

**OmicsArray™ Paraneoplastic Neurological Syndrome (PNS) Antigen Array, 34-Plex - Catalog No. PA024**

Seial#	Antigen ID	Antigen Name	UniProt ID	Symbol	Associated Diseases
1	AMPH	Amphiphysin (AMPH)	<a href="#">P49418</a>	AMPH	Paraneoplastic neurological syndrome (PNS); Stiff-person syndrome (SPS); Limbic encephalitis; Brainstem encephalitis; Myelopathy; Peripheral neuropathy
2	AQP4	Aquaporin-4	<a href="#">P55087</a>	AQP4	Neuromyelitis optica spectrum disorder (NMOSD)
3	CASPR2	Contactin-associated protein 2 (CASPR2)	<a href="#">Q9JHC6</a>	CNTNAP2	Autoimmune encephalitis; Neuroomyotonia (Morvan's syndrome); Neuropathic pain
4	CDR2	Cerebellar degeneration-related protein 2 (CDR2)	<a href="#">Q01850</a>	CDR2	Paraneoplastic cerebellar degeneration (PCD)
5	CRMP5/DPYSL5/CV2	Collapsin response mediator protein 5 (CRMP5)/Dihydropyrimidinase-related protein 5	<a href="#">Q98PU6</a>	DPYSL5	Paraneoplastic neurological syndrome (PNS)
6	CSPG4/NG2/MCSP	Chondroitin sulfate proteoglycan 4 (CSPG4)/Chondroitin sulfate proteoglycan NG2/Melanoma chondroitin sulfate proteoglycan (MCSP)	<a href="#">Q6LJK1</a>	CSPG4	Multiple Sclerosis (MS)
7	ELAVL1/HuR (Hu A)	ELAV-like protein 1/ELAVL1	<a href="#">Q15717</a>	ELAVL1	Paraneoplastic neurological syndromes
8	ELAVL2/HuB	ELAV-LIKE Protein 2 (ELAVL2)/Hu-antigen B (HuB)	<a href="#">Q12926</a>	ELAVL2	Paraneoplastic neurological syndromes (PNS), particularly those associated with small-cell lung cancer (SCLC).
9	ELAVL3/HuC	ELAV-Like Protein 3 (ELAVL3)/Hu-antigen C (HuC)	<a href="#">Q14576</a>	ELAVL3	Paraneoplastic encephalomyelitis; sensory neuronopathy (PEM/PSN), amyotrophic lateral sclerosis (ALS).
10	ELAVL4/HuD	ELAV-like protein 4 (ELAVL4)/Hu-antigen D (HuD)	<a href="#">P26378</a>	ELAVL4	Paraneoplastic neurological syndromes (PNS) with small cell lung cancer (SCLC)
11	GAD2/GAD65	Glutamate decarboxylase 2 (GAD2/65 kDa glutamic acid decarboxylase (GAD-65)	<a href="#">Q05329</a>	GAD2	Type 1 Diabetes; Stiff-Person Syndrome; autoimmune encephalitis, cerebellar ataxia, and certain types of epilepsy.
12	GJA1/CX43	Gap junction alpha-1 (GJA1) / connexin 43 (Cx43)	<a href="#">P17302</a>	GJA1	Neuromyelitis Optica Spectrum Disorder (NMOSD); general immune-mediated diseases
13	GLUD2	Glutamate dehydrogenase 2, mitochondrial	<a href="#">P49448</a>	GLUD2	Opsoclonus-Myoclonus Syndrome (OMS)
14	GluN1/NMDAR1	Glutamate receptor ionotropic, NMDA 1/N-methyl-D-aspartate receptor subunit NR1 (NMD-R1)	<a href="#">Q05586</a>	GRIN1	anti-NMDAR encephalitis; NPSLE
15	GluR-2	Glutamate receptor 2/AMPA-selective glutamate receptor 2	<a href="#">P42262</a>	GRIA2	Limbic Encephalitis; Rasmussen's Syndrome; Progressive Degenerative Diseases; Epilepsy
16	GRID2	Glutamate receptor ionotropic, delta-2/Orphan glutamate receptor delta2 (GRID2)	<a href="#">Q43424</a>	GRID2	Neurological disorders such as cerebellar ataxia and schizophrenia.
17	GRIN1	G protein-regulated inducer of neurite outgrowth 1/Glutamate receptor, ionotropic, N-methyl D-aspartate 1 (GRIN1, transcript variant NR1-1	<a href="#">Q772K8</a>	GPRIN1	Autoimmune encephalitis
18	IgLON5	IgLON family member 5	<a href="#">A6NGN9</a>	IGLON5	Anti-IgLON5 disease, an autoimmune and neurodegenerative disorder typically presents with a combination of sleep disorders, gait instability, bulbar dysfunction (difficulty swallowing or speaking), and cognitive issues.
