

OmicsArray™ Brain and Neuronal-associated Antigen Array Antigen List - Catalog No. PA002

Antigen Number	Antigen Name	Antigen Description
1	AQP-4	Aquaporin-4 (AQP4)
2	HSPD1/Hsp60	Heat shock 60kDa protein 1 (chaperonin) (HSPD1)
3	GalC	Galactocerebroside
4	AchR3	Acetylcholine Receptor Subunit Gamma (CHRNA3)
5	GAD1/GAD67	Glutamate decarboxylase 1 (GAD1)
6	CHRM3	Cholinergic Receptor Muscarinic 3 (CHRM3)
7	MAG	Myelin Associated Glycoprotein (MAG)/Siglec-4a
8	ADGRB3/BAI3	Adhesion G protein-coupled receptor B3 (ADGRB3)
9	AMPH	Amphiphysin (AMPH)
10	GJB1	Gap Junction Protein Beta 1 (GJB1)
11	GJC2/CX47	Gap Junction Protein, gamma 2 (GJC2)/Connexin-47
12	ELAVL4/HuD	ELAV-like protein 4 (ELAVL4, HuD)
13	MMP-14	Matrix metalloproteinase-14 (MMP-14)
14	YWHAE	14-3-3 protein epsilon (YWHAE)
15	APLP-1	Amyloid beta precursor like protein 1
16	mGluR5	Metabotropic glutamate receptor 5 (mGluR5)
17	mGluR3	Metabotropic glutamate receptor 3 (mGluR3)
18	CRYAB	Alpha beta-crystallin (CRYAB)
19	GFAP	Glial Fibrillary Acidic Protein (GFAP)
20	GRID2	Orphan glutamate receptor delta2 (GRID2)
21	LGI-2	leucine-rich glioma-inactivated 2 (LGI2)
22	PNMA2	Paraneoplastic antigen Ma2 (PNMA2)
23	MAP2	Microtubule Associated Protein 2 (MAP2)
24	mGluR1/GRM1	Metabotropic glutamate receptor type 1 (mGluR1)/Glutamate receptor metabotropic 1 (GRM1)
25	PCDHA1	Protocadherin Alpha 1 (PCDHA1)
26	ITPR1	Inositol 1,4,5-trisphosphate receptor type 1 (ITPR1)
27	CADM2	Cell Adhesion Molecule 2 /CADM2
28	CHI3L1/YKL40	Chitinase-3 like-protein-1 (CHI3L1)/YKL40
29	MBP	Myelin Basic Protein

30	GRIA2	Glutamate Ionotropic Receptor AMPA Type Subunit 2 (GRIA2)
31	KCNJ10/Kir4.1	Potassium Inwardly Rectifying Channel Subfamily J Member 10 (KCNJ10)
32	OLIG2	Oligodendrocyte transcription factor (OLIG2)
33	MOBP	Myelin-associated oligodendrocyte basic protein (MOBP)
34	NEFL	Neurofilament light polypeptide (NEFL)
35	GRIN1	Glutamate receptor, ionotropic, N-methyl D-aspartate 1 (GRIN1), transcript variant NR1-1
36	S100B	S100 calcium binding protein B (S100B)
37	SYN1	Synapsin I (SYN1), transcript variant Ia
38	GAPDH	Glyceraldehyde-3-phosphate dehydrogenase (GAPDH)
39	GLUD2	Glutamate dehydrogenase 2 (GLUD2)
40	PKC-r	Protein kinase C, gamma (PRKCG)
41	SPAG16	Sperm associated antigen 16 (SPAG16)
42	EAAT2	Excitatory Amino Acid Transporter 2 (EAAT2)
43	DPP6	Dipeptidyl Peptidase 6 (DPP6)
44	PLP1	Myelin proteolipid protein 1 (PLP1)
45	HOMER3	Homer Homolog 3 (Drosophila) (HOMER3)
46	RBPJ	Recombination Signal Binding Protein For Immunoglobulin Kappa J Region (RBPJ)
47	REEP2	Receptor Accessory Protein 2 (REEP2)
48	UCH-L1	Ubiquitin Carboxyl-Terminal Hydrolase Isozyme L1 (UCH-L1)/UCH-L1
49	PMP2	Myelin P2 Protein/Peripheral myelin protein 2 (PMP2)
50	VILIP/VSNL1	Visinin-Like Protein 1 (VILIP)/VSNL1
51	NPTX1	Neuronal Pentraxin-1 (NPTX1)
52	HepaCAM	Hepatocyte Cell Adhesion Molecule/HepaCAM
53	LY6H	Lymphocyte Antigen 6H/LY6H
54	BCAN/BEHAB	Brevican Core Protein/BCAN/BEHAB
55	SULT4A1	Sulfotransferase 4A1 (SULT4A1)
56	DLG4/PSD95/SAP90	Disks Large Homolog 4 (DLG4)/SAP90//PSD95
57	SNAP-a/NAPA	Alpha-soluble NSF attachment protein 1 (ASnap1)/SNAP-a/NAPA
58	TALDO1	Transaldolase 1 (TALDO1)
59	LRRN2/LRRTM2	Leucine-rich repeat-containing protein 2 (LRRN2)
60	Contactin-2 (CNTN2)/TAG-1	Contactin-2 (CNTN2)/TAG-1
61	NCAM-1/CD56	Neural Cell Adhesion Molecule 1 (NCAM-1)/CD56
62	SLC17A6	Solute Carrier Family 17 (Vesicular Glutamate Transporter)
63	Lactoferrin	Lactoferrin Native Protein

