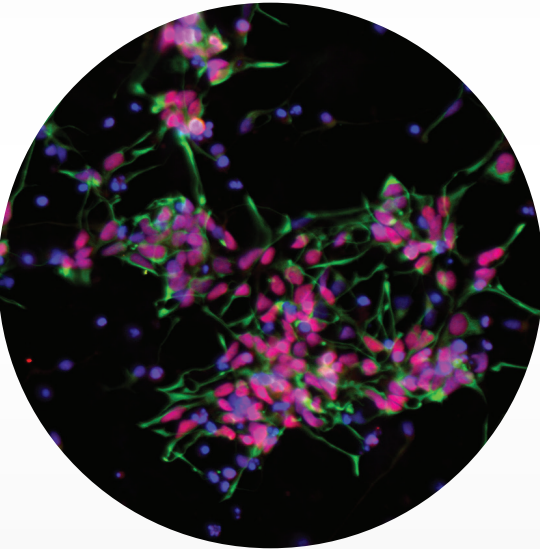
Oligodendrocyte progenitor cell and oligodendrocyte markers

Target	Cat. No. (clone number)
A2B5	43-3110 (105)
APC	710859, PA5-16883
GALC	PA5-42652, PA5-27352
MBP	PA5-16413, MA5-15922 (2H9)
MOG	PA5-19602
NG2	37-2300 (N143.8), 37-2700 (D120.43/D4.11/N143.8/N109.6), 41-2000 (LHM2)
Nogo-A	36-6600
OLIG1	PA5-21613, MA5-23954, MA5-22956
OLIG2	P21954, MA5-15810 (1G11)
PDGFRA	PA5-16571, PA5-16742, 701142 (7H13L1), 710169 (7HCLC)



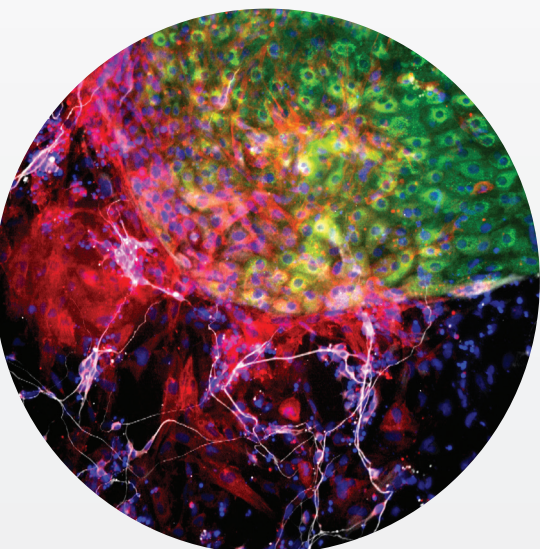
Neural stem cell markers

Target	Cat. No. (clone number)
FOXG1	PA5-26794, PA1-31654
FZD4	710731 (22HCLC), 720062
Nestin	MA1-110 (10C2)
PAX3	701147, 710178, 38-1801, PA1-107
PAX6	42-6600, MA1-109 (13B10-1A10), MA1-109D488
SOX1	PA5-23351, PA5-23370
SOX2	48-1400 (20G5), MA1-014 (20G5), MA1-014D488, MA1-014D550, MA1-014D650, MA1-014HRP, PA1-094



