

PA001-V2 Systemic Autoimmune-associated Antigen Array (120-Plex)					
Antigen Number	Antigen ID	Protein Name	Alternative name / Synonyms	Gene symbol	UniProt ID
1	Aggrecan	Aggrecan	cartilage-specific proteoglycan core protein (CSPCP) or chondroitin sulfate proteoglycan 1	ACAN	P16112
2	AGTR	Type-1 angiotensin II receptor	Type-1 angiotensin II receptor	AGTR1	P30556
3	Alpha Fodrin	Alpha Fodrin	Spectrin alpha chain, non-erythrocytic 1	SPTAN1	Q13813
4	Alpha-actinine	Alpha-actinine	Alpha-actinin-1; F-actin cross-linking protein	ACTN1	P12814
5	Amyloid	Amyloid	Amyloid-beta precursor; proteinAlzheimer disease amyloid protein; Protease nexin-II	APP	P05067
6	AQP4	Aquaporin-4	Aquaporin-4	AQP4	P55087
7	B2 glycoprotein 1	B2-glycoprotein 1 (B2GPI)	Apolipoprotein H (APO-H)	APOH	P02749
8	B2-microglobulin	B2-microglobulin	Beta-2-microglobulin	B2M	P61769
9	B-actin	B-actin	Actin, cytoplasmic 1	ACTB	P60709
10	BPI	Bactericidal permeability-increasing protein	Bactericidal permeability-increasing protein	BPI	P17213
11	Calprotectin	Calprotectin	Protein S100, calgranulin, leucocyte L1 complex, migration inhibitory factor-related protein, MRP8/14, S100 calcium-binding protein	S100A8	P05109
				S100A9	P06702
12	Cardiolipin	Cardiolipin	Diphosphatidylglycerol	Cardiolipin sodium salt from bovine heart	C0563
13	CD4	T-cell surface glycoprotein CD4	T-cell surface glycoprotein CD4	CD4	P01730
14	CD8	T-cell surface glycoprotein CD8 beta chain	T-cell surface glycoprotein CD8 beta chain	CD8B	P10966
15	CENP-A	Centromere protein A	Histone H3-like centromeric protein A, Centromere protein A	CENPA	P49450
16	CENP-B	Centromere protein B	Centromere protein B	CENPB	P07199
17	Chondroitin Sulfate C	Chondroitin Sulfate C	Chondroitin Sulfate Proteoglycan (CSPG), sulfated glycosaminoglycan polysaccharide	Chondroitin sulfate sodium salt from shark cartilage	C4384
18	COL17A1	Collagen alpha-1(XVII) chain	COL17A1; collagen, type XVII, alpha 1; BPAG2	COL17A1	Q9UMD9
19	Collagen I	Collagen type I	Collagen alpha-1(I) chain	COL1A1	P02452
20	Collagen II	Collagen Type II	Collagen alpha-1(II) chain	COL2A1	P02458
21	Collagen III	Collagen type III	Collagen alpha-1(III) chain/COL3A1	COL3A1	P02461
22	Collagen IV	Collagen type IV	Collagen alpha-1(IV) chain	COL4A1	P02462
23	Collagen V	Collagen type V	Collagen alpha-1(V) chain	COL5A1	P20908
24	Collagen VI	Collagen type VI	Collagen alpha-1(VI) chain	COL6A1	P12109
25	complement C1q	complement C1q	Complement C1q subcomponent subunit A	C1QA	P02745
26	complement C3/C3a	complement C3/C3a	Complement C3	C3	P01024
27	Core Histone	Core Histone	H1, H2A, H2B, H3 and H4	NA	
28	CRP	C-reactive protein	C-reactive protein	CRP	P02741
29	dsDNA	double strand DNA	double strand DNA	NA	
30	EBNA1	Epstein-Barr nuclear antigen 1	Epstein-Barr nuclear antigen 1	EBNA1	Q3KSS4
31	Elastin	Elastin	Tropoelastin	ELN	P15502
32	Factor B	Complement factor B	Complement factor B	CFB	P00751
