

OmicsArray™ Brain and Neuronal-associated Antigen Array Antigen List - Catalog No. PA002			
Antigen Number	Antigen Name	Antigen Description	UniProt ID
1	AQP-4	Aquaporin-4 (AQP4)	<a href="#">P55087</a>
2	HSPD1/Hsp60	Heat shock 60kDa protein 1 (chaperonin) (HSPD1)	<a href="#">P10809</a>
3	GALC	Galactocerebrosidase	<a href="#">P54803</a>
4	CHRNA3/AChR3	Acetylcholine Receptor Subunit Gamma (CHRNA3)	<a href="#">P07510</a>
5	GAD1/GAD67	Glutamate decarboxylase 1 (GAD1)	<a href="#">Q99259</a>
6	CHRM3	Cholinergic Receptor Muscarinic 3 (CHRM3)	<a href="#">P20309</a>
7	MAG	Myelin Associated Glycoprotein (MAG)/Siglec-4a	<a href="#">P20916</a>
8	ADGRB3/BAI3	Adhesion G protein-coupled receptor B3 (ADGRB3)	<a href="#">O60242</a>
9	AMPH	Amphiphysin (AMPH)	<a href="#">P49418</a>
10	GJB1	Gap Junction Protein Beta 1 (GJB1)	<a href="#">P08034</a>
11	GJC2/CX47	Gap Junction Protein, gamma 2 (GJC2)/Connexin-47	<a href="#">Q5T442</a>
12	ELAVL4/HuD	ELAV-like protein 4 (ELAVL4, HuD)	<a href="#">P26378</a>
13	MMP-14	Matrix metalloproteinase-14 (MMP-14)	<a href="#">P50281</a>
14	YWHAE	14-3-3 protein epsilon (YWHAE)	<a href="#">P62258</a>
15	APLP-1	Amyloid beta precursor like protein 1	<a href="#">P51693</a>
16	mGluR5	Metabotropic glutamate receptor 5 (mGluR5)	<a href="#">P41594</a>
17	mGluR3	Metabotropic glutamate receptor 3 (mGluR3)	<a href="#">Q14832</a>
18	CRYAB	Alpha beta-crystallin (CRYAB)	<a href="#">P02511</a>
19	GFAP	Glial Fibrillary Acidic Protein (GFAP)	<a href="#">P14136</a>
20	GRID2	Orphan glutamate receptor delta2 (GRID2)	<a href="#">O43424</a>
21	LGI-2	leucine-rich glioma-inactivated 2 (LGI2)	<a href="#">Q8NOV4</a>
22	PNMA2	Paraneoplastic antigen Ma2 (PNMA2)	<a href="#">Q9UL42</a>
23	MAP2	Microtubule Associated Protein 2 (MAP2)	<a href="#">P11137</a>
24	mGluR1/GRM1	Metabotropic glutamate receptor type 1 (mGluR1)/Glutamate receptor metabotropic 1 (GRM1)	<a href="#">Q13255</a>
25	PCDHA1	Protocadherin Alpha 1 (PCDHA1)	<a href="#">Q9Y5I3</a>
26	ITPR1	Inositol 1,4,5-trisphosphate receptor type 1 (ITPR1)	<a href="#">Q14643</a>
27	CADM2	Cell Adhesion Molecule 2 /CADM2	<a href="#">Q8N3J6</a>
28	CHI3L1/YKL40	Chitinase-3 like-protein-1 (CHI3L1)/YKL40	<a href="#">P36222</a>
29	MBP	Myelin Basic Protein	<a href="#">P02686</a>
30	GRIA2	Glutamate Ionotropic Receptor AMPA Type Subunit 2 (GRIA2)	<a href="#">P42262</a>
31	KCNJ10/Kir4.1	Potassium Inwardly Rectifying Channel Subfamily J Member 10 (KCNJ10)	<a href="#">P78508</a>
32	OLIG2	Oligodendrocyte transcription factor (OLIG2)	<a href="#">Q13516</a>
33	MOBP	Myelin-associated oligodendrocyte basic protein (MOBP)	<a href="#">Q13875</a>
34	NEFL	Neurofilament light polypeptide (NEFL)	<a href="#">P07196</a>
35	GRIN1	Glutamate receptor, ionotropic, N-methyl D-aspartate 1 (GRIN1), transcript variant NR1-1	<a href="#">Q05586</a>
36	S100B	S100 calcium binding protein B (S100B)	<a href="#">P04271</a>
37	SYN1	Synapsin I (SYN1), transcript variant Ia	<a href="#">P17600</a>
38	GAPDH	Glyceraldehyde-3-phosphate dehydrogenase (GAPDH)	<a href="#">P04406</a>
39	GLUD2	Glutamate dehydrogenase 2 (GLUD2)	<a href="#">P49448</a>
40	PKC-r	Protein kinase C, gamma (PRKCG)	<a href="#">P05129</a>