19	LGI1	Leucine-rich glioma-inactivated protein 1	<a href="#">Q95970</a>	LGI1	LGI1 autoantibody encephalitis
20	mGluR1	Metabotropic glutamate receptor type 1 (mGluR1)/Glutamate receptor metabotropic 1 (GRM1)	<a href="#">Q13255</a>	GRM1	Autoimmune Encephalitis; Cerebellar Ataxia; Paraneoplastic syndrome
21	mGluR2	Metabotropic glutamate receptor type 2 (mGluR2)	<a href="#">Q14416</a>	GRM2	Paraneoplastic cerebellar ataxia; Seizures; altered consciousness, and brainstem atrophy
22	mGluR3	Metabotropic glutamate receptor 3 (mGlu3)	<a href="#">Q14832</a>	GRM3	Rasmussen's encephalitis; Epilepsies; Systemic lupus erythematosus (SLE), Autoimmune encephalitis, Central nervous system ischemia
23	mGluR5	Metabotropic glutamate receptor 5 (mGluR5)	<a href="#">P41594</a>	GRM5	Neurological and psychiatric symptoms (mood changes, anxiety, agitation, psychosis, hallucinations, and personality shifts); Cognitive impairment; Seizures; Movement disorders; Sleep dysfunction; Altered consciousness
24	MNDA	Myeloid cell nuclear differentiation antigen (MNDA)	<a href="#">P41218</a>	MNDA	anti-NMDA receptor encephalitis (brain inflammation with symptoms including psychiatric changes, seizures, movement disorders, and cognitive dysfunction)
25	MOG	Myelin-oligodendrocyte glycoprotein	<a href="#">Q16653</a>	MOG	MOG antibody-associated disease (MOGAD)(Symptoms can include vision loss, weakness, numbness, tingling, and loss of bowel or bladder control)
26	NCAM-1/CD56	Neural Cell Adhesion Molecule 1 (NCAM-1)/CD56	<a href="#">P13591</a>	NCAM1	schizophrenia; membranous lupus nephritis; long COVID.
27	NOVA1	RNA-binding protein Nova-1/Onconeural antigen Nova-1/Paraneoplastic Ri antigen	<a href="#">P51513</a>	NOVA1	Paraneoplastic opsoclonus-myoclonus ataxia (POMA); Paraneoplastic syndromes; Neurological disorders; hepatocellular carcinoma
28	NOVA2	Neuro-oncological ventral antigen 2 (NOVA2) /RNA-binding protein Nova-2	<a href="#">Q9JUNW9</a>	NOVA2	Paraneoplastic opsoclonus myoclonus ataxia (POMA); Dementia; Cancer
29	PKC-gamma (PKCG)	Protein kinase C gamma type	<a href="#">P05129</a>	PRKCG	Paraneoplastic neurological syndromes; Paraneoplastic cerebellar degeneration (PCD)
30	PNMA1	Paraneoplastic antigen MA1 (PNMA1)/37 kDa neuronal protein/Neuron- and testis-specific protein 1	<a href="#">Q8ND90</a>	PNMA1	Paraneoplastic neurological syndromes (PNS)
31	PNMA2	Paraneoplastic antigen Ma2 (PNMA2)/40 kDa neuronal protein/Onconeural antigen Ma2	<a href="#">Q9JL42</a>		Paraneoplastic Neurological Syndromes (PNS); Neurological symptoms (limbic encephalitis, cerebellar degeneration, memory problems, and learning deficits)
32	RCV1/RCVRN	Recoverin (RCVRN)/Cancer-associated retinopathy protein (Protein CAR)	<a href="#">P35243</a>	RCVRN	Cancer-associated retinopathy (CAR), Paraneoplastic syndrome; Vision loss
33	SOX1	Transcription factor SOX-1	<a href="#">Q00570</a>	SOX1	Paraneoplastic neurological syndromes (PNS), Small-cell lung cancer (SCLC); Lambert-Eaton myasthenic syndrome (LEMS), paraneoplastic cerebellar degeneration (PCD), and paraneoplastic
34	ZIC4	Zinc finger protein ZIC 4/Zinc finger protein of the cerebellum 4	<a href="#">Q8N9L1</a>	ZIC4	Paraneoplastic cerebellar degeneration (PCD), a neurological disorder linked to small cell lung carcinoma (SCLC).
35	Ig control 1	control			
36	Ig control 2	control			
37	Ig control 3	control			
38	Ig control 4	control			
39	anti-Ig control 1	control			
40	anti-Ig control 2	control			
41	anti-Ig control 3	control			
42	anti-Ig control 4	control			