

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

Antigen	Antigen ID	Name	UniProtID
1	CHRNA3	Acetylcholine receptor subunit gamma (AChR)	P07510
2	AFP	Alpha-fetoprotein	P02771
3	A2ML1	Alpha-2-macroglobulin-like protein 1	A8K2U0
4	SPTAN1	alpha-fodrin/Spectrin alpha chain, non-erythrocytic 1	Q13813
5	AGTR1	Type-1 angiotensin II receptor	P30556
6	ANXA2	Annexin A2	P07355
7	APEX1	DNA-(apurinic or apyrimidinic site) endonuclease	P27695
8	APOB/APOE	Apolipoprotein B/Apolipoprotein E	P04114, P02649
9	AURKA	Aurora Kinase A	O14965
10	AZU1	Azurocidin	P20160
11	BAFF/TNFSF13B	B-cell-activating factor/Tumor necrosis factor ligand superfamily member 13B	Q9Y275
12	BAG3	BAG family molecular chaperone regulator 3	O95817
13	BGLR	Beta-glucuronidase	P08236
14	BRCA1	Breast cancer type 1 susceptibility protein	P38398
15	CA125/MUC16	Ovarian carcinoma antigen CA125/Mucin-16	Q8WXI7
16	CA15-3/MAGEA3/Mucin-1	Breast carcinoma-associated antigen DF3, Melanoma-associated antigen 3	P43357
17	CA 19-9	CA19-9 (carbohydrate antigen 19-9, also called cancer antigen 19-9 or sialylated Lewis (a) antigen). Native Human CA 19-9 was purified from Human Metastatic Liver Carcinoma	
18	CA 242	CA 242 is a tumor marker for sialylated Lewis carbohydrates associated with adenocarcinomas and e-selectin-mediated metastatic risk. Native Human CA-242 was purified from Human Tissue Culture	
19	CA 27-29	MUC-1, CA 27-29, Breast Tumor protein	P15941
20	CA50	CA-50, Cancer Antigen 50, carbohydrate antigen 50	
21	CA72-4	Cancer Antigen 72-4, tumor-associated glycoprotein-72 (TAG72-4)	
22	CALM	Calmodulin-1	P0DP23
23	CA6	Carbonic Anhydrase 6	P23280
24	CAT	Catalase	P04040

25	CTSG	Cathepsin G	P08311
26	CCL11	Eotaxin	P51671
27	CCL2	C-C motif chemokine 2	P13500
28	CCL3	C-C motif chemokine 3	Q68A92
29	CCL5	C-C motif chemokine 5	P13501
30	CD74	HLA class II histocompatibility antigen gamma chain	P04233
31	CDK2	Cyclin-dependent kinase 2	P24941
32	CEA	Carcinoembryonic antigen	Q13985
33	CHRM3	Muscarinic acetylcholine receptor M3	P20309
34	CMV	Cytomegalovirus (CMV)	
35	C-MYC	Myc proto-oncogene protein	P01106
36	CTLA4	Cytotoxic T-lymphocyte protein 4	P16410
37	CXCL10	C-X-C motif chemokine 10	P02778
38	KRT19	Cytokeratin 19, Keratin, type I cytoskeletal 19	P08727
39	DDX53	Probable ATP-dependent RNA helicase DDX53	Q86TM3
40	dsDNA	double strand DNA	
41	EBNA1	Epstein-Barr nuclear antigen 1	P03211
42	Elastase	Neutrophil elastase	P08246
43	ENO1	Alpha-enolase	P06733
44	ErbB2	Receptor tyrosine-protein kinase erbB-2	P04626
45	ErbB3	Receptor tyrosine-protein kinase erbB-3	P21860
46	ErbB4	Receptor tyrosine-protein kinase erbB-4	Q15303
47	ERP29	Endoplasmic reticulum resident protein 29	P30040
48	Ferritin	Ferritin	Q8TD27
49	FUCA1	Tissue alpha-L-fucosidase	P04066
50	GAD65/GAD2	Glutamate decarboxylase 2	Q04329
51	GALC	Galactocerebrosidase	P54803
52	GBM	Glomerular Basement Membrane, Collagen alpha-4(IV) chain	P53420
53	GBU4-5/TDRD-12	cancer-associated gene protein (CAGE), Putative ATP-dependent RNA helicase TDRD12	Q587J7
54	GNAL	Guanine nucleotide-binding protein G(olf) subunit alpha	P38405
55	H. Pylori	Helicobacter pylori	85962
56	HBsAg	Hepatitis B virus surface antigen	10407
57	HMGB1	High mobility group protein B1	P09429

58	HPV11	Human papillomavirus 11 E1, E2, E4, E5B, E6, E7, L1	10580
59	HPV16	Human papillomavirus type 16 (E6, E7)	333760
60	HPV18	Human papillomavirus type 18	333761
61	HPV-6 (mix 6)	Human papillomavirus type 6 (E2, E4, E5, E6, E7, L1, L2)	37122
62	IFNA2	Interferon alpha-2	P01563
63	IFNB1	Interferon beta	P01574
64	IFNG	Interferon gamma	P01579
65	IL-12 (p70)	Interleukin-12 subunit beta/alpha (p40/p35)	P29460, P29459
66	IL17A	Interleukin-17A	Q16552
67	IL-1A	Interleukin-1 alpha	P01583
68	IL-1B	Interleukin-1 beta	P01584
69	IL-2	Interleukin-2	P60568
70	IL-6	Interleukin-6	P05231
71	IL-8	Interleukin-8	P10145
72	ICA 512/PTPRN	Insulinoma-associated protein IA-2 (ICA 512)/Receptor-type tyrosine-protein phosphatase-like N	Q16849
73	ITM2B	Integral membrane protein 2B	Q9Y287
74	Jo-1/HARS1	Histidyl-tRNA Synthetase (Jo-1)/Histidine--tRNA ligase	P12081
75	Lactoferrin	Lactotransferrin	P02788
76	LC1/FTCD	liver cytosol type 1 antigen (LC1)/Formiminotransferase Cyclodeaminase (FTCD)	O95954
77	LKM 1/CYP2D6	Cytochrome P450 2D6 monooxygenase	P10635
78	LAMP2	Lysosome-associated membrane glycoprotein 2	P13473
79	MAG	Myelin-associated glycoprotein	P20916
80	MAGEA3	Melanoma-associated antigen 3	P43357
81	Matrigel		
82	MBP	Myelin basic protein	P02686
83	MMP2/GelA/Clg4a	Matrix metalloproteinase 2, 72 kDa type IV collagenase, Gelatinase A	P08253
84	PLP1	Myelin proteolipid protein	P60201
85	Myosin	Myosin heavy and light chain from bovine muscle	M6643
86	NSE/ENO2	Neuron Specific Enolase (NSE), Gamma-enolase	P09104
87	Nucleosome	Nucleosome from Bovin, non-recombinant	
88	NY-ESO-1	Cancer/testis antigen 1, Autoimmunogenic cancer/testis antigen NY-ESO-1	P78358
89	OX40L/TNFSF4	OX40 ligand (OX40L), Tumor necrosis factor ligand superfamily member 4	P23510

90	P53	Cellular tumor antigen p53	P04637
91	PTH	Parathyroid Hormone (PTH)	P01270
92	Parotid Secretory Protein (PSP)	BPI fold-containing family A member 2	Q96DR5
93	PD-1	Programmed cell death protein 1	Q15116
94	PD-L1	Programmed cell death 1 ligand 1	Q9NZQ7
95	Peptidyl Arginine Deiminases	Peptidyl Arginine Deiminase from rabbit skeletal muscle	P1584
96	Peroxiredoxin 1 (PRDX1)	Peroxiredoxin-1	Q06830
97	Phosphatidyl-L-serine	3-sn-Phosphatidyl-L-serine sodium salt from bovine brain	P5660
98	PKM2	Pyruvate kinase PKM	P14618
99	Prostate Specific Antigen (PSA)	Prostate-specific antigen	P07288
100	Prostate Specific Membrane	Glutamate carboxypeptidase 2	Q04609
101	Prostatic Acid Phosphatase	Prostatic acid phosphatase	P15309
102	PSMA	Glutamate carboxypeptidase 2	Q04609
103	HSPG2	Human Heparan Sulfate Proteoglycan 2/Perlecan protein	P98160
104	Ribosomal phosphoproteins	60S acidic ribosomal protein P0, P1, P2	P05388, P05386, P05387
105	SCCA/SERPINA4	Kallistatin, Serpin A4	P29622
106	ScL-70	DNA Topoisomerase I (ScL-70; full length)	P11387
107	Sm/RNP	Bovin RNP/Sm ribonuclear protein complex (U-snRNPs)	
108	SOX2	Transcription factor SOX-2	P48431
109	Sphingomyelin	N-Acyl-4-sphingenyl-1-O-phosphorylcholine, N-Acyl-D-sphingosine-1-phosphocholine	S0756
110	SSA/SSB	SSA/Ro and SSB/La	P19474, P10155, P05455
111	TGFB1	Transforming growth factor beta-1 proprotein	P01137
112	Thrombopoietin (TPO)	Thrombopoietin	P40225
113	Thyroglobulin	Thyroglobulin (TG)	P01266
114	Tissue Transglutaminase (TTG)	Tissue transglutaminase (tTG; syn. TGase-2)	P21980
115	TNF-alpha	Tumor necrosis factor	P01375
116	TNF-beta	Lymphotoxin-alpha	P01374
117	Troponin I-T-C	Troponin C, Troponin I, and Troponin T	P19429, P02585, P45379

118	UBE2S	Ubiquitin-conjugating enzyme E2 S	Q16763
119	uPA	Urokinase-type plasminogen activator	P00749
120	VEGF-165/VEGF	Vascular endothelial growth factor A	P15692
121	Ig control 1	Ig control dilution 1	
122	Ig control 2	Ig control dilution 2	
123	Ig control 3	Ig control dilution 3	
124	Ig control 4	Ig control dilution 4	
125	anti-Ig control 1	Anti-Ig control dilution 1	
126	anti-Ig control 2	Anti-Ig control dilution 2	
127	anti-Ig control 3	Anti-Ig control dilution 3	
128	anti-Ig control 4	Anti-Ig control dilution 4	