64	Kappa Light Chains	Kappa Light Chains (Free) from human myeloma plasma
65	GJA1/CX43	Gap junction alpha-1 (GJA1) / connexin 43 (Cx43)
66	ELAVL3/HuC	ELAV Like RNA Binding Protein 3 (ELAVL3)/HuC
67	NOVA2	Neuro-oncological ventral antigen 2 (NOVA2)
68	OSP	Oligodendrocyte specific protein (OSP)
69	LTBP3	Latent-transforming growth factor beta-binding protein 3 (LTBP3)
70	AchRm2	M2 Muscarinic Acetylcholine Receptor
71	Amyloid B-peptide	Amyloid b-peptide (1-42)
72	PGLYRP1/PGRP-S	Peptidoglycan recognition protein 1 (PGLYRP1)/PGRP-S
73	Lipocalin-2/NGAL	Lipocalin-2/NGAL Protein
74	Profilin 1	Profilin 1 Protein
75	Transketolase	Transketolase Protein
76	S100A8/A9	Alarmin S100A8/A9 Heterodimer
77	Cytokeratin 10 (CK10)	Cytokeratin 10 (CK10)
78	Coronin-1	Coronin-1
79	Plastin-2/LCP1	Plastin-2/LCP1
80	Myosin-9 (MYH9)	Myosin-9 (MYH9)
81	S100A12	S100 calcium-binding protein A12 (S100A12)
82	Resistin	Resistin
83	CAP1/CAP2	Adenylyl cyclase-associated protein 1 and 2 (CAP1/CAP2)
84	RCV1/RCVRN	Recoverin (RCVRN)
85	CRMP5/CV-2	Collapsin response mediator protein 5 (CRMP5)/CV-2
86	CDR2	Cerebellar degeneration-related protein 2 (CDR2)
87	GAD2/GAD65	Glutamate decarboxylase (GAD2/GAD65)
88	NOVA1	Onconeural antigen Nova-1
89	PNMA1	Paraneoplastic Ma proteins (PNMA1)
90	ELAVL1/HuR (Hu A)	ELAV-like protein 1/ELAVL1
91	ERP29	Endoplasmic reticulum protein 29 (EPR29)
92	HSPA5 (GRP78)	Heat Shock Protein Family A (Hsp70) Member 5 · GRP78
93	SOX1	Sry-like high mobility group box 1 (SOX1)
94	Titin (TTN)	Titin (TTN)
95	EBNA1BP2	EBNA1 binding protein (Ebna1bp2)
96	ADCY	Adenylate Cyclase (ADCY 1-2-4-5)
97	ELAVL2/HuB	ELAV Like RNA Binding Protein 2 (ELAVL2)/HuB

98	Neurofascin	Neurofascin Protein
99	mGluR2	Metabotropic glutamate receptor type 2 (mGluR2)
100	NG2/MCSP/CSPG4	Neural/glial antigen 2 (NG2)/NG2 proteoglycan
101	Caspr2	Contactin-associated protein 2 (CASPR2)
102	MNDA	Myeloid cell nuclear differentiation antigen (MNDA)
103	ANXA1	Annexin A1(ANXA1)
104	TSHR	Thyroid Stimulating Hormone Receptor (TSHR)
105	ZIC4	Zic Family Member 4 (ZIC4)
106	MOG	Myelin Oligodendrocyte Glycoprotein
107	POLR3D	POLR3D (RNA Polymerase III Subunit D)
108	SepSecS/SLA/LP	SEPSECS (O-phosphoseryl-tRNA(Sec) selenium transferase)
109	FARS2	phenylalanyl-tRNA synthetase 2 (FARS2)
110	RPP25	RPP25 (Th/To complex)
111	OLFM4	Olfactomedin-4
112	NRP1	Neuropilin-1 (NRP1)
113	MARCO	Macrophage receptor with collagenous structure
114	EGFR	Human EGFR/HER1/ErbB1 protein
115	MuSk (ECD)	MuSK (extra-cellular domain)
116	MuSk (ICD)	MuSK (intra-cellular domain)
117	NMDAR1-M1	NMDAR1 -M1
118	NMDAR1-S1S2	NMDAR1-S1-GT-S2
119	Ganglioside	Ganglioside mixture
120	Sphingomyelin	Sphingomyelin
121	Ig control	Ig control 1
122	Ig control	Ig control 2
123	Ig control	Ig control 3
124	Ig control	Ig control 4
125	anti-Ig control	anti-Ig control 1
126	anti-Ig control	anti-Ig control 2
127	anti-Ig control	anti-Ig control 3
128	anti-Ig control	anti-Ig control 4