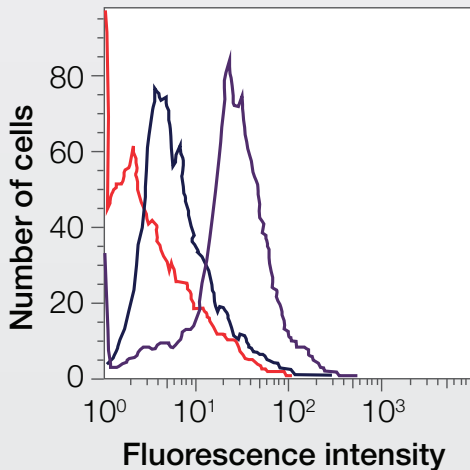
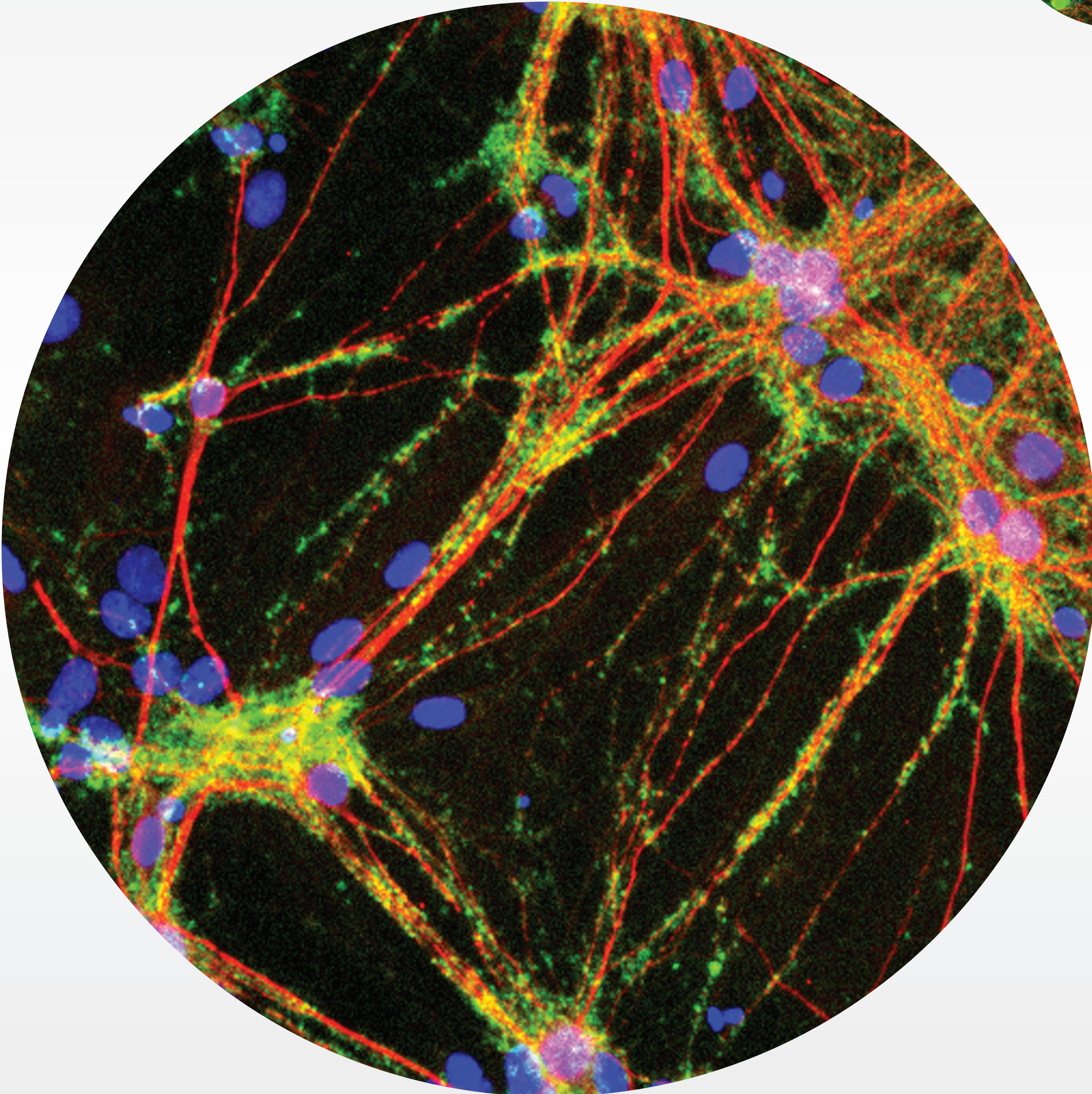
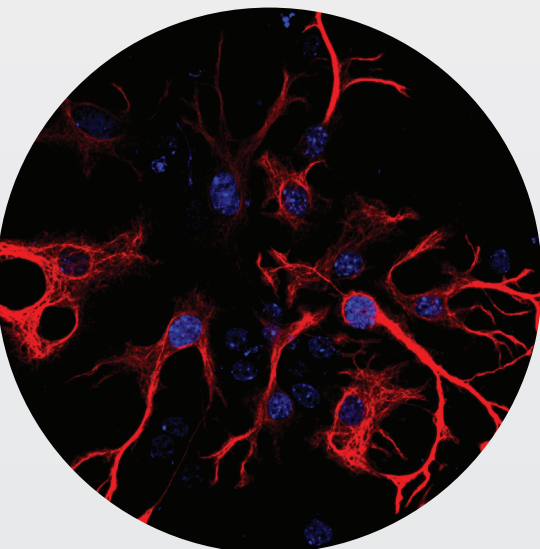
Neuroectoderm markers

