

OmicsArray™ Coronavirus Infection-associated Autoimmune Antigen Array Antigen List - Catalog No. PA012						
No.	Antigen Name	Alternative name / Synonyms	Category	Uniprot ID	Associate disease	Full name of Disease
1	TIF1-r	Transcription intermediary factor 1 gamma	autoAg	<a href="#">Q9UPN9</a>	Cancer-associated DM	Cancer-associated DM
2	B2-GP1 (Native)	Beta-2 glycoprotein	autoAg	<a href="#">P20231</a>	APS	Antiphospholipid syndrome
3	RNP/Sm (Native)	Small nuclear ribonucleoprotein Sm	autoAg	<a href="#">P62314</a>	SLE, RA	Systemic lupus erythematosus,rheumatoid arthritis
4	SmD (Recombinant)	Small nuclear ribonucleoprotein Sm	autoAg		SLE	Systemic lupus erythematosus
5	TPO	Thyroid Peroxidase	autoAg	<a href="#">P07202</a>	Chronic thyroiditis	Chronic thyroiditis
6	Thyroglobulin	Thyroglobulin	autoAg	<a href="#">P01266</a>	Thyroid cancer	Thyroid cancer
7	dsDNA	double stranded DNA	autoAg		SLE	Systemic lupus erythematosus
8	ScI-70/Topo Isomerase I	topoisomerase I	autoAg	<a href="#">P11387</a>	SLE, SSc	Systemic lupus erythematosus, Systemic sclerosis
9	CENP-A	Centromere protein A	autoAg	<a href="#">P49450</a>	SSc	Systemic sclerosis
10	CENP-B	Centromere protein B	autoAg	<a href="#">P07199</a>	SS, Breast cancer	Sjögren's syndrome, Breast cancer
11	Ro-52	52 kDa ribonucleoprotein autoantigen Ro/SS-A	autoAg	<a href="#">P19474</a>	SS	Secondary Sjögren's syndrome
12	Ro-60 (native)	RNA-binding protein RO60	autoAg	<a href="#">P10155</a>	SLE,SS	Systemic lupus erythematosus and Sjogren's
13	La/SS-B	Lupus La protein/Sjogren syndrome type B antigen	autoAg	<a href="#">P05455</a>	SLE,SS	Systemic lupus erythematosus and Sjogren's syndrome
14	Jo-1	Histidyl-tRNA synthetase	autoAg	<a href="#">P12081</a>	Antisynthetase syndrome	Antisynthetase syndrome
15	U1-snRNP-68	U1 small nuclear ribonucleoprotein 70 kDa	autoAg	<a href="#">P08621</a>	SLE	Systemic lupus erythematosus
16	U1-snRNP-A	U1 small nuclear ribonucleoprotein A	autoAg	<a href="#">P09012</a>	SLE	Systemic lupus erythematosus
17	U1-snRNP-C	U1 small nuclear ribonucleoprotein C	autoAg	<a href="#">P09234</a>	SLE	Systemic lupus erythematosus
18	U1-snRNP-BB'	Small nuclear ribonucleoprotein-associated proteins B and B'	autoAg	<a href="#">P14678</a>	SLE	Systemic lupus erythematosus
19	LC1	Liver cytosol type 1	autoAg	<a href="#">Q95954</a>	AIH	Autoimmune hepatitis
20	GAD65	Glutamate decarboxylase 65 kDa isoform	autoAg	<a href="#">Q05329</a>	T1D, SPS	Type 1 diabetes, Stiff-person syndrome
21	RPLP0	Ribosomal protein P0	autoAg	<a href="#">P05388</a>	SLE, Gynecologic tumors	Systemic lupus erythematosus, Gynecologic tumors
22	RPLP1	Ribosomal protein P1	autoAg	<a href="#">P05386</a>	Gynecologic tumors	Gynecologic tumors
23	RPLP2	Ribosomal protein P2	autoAg	<a href="#">P05387</a>	Gynecologic tumors	Gynecologic tumors
