

OmicsArray™ APS-NETs antigen Array Antigen List - Catalog No. PA004

Antigen No.	Antigen ID	Antigen Name
1	a-actin	Alpha-actinine
2	Apo B	Apolipoprotein B
3	Apo E	Apolipoprotein E
4	Azurocidin	Azurocidin
5	B -actin	β -Actin
6	B2GPI	B2 Glycoprotein I
7	B2-microglobulin	β 2-microglobulin
8	B4-integrin	β 4-integrin
9	B-glucuronidase	β -glucuronidase
10	BPI	Bactericidal/permeability-increasing protein
11	C1q	complement C1q
12	Calmodulin	Calmodulin
13	Calreticulin	Calreticulin
14	CAP1	Adenylate Cyclase-Associated Protein 1
15	CAP2	Adenylate Cyclase-Associated Protein 2
16	Cardiolipin	Cardiolipin
17	Catalase	Catalase
18	Cathepsin G	Cathepsin G
19	CCL11	Eotaxin/C-C motif chemokine 11
20	CCL2	C-C motif chemokine 2
21	CCL3	C-C motif chemokine 3
22	CCP	Cyclic citrullinated peptide
23	CENP-A	Centromere protein A (CENP-A)
24	CENP-B	Centromere protein B (CENP-B)
25	Chromatin	Chromatin
26	Collagen I	Collagen I
27	Collagen II	Collagen II
28	Collagen III	Collagen III
29	Collagen IV	Collagen IV
30	Collagen V	Collagen V
31	Collagen VI	Collagen VI
32	Coronin-1	Coronin-1
33	CTLA4	Cytotoxic T-lymphocyte protein 4
34	CXCL10	C-X-C motif chemokine 10
35	Cytokeratin 10	Cytokeratin 10
36	Cytokeratin 19	Cytokeratin 19
37	DEFA1	Defensin α 1
38	DEFA3	Defensin α 3
39	DNase I	DNase I
40	dsDNA	double strand DNA
41	EGFR	EGF receptor
42	Elastase	Elastase
43	Enolase	Enolase
44	Fibrillarin	Fibrillarin

45	Fibrinogen	Fibrinogen
46	Fibrinogen IV	Fibrinogen IV
47	Fibrinogen-cit	Fibrinogen-citrullinated
48	Fibronectin	Fibronectin
49	Filaggrin	Filaggrin
50	Filamin-A	Filamin-A
51	Flagellin	Flagellin
52	Genomic DNA	Genomic DNA
53	GM2	Monosialoganglioside GM2
54	GSTP1	Glutathione S-transferase P
55	Heparan	Heparan
56	Heparan sulfate	Heparan sulfate
57	Histone H1	Histone H1
58	Histone H2A	Histone H2A
59	Histone H2B	Histone H2B
60	Histone H3	Histone H3
61	Histone H3-cit	citrullinated Histone H3
62	Histone H4	Histone H4
63	HMGB1	High mobility group protein B1
64	HMGB2	High mobility group protein B2
65	Hsp70	Heat shock 70 kDa protein
66	HSPG	Heparan sulfate proteoglycan
67	IFN-A2	Interferon alpha-2
68	IFN-r	Interferon gamma
69	IL 1b	Interleukin-1 beta
70	IL 4	Interleukin-4
71	IL 6	Interleukin-6
72	Ku (p70/p80)	Ku(p70/p80)
73	La/SSB	Lupus La protein/Sjogren syndrome type B antigen (SS-B)
74	lactoferrin	Lactoferrin
75	Laminin	Laminin
76	Laminin r1	Laminin subunit gamma-1
77	LAMP2	Lysosome-associated membrane glycoprotein 2
78	Lipocalin-2/NGAL	Lipocalin (neutrophil gelatinase-associated)
79	LL-37	Cathelicidin antimicrobial peptide
80	Lysozyme	Lysozyme
81	MAGEA4	Melanoma-associated antigen 4
82	MDA5	Melanoma differentiation-associated protein 5 (MDA-5)
83	Mi-2	ATP-dependent helicase Mi-2
84	MMP9	Matrix metalloproteinase-9
85	MNDA	Myeloid cell nuclear differentiation antigen
86	MPO/DNA	MPO/DNA complex
87	MPO	Myeloperoxidase
88	Myosin	Myosin
89	Myosin-9	Myosin-9

90	Nucleolin	Nucleolin
91	Nucleosome	Nucleosome
92	Nup62	Nuclear pore glycoprotein p62
93	Olfactomedin-4	Olfactomedin-4
94	PAD1	Protein-arginine deiminase type-1
95	PAD2	Protein-arginine deiminase type-2
96	PAD3	Protein-arginine deiminase type-3
97	PAD4	Protein-arginine deiminase type-4
98	PCNA	Proliferating cell nuclear antigen
99	Peroxiredoxin 1	Peroxiredoxin 1
100	PGLYRP1	Peptidoglycan recognition protein 1
101	Phosphatidylinositol	Phosphatidylinositol
102	Plastin-2	Plastin-2
103	PR3	Proteinase 3
104	Profilin-1	Profilin-1
105	Resistin	Resistin
106	Ribo P0	Phosphoprotein P0
107	Ribo P1	Phosphoprotein P1
108	Ribo P1	Phosphoprotein P2
109	Ro/SSA	Ribonucleoprotein autoantigen Ro/SS-A
110	S100-A12	Protein S100-A12
111	S100-A8/A9	Protein S100-A8 and A9 complex
112	Sm/RNP	U-small nuclear riboprotein complexes
113	SmD	Small nuclear ribonucleoprotein Sm
114	SOX2	Transcription factor SOX-2
115	ssDNA	Single strand DNA
116	Total histone	Total histone
117	Transketolase	Transketolase
118	TTG	Tissue transglutaminase
119	Vimentin	Vimentin
120	Vitronectin	Vitronectin
121	Ig control_1	Ig control
122	Ig control_2	Ig control
123	Ig control_3	Ig control
124	Ig control_4	Ig control
125	anti-Ig control_1	anti-Ig control
126	anti-Ig control_2	anti-Ig control
127	anti-Ig control_3	anti-Ig control
128	anti-Ig control_4	anti-Ig control