

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

Antigen number	Antigen name	Description
1	14-3-3 protein	14-3-3 protein alpha/beta/epsilon
2	AchR3	Acetylcholine Receptor Subunit Gamma(CHRNG)
3	AchRm2	M2 Muscarinic Acetylcholine Receptor
4	ADCY (mix)	Adenylate Cyclase (ADCY 1,2,4,5 mix)
5	ADGRB3/BAI3	Adhesion G protein-coupled receptor B3 (ADGRB3)
6	AMPH	Amphiphysin (AMPH)
7	Amyloid B-peptide	Amyloid b-peptide (1-42)
8	ANXA1	Annexin A1(ANXA1), 20ug
9	APLP-1	Amyloid beta precursor like protein 1
10	AQP-4	Aquaporin-4 (AQP4)
11	BCAN/BEHAB	Brevican Core Protein/BCAN/BEHAB
12	CADM2	Cell Adhesion Molecule 2 /CADM2
13	CAP1/CAP2	Adenylyl cyclase-associated protein 1 and 2 (CAP1/CAP2)
14	Caspr2	contactin-associated protein-like 2 (Caspr2)
15	CDR2	Cerebellar degeneration-related protein 2 (CDR2)
16	CHI3L1/YKL40	Chitinase-3 like-protein-1 (CHI3L1)/YKL40
17	CHRM3	Cholinergic Receptor Muscarinic 3 (CHRM3)
18	Contactin-2/CNTN2/TAG-	Contactin-2 (CNTN2)/TAG-1
19	Coronin-1	Coronin-1
20	CRMP5/CV-2	collapsin response mediator protein 5 (CRMP5)/CV-2
21	CRYAB	Alpha beta-crystallin (CRYAB)
22	Cytokeratin 10 (CK10)	Cytokeratin 10
23	DLG4/PSD95/SAP90	Disks Large Homolog 4 (DLG4)/SAP90//PSD95
24	DPP6	Dipeptidyl Peptidase 6 (DPP6)
25	EAAT2	Excitatory Amino Acid Transporter 2 (EAAT2)
26	EBNA1 binding protein 2	EBNA1 binding protein (Ebna1bp2)
27	ELAVL2/HuB	ELAV Like RNA Binding Protein 2 (ELAVL2)/HuB
28	ELAVL3/HuC	ELAV Like RNA Binding Protein 3 (ELAVL3)/HuC
29	ELAVL4/HuD	ELAV-like protein 4 (ELAVL4, HuD)

30	ERP29	Endoplasmic reticulum protein 29 (EPR29)
31	FARS2	phenylalanyl-tRNA synthetase 2 (FARS2)
32	GAD1/GAD67	Glutamate decarboxylase 1 (GAD1)
33	GAD2/GAD65	Glutamate decarboxylase (GAD2/GAD65)
34	GalC	galactocerebroside
35	GAPDH	Glyceraldehyde-3-phosphate dehydrogenase (GAPDH)
36	GFAP	Glial Fibrillary Acidic Protein (GFAP)
37	GJA1/CX43	Gap junction alpha-1 (GJA1) / connexin 43 (Cx43)
38	GJB1	Gap Junction Protein Beta 1 (Gjb1)
39	GJB6/CX30	Gap junction beta-6 (GJB6), connexin 30 (Cx30)
40	GJC2/CX47	Gap Junction Protein, gamma 2 (GJC2)/Connexin-47
41	GJC3/GJE1/CX29	Gap junction gamma-3 (GJC3), connexin-29 (Cx29) or gap junction epsilon-1 (GJE1)
42	GLUD2	Glutamate dehydrogenase 2 (GLUD2)
43	GM1, GD1a, GD1b, GT1b	Ganglioside mixture
44	GM2 ganglioside	Monosialoganglioside GM2
45	GRIA2	Glutamate Ionotropic Receptor AMPA Type Subunit 2 (GRIA2)
46	GRID2	Orphan glutamate receptor delta2 (GRID2)
47	GRIN1	Glutamate receptor, ionotropic, N-methyl D-aspartate 1 (GRIN1), transcript variant NR1-1
48	GRM1	Glutamate receptor metabotropic 1 (GRM1)
49	HepaCAM	Hepatocyte Cell Adhesion Molecule/HepaCAM
50	HOMER3	Homer Homolog 3 (Drosophila) (HOMER3) (AA 1-361) protein
51	HSPA5	HSPA5/GRP78
52	HSPD1/Hsp60	Heat shock 60kDa protein 1 (chaperonin) (HSPD1)
53	HuR (Hu A)	ELAV-like protein 1/ELAVL1
54	ITPR1	Inositol 1,4,5-trisphosphate receptor type 1 (ITPR1)
55	Kappa Light Chains	Kappa Light Chains (Free) from human myeloma plasma
56	KCNJ10/Kir4.1	Potassium Inwardly Rectifying Channel Subfamily J Member 10 (KCNJ10)
57	Lactoferrin	Lactoferrin Native Protein
58	LGI-2	leucine-rich glioma-inactivated 2 (LGI2)
59	Lipocalin-2/NGAL	Lipocalin-2/NGAL Protein
60	LRRN2/LRRTM2	Leucine-rich repeat-containing protein 2 (LRRN2)
61	LTBP3	Latent-transforming growth factor beta-binding protein 3 (LTBP3)
62	LY6H	Lymphocyte Antigen 6H/LY6H
63	MAG	Myelin Associated Glycoprotein (MAG)/Siglec-4a