24	B2GP1 (Recombinant)	Beta 2 glycoprotein 1	autoAg	<a href="#">P02749</a>	APS	Antiphospholipid syndrome
25	TTG	Tissue transglutaminase	autoAg	<a href="#">P21980</a>	CD	celiac disease
26	PCNA	Proliferating cell nuclear antigen	autoAg	<a href="#">P12004</a>	SLE	Systemic lupus erythematosus
27	PL-7	Threonyl-tRNA synthetase	autoAg	<a href="#">Q9BW92</a>	IIM	Idiopathic inflammatory myopathies
28	PL-12	Alanyl-tRNA synthetase	autoAg	<a href="#">Q5JT29</a>	IIM	Idiopathic inflammatory myopathies
29	PM/ScI100	Polymyositis/scleroderma autoantigen 100 kDa	autoAg	<a href="#">Q01780</a>	SSc	Systemic sclerosis
30	IF	Intrinsic factor	autoAg	<a href="#">P27352</a>	Autoimmune atrophic gastritis	Autoimmune atrophic gastritis
31	GBM	glomerular basement membrane	autoAg		Anti-Glomerular Basement Membrane Disease	Anti-Glomerular Basement Membrane Disease
32	Neuropilin-1	Neuropilin-1	autoAg	<a href="#">Q14786</a>	LN	Lupus Nephritis
33	PM/ScI75	Polymyositis/scleroderma autoantigen 75 kDa	autoAg	<a href="#">Q06265</a>	SSc	Systemic sclerosis
34	Ku (p70/p80)	ATP-dependent DNA helicase 2	autoAg	<a href="#">P12956</a>	SLE	Systemic lupus erythematosus
				<a href="#">P13010</a>		
35	Sm (Native)	Small nuclear ribonucleoprotein	autoAg		SLE	Systemic lupus erythematosus
36	BCOADC-E2	Branched chain 2-oxo-acid dehydrogenase complex component	autoAg	<a href="#">P11182</a>	Primary biliary cirrhosis	Primary biliary cirrhosis
37	OGDC-E2	2-oxoglutarate dehydrogenase complex component E2	autoAg	<a href="#">P36957</a>	Primary biliary cirrhosis	Primary biliary cirrhosis
38	PDC-E2	Pyruvate dehydrogenase complex component E2	autoAg	<a href="#">P10515</a>	Primary biliary cirrhosis	Primary biliary cirrhosis
39	M2 (AMA-M2)	mitochondrial subtype 2	autoAg	<a href="#">P11182</a>	Myopathies	Myopathies
40	Mi-2	Mi-2 autoantigen 218 kDa protein	autoAg	<a href="#">Q14839</a>	Dermatomyositis	Dermatomyositis
41	SRP54	Signal recognition particle 54 kDa protein	autoAg	<a href="#">P61011</a>	Myopathies	Myopathies
42	MPO	Myeloperoxidase	autoAg	<a href="#">P05164</a>	Glomerulonephritis	Glomerulonephritis
43	PR3	Leukocyte proteinase 3	autoAg	<a href="#">P24158</a>	Vasculiti	Vasculiti
44	SP100	Nuclear autoantigen Sp-100	autoAg	<a href="#">P23497</a>	PBC	Primary biliary cholangitis
45	Gp210	Nuclear pore membrane glycoprotein 210	autoAg	<a href="#">Q8TEM1</a>	PBC	Primary biliary cholangitis
46	BPI	bactericidal/permeability-increasing protein (BPI)	autoAg	<a href="#">P17213</a>	Cystic fibrosis, inflammatory bowel disease, vasculitis, and primary sclerosing cholangitis	Cystic fibrosis, inflammatory bowel disease, vasculitis, and primary sclerosing cholangitis
47	Nup62	Nuclear pore glycoprotein p62	autoAg	<a href="#">P37198</a>	AGS	Aicardi-Goutières syndrome