33	Factor D	Complement factor D	Complement factor D	CFD	P00746
34	Factor H	Complement factor H	Complement factor H	CFH	P08603
35	Factor I	Complement factor I	Complement factor I	CFI	P05156

36	Factor P	Complement factor P	Complement factor P/Properdin	CFP	P27918
37	Fibrinogen IV	Fibrinogen IV	Fibrinogen alpha chain/Fibrinogen beta chain	FGA	P02672
38	Fibrinogen S	Fibrinogen Type I-S	Fibrinogen gamma-B chain	FGB	P02676
					P12799
39	Fibronectin	Fibronectin	Fibronectin	FN1	P02751
40	GAD65	Glutamate decarboxylase 2	Glutamate decarboxylase 2	GAD2/GAD65	Q05329
41	GBM	Glomerular Basement Membrane	Glomerular Basement Membrane		
42	GDF-15	Growth/differentiation factor 15	Growth/differentiation factor 15	GDF15	Q99988
43	GM-CSF	Granulocyte-macrophage colony-stimulating factor	Granulocyte-macrophage colony-stimulating factor	CSF2/GM-CSF	P04141
44	GP2	Pancreatic secretory granule membrane major glycoprotein GP2	Pancreatic secretory granule membrane major glycoprotein GP2	GP2	P55259
45	GP210	Nuclear pore membrane glycoprotein 210	Nuclear pore membrane glycoprotein 210	NUP210/gp210	Q8TEM1
46	Hemocyanin	Hemocyanin	Hemocyanin		
47	Heparan Sulphate	Heparan Sulphate	Basement membrane-specific heparan sulfate proteoglycan core protein	HSPG2	P98160
48	Heparin	Heparin	Heparin		
49	Histone H1	Histone H1	Histone H1 human	H1-0	P07305
50	Histone H2A	Histone H2A	Histone H2A type 1	H2AC11	P0C0S8
51	Histone H2B	Histone H2B	Histone H2B	Histone H2B type 1-C/E/F/G/I	P62807
52	Histone H3	Histone H3	Histone H3.1	H3C1	P68431
53	Histone H4	Histone H4	Histone H4	H4C1	P62805
54	IA2	Insulinoma-associated protein (IA-2)	islet cell antigen 512 (ICA 512); Receptor-type tyrosine-protein phosphatase-like N	PTPRN	Q16849
55	IFN-A1	Interferon alpha-1/13	Interferon alpha-1/13	IFNA1	P01562
56	IFN-B	Interferon beta	Interferon beta	IFNB1	P01574
57	IFN-E	Interferon epsilon	Interferon epsilon	IFNE	Q86WN2
58	IFN-L2	Interferon lambda-2	Interferon lambda-2	IFNL2	Q8IZJ0
59	IFN-r	Interferon gamma	Interferon gamma	IFNG	P01579
60	IL-12	Interleukin-12 subunit alpha/beta	Interleukin-12 subunit alpha/beta	IL12A	P29459
				IL12B	P29460
61	IL-15	Interleukin-15	Interleukin-15	IL15	P40933
62	Insulin	Insulin	Insulin	INS	P01308
63	Jo-1	Histidyl-tRNA Synthetase (Jo-1)	Histidine--tRNA ligase, cytoplasmic	HARS1	P12081
64	KU (P70/P80)	Human KU protein, composed of P70 and P80 subunits	XRCC5 and XRCC6	XRCC5	P12956
				XRCC6	P13010
65	La/SSB	Lupus La protein homolog	Lupus La protein homolog	SSB	P10881
66	Laminin	Laminin	Laminin subunit gamma-1	LAMC1	P11047
67	LC1 (liver cytosol)	LC1 (liver cytosol)	Formimidoyltransferase-cyclodeaminase	FTCD	Q95954
68	LPS	Lipopolysaccharide	Lipopolysaccharide		
69	M2	Mitochondrial 2-oxo acid dehydrogenase multienzyme complexes	PDC-E2, OGDC-E2, and BCOADC-E2	DLAT	P10515
70	MDA5	melanoma differentiation-associated protein 5 (MDA5; CADM 140)	Interferon-induced helicase C domain-containing protein 1	IFIH1	Q9BYX4

71	Mi-2	nucleosome remodeling deacetylase (NuRD) complex, subunit	Chromodomain-helicase-DNA-binding protein	CHD3	Q12873
				CHD4	Q14839
72	Mitochondrial antigen	Mitochondrial antigen	Pyruvate dehydrogenase complex, mitochondrial	PDH	P11966
					Q9N1X8
73	MPO	Myeloperoxidase	Myeloperoxidase	MPO	P05164
74	Muscarinic receptor	Muscarinic receptor	Muscarinic acetylcholine receptor M5	CHRM5	P08912
75	Myelin basic protein (MBP)	Myelin basic protein (MBP)	Myelin basic protein	MBP	P02686
76	Myosin	Myosin	Myosin Calcium activated from bovine muscle	MYH	
77	Myosin heavy & light chain	Myosin heavy and light chain	Myosin Heavy and Light Chain from human heart	MYH1	P12882
					P05976
					P08590
					P10916
78	Neuropilin-1	Neuropilin-1	Neuropilin-1	NRP1	Q14786
79	Nucleolin	Nucleolin	Nucleolin	NCL	P19338
80	Nucleosome antigen	Nucleosome antigen	Nucleosome (Native. Purified from bovine tissues)		
81	Nup 62	Nucleoporin Nuclear pore glycoprotein p62	Nucleoporin Nuclear pore glycoprotein p62	Nup62	P37198
82	PCNA	Proliferating cell nuclear antigen	Proliferating cell nuclear antigen	PCNA	P12004
83	Phosphatidyl-serine	1,2-Diacyl-sn-glycero-3-phospho-L-serine	1,2-Diacyl-sn-glycero-3-phospho-L-serine	DGPS	
84	PL-12	Alanyl-tRNA Synthetase (PL-12)	Alanyl-tRNA Synthetase (PL-12)	AARS1	P49588
85	PL-7	Threonyl-tRNA Synthetase (PL-7)	Threonyl-tRNA Synthetase (PL-7)	TARS1	P26639
86	PM/Scf 100	Exosome component 10	Exosome component 10	EXOSC10	Q01780
87	PM/Scf-75	Exosome complex component RRP45	Exosome complex component RRP45	EXOSC9	Q06265
88	PR3	Proteinase 3/Myeloblastin	Proteinase 3/Myeloblastin	PRTN3	P24158
89	Proteoglycan	Proteoglycan 4	Proteoglycan 4	PRG4	F1MDS0
90	Prothrombin protein	Prothrombin	Prothrombin	F2	P00734
91	Rheumatoid factor (RF)	Rheumatoid factor (RF)	Human IgG Fc fragment		
92	Ribo Phosphoprotein P0	Ribo Phosphoprotein P0	Ribosomal Phosphoprotein P0	RPLP0	P05388
93	Ribo Phosphoprotein P1	Ribo Phosphoprotein P1	Ribosomal Phosphoprotein P1	RPLP1	P05386
94	Ribo Phosphoprotein P2	Ribo Phosphoprotein P2	Ribosomal Phosphoprotein P2	RPLP2	P05387
95	Ro/SSA (52 Kda)	E3 ubiquitin-protein ligase TRIM21	E3 ubiquitin-protein ligase TRIM21	TRIM21	P19474
96	Ro/SSA (60 Kda)	RNA-binding protein RO60	RNA-binding protein RO60	RO60	P10155
97	S100	S100 Calcium Binding Protein B	S100 Calcium Binding Protein B	S100B	P04271
98	Scf-70 /Topoisomerase I	DNA Topoisomerase I (Scf-70; full length)	DNA Topoisomerase I (Scf-70; full length)	TOP1	P11387
99	Sm	Small nuclear ribonucleoprotein Sm from Bovine tissue	Small nuclear ribonucleoprotein Sm from Bovine tissue	Sm	
100	Sm/RNP	RNP/Sm ribonucleoprotein complex native protein from bovine tissue	RNP/Sm ribonucleoprotein complex native protein from bovine tissue	RNP/Sm	
101	SmD	Small nuclear ribonucleoprotein Sm, recombinant D1,D2,D3 mixture	Small nuclear ribonucleoprotein Sm, recombinant D1,D2,D3 mixture	SNRPD1	P62314