Target	Cat. No. (clone number)
Beta-3 tubulin	MA1-118 (2G10)
GBX2	PA5-18561, PA5-40441
NCAM	701379 (3H15L12), 710388 (3HCLC), MA5-11563 (56C04), MA5-12685 (123C3.D5), MHCD5600 (MEM-188)
Nestin	MA1-110 (10C2)
OTX2	MA5-15854 (1H12C4B5), MA5-15855 (1H12G8B2), PA5-23406, PA5-29914
PAX3	701147, 710178, 38-1801
PAX6	42-6600, MA1-109 (13B10-1A10), MA1-109D488
SOX1	PA5-23351, PA5-23370
SOX2	48-1400 (20G5), MA1-014 (20G5), MA1-014D488, MA1-014D550, MA1-014D650, MA1-014HRP, PA1-094



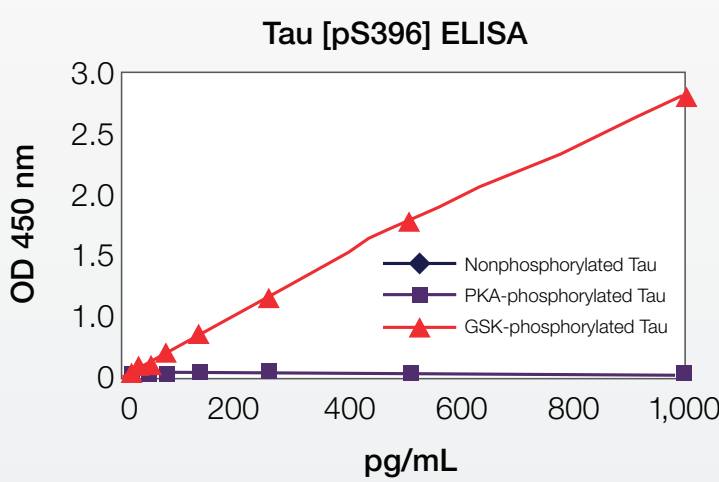
Astrocyte markers

Target	Cat. No. (clone number)
ALDH1L1	720169
EAAT2 (GLT-1)	711020, PA3-040A
GFAP	13-0300 (2.2B10), MA5-12023 (ASTRO6), PA5-16291, A21282 (131-1771), A21294, A21295
Glutamine synthetase	710963 (7HCLC)
S100beta	701340 (16H24L21), 710363 (16HCLC)



Microglia markers

Target	Cat. No. (clone number)
CD11b	CD11b00 (VIM12), RM2804 (M1/70.15)
CD68	MA5-12407 (PG-M1), MA5-13324 (KP1)
CX3CR1	PA5-19910, PA5-19935
HLA-DR	MA5-11966 (LN3), H11212 (Tu36)
IBA1	MA1-40227 (1022-5), PA5-21274, PA5-27436
TREM2	PA5-46980



Neurodegenerative disease research targets

Disease type	Target	Cat. No. (clone number)
Alzheimer's disease	Amyloid precursor protein	OMA103132 (mAbP2-1)
	APOE	701241 (16H22L18)
	Beta amyloid	MA5-11617 (BAM01 (6F/3D)), OPA1-03131
	PSEN1	MA1-752 (APS 18)
	PSEN2	MA1-754 (APS 26)
	Tau	44-738G
Amyotrophic lateral sclerosis	C9orf72	PA5-31565
	FUS	PA5-27531
	SOD1	LF-PA0013
	SQSTM1	PA5-27247
	TARDBP/TDP43	PA5-29949
Huntington's disease	HTT	710695 (3HCLC), MA1-115 (3-19), MA1-058 (4-13), 720029, PA1-003, PA1-002
Multiple sclerosis	ASIC1	PA5-28241
	IL-17A	MA1068 (3-36.2.7.3), MA1069 (3.41.2.6.1), PA1-24811
	TREM2	PA5-18763
	TRPM4	PA5-34283
Parkinson's disease	Alpha synuclein	701085 (14H2L1), MA5-15290 (2B2D1), PA5-16738
	GBA	PA5-21347, PA5-22326
	LRRK2	PA1-16770
	PARKIN	39-0900 (PRK109), PA1-751
	PINK1	MA5-11153 (38CT20.8.5), PA5-23072