

OmicsArray™ General Autoimmune Antigen Array 82-plex Antigen List			Catalog No. PA018	
Antigen Number	Antigen ID	Antigen Name	Gene symbol	UniProt ID
1	CHRNA3	Acetylcholine receptor subunit gamma	CHRNA3	P07510
2	AQP-4 (partial))	Aquaporin-4 (aa208-323)	AQP4	P55087
3	AQP-4 (full length)	Aquaporin-4 (aa208-323)	AQP4	P55087
4	CHRM2	M2 Muscarinic Acetylcholine Receptor	CHRM2	P08172
5	ADCY1	Adenylate cyclase type 1	ADCY1	Q08828
6	ADCY2	Adenylate cyclase type 2	ADCY2	Q08462
7	ADCY3	Adenylate cyclase type 3	ADCY3	Q60266
8	ADCY4	Adenylate cyclase type 4	ADCY4	Q8NFM4
9	ADGRB3_BAI3	Adhesion G protein-coupled receptor B3	ADGRB3	Q60242
10	ARS	Aminoacyl-tRNA Synthetase from Escherichia coli	ARS	P0A8L1
11	BAFF	B cell activating factor belonging to the TNF family (BAFF)	TNFSF13B	Q9Y275
12	BPI	Bactericidal permeability-increasing protein	BPI	P17213
13	CD8B	T-cell surface glycoprotein CD8 beta chain	CD8B	P10966
14	CENP-A	Histone H3-like centromeric protein A	CENPA	P49450
15	Collagen V	Human Collagen Type V	COL5A1	P20908
16	EBNA1	Epstein-Barr nuclear antigen 1	EBNA1	EBNA1
17	EBNA1BP2	EBNA1-binding protein 2	EBNA1BP2	Q99848
18	ELAVL2 (HuB)	ELAV-like protein 2/Hu-antigen B(HuB)	ELAVL2	Q12926
19	ELAVL3 (HuC)	ELAV-like protein 3/Hu-antigen C (HuC)	ELAVL3	Q14576
20	FARS2	Phenylalanine--tRNA ligase, mitochondrial	FARS2	Q95363
21	GBM	Glomerular Basement Membrane (GBM; dissociated)		
22	GDF15	Growth/differentiation factor 15	GDF15	Q99988
23	Gp210	Nuclear pore protein gp210	NUP210	Q8TEM1
24	GRIA2	Glutamate receptor 2	GRIA2	P42262
25	HCOV-229E_S	Human coronavirus (HCoV-229E) Spike Protein	S	P15423
26	HCOV-NL63_S	Human coronavirus (HCoV-NL63) Spike Protein	S	Q6Q1S2
27	HCOV-HKU1_S	Human coronavirus HKU1 (HCoV-HKU1) Spike Protein	S	Q5MQD0
28	Histone	Bovine Histone (natural antigen)		
29	IAV_H1N1	Influenza A (H1N1) Antigen		
30	IBV	Influenza B Antigen		
31	IFNA	Mouse interferon alpha recombinant	ifna1	P01572
32	IFNB1	Mouse interferon beta recombinant	ifnb1	P01575
33	IFN epsilon	Human IFN-epsilon Protein	IFNE	Q86WN2
34	IFN lambda1	human Interferon lambda-1	IFNL1	Q8IU54
35	IL-10	Interleukin-10	IL10	P22301
36	IL-17A	Interleukin-17A	IL17A	Q16552
37	IL-17F	Interleukin-17F	IL17F	Q96PD4

