

OmicsArray™ Infection and Vaccination-associated Autoimmune Antigen Array Antigen List - Catalog No. PA010

Antigen number	Antigen name	Category
1	Arrestin, Beta 1/(Human ARRB1)	autoantigen
2	AChR3 (muscle)	autoantigen
3	Adenovirus (ADV type 5) natural antigen	Virus
4	Alpha enolase	autoantigen
5	Alpha Fodrin	autoantigen
6	Aquaporin 4 (AQP4)	autoantigen
7	B cell activating factor (BAFF)	autoantigen
8	Beef_Bos taurus	Allergen
9	Bermuda_Cynodon dactylon	Allergen
10	CANDIDA ALBICANS	Allergen
11	Carboxypeptidase H	autoantigen
12	Casein (milk protein)	autoantigen
13	Cat dander	Allergen
14	Cedar,Red_Juniper rus virginiana	Allergen
15	CENP-A	autoantigen
16	Chitinase 3-Like 1 (Cartilage Glycoprotein-39)	autoantigen
17	Cholinergic Receptor Muscarinic 3 (CHRM3)	autoantigen
18	CMV Ext-2, enriched for cell surface glycoprotein antigens	Virus
19	CMV III Antigen, enriched for pp65	Virus
20	CMV-G, whole cell extract	Virus
21	CMV-M Concentrate, nuclear extract and ER antigens	Virus
22	Cockroach	Allergen
23	Complement C1q	autoantigen
24	Core Histone_Citrullinated	autoantigen
25	CoV SARS SPIKE /RBD(SB)	Virus
26	CoV -HKU1 (N5) SPIKE S 1/2 (SB)	Virus
27	CoV HKU1 SPIKE S1 (CD)	Virus
28	CoV MERS SPIKE/ECD (SB)	Virus
29	CoV -NL63 SPIKE S 1/2 (SB)	Virus

30	CoV-229 E SPIKE S 1/2 (SB)	Virus
31	COVID 19 NCP (CD)	Virus
32	COVID 19 SPIKE S1 (CD)	Virus
33	COVID 19 SPIKE S1(SB)	Virus
34	COVID 19 SPIKE S1/2(SB)	Virus
35	COVID 19 SPIKE S2/ECD(SB)	Virus
36	CoV-oc43 SPIKE S 1/2 (SB)	Virus
37	Coxsackievirus (B1) Native Antigen	Virus
38	Dengue virus (DENV)	Virus
39	Diphtheria Toxin	Bacterial
40	Dog Dander, Mixed-Breed_Canis familiaris	Allergen
41	Dog Epithelia, Mixed-Breed_Canis familiaris	Allergen
42	dsDNA/ssDNA	autoantigen
43	EBV EBNA1 ANTIGEN	Virus
44	Echovirus type 9	Virus
45	Egg-white Lysozyme (HEL)	autoantigen
46	Endothelial cells antigen	autoantigen
47	Enterovirus 71 VP1 (EV71)	Virus
48	Esterase Pseudomonas fluorescens	Pseudomonas
49	Fibrinogen IV	autoantigen
50	GBM (kidney)	autoantigen
51	Glucose-6-Phosphatase, Catalytic, 2 /(G6PC2)	autoantigen
52	Glutamate decarboxylase 1 (GAD1)	autoantigen
53	Glutamate decarboxylase 2 (GAD2/GAD65)	autoantigen
54	GPIb Alpha/GPIIbIIIa/GPIV	autoantigen
55	Haemophilus influenzae OmpP2	Bacterial
56	Hepatitis A Virus (HAV) Antigen	Virus
57	Hepatitis A Virus (HAV) Antigen, concentrate	Virus
58	Hepatitis B surface antigen	Virus
59	Herpes simplex virus-1 (HSV-1)	Virus
60	Herpes simplex virus-2 (HSV-2)	Virus
61	Histone H1/H2A/H2B/H3/H4	autoantigen
62	Honey Bee_Apis mellifera	Allergen
63	House Dust	Allergen

64	Influenza A antigen	Virus
65	Influenza B antigen	Virus
66	Influenza H1N1	Virus
67	Jo-1	autoantigen
68	KU -P70/P80	autoantigen
69	La/SSB	autoantigen
70	Liver Cytosol Type 1 (LC1)	autoantigen
71	Liver Kidney Microsome(LKM1)	autoantigen
72	M. Tuberculosis CFP10	Virus
73	M. Tuberculosis ESAT-6 Antigen	Virus
74	MDA5	autoantigen
75	MERS-CoV NCP	Virus
76	Mite, House Dust_ Blomia tropicalis	Allergen
77	Mitochondrial Ribosomal Protein S31	autoantigen
78	Mumps virus antigen Inactivated	Virus
79	Mycoplasma Pneumoniae Antigen	Virus
80	Myelin Basic Protein (MBP)	autoantigen
81	Myelin oligodendrocytic glucoprotein (MOG)	autoantigen
82	Myelin-associated Glycoprotein (MAG)	autoantigen
83	Myosin	autoantigen
84	Native cytomegalovirus	Virus
85	Native Influenza A Virus H3N2	Virus
86	Native Influenza B Virus Antigen	Virus
87	Norovirus vp1	Virus
88	Nucleosome	autoantigen
89	Ovalbumin (OVA)	autoantigen
90	Parvovirus VP1 (aa 1-781)	Virus
91	Poliorirus tape 1,2,3	Virus
92	Proteolipid protein(PLP)	autoantigen
93	Respiratory Syncytial Virus (RSV) long strain	Virus
94	Ro/SSA (52kd and 60kd)	autoantigen
95	ROTAVIRUS SA-11 ANTIGEN	Virus
96	RSV(DAG211)	Virus
97	RUBELLA GRADE IV ANTIGEN PBS	Virus

98	Rubella RSVP Antigen	Virus
99	RUBELLA VIRUS GRADE IV ANTIGEN	Virus
100	RUBEOLA ANTIGEN	Virus
101	SARS1 S1	Virus
102	SARS2_S1_RBD (F342L)	Virus
103	SARS-CoV-1 NCP	Virus
104	SARS-CoV-2 NCP	Virus
105	SARS-CoV-2 S1 RBD	Virus
106	Scl-70	autoantigen
107	Scl-75/Scl-100	autoantigen
108	Shrimp_Penacidae	Allergen
109	Sm/SmD	autoantigen
110	Streptococcus pneumoniae	Bacterial
111	TETANUS TOXIN	Bacterial
112	Thyroglobulin (TG)	autoantigen
113	Tissue Transglutaminase (TTG)	autoantigen
114	Toxoplasma Antigen (Toxo) grade II antigen	Virus
115	U1-sn-RNP (A, BB', C, 68)	autoantigen
116	Varicella Zoster Virus (VZV) Antigen	Virus
117	Varicella Zoster Virus (VZV) II Antigen	Virus
118	Vitronectin	autoantigen
119	West Nile Virus (WNV) Envelope protein	Virus
120	Wheat, Whole_Triticum aestivum	Allergen
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