

OmicsArray™ Systemic Autoimmune-associated Antigen Array Antigen List - Catalog No. PA001

Antigen number	Antigen Name	Category	Associated Disease
1	Aggrecan	extracellular matrix	Rheumatoid arthritis
2	AGTR	receptor	Essential hypertension
3	Alpha Fodrin	cytoskeleton	Sjogren's syndrome, Systemic lupus erythematosus
4	Alpha-actinine	cytoskeletal	Systemic lupus erythematosus, Multiple sclerosis
5	Amyloid	extracellular	Alzheimer's disease
6	AQP4	membrane	Neuromyelitis
7	B2 glycoprotein 1	extracellular	Anti phospholipid syndrome
8	B2-microglobulin	membrane	Cancer
9	B-actin	cytoskeletal	Autoimmune hepatitis
10	BPI	cytoplasmic	Irritable bowel syndrome, Systemic lupus erythematosus
11	Calprotectin	extracellular	Irritable bowel syndrome/Irritable bowel disease
12	Cardolipin	Phaspholipid	Anti phospholipid syndrome
13	CENP-A	extracellular matrix	Systemic sclerosis, Systemic lupus erythematosus, Rheumatoid arthritis
14	CENP-B	extracellular matrix	Systemic sclerosis, Systemic lupus erythematosus, Rheumatoid arthritis
15	Chondroitin Sulfate C	extracellular matrix	Rheumatoid arthritis, Lambert-Eaton syndrome
16	COL17A1	extracellular matrix	Irritable bowel syndrome
17	Collagen I	extracellular matrix	Rheumatoid arthritis, Systemic lupus erythematosus
18	Collagen II	extracellular matrix	Rheumatoid arthritis, Systemic lupus erythematosus
19	Collagen III	extracellular matrix	Rheumatoid arthritis, Systemic lupus erythematosus
20	Collagen IV	extracellular matrix	Rheumatoid arthritis, Systemic lupus erythematosus
21	Collagen V	extracellular matrix	Rheumatoid arthritis, Systemic lupus erythematosus
22	Collagen VI	extracellular matrix	Rheumatoid arthritis, Systemic lupus erythematosus
23	complement C1q	complement	Systemic lupus erythematosus, Lupus nephritis
24	complement C3/C3a	complement	Systemic lupus erythematosus, Lupus nephritis
25	Core Histone	nuclear antigen	Systemic lupus erythematosus, Lupus nephritis
26	Crohn's disease4	cytokine	Rheumatoid arthritis
27	Crohn's disease8	cytokine	Rheumatoid arthritis
28	CRP	blood factor	Rheumatoid arthritis, Systemic lupus erythematosus
29	dsDNA	nuclear antigen	Systemic lupus erythematosus

30	EBNA1	Viral antigen	Systemic lupus erythematosus
31	Elastin	matrix	Chronic obstructive pulmonary disease
32	Factor B	blood factor	Dense deposit disease
33	Factor D	blood factor	Dense deposit disease
34	Factor H	blood factor	Membranous nephropathy
35	Factor I	blood factor	Membranous nephropathy
36	Factor P	blood factor	Membranous nephropathy
37	Fibrinogen IV	extracellular	Systemic lupus erythematosus, Rheumatoid arthritis
38	Fibrinogen S	extracellular	Systemic lupus erythematosus, Rheumatoid arthritis
39	Fibronectin	extracellular	Systemic lupus erythematosus, Rheumatoid arthritis
40	GAD65	Enzyme	Type 1 diabetes, Central Nervous System
41	GBM	membrane	Lupus nephritis, Goodpasture disease
42	GDF-15	cytokine	Cardiovascular disease, Anti phospholipid syndrome
43	GM-CSF	cytokine	Multiple sclerosis, Rheumatoid arthritis
44	GP2	Glycoprotein	Anti phospholipid syndrome, Crohn's disease, Irritable bowel syndrome
45	GP210	Glycoprotein	Primary biliary cirrhosis
46	Hemocyanin	extracellular	Autoimmune hemolytic anemia
47	Heparan Sulphate	cell matrix	Systemic lupus erythematosus
48	Heparin	cell matrix	Systemic lupus erythematosus
49	Histone H1	nuclear antigen	Systemic lupus erythematosus, Connective tissue disease
50	Histone H2A	nuclear antigen	Systemic lupus erythematosus, Connective tissue disease
51	Histone H2B	nuclear antigen	Systemic lupus erythematosus, Connective tissue disease
52	Histone H3	nuclear antigen	Systemic lupus erythematosus, Connective tissue disease
53	Histone H4	nuclear antigen	Systemic lupus erythematosus, Connective tissue disease
54	IA2	extracellular	Type 1 diabetes, Cancer
55	IFN-A1	cytokine	Systemic lupus erythematosus, COVID
56	IFN-B	cytokine	Systemic lupus erythematosus, COVID
57	IFN-E	cytokine	Systemic lupus erythematosus, COVID
58	IFN-L2	cytokine	Systemic lupus erythematosus, COVID
59	IFN-r	cytokine	Systemic lupus erythematosus, COVID
60	IL-12	cytokine	Systemic lupus erythematosus, COVID
61	IL-15	cytokine	Systemic lupus erythematosus, COVID
62	Insulin	hormone	Type 1 diabetes
63	Jo-1	extracellular	Infectious mononucleosis

