

OmicsArray™ Coronavirus infection-associated Autoimmune Antigen Array Antigen List - Catalog No. PA012

Antigen#	ID	Antigen name	Uniprot	Category
1	ACE2	Angiotensin-converting enzyme 2	Q9BYF1	autoAg
2	Aldolase C	Fructose-bisphosphate aldolase C	P09972	autoAg
3	ARS	Aminoacyl-tRNA Synthetase	P49588	autoAg
4	B2GP1 (Native)	Beta-2 glycoprotein	P20231	autoAg
5	B2GP1 (Recombinant)	Beta 2 glycoprotein 1	P02749	autoAg
6	BAFF	B cell-activating factor (BAFF)	Q9Y275	cytokines
7	BCOADC-E2	Branched chain 2-oxo-acid dehydrogenase complex component E2	P11182	autoAg
8	BPI	Bactericidal/permeability-increasing protein (BPI)	P17213	autoAg
9	Calprotectin	Calprotectin	P05109 P06702	autoAg
10	CD4	T-cell surface glycoprotein CD4	P01730	cytokines
11	CD8	T-cell surface glycoprotein CD8	P01732 P10966	cytokines
12	CENP-A	Centromere protein A	P49450	autoAg
13	CENP-B	Centromere protein B	P07199	autoAg
14	CMV	NATIVE CYTOMEGALOVIRUS Inactivated		viral Ags
15	Collagen V	Collagen V	P20908	autoAg
16	Collagen VI	Collagen VI	P12109	autoAg

17	Complement C1q	Complement C1q	P02745 P02746	autoAg
18	DFS70	Dense fine speckles 70 kDa protein	O75475	autoAg
19	dsDNA	double stranded DNA		autoAg
20	EBV	Epstein-Barr virus nuclear antigen 1 (EBNA1)	P03211	viral Ags
21	EJ	Glycyl-tRNA Synthetase	P41250	autoAg
22	Fibrinogen	Fibrinogen	P02671 P02675	autoAg
23	GAD65	Glutamate decarboxylase 65 kDa isoform	Q05329	autoAg
24	GBM	Glomerular basement membrane/Collagen alpha-4(IV) chain	P53420	autoAg
25	GDF-15	Growth/differentiation factor 15	Q99988	autoAg
26	GM-CSF	Granulocyte-macrophage colony-stimulating factor	P04141	cytokines
27	GP2	Glycoprotein 2	P55259	autoAg
28	Gp210	Nuclear pore membrane glycoprotein 210	Q8TEM1	autoAg
29	HCOV-229E S	HCOV-229E Spike glycoprotein	P15423	viral Ags
30	HCOV-HKU1 S	HCOV-HKU1 Spike glycoprotein	Q5MQD0	viral Ags
31	HCOV-NL63 S	HCOV-NL63 Spike glycoprotein	P0C6U6	viral Ags
32	HCOV-OC43 S	HCOV-OC43 Spike glycoprotein	P36334	viral Ags
33	Histone	Core Histone	P16403	autoAg
34	HSPG	Heparan sulfate proteoglycan	P34741	autoAg

35	IAV (H1N1)	Inactivated Influenza A virus subtype H1N1		viral Ags
36	IAV (H3N2)	Inactivated Influenza A virus subtype H3N2		viral Ags
37	IBV	Inactivated Influenza B virus		viral Ags
38	IF	Intrinsic factor	P27352	autoAg
39	IFN-a	Mouse IFN- α	Q61718	cytokines
40	IFN-a1	Interferon alpha-1/13	P01562	cytokines
41	IFN-a2	Interferon alpha-2	P01563	cytokines
42	IFN-B	Interferon beta	P01574	cytokines
43	IFN-B1	Interferon beta	P01574	cytokines
44	IFN-epsilon	Interferon epsilon	Q86WN2	cytokines
45	IFN-gamma	Interferon gamma	P01579	cytokines
46	IFN-lambda1	Interferon lambda-1	Q8IU54	cytokines
47	IFN-lambda2/IL28A	Interferon lambda-2	Q8IZJ0	cytokines
48	IFN-lambda3	Interferon lambda-3	Q8IZI9	cytokines
49	IFN-omega	Interferon omega-1	P05000	cytokines
50	IL-10	Interleukin-10	P22301	cytokines
51	IL-12 (p70)	Interleukin-12	P29459 P29460	cytokines
52	IL-15	Interleukin-15	P40933	cytokines

53	IL-17A	Interleukin-17A	Q16552	cytokines
54	IL-17F	Interleukin-17F	Q96PD4	cytokines
55	IL-1a	Interleukin-1 alpha	P01583	cytokines
56	IL-1b	Interleukin-1 beta	P01584	cytokines
57	IL-2	Interleukin-2	P60568	cytokines
58	IL-22	Interleukin-22	Q9GZX6	cytokines
59	IL-4	Interleukin-4	P05112	cytokines
60	IL-6	Interleukin-6	P05231	cytokines
61	IL-8	Interleukin-8	P10145	cytokines
62	IA-2	Insulinoma-associated protein 2	Q96T92	autoAg
63	Jo-1	Histidyl-tRNA Synthetase (Jo-1)	P12081	autoAg
64	KS/AsnRS	Asparaginyl-tRNA Synthetase	O43776	autoAg
65	Ku (p70/p80)	ATP-dependent DNA helicase 2	P12956 P13010	autoAg
66	La/SS-B	Lupus La protein/Sjogren syndrome type B antigen	P05455	autoAg
67	Laminin	Laminin	P07942	autoAg
68	LC1	Liver cytosol type 1	O95954	autoAg
69	LKM-1/CYP2D6	Cytochrome P450 IID6	P10635	autoAg

