

OmicsArray™ Common Pathogen Antigen Array, 64-plex, Antigen List - Catalog No. PA021

Antigen#	Antigene ID	Antigen Name
1	IAV H5N1	Native IAV H5N1 Antigen
2	IAV H3N2	Native IAV H3N2 Antigen
3	IAV H1N1	Inactivated IAV H1N1 Antigen
4	IBV	Native Influenza B Virus Antigen
5	RSV	Native RSV Antigen
6	RSV	RSV Grade II antigen
7	SARS2 S	SARS-CoV-2 (2019-nCoV) Spike S1+S2 ECD
8	SARS2 RBD	SARS-CoV-2 (2019-nCoV) Spike RBD
9	HAV	Hepatitis A virus (HAV) Antigen
10	HBV sAg	Hepatitis B (HBV) surface antigen
11	CMV	Native cytomegalovirus Antigen
12	Coxsackie virus	Coxsackievirus (B1) Native Antigen
13	Echovirus	Echovirus type 9 Native Antigen
14	Tuberculosis	M. Tuberculosis ESAT-6 antigen (a.a.1-99) and CFP10 + MTB81
15	S. Pneumoniae	Streptococcus pneumoniae native antigen
16	Parainfluenza type-1	Human Para influenza virus Type-1 (HPIV-1)
17	Haemophilus influenzae	Haemophilus influenzae OmpP2 (HIV OmpP2)
18	HSV-1	HERPES (HSV)-1 Antigen
19	HSV-2	HERPES (HSV) 2 Antigen
20	IBV Hemagglutinin	Influenza B virus (IBV) Hemagglutinin
21	Dengue virus (DENV)	Recombinant DENV Polyvalent Antigen from E. coli
22	Enterovirus (EV71)	Enterovirus 71 VP1, recombinant protein from E. coli
23	Parvovirus B19 VP1	Recombinant Parvovirus B19 VP1 Protein (aa 1-781)
24	West Nile Virus	West Nile Virus (WNV) Envelope protein recombinant protein from E. coli
25	TOXO Gon	Toxoplasma gondii Antigen
26	VZV	Varicella Zoster Virus antigen
27	Rotavirus	Rotavirus SA-11 antigen
28	Parvovirus B19 VP2	Parvovirus B19 VLP VP2 recombinant protein
29	Rubeola	Rubeola Antigen Partially Purified
30	HBV cAg	HBCAG (HBV CORE δ ANTIGEN) recombinant protein in E Coli
31	M. Pneumoniae	Mycoplasma Pneumoniae native antigen
32	HCoV NL63 Spike protein	Human coronavirus (HCoV-NL63) Spike Protein (S1+S2 ECD)
33	HCoV 229E Spike protein	Human coronavirus (HCoV-229E) Spike Protein (S1+S2 ECD)
34	HCoV HKU1 Spike protein	Human coronavirus HKU1 (HCoV-HKU1) Spike Protein (S1+S2 ECD)
35	HCoV OC43 Spike protein	Human coronavirus (HCoV-OC43) Spike Protein (S1+S2 ECD)
36	MERS Spike protein	MERS-CoV Spike Protein (S1+S2 ECD, aa 1-1297)
37	SARS1 Spike/S1 Protein	SARS-CoV Spike/S1 Protein (S1 Subunit)
38	H.Pylori Antigen	Helicobacter pylori Antigen
39	Tetanus toxin	Tetanus Txon from clostridium tetani
40	HIV	HIV Chimeric Recombinant
41	E.coli	E. coli O157

42	Poliovirus type 1	Salk Inactivated poliovirus type 1 (sIPV-1), inactivated native antigen
43	Poliovirus type 2	Salk Inactivated poliovirus type 2 (sIPV-2), inactivated native antigen
44	Poliovirus type 3	Salk Inactivated poliovirus type 3 (sIPV-3), inactivated native antigen
45	Norovirus GII.4 VP1 Antigen	Norovirus GII.4 VP1 recombinant antigen produced in mammalian HEK293 cells
46	Diphtheria Toxin	Diphtheria Toxin from <i>Corynebacterium diphtheriae</i>
47	RUBELLA RSV	Native Rubella Virus RSV Antigen
48	Epstein-Barr virus (EBV)	Epstein-Barr Virus (HHV-4) EBNA1 Mosaic Recombinant
49	HPV-6 L1	Human HPV-6 L1 protein (CDK37669) expressed in Yeast
50	HPV-6 L2	Human HPV-6 L2 protein(CDK37668) expressed in Yeast
51	HPV-6 E2	Human HPV-6 E2 Protein expressed in <i>E. coli</i>
52	HPV-6a E4 E5	Human HPV-6a E4 and E5 Protein expressed in <i>E. coli</i>
53	HPV-6 E6	Human HPV-6 E6 Protein expressed in <i>E. coli</i>
54	HPV-6 E7	Human HPV-6 E7 Protein expressed in <i>E. coli</i>
55	HPV-11 E1	Human HPV-11 E1 protein
56	HPV-11 E2	Human HPV-11 E2 (aa 1-367) Protein
57	HPV-11 E4	Human HPV-11 E4 & E5B Protein
58	HPV-11 E6	Human HPV-11 E6 (aa 1-150) Protein
59	HPV-11 E7	Human HPV-11 E7 (aa 1-98) Protein
60	HPV-11 L1	Human HPV-11 L1 Protein (aa 1- 501) from <i>E.coli</i>
61	HPV-16	Human HPV-16 full length protein expressed in <i>E. coli</i>
62	HPV16 E6E7	Human HPV16 E6 and E7 full length protein
63	HPV-18	Human HPV-18 full length large capsid protein
64	HPV18 E7	Human HPV 18 E7 protein
65	Ig control 1	Ig control 1
66	Ig control 2	Ig control 2
67	Ig control 3	Ig control 3
68	Ig control 4	Ig control 4
69	alg control 1	alg control 1
70	alg control 2	alg control 2
71	alg control 3	alg control 3
72	alg control 4	alg control 4