64	KU (P70/P80)	nuclear antigen	Systemic lupus erythematosus
65	La/SSB	nuclear antigen	Sjogren's syndrome, Systemic sclerosis
66	Laminin	extracellular matrix	Pemphigoid
67	LC1 (liver cytosol)	cytosol	Autoimmune chronic active hepatitis type 2
68	LPS	bacterial	Autoimmune arthritis
69	M2	mitochondrial	Primary biliary cirrhosis
70	MDA5	enzyme	Dermatomyositis/polyMyositis
71	Mi-2	nuclear antigen	Dermatomyositis
72	Mitochondrial antigen	mitochondrial	Primary biliary cholangitis
73	MPO	enzyme	Systemic vasculitis
74	Muscarinic receptor	receptor	Primary biliary cirrhosis, Sjogren's syndrome
75	Myelin basic protein (MBP)	membrane	Multiple sclerosis
76	Myosin	Muscle cell structure	Cardiac dysfunction
77	Myosin heavy & light chain	Muscle cell structure	Cardiac dysfunction
78	Neuropilin-1	transmembrane	cardiovascular and CNS
79	Nucleolin	nuclear antigen	Systemic lupus erythematosus
80	Nucleosome antigen	nuclear antigen	Systemic lupus erythematosus
81	Nup 62	nuclear antigen	Primary biliary cirrhosis
82	PCNA	nuclear antigen	Systemic lupus erythematosus
83	Phosphatidyl-serine	Phospholipid	Anti phospholipid syndrome
84	PL-12	tRNA Synthetase	Interstitial lung disease
85	PL-7	tRNA Synthetase	Interstitial lung disease
86	PM/Scl 100	nuclear antigen	Myositis
87	PM/Scl-75	nuclear antigen	PolyMyositis/Scleroderma
88	PR3	enzyme	Systemic vasculitis
89	Proteoglycan	extracellular matrices	Lupus nephritis
90	Prothrombin protein	blood factor	Anti phospholipid syndrome
91	Rheumatoid factor (RF)	blood factor	Systemic lupus erythematosus, Rheumatoid arthritis
92	Ribo Phosphoprotein P0	nuclear antigen	Systemic lupus erythematosus, Anti phospholipid syndrome, Rheumatoid arthritis
93	Ribo Phosphoprotein P1	nuclear antigen	Systemic lupus erythematosus, Anti phospholipid syndrome, Rheumatoid arthritis
94	Ribo Phosphoprotein P2	nuclear antigen	Systemic lupus erythematosus, Anti phospholipid syndrome, Rheumatoid arthritis
95	Ro/SSA (52 Kda)	ribonucleic	Sjogren's syndrome, Systemic sclerosis
96	Ro/SSA (60 Kda)	ribonucleic	Sjogren's syndrome, Systemic sclerosis
97	S100	cytosolic	Vascular disease, Auto-Inflammatory diseases

98	Scl-70/Topoisomerase I	cytosolic	Sjogren's syndrome
99	Sm	nuclear antigen	Systemic lupus erythematosus
100	Sm/RNP	nuclear antigen	Systemic lupus erythematosus
101	SmD	nuclear antigen	Systemic lupus erythematosus
102	SP100	Nuclear Antigen	Primary biliary cirrhosis
103	Sphingomyelin	membranes	Multiple sclerosis, Experimental autoimmune encephalitis
104	SRP54	cytosolic	Myopathy
105	ssDNA	nuclear antigen	Systemic lupus erythematosus, Systemic sclerosis
106	ssRNA	nuclear antigen	Systemic lupus erythematosus, Systemic sclerosis
107	TGF-B1	secreted protein	Systemic lupus erythematosus, Autoimmune hepatitis
108	Thyroglobulin	glycoprotein	Autoimmune thyroiditis
109	TIF1 gama	ubiquitin ligase	Myositis, Systemic lupus erythematosus, Tumor
110	TNF-a	cytokines	Vasculitis, Systemic lupus erythematosus, Interstitial lung disease
111	TPO	enzyme	Autoimmune thyroiditis
112	Troponin-I	cardiac muscle fibers	Myocardial dysfunction
113	TTG	enzyme	Celiac disease
114	Tweek	cytokines	Systemic lupus erythematosus
115	U1-snRNP 68/70	Nuclear Antigen	Connective tissue disease, Systemic lupus erythematosus
116	U1-snRNP A	Nuclear Antigen	Connective tissue disease, Systemic lupus erythematosus
117	U1-snRNP B/B'	Nuclear Antigen	Connective tissue disease, Systemic lupus erythematosus
118	U1-snRNP C	Nuclear Antigen	Connective tissue disease, Systemic lupus erythematosus
119	Vimentin	intermediate filament	Rheumatoid arthritis Sjogren's syndrome, Systemic lupus erythematosus
120	Vitronectin	extracellular matrix	Myasthenia gravis, Systemic lupus erythematosus
121	Ig control	Ig control 1	N/A
122	Ig control	Ig control 2	N/A
123	Ig control	Ig control 3	N/A
124	Ig control	Ig control 4	N/A
125	anti-Ig control	anti-Ig control 1	N/A
126	anti-Ig control	anti-Ig control 2	N/A
127	anti-Ig control	anti-Ig control 3	N/A
128	anti-Ig control	anti-Ig control 4	N/A