48	GDF-15	Growth/differentiation factor 15	autoAg	<a href="#">Q99988</a>	SLE, myositis	Systemic lupus erythematosus, myositis
49	GP2	Glycoprotein 2	autoAg	<a href="#">P55259</a>	Primary sclerosing cholangitis	Primary sclerosing cholangitis
50	LKM-1/CYP2D6	Cytochrome P450 IID6	autoAg	<a href="#">P10635</a>	Autoimmune hepatitis	Autoimmune hepatitis
51	MDA5	Melanoma differentiation-associated protein 5	autoAg	<a href="#">Q9BYX4</a>	Dermatomyositis	Dermatomyositis
52	DFS70	Dense fine speckles 70 kDa protein	autoAg	<a href="#">Q75475</a>	AARD	ANA associated rheumatic diseases

53	Insulinoma-associated (IA-2)	Insulinoma-associated protein 2	autoAg	<a href="#">Q16849</a>	T1D	Type 1 diabetes
54	Nucleosome		autoAg		SLE	Systemic lupus erythematosus
55	Histone	Histone	autoAg	<a href="#">P16403</a>	SLE, RA	Systemic lupus erythematosus, Rheumatoid arthritis
56	Laminin	Laminin	autoAg		AIP	Autoimmune pancreatitis
57	Collagen V	Collagen V	autoAg	<a href="#">P20908</a>	SSc	Systemic sclerosis
58	Collagen VI	Collagen VI	autoAg	<a href="#">P12109</a>	Vasculitis, SLE	Vasculitis, Systemic lupus erythematosus
59	Vimentin	Vimentin	autoAg	<a href="#">P08670</a>	Secondary Sjögren's syndrome (SSS), RA	Secondary Sjögren's syndrome, Rheumatoid arthritis
60	Troponin I	Troponin I	autoAg	<a href="#">P19429</a>	DCM	Dilated cardiomyopathy
61	Aldolase C		autoAg	<a href="#">P09972</a>	Diabetic retinopathy	Diabetic retinopathy
62	ARS	Aminoacyl-tRNA Synthetase	autoAg		Myositis	Myositis
63	Complement C1q	Complement C1q	autoAg	<a href="#">P02745</a>	SLE	Systemic lupus erythematosus
				<a href="#">P02746</a>		
				<a href="#">P02747</a>		
64	HSPG	Heparan sulfate proteoglycan	autoAg	<a href="#">P34741</a>	SLE	Systemic lupus erythematosus
65	Vitronectin	Vitronectin	autoAg	<a href="#">P04004</a>	PBC, SLE	Primary biliary cirrhosis, Systemic lupus
66	Fibrinogen	Fibrinogen	autoAg	<a href="#">P02671</a>	SLE	Systemic lupus erythematosus
				<a href="#">P02675</a>		
67	rHuPTH	Parathyroid Hormone	autoAg	<a href="#">P01270</a>	IHP	Idiopathic hypoparathyroidism
68	EJ	Glycyl-tRNA Synthetase	autoAg	<a href="#">P41250</a>	Myositis	Myositis
69	KS/AsnRS	Asparaginyl-tRNA Synthetase	autoAg	<a href="#">O43776</a>	ILD	Myositis and interstitial lung disease
70	Calprotectin	Calprotectin	autoAg	<a href="#">P05109</a>	Myasthenia gravis, RA	Myasthenia gravis, Rheumatoid arthritis
				<a href="#">P06702</a>		
71	Rheumatoid Factor (RF)	Rheumatoid Factor (RF)	autoAg		RA	Rheumatoid arthritis
72	ACE2	angiotensin-converting enzyme 2	autoAg	<a href="#">Q9BYF1</a>	COVID-19, Parkinson's disease	COVID-19, Parkinson's disease
73	BAFF	B cell-activating factor (BAFF)	cytokines	<a href="#">Q9Y275</a>	SLE,	Systemic lupus erythematosus