64	MAP2	Microtubule Associated Protein 2 (MAP2)
65	mGluR1/GRM1	Metabotropic glutamate receptor type 1 (mGluR1)
66	mGluR2	Metabotropic glutamate receptor type 2 (mGluR2)
67	mGluR3	Metabotropic glutamate receptor 3 (mGlu3)
68	mGluR5	Metabotropic glutamate receptor 5 (mGluR5)
69	MMP-14	Matrix metalloproteinase-14 (MMP-14)
70	MNDA	Myeloid cell nuclear differentiation antigen (MNDA)
71	MOBP	Myelin-associated oligodendrocyte basic protein (MOBP)
72	MOG	Myelin Oligodendrocyte Glycoprotein
73	Myosin-9 (MYH9)	Myosin-9 (MYH9)
74	NCAM-1/CD56	Neural Cell Adhesion Molecule 1 (NCAM-1)/CD56
75	NEFL	Neurofilament light polypeptide (NEFL)
76	Neurofascin	Neurofascin Protein
77	NG2/MCSP/CSPG4	Neural/glial antigen 2 (NG2)/NG2 proteoglycan
78	NOVA1	Onconeural antigen Nova-1
79	NOVA2	neuro-oncological ventral antigen 2 (NOVA2)
80	NPTX1	Neuronal Pentraxin-1 (NPTX1)
81	NRP1	Neuropilin-1 (NRP1)
82	OLFM4	Olfactomedin-4
83	OLIG2	Oligodendrocyte transcription factor (OLIG2)
84	OSP	Oligodendrocyte specific protein (OSP)
85	PCDHA1	Protocadherin Alpha 1 (PCDHA1)
86	PGLYRP1/PGRP-S	Peptidoglycan recognition protein 1 (PGLYRP1)/PGRP-S
87	Phosphatidylinositol (PL)	Phosphatidylinositol
88	Plastin-2/LCP1	Plastin-2/LCP1
89	PLP1	Myelin proteolipid protein 1 (PLP1)
90	PMP2	Myelin P2 Protein/Peripheral myelin protein 2 (PMP2)
91	PNMA1 (anti-Ma)	Paraneoplastic Ma proteins (PNMA1)
92	PNMA2 (Anti-Ta)	Paraneoplastic antigen Ma2 (PNMA2)
93	POLR3D	POLR3D (RNA Polymerase III Subunit D)
94	PRKCG	Protein kinase C, gamma (PRKCG)
95	Profilin 1	Profilin 1 Protein
96	RBPJ	Recombination Signal Binding Protein For Immunoglobulin Kappa J Region (RBPJ) (Transcript Variant 3)
97	RCV1/RCVRN	Recoverin (RCVRN)

98	REEP2	Receptor Accessory Protein 2 (REEP2) (Myc-DYKDDDDK tag)
99	Resistin	Resistin
100	RPP25	RPP25 (Th/To complex)
101	S100A12	S100 calcium-binding protein A12 (S100A12)
102	S100A8/A9	Alarmin S100A8/A9 Heterodimer
103	S100B	S100 calcium binding protein B (S100B)
104	SepSecS or SLA/LP	SEPSECS (O-phosphoserine-tRNA(Sec) selenium transferase)
105	SLC17A6	Solute Carrier Family 17 (Vesicular Glutamate Transporter), Member 6 (SLC17A6) (AA 1-582)
106	SNAP-a/NAPA	Alpha-soluble NSF attachment protein 1 (ASNA1)/SNAP-a/NAPA
107	SOX1	Sry-like high mobility group box 1 (SOX1)
108	SPAG16	Sperm associated antigen 16 (SPAG16), transcript variant 1
109	Sphingomyelin	Sphingomyelin
110	Sulfatide	Sulfated galactocerebroside
111	SULT4A1	Sulfotransferase 4A1 (SULT4A1)
112	SYN1	Synapsin I (SYN1), transcript variant Ia
113	SYP	Synaptophysin (SYP)
114	TALDO1	Transaldolase 1 (TALDO1)
115	Titin (TTN)	Titin (TTN)
116	Transketolase	Transketolase Protein
117	TSHR	Thyroid Stimulating Hormone Receptor (TSHR)
118	UCH-L1	Ubiquitin Carboxyl-Terminal Hydrolase Isozyme L1 (UCH-L1)/UCH-L1
119	VILIP/VSNL1	Visinin-Like Protein 1 (VILIP)/VSNL1
120	ZIC4	Zic Family Member 4 (ZIC4)
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