

PA001-V3 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	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	AlaRS	Aminoacyl tRNA synthetase	Aminoacyl-tRNA Synthetase from Escherichia coli	alaS	P00957
8	AsnRS/KS	Asparaginyl-tRNA Synthetase	Human Asparaginyl-tRNA Synthetase (KS)	NARS1	Q43776
9	B2 glycoprotein 1	B2-glycoprotein 1 (B2GPI)	Apolipoprotein H (APO-H)	APOH	P02749
10	B2-microglobulin	B2-microglobulin	Beta-2-microglobulin	B2M	P61769
11	B-actin	B-actin	Actin, cytoplasmic 1	ACTB	P60709
12	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
14	Cardiolipin	Cardiolipin	Diphosphatidylglycerol	Cardiolipin sodium salt from bovine heart	C0563
15	CCP	Cyclic citrullinated peptide			
16	CD4	T-cell surface glycoprotein CD4	T-cell surface glycoprotein CD4	CD4	P01730
17	CD8	T-cell surface glycoprotein CD8 beta chain	T-cell surface glycoprotein CD8 beta chain	CD8B	P10966
18	CENP-A	Centromere protein A	Histone H3-like centromeric protein A, Centromere protein A	CENPA	P49450
19	CENP-B	Centromere protein B	Centromere protein B	CENPB	P07199
20	Chondroitin Sulfate C	Chondroitin Sulfate C	Chondroitin Sulfate Proteoglycan (CSPG), sulfated glycosaminoglycan polysaccharide	Chondroitin sulfate sodium salt from shark	C4384
21	COL17A1	Collagen alpha-1(XVII) chain	COL17A1; collagen, type XVII, alpha 1; BPAG2	COL17A1	Q9UMD9
22	Collagen I	Collagen type I	Collagen alpha-1(I) chain	COL1A1	P02452
23	Collagen II	Collagen Type II	Collagen alpha-1(II) chain	COL2A1	P02458
24	Collagen III	Collagen type III	Collagen alpha-1(III) chain/COL3A1	COL3A1	P02461
25	Collagen IV	Collagen type IV	Collagen alpha-1(IV) chain	COL4A1	P02462
26	Collagen V	Collagen type V	Collagen alpha-1(V) chain	COL5A1	P20908
27	Collagen VI	Collagen type VI	Collagen alpha-1(VI) chain	COL6A1	P12109
28	complement C1q	complement C1q	Complement C1q subcomponent subunit A	C1QA	P02745
29	complement C3/C3a	complement C3/C3a	Complement C3	C3	P01024
30	Core Histone	Core Histone	H2A, H2B, H3 and H4		
31	CRP	C-reactive protein	C-reactive protein	CRP	P02741
32	DFS70	Dense fine speckles 70 kDa protein	Transcriptional coactivator p75/p52	PSIP1	Q75457
33	dsDNA	double strand DNA	double strand DNA		
34	Elastin	Elastin	Tropoelastin	ELN	P15502
35	Factor B	Complement factor B	Complement factor B	CFB	P00751
36	Factor D	Complement factor D	Complement factor D	CFD	P00746
37	Factor H	Complement factor H	Complement factor H	CFH	P08603
38	Factor I	Complement factor I	Complement factor I	CFI	P05156
39	Factor P	Complement factor P	Complement factor P/Properdin	CFP	P27918
40	FARS2	Phenylalanine--tRNA ligase, mitochondrial	Phenylalanyl-tRNA synthetase (PheRS)	FARS2	Q95363
41	Fibrinogen IV	Fibrinogen IV	Fibrinogen alpha chain/Fibrinogen beta chain	FGA	P02672
42	Fibrinogen S	Fibrinogen Type I-S	Fibrinogen gamma-B chain	FGB	P02676
					P12799
43	Fibronectin	Fibronectin	Fibronectin	FN1	P02751
44	GAD65	Glutamate decarboