

OmicsArray™ Interferon, Infection and autoimmune panel, 64-plex antigen list - PA011

Antigen Number	Antigen Name
1	Aldolase C
2	Aminoacyl-tRNA Synthetase (ARS)
3	Beta-2 Glycoprotein I (B2GP1)
4	CD4
5	CD8
6	Collagen V
7	Collagen VI
8	Complement C1q
9	dsDNA
10	GAD65
11	Histone
12	IFN-a 4B
13	IFN-a A
14	IFN-a B2
15	IFN-a C
16	IFN-a D
17	IFN-a F
18	IFN-a G
19	IFN-a H2
20	IFN-a I
21	IFN-a J1
22	IFN-a K
23	IFN-a Wa
24	IFN-a/IFN-a1
25	IFN-a2
26	IFN-b/IFN-b1
27	IFN-r
28	IFN-e
29	IFN-L1
30	IFN-L2/IL28A
31	IFN-L3
32	IFN-Omega
33	CXCR4/PF4
34	Ku (p70/p80)
35	La/SS-B
36	MPO
37	Neuropilin-1
38	Nucleosome
39	Nup62
40	Parathyroid Hormone (rHuPTH)

41	PCNA
42	RNP/Sm
43	Ro/SS-A (52 + 60 kDa)
44	CMV
45	EBV
46	Influenza A H1N1
47	Influenza A H3N2
48	Influenza B
49	RSV
50	HCOV-229E S
51	HCOV-HKU1 S
52	HCOV-NL63 S
53	HCOV-OC43 S
54	MERS-CoV S
55	SARS-CoV-1 S1
56	SARS-CoV-2 NCP
57	SARS-CoV-2 S
58	SARS-CoV-2 S1
59	SARS-CoV-2 S1RBD
60	SARS-CoV-2 S2
61	Sm
62	SmD
63	Thyroglobulin
64	Thyroid Peroxidase (TPO)
65	Ig control 1
66	Ig control 2
67	Ig control 3
68	Ig control 4
69	anti-Ig control 1
70	anti-Ig control 2
71	anti-Ig control 3
72	anti-Ig control 4