

OmicsArray™ SARS-CoV-2 and Variants Antigen Array Antigen List - Catalog No. PA013

Antigen number	Antigen name	Variant strain	Origin
1	MERS-COV NCP		
2	MERS-CoV S1		
3	SARS-CoV-1 NCP		
4	SARS-CoV-1 S1		
5	SARS-CoV-1 S1 RBD		
6	SARS-CoV-1 Plpro		
7	SARS-CoV-2 NCP		
8	SARS-CoV-2-NCP (EC)		
9	SARS-CoV-2 NCP (094)		
10	SARS-CoV-2 S		
11	SARS-CoV-2 S2		
12	SARS-CoV-2 S1		
13	SARS-CoV-2 S1 (091)		
14	SARS-CoV-2 S1RBD		
15	SARS-CoV-2 S1RBD (No tag)		
16	SARS-CoV-2 Plpro		
17	SARS-CoV-2 ME		
18	HCoV-NL63 S		
19	HCoV-229E S		
20	HCoV-HKU1 S		
21	HCoV-HKU1 S1		
22	HCoV-OC43 S		
23	SARS-CoV-2 E		
24	SARS-CoV-2 M		
25	SARS-CoV-2 3CL		
26	SARS-CoV-2 ORF3A		
27	SARS-CoV-2 ORF7A		
28	SARS-CoV-2 ORF8		
29	SARS-CoV-2 ORF9		

30	SARS-CoV-2 NSP5		
31	SARS-CoV-2 S1RBD (F342L)		
32	SARS-CoV-2 S1RBD (N501Y)	B.1.1.7	UK
33	SARS-CoV-2 S1 (HV69-70 Del, Y144 Del, N501Y, A570D, D614G, P681H)		
34	SARS-CoV-2 S1 (HV69-70 Del, N501Y, D614G)		
35	SARS-CoV-2 S1 (K417N, E484K, N501Y, D614G)	B.1.351 (Beta)	South Africa
36	SARS-CoV-2 NCP (D3L, R203K, G204R, S235F)		
37	SARS-CoV-2 S1 (HV69-70del, Y453F, D614G)		
38	SARS-CoV-2 NCP (R203K, G204R)		
39	SARS-CoV-2 S (HV69-70del, Y144del, N501Y, A570D, D614G, P681H, T716I, S982A, D1118H)		
40	SARS-CoV-2 S1 (D614G)	B.1.525	Europe
41	SARS-CoV-2 S1 (P618H)		
42	SARS-CoV-2 S1 (E484K)	B.1.525	Europe
43	SARS-CoV-2 S1 (K417N)	B.1.351 (Beta)	South Africa
44	SARS-CoV-2 S1 (L452R)	B.1.526.1, B.1.427 (Epsilon)	Europe
45	SARS-CoV-2 S (W152C, L452R, D614G)	B.1.429	U.S.
46	SARS-CoV-2 S1RBD (L452R, E484Q)	B.1.617	India
47	SARS-CoV-2 S1RBD (K417N, E484K, N501Y)	B.1.351	South Africa
48	SARS-CoV-2 S1RBD (K417T, E484K, N501Y)	P.1	Brazil
49	SARS-CoV-2 S1 (E154K, L452R, E484Q, D614G, P681R)	B.1.617	India
50	ACE2		
51	Influenza A IAV H1N1		
52	Influenza A IAV H3N2		
53	Influenza B		
54	CMV Native		
55	RSV Inactivate		
56	Adenovirus (ADV type 5) natural antigen		
57	Coxsackievirus (B1) Native Antigen		
58	Echovirus type 9		
59	Ig control		
60	Ig control		
61	Ig control		

62	Ig control		
63	anti-Ig control		
64	anti-Ig control		
65	anti-Ig control		
66	anti-Ig control		