

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

Antigen number	Antigen name	Category
1	ACE2	Autoantigen
2	Aldolase C	Autoantigen
3	Aminoacyl-tRNA Synthetase (ARS)	Autoantigen
4	Asparaginyl-tRNA Synthetase (KS)	Autoantigen
5	B2-GP1 (Native)	Autoantigen
6	B2GP1 (Recombinant)	Autoantigen
7	BAFF	Cytokine
8	BCOADC-E2	Autoantigen
9	BPI	Autoantigen
10	Calprotectin	Autoantigen
11	CD4	Cytokine
12	CD8	Cytokine
13	CENP-A	Autoantigen
14	CENP-B	Autoantigen
15	CMV	Viral antigen
16	Collagen V	Autoantigen
17	Collagen VI	Autoantigen
18	Complement C1q	Autoantigen
19	DFS70	Autoantigen
20	dsDNA	Autoantigen
21	EBV	Viral antigen
22	Fibrinogen	Autoantigen
23	GAD65	Autoantigen
24	GBM	Autoantigen
25	GDF-15	Autoantigen
26	Glycyl-tRNA Synthetase (EJ)	Autoantigen
27	GM-CSF	Cytokine
28	GP2	Autoantigen
29	Gp210	Autoantigen

30	HCOV-229E S	Viral antigen
31	HCOV-HKU1 S	Viral antigen
32	HCOV-NL63 S	Viral antigen
33	HCOV-OC43 S	Viral antigen
34	Heparan sulfate proteoglycan (HSPG)	Autoantigen
35	Histone	Autoantigen
36	IAV (H1N1)	Viral antigen
37	IAV (H3N2)	Viral antigen
38	IBV (Influenza B)	Viral antigen
39	IFN-a	Cytokine
40	IFN-a1	Cytokine
41	IFN-a2	Cytokine
42	IFN-B	Cytokine
43	IFN-B1	Cytokine
44	IFN-epsilon	Cytokine
45	IFN-gamma	Cytokine
46	IFN-lambda1	Cytokine
47	IFN-lambda2/IL28A	Cytokine
48	IFN-lambda3	Cytokine
49	IFN-omega	Cytokine
50	IL-10	Cytokine
51	IL-12 (p70)	Cytokine
52	IL-15	Cytokine
53	IL-17A	Cytokine
54	IL-17F	Cytokine
55	IL-1a	Cytokine
56	IL-1b	Cytokine
57	IL-2	Cytokine
58	IL-22	Cytokine
59	IL-4	Cytokine
60	IL-6	Cytokine
61	IL-8	Cytokine
62	Insulinoma-associated (IA-2)	Autoantigen
63	Intrinsic factor (IF)	Autoantigen

64	Jo-1	Autoantigen
65	Ku (p70/p80)	Autoantigen
66	La/SS-B	Autoantigen
67	Laminin	Autoantigen
68	Liver cytosol type 1 (LC1)	Autoantigen
69	LKM-1	Autoantigen
70	M2	Autoantigen
71	MDA5	Autoantigen
72	MERS S	Viral antigen
73	Mi-2	Autoantigen
74	MPO	Autoantigen
75	Neuropilin-1	Autoantigen
76	Nucleosome	Autoantigen
77	Nup62	Autoantigen
78	OGDC-E2	Autoantigen
79	Parathyroid Hormone (rHuPTH)	Autoantigen
80	PCNA	Autoantigen
81	PDC-E2	Autoantigen
82	PL-12	Autoantigen
83	PL-7	Autoantigen
84	PM/Scl100	Autoantigen
85	PM/Scl75	Autoantigen
86	PR3	Autoantigen
87	Rheumatoid Factor (RF)	Autoantigen
88	ribosomal protein P0 (RPLP0)	Autoantigen
89	ribosomal protein P1 (RPLP1)	Autoantigen
90	ribosomal protein P2 (RPLP2)	Autoantigen
91	RNP/Sm (Native)	Autoantigen
92	Ro-52	Autoantigen
93	Ro-60 (native)	Autoantigen
94	RSV	Viral antigen
95	SARS S2	Viral antigen
96	SARS1 S1	Viral antigen
97	SARS2 NCP	Viral antigen

98	SARS2 S	Viral antigen
99	SARS2 S (B.1.1.529)	Viral antigen
100	SARS2 S1	Viral antigen
101	SARS2 S1RBD	Viral antigen
102	Scl-70/Topo Isomerase I	Autoantigen
103	Sm (Native)	Autoantigen
104	SmD (Recombinant)	Autoantigen
105	SP100	Autoantigen
106	SRP54	Autoantigen
107	TGFB1	Cytokine
108	Thyroglobulin	Autoantigen
109	Thyroid Peroxidase (TPO)	Autoantigen
110	TIF1-r	Autoantigen
111	Tissue transglutaminase (TTG)	Autoantigen
112	TNF-B (LT-a)	Cytokine
113	Troponin I	Autoantigen
114	TWEAK (CD255)	Cytokine
115	U1-snRNP-68	Autoantigen
116	U1-snRNP-A	Autoantigen
117	U1-snRNP-BB'	Autoantigen
118	U1-snRNP-C	Autoantigen
119	Vimentin	Autoantigen
120	Vitronectin	Autoantigen
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