70	M2	Mitochondrial 2-oxo acid dehydrogenase multienzyme complexes		autoAg
71	MDA5	Melanoma differentiation-associated protein 5	Q9BYX4	autoAg
72	MERS S	Spike glycoprotein	K9N5Q8	viral Ags
73	Mi-2	Mi-2 autoantigen 218 kDa protein	Q14839	autoAg
74	MPO	Myeloperoxidase	P05164	autoAg
75	Neuropilin-1	Neuropilin-1	O14786	autoAg
76	Nucleosome	Nucleosome cores from Bovine		autoAg
77	Nup62	Nuclear pore glycoprotein p62	P37198	autoAg
78	OGDC-E2	2-oxoglutarate dehydrogenase complex component E2	P36957	autoAg
79	PCNA	Proliferating cell nuclear antigen	P12004	autoAg
80	PDC-E2	Pyruvate dehydrogenase complex component E2	P10515	autoAg
81	PL-12	Alanyl-tRNA synthetase	Q5JTZ9	autoAg
82	PL-7	Threonyl-tRNA synthetase	Q9BW92	autoAg
83	PM/Scl100	Polymyositis/scleroderma autoantigen 100 kDa	Q01780	autoAg
84	PM/Scl75	Polymyositis/scleroderma autoantigen 75 kDa	Q06265	autoAg
85	PR3	Leukocyte proteinase 3	P24158	autoAg
86	RF	Rheumatoid Factor (RF)		autoAg

87	PTH	Parathyroid Hormone	P01270	autoAg
88	RNP/Sm (Native)	U-small nuclear riboprotein complexes	P62314	autoAg
89	Ro-52	52 kDa ribonucleoprotein autoantigen Ro/SS-A	P19474	autoAg
90	Ro-60 (native)	RNA-binding protein RO60	P10155	autoAg
91	RPLP0	Ribosomal Phosphoprotein P0	P05388	autoAg
92	RPLP1	Ribosomal Phosphoprotein P1	P05386	autoAg
93	RPLP2	Ribosomal Phosphoprotein P2	P05387	autoAg
94	RSV	Inactivated Respiratory Syncytial Virus Antigen		viral Ags
95	SARS2 S2	SARS-CoV-2 Spike protein S2	P0DTC2	viral Ags
96	SARS1 S1	SARS-CoV-1 Spike S1 Protein	P59594	viral Ags
97	SARS2 NCP	SARS-CoV-2 Nucleocapsid protein	P0DTC2	viral Ags
98	SARS2 S	SARS-CoV-2 Spike protein	P0DTC2	viral Ags
99	SARS2 S (B.1.1.529)	SARS-CoV-2 Spike protein (Omicron)	P0DTC2	viral Ags
100	SARS2 S1	SARS-CoV-2 Spike S1 protein	P0DTC2	viral Ags
101	SARS2 S1RBD	SARS-CoV-2 Spike S1 receptor binding domain	P0DTC2	viral Ags
102	Scl-70	Scl-70/Topo Isomerase I	P11387	autoAg
103	Sm (Native)	Small nuclear ribonucleoprotein		autoAg
104	SmD (Recombinant)	SmD comprises equal masses of SmD1, SmD2, and SmD3	P62314	autoAg

105	SP100	Nuclear autoantigen Sp-100	P23497	autoAg
106	SRP54	Signal recognition particle 54 kDa protein	P61011	autoAg
107	TGFB1	Transforming growth factor beta-1 proprotein	P01137	cytokines
108	Thyroglobulin	Thyroglobulin	P01266	autoAg
109	TPO	Thyroid Peroxidase	P07202	autoAg
110	TIF1-r	Transcription intermediary factor 1 gamma	Q9UPN9	autoAg
111	TTG	Tissue transglutaminase	Q9BXQ0	autoAg
112	TNF-B (LT-a)	TNF-beta (Lymphotoxin-alpha)	P01374	cytokines
113	Troponin I	Troponin I	P19429	autoAg
114	TWEAK (CD255)	TNF weakly inducer of apoptosis	O43508	cytokines
115	U1-snRNP-68/70	U1 small nuclear ribonucleoprotein 70 kDa	P08621	autoAg
116	U1-snRNP-A	U1 small nuclear ribonucleoprotein A	P09012	autoAg
117	U1-snRNP-BB'	U-snRNP B/B'	P14678	autoAg
118	U1-snRNP-C	U1 small nuclear ribonucleoprotein C	P09234	autoAg
119	Vimentin	Vimentin	P08670	autoAg
120	Vitronectin	Vitronectin	P04004	autoAg
121	Ig control 1			control
122	Ig control 2			control

123	Ig control 3			control
124	Ig control 4			control
125	anti-Ig control 1			control
126	anti-Ig control 2			control
127	anti-Ig control 3			control
128	anti-Ig control 4			control