74	CD4	T-cell surface glycoprotein CD4	cytokines	<a href="#">P01730</a>	SLE, HIV infection	Systemic lupus erythematosus, HIV infection
75	CD8	T-cell surface glycoprotein CD8	cytokines	<a href="#">P01732</a>	SLE, HIV infection	Systemic lupus erythematosus, HIV infection
				<a href="#">P10966</a>		
76	GM-CSF	Granulocyte-macrophage colony-stimulating factor	cytokines	<a href="#">P04141</a>	PAP	Pulmonary alveolar proteinosis
77	IFN-a		cytokines		SLE	Systemic lupus erythematosus
78	IFN-a1	Interferon alpha-1/13	cytokines	<a href="#">P01562</a>	SLE	Systemic lupus erythematosus
79	IFN-a2	Interferon alpha-2	cytokines	<a href="#">P01563</a>	SLE, Esophageal cancer	Systemic lupus erythematosus, Esophageal cancer
80	IFN-B	Interferon beta	cytokines	<a href="#">P01574</a>	Multiple sclerosis	Multiple sclerosis
81	IFN-B1	Interferon beta	cytokines	<a href="#">P01574</a>	SLE	Systemic lupus erythematosus
82	IFN-gamma	Interferon gamma	cytokines	<a href="#">P01579</a>	Nontuberculous mycobacterial infection	Nontuberculous mycobacterial infection
83	IFN-lambda1	Interferon lambda-1	cytokines	<a href="#">Q8IU54</a>	RA, CLE	Rheumatoid arthritis, Cutaneous lupus erythematosus (CLE)
84	IFN-lambda2/IL28A	Interferon lambda-2	cytokines	<a href="#">Q8IZJ0</a>	Cancer and autoimmunity	Cancer and autoimmunity
85	IFN-lambda3	Interferon lambda-3	cytokines	<a href="#">Q8IZI9</a>	Cancer and autoimmunity	Cancer and autoimmunity
86	IFN-epsilon	Interferon epsilon	cytokines	<a href="#">Q86WN2</a>	COVID-19	COVID-19
87	IFN-omega	Interferon omega-1	cytokines	<a href="#">P05000</a>	COVID-19	COVID-19
88	IL-10	Interleukin-10	cytokines	<a href="#">P22301</a>	SLE	Systemic lupus erythematosus
89	IL-12 (p70)	Interleukin-12	cytokines	<a href="#">P29459</a>	SLE	Systemic lupus erythematosus
				<a href="#">P29460</a>		
90	IL-15	Interleukin-15	cytokines	<a href="#">P40933</a>	RA	Rheumatoid arthritis
91	IL-17A	Interleukin-17A	cytokines	<a href="#">Q16552</a>	CMC	Chronic mucocutaneous candidiasis
92	IL-17F	Interleukin-17F	cytokines	<a href="#">Q96PD4</a>	CMC	Chronic mucocutaneous candidiasis
93	IL-1a	Interleukin-1 alpha	cytokines	<a href="#">P01583</a>	RA	Rheumatoid arthritis
94	IL-1b	Interleukin-1 beta	cytokines	<a href="#">P01584</a>	RA	Rheumatoid arthritis
95	IL-2	Interleukin-2	cytokines	<a href="#">P60568</a>	T1D	Type 1 diabetes
96	IL-22	Interleukin-22	cytokines	<a href="#">Q9GZX6</a>	APECED	Autoimmune polyendocrinopathy-candidiasis-ectodermal dystrophy
97	IL-4	Interleukin-4	cytokines	<a href="#">P05112</a>	Myasthenia gravis	Myasthenia gravis
98	IL-6	Interleukin-6	cytokines	<a href="#">P05231</a>	APECED, SSc	autoimmune polyendocrinopathy-candidiasis-ectodermal dystrophy (APECED) ;systemic sclerosis
99	IL-8	Interleukin-8	cytokines	<a href="#">P10145</a>	ARDS, Helicobacter pylori infection	acute respiratory distress syndrome (ARDS) ; Helicobacter pylori infection