38	IA-2	Insulinoma-associated protein 2	INSM2	Q96T92
39	Intrinsic factor	Cobalamin binding intrinsic factor	CBLIF	P27352
40	Kappa Light Chain	Immunoglobulin kappa	IGKC	P01834
41	KCNJ10_Kir4	ATP-sensitive inward rectifier potassium channel 10	KCNJ10	P78508
42	LKM 1	Cytochrome p450 2D6/liver-kidney microsome 1	CYP2D6	Q2XND0
43	M2	Mitochondrial 2-oxo acid dehydrogenase multienzyme		
44	MBP	Myelin basic protein	MBP	P02686
45	MERS-S	MERS-CoV Spike Protein	S	K9N5Q8
46	mGluR5	Metabotropic glutamate receptor 5	GRM5	P41594
47	Mi-2	Nucleosome remodeling deacetylase (NuRD) complex		
48	NMDAR1_M1 (aa19-393)	Glutamate receptor ionotropic, NMDA 1	GRIN1	Q05586
49	NMDAR1_S1S2 (aa384-544, aa663-	Glutamate receptor ionotropic, NMDA 1	GRIN1	Q05586
50	NMDAR1 (aa21-120)	Glutamate receptor ionotropic, NMDA 1	GRIN1	Q05586
51	MMP14	Matrix metalloproteinase-14	MMP14	Q5RES1
52	Nucleosome	Nucleosome (non-recombinant, bovine)		
53	OLIG2	Oligodendrocyte transcription factor 2	OLIG2	Q13516
54	PCDHA1	Protocadherin alpha-1	PCDHA1	Q9Y5I3
55	PCNA	Proliferating cell nuclear antigen	PCNA	P12004
56	Plastin-2 (LCP1)	Plastin-2	LCP1	P13796
57	PLP1	Myelin proteolipid protein	PLP1	P60201
58	PM/Scl100	Human PM/Scl 100 protein component of nuclear exosome		
59	PM/Scl75	Human PM/Scl 75 protein component of nuclear exosome		
60	POLR3D	DNA-directed RNA polymerase III subunit RPC4	POLR3D	P05423
61	PR3	Leukocyte proteinase 3	PRTN3	P24158
62	RCVRN/RCV1	Recoverin	RCVRN	P35243
63	RCVRN (Abcam)	Recoverin	RCVRN	P35243
64	Ribo P2	Ribosomal Phosphoprotein P2	RPLP2	Q60841
65	RPP25	Ribonuclease P protein subunit p25	RPP25	Q9BUL9
66	RSV	Respiratory Syncytial virus antigen	G	P20895
67	SARS2_NCP	SARS-CoV-2 (2019-nCoV) Nucleocapsid protein	N	P0DTC9
68	SARS2_S1	SARS-CoV-2 (2019-nCoV) Spike S1	S	P0DTC2
69	SARS2_S2	SARS-CoV-2 (2019-nCoV) Spike S2 ECD	S	P0DTC2
70	SARS2-S	ARS-CoV-2 (2019-nCoV) Spike S1+S2 ECD	S	P0DTC2
71	Scl70-Topo I	DNA topoisomerase 1	TOP1	P11387
72	SOX1	ranscription factor SOX-1	SOX1	O00570
73	SP100	Nuclear autoantigen Sp-100	SP100	P23497
74	TSHR	Thyrotropin receptor	TSHR	P16473
75	U1-snRNP BB	Small nuclear ribonucleoprotein-associated proteins B and B'	SNRPB	P14678

76	U1-snRNP C	U1 small nuclear ribonucleoprotein C	SNRPC	P09234
77	CHRM3	Muscarinic acetylcholine receptor M3	CHRM3	P20309
78	HSP27/HSPB1	Heat shock 27 kDa protein (HSP 27)	HSPB1	P04792
79	Col I	Human Collagen Type I		P02666
80	Histone H2B/H2F	Histone H2B type 2-F	H2BC18	Q5QNW6
81	COLQ	Acetylcholinesterase collagenic tail peptide	COLQ	Q9Y215
82	Collagen X	Collagen alpha-1(X) chain	COL10A1	Q03692
83	Buffer	negative control		
84	Ig control 1	Control		
85	Ig control 2	Control		
86	Ig control 3	Control		
87	Ig control 4	Control		
88	anti-Ig control 1	Control		
89	anti-Ig control 2	Control		
90	anti-Ig control 3	Control		
91	anti-Ig control 4	Control		