102	SP100	Nuclear dot-associated Sp100 protein; Speckled 100 kDa	Nuclear dot-associated Sp100 protein; Speckled 100 kDa	SP100	P23497
103	Sphingomyelin	N-Acyl-D-sphingosine-1-phosphocholine, from egg yolk	N-Acyl-D-sphingosine-1-phosphocholine, from egg yolk		

104	SRP54	Signal recognition particle 54 kDa protein, SRP54	Signal recognition particle 54 kDa protein, SRP54	SRP54	P61011
105	ssDNA	Single strand DNA	Single strand DNA		
106	ssRNA	Single strand RNA	Single strand RNA		
107	TGF-B1	Transforming growth factor beta-1 proprotein	Transforming growth factor beta-1 proprotein,	TGFB1; TGFB	P01137
108	Thyroglobulin	Native Thyroglobulin Isolated from human thyroid glands	Native Thyroglobulin Isolated from human thyroid glands	TG	P01266
109	TIF1 gama	E3 ubiquitin-protein ligase TRIM33	E3 ubiquitin-protein ligase TRIM33	TRIM33; KIAA1113, RFG7, TIF1G	Q9UPN9
110	TNF-a	Tumor necrosis factor	Tumor necrosis factor	TNF TNFA, TNFSF2	P01375
111	TPO	Thyroid peroxidase; TPO	Thyroid peroxidase; TPO	TPO	P07202
112	Troponin-I	Troponin I, cardiac muscle	Troponin I, cardiac muscle	TNNI3 TNNC1	P19429
113	TTG	Tissue transglutaminase (Ttg)/Protein-glutamine gamma-glutamyltransferase	Tissue transglutaminase (Ttg)/Protein-glutamine gamma-glutamyltransferase 2 (TGase-2)	TGM2	P21980
114	Tweek	Tumor necrosis factor ligand superfamily, member 12, TNFSF12, YNF-	Tumor necrosis factor ligand superfamily, member 12, TNFSF12, YNF-related weak inducer of apoptosis, Apo-3 ligand (Apo3L).	TNFSF12	O43508
115	U1-snRNP 68/70	U1 small nuclear ribonucleoprotein 70 kD, U1 snRNP 70 kDa, U1-70K, snRNP70	U1 small nuclear ribonucleoprotein 70 kD, U1 snRNP 70 kDa, U1-70K, snRNP70	SNRNP70 RNPU1Z, RPU1, SNRP70, U1AP1	P08621
116	U1-snRNP A	U1 small nuclear ribonucleoprotein A, U1 snRNP A, U1-A, U1A	U1 small nuclear ribonucleoprotein A, U1 snRNP A, U1-A, U1A	SNRPA	P09012
117	U1-snRNP B/B'	Small nuclear ribonucleoprotein-associated proteins B and B', snRNP-B,	Small nuclear ribonucleoprotein-associated proteins B and B', snRNP-B, Sm protein B/B', Sm-B/B', SmB/B'	SNRPB, COD, SNRPB1	P14678
118	U1-snRNP C	U1 small nuclear ribonucleoprotein C, U1 snRNP C, U1-C, U1C	U1 small nuclear ribonucleoprotein C, U1 snRNP C, U1-C, U1C	SNRPC	P09234
119	Vimentin	Vimentin	Vimentin	VIM	P08670
120	Vitronectin	Vitronectin	Vitronectin, VN, S-protein, Serum-spreading factor, V75	VTN	P04004
121	Ig control 1	Ig control 1	Human and mouse Ig control mix		
122	Ig control 2	Ig control 2	Human and mouse Ig control mix		
123	Ig control 3	Ig control 3	Human and mouse Ig control mix		
124	Ig control 4	Ig control 4	Human and mouse Ig control mix		
125	anti-Ig control 1	anti-Ig control 1	Anti-human and anti-mouse Ig control mix		
126	anti-Ig control 2	anti-Ig control 2	Anti-human and anti-mouse Ig control mix		
127	anti-Ig control 3	anti-Ig control 3	Anti-human and anti-mouse Ig control mix		
128	anti-Ig control 4	anti-Ig control 4	Anti-human and anti-mouse Ig control mix		