

## OmicsArray™ Cancer and Neoplasm-associated Antigen Array Antigen List - Catalog No. PA003

Antigen number	Antigen name
1	ALDOC
2	AChR3
3	AFP
4	Aldolase (muscle)
5	Alpha-2-macroglobulin-like 1
6	alpha-fodrin
7	Angiotensin II type 1 Receptor
8	ANXA1
9	Apolipoprotein B/E
10	Azurocidin
11	BAFF
12	beta-glucuronidase
13	BRCA2/BRCA1
14	CA125
15	CA15-3
16	CA19-9
17	CA242
18	CA27-29
19	CA50
20	CA72-4
21	Calmodulin (brain)
22	Calreticulin(rabbit calreticulin)
23	Carbonic Anhydrase 6 (CA6)
24	Catalase
25	Cathepsin G
26	CCL11
27	CCL2
28	CCL3
29	CCL5

30	CD74 Protein
31	CDK2
32	CEA
33	Cholinergic Receptor Muscarinic 3 (CHRM3)
34	CMV native protein
35	C-MYC
36	CTLA4
37	CXCL10
38	Cytokeratin 19 Ag
39	DDX53
40	dsDNA
41	EBV EBNA1
42	EGF Receptor
43	Elastase
44	Endothelial Cell Extract
45	Enolase
46	ErbB2
47	ErbB3 + ErbB4
48	ERP29
49	Ferritin
50	FUCA1
51	Galactocerebroside
52	GBM
53	GBU4-5/TDRD-12
54	Glutamate decarboxylase-65/GAD2
55	GNAL Protein
56	Hepatitis B antigen
57	HPV-11 E2 (aa 1-367)
58	HPV-11 E4 (aa 1-108)
59	HPV-11 E5B (aa 1-74)
60	HPV-11 E6 (aa 1-150)
61	HPV-11 E7 (aa 1-98)
62	HPV-16 E6
63	HPV-16 E7

64	HPV-18 E7
65	HPV-6 E2
66	HPV-6 E6
67	HPV-6 E7
68	HPV-6 L1
69	HPV-6 L2
70	HPV-6a E4 (aa 1-99)
71	HuD/ELAVL4
72	IFN-?
73	IFN-b/IFN-a2
74	IL-12 (p70)
75	IL-17A
76	IL-1a/IL-1b
77	IL-2
78	IL-6
79	IL-8
80	ITM2B Protein
81	Lactoferrin
82	LC1
83	Lysosomal Associated Protein 2 (LAMP-2)
84	MAG
85	MAGEA3
86	MAGEA4
87	MBP
88	MMP-2
89	NSE
90	Nucleosome
91	NY-ESO-1
92	Ox40L
93	P53
94	Parotid Secretory Protein (PSP)
95	PD-1
96	PD-L1
97	Peptidyl Arginine Deiminases 1&2&3&4

98	Phosphatidyl-l-serine
99	PKM2
100	Prostate Specific Antigen (PSA)
101	Prostate Specific Membrane Antigen
102	Prostatic Acid Phosphatase
103	PSMA
104	rhHSPG2
105	Ribo_PO.P1.P2 mix
106	Ro-SSA (52+60)
107	SCCA
108	Sm/RNP
109	SOX2
110	Sphingomyelin
111	TGF- $\beta$ 1
112	Thrombopoietin (TPO)
113	Thyroglobulin
114	Tissue Transglutaminase (TTG)
115	TNF-a/TNF-b
116	Troponin I
117	Troponin I-T-C ternary complex mixture
118	TWEAK (CD255)
119	uPA
120	VEGF-165
121	Ig control
122	Ig control
123	Ig control
124	Ig control
125	anti-Ig control
126	anti-Ig control
127	anti-Ig control
128	anti-Ig control