41	SPAG16	Sperm associated antigen 16 (SPAG16)	<a href="#">Q8N0X2</a>
42	EAAT2	Excitatory Amino Acid Transporter 2 (EAAT2)	<a href="#">P43004</a>
43	DPP6	Dipeptidyl Peptidase 6 (DPP6)	<a href="#">P42658</a>
44	PLP1	Myelin proteolipid protein 1 (PLP1)	<a href="#">P60201</a>
45	HOMER3	Homer Homolog 3 (Drosophila) (HOMER3)	<a href="#">Q9NSC5</a>
46	RBPJ	Recombination Signal Binding Protein For Immunoglobulin Kappa J Region (RBPJ)	<a href="#">Q06330</a>
47	REEP2	Receptor Accessory Protein 2 (REEP2)	<a href="#">Q9BRK0</a>
48	UCH-L1	Ubiquitin Carboxyl-Terminal Hydrolase Isozyme L1 (UCH-L1)/UCH-L1	<a href="#">P09936</a>
49	PMP2	Myelin P2 Protein/Peripheral myelin protein 2 (PMP2)	<a href="#">P02689</a>
50	VSNL1/VILIP	Visinin-Like Protein 1 (VILIP)/VSNL1	<a href="#">P62760</a>
51	NPTX1	Neuronal Pentraxin-1 (NPTX1)	<a href="#">Q15818</a>
52	HepaCAM	Hepatocyte Cell Adhesion Molecule/HepaCAM	<a href="#">Q14CZ8</a>
53	LY6H	Lymphocyte Antigen 6H/LY6H	<a href="#">O94772</a>
54	BCAN/BEHAB	Brevican Core Protein/BCAN/BEHAB	<a href="#">Q96GW7</a>
55	SULT4A1	Sulfotransferase 4A1 (SULT4A1)	<a href="#">Q9BR01</a>
56	DLG4/PSD95/SAP90	Disks Large Homolog 4 (DLG4)/SAP90//PSD95	<a href="#">P78352</a>
57	SNAP-a/NAPA	Alpha-soluble NSF attachment protein 1 (ASNA1)/SNAP-a/NAPA	<a href="#">P54920</a>
58	TALDO1	Transaldolase 1 (TALDO1)	<a href="#">P37837</a>
59	LRRN2/LRRTM2	Leucine-rich repeat-containing protein 2 (LRRN2)	<a href="#">Q9BYS8</a>
60	Contactin-2 (CNTN2)/TAG-1	Contactin-2 (CNTN2)/TAG-1	<a href="#">Q02246</a>
61	NCAM-1/CD56	Neural Cell Adhesion Molecule 1 (NCAM-1)/CD56	<a href="#">P13591</a>
62	SLC17A6	Solute Carrier Family 17 (Vesicular Glutamate Transporter)	<a href="#">Q9P2U8</a>
63	Lactoferrin	Lactoferrin Native Protein	<a href="#">P02788</a>
64	Kappa Light Chains	Kappa Light Chains (Free) from human myeloma plasma	<a href="#">P0DOX7</a>
65	GJA1/CX43	Gap junction alpha-1 (GJA1) / connexin 43 (Cx43)	<a href="#">P17302</a>
66	ELAVL3/HuC	ELAV Like RNA Binding Protein 3 (ELAVL3)/HuC	<a href="#">Q14576</a>
67	NOVA2	Neuro-oncological ventral antigen 2 (NOVA2 ) /RNA-binding protein Nova-2	<a href="#">Q9UNW9</a>
68	Claudin-11/OSP	Claudin-11/Oligodendrocyte specific protein (OSP)	<a href="#">O75508</a>
69	LTBP3	Latent-transforming growth factor beta-binding protein 3 (LTBP3)	<a href="#">Q9NS15</a>
70	CHRM2/AchRm2	Muscarinic acetylcholine receptor M2	<a href="#">P08172</a>
71	Amyloid B-peptide	Amyloid b-peptide (1-42)	<a href="#">P86906</a>
72	PGLYRP1/PGRP-S	Peptidoglycan recognition protein 1 (PGLYRP1)/PGRP-S	<a href="#">O75594</a>
73	Lipocalin-2/NGAL	Lipocalin-2/NGAL Protein	<a href="#">P80188</a>
74	Profilin 1	Profilin 1 Protein	<a href="#">P07737</a>
75	TKT	Transketolase Protein	<a href="#">P29401</a>
76	S100A8/A9	Alarmin S100A8/A9 Heterodimer	<a href="#">P05109,</a> <a href="#">P06702</a>
77	Cytokeratin 10 (CK10)	Cytokeratin 10 (CK10)	<a href="#">P13645</a>
78	Coronin-1	Coronin-1	<a href="#">P31146</a>
79	Plastin-2/LCP1	Plastin-2/LCP1	<a href="#">P13796</a>
80	Myosin-9 (MYH9)	Myosin-9 (MYH9)	<a href="#">P35579</a>
81	S100A12	S100 calcium-binding protein A12 (S100A12)	<a href="#">P80511</a>