100	TGFB1	Transforming growth factor beta-1 proprotein	cytokines	<a href="#">P01137</a>	SLE	Systemic lupus erythematosus
101	TNF-B (LT-a)	TNF-beta(Lymphotoxin-alpha)	cytokines	<a href="#">P01374</a>	SLE	Systemic lupus erythematosus
102	TWEAK (CD255)	TNF weakly inducer of apoptosis	cytokines	<a href="#">Q43508</a>	Psoriatic arthritis (PsoA); systemic lupus erythematosus (SLE) and primary Sjögren's syndrome	Psoriatic arthritis (PsoA); systemic lupus erythematosus (SLE) and primary Sjögren's syndrome
103	SARS1 S1	Spike glycoprotein	viral Ags	<a href="#">P59594</a>	SARS	Severe acute respiratory syndrome coronavirus
104	SARS2 S	Spike glycoprotein	viral Ags	<a href="#">P0DTC2</a>	COVID-19	COVID-19
105	SARS S2	Spike protein S2	viral Ags	<a href="#">P59594</a>		Severe acute respiratory syndrome coronavirus
106	SARS2 S1RBD	SARS-CoV-2 spike receptor binding domain	viral Ags	<a href="#">P0DTC2</a>	COVID-19	COVID-19

107	SARS2 S1	Spike glycoprotein	viral Ags	<a href="#">P0DTC2</a>	COVID-19	COVID-19
108	SARS2 NCP	Spike glycoprotein	viral Ags	<a href="#">P0DTC2</a>	COVID-19	COVID-19
109	MERS S	Spike glycoprotein	viral Ags	<a href="#">K9N5Q8</a>	MERS-CoV	Middle East respiratory syndrome coronavirus
110	HCOV-NL63 S	Spike glycoprotein	viral Ags	<a href="#">P0C6U6</a>	SARS	Severe acute respiratory syndrome coronavirus
111	HCOV-229E S	Spike glycoprotein	viral Ags	<a href="#">P15423</a>	SARS	Severe acute respiratory syndrome coronavirus
112	HCOV-HKU1 S	Spike glycoprotein	viral Ags	<a href="#">Q5MQD0</a>	SARS	Severe acute respiratory syndrome coronavirus
113	HCOV-OC43 S	Spike glycoprotein	viral Ags	<a href="#">P36334</a>	SARS	Severe acute respiratory syndrome coronavirus
114	SARS2 S (B.1.1.529)	Spike glycoprotein	viral Ags	<a href="#">P0DTC2</a>	COVID-19	COVID-19
115	IAV (H1N1)	Influenza A virus	viral Ags		Influenza A	Influenza A
116	IAV (H3N2)	Influenza A virus	viral Ags		Influenza A	Influenza A
117	IBV (Influenza B)	Influenza B virus	viral Ags		Influenza B	Influenza B
118	RSV	Respiratory syncytial virus (RSV)	viral Ags	<a href="#">Q85431</a>	Respiratory illness among children and infants	Respiratory illness among children and infants
119	CMV	Cucumber mosaic virus	viral Ags	<a href="#">P17769</a> <a href="#">P0C783</a> <a href="#">P16490</a>	SLE, Guillain-Barré syndrome	Systemic lupus erythematosus, Guillain-Barre syndrome
120	EBV	Epstein-Barr virus	viral Ags		RA, SLE	Rheumatoid arthritis, Systemic lupus erythematosus
121	Ig control 1		control			
122	Ig control 2		control			
123	Ig control 3		control			
124	Ig control 4		control			
125	anti-Ig control 1		control			
126	anti-Ig control 2		control			
127	anti-Ig control 3		control			
128	anti-Ig control 4		control			