82	Resistin	Resistin	<a href="#">Q9HD89</a>

83	CAP1/CAP2	Adenylyl cyclase-associated protein 1 and 2 (CAP1/CAP2)	<a href="#">Q01518</a> , <a href="#">P40123</a>
84	RCV1/RCVRN	Recoverin (RCVRN)	<a href="#">P35243</a>
85	CRMP5/CV-2	Collapsin response mediator protein 5 (CRMP5)/CV-2	<a href="#">Q9BPU6</a>
86	CDR2	Cerebellar degeneration-related protein 2 (CDR2)	<a href="#">Q01850</a>
87	GAD2/GAD65	Glutamate decarboxylase (GAD2/GAD65)	<a href="#">Q04329</a>
88	NOVA1	Onconeural antigen Nova-1/RNA-binding protein Nova-1	<a href="#">P51513</a>
89	PNMA1	Paraneoplastic Ma proteins (PNMA1)	<a href="#">Q8ND90</a>
90	ELAVL1/HuR (Hu A)	ELAV-like protein 1/ELAVL1	<a href="#">Q15717</a>
91	ERP29	Endoplasmic reticulum protein 29 (EPR29)	<a href="#">P30040</a>
92	HSPA5 (GRP78)	Heat Shock Protein Family A (Hsp70) Member 5 - GRP78	<a href="#">P11021</a>
93	SOX1	Sry-like high mobility group box 1 (SOX1)	<a href="#">O00570</a>
94	Titin (TTN)	Titin (TTN)	<a href="#">Q8WZ42</a>
95	EBNA1BP2	EBNA1 binding protein (Ebna1bp2)	<a href="#">Q99848</a>
96	ADCY	Adenylate Cyclase 1, 2, 4, and 5	<a href="#">Q08828</a> ,
			<a href="#">Q08462</a> ,
			<a href="#">Q8NFM4</a> ,
			<a href="#">Q95622</a>
97	ELAVL2/HuB	ELAV Like RNA Binding Protein 2 (ELAVL2)/HuB	<a href="#">Q12926</a>
98	Neurofascin	Neurofascin Protein	<a href="#">Q94856</a>
99	mGluR2	Metabotropic glutamate receptor type 2 (mGluR2)	<a href="#">Q14416</a>
100	NG2/MCSP/CSPG4	Neural/glial antigen 2 (NG2)/NG2 proteoglycan	<a href="#">Q6UVK1</a>
101	Caspr2	Contactin-associated protein 2 (CASPR2)	<a href="#">Q9UHC6</a>
102	MNDA	Myeloid cell nuclear differentiation antigen (MNDA)	<a href="#">P41218</a>
103	ANXA1	Annexin A1(ANXA1)	<a href="#">P04083</a>
104	TSHR	Thyroid Stimulating Hormone Receptor (TSHR)	<a href="#">P16473</a>
105	ZIC4	Zic Family Member 4 (ZIC4)	<a href="#">Q8N9L1</a>
106	MOG	Myelin Oligodendrocyte Glycoprotein	<a href="#">Q16653</a>
107	POLR3D	POLR3D (RNA Polymerase III Subunit D)	<a href="#">P05423</a>
108	SepSecS/SLA/LP	SEPSECS (O-phosphoseryl-tRNA(Sec) selenium transferase)	<a href="#">Q9HD40</a>
109	FARS2	phenylalanyl-tRNA synthetase 2 (FARS2)/Phenylalanine --tRNA ligase, mitochondrial	<a href="#">Q95363</a>
110	RPP25	Ribonuclease P protein subunit p25 (Th/To complex)	<a href="#">Q9BUL9</a>
111	OLFM4	Olfactomedin-4	<a href="#">Q6UX06</a>
112	NRP1	Neuropilin-1 (NRP1)	<a href="#">O14786</a>
113	MARCO	Macrophage receptor with collagenous structure	<a href="#">Q9UEW3</a>
114	EGFR	Human EGFR/HER1/ErbB1 protein	<a href="#">P00533</a>
115	MuSk (ECD)	MuSK (extra-cellular domain)	<a href="#">Q15146</a>
116	MuSK (ICD)	MuSK (intra-cellular domain)	<a href="#">Q15146</a>
117	NMDAR1-M1	Glutamate receptor ionotropic, NMDA 1/N-methyl-D-aspartate receptor subunit NR1 (NMD-R1)	<a href="#">Q05586</a>
118	NMDAR1-S1S2	Glutamate receptor ionotropic, NMDA 1/N-methyl-D-aspartate receptor subunit NR1 (NMD-R1)	<a href="#">Q05586</a>
119	Ganglioside	Ganglioside Mixture, Bovine Brain	<a href="#">345717-M</a>
120	Sphingomyelin	Sphingomyelin, Bovine Spinal Cord	<a href="#">SKU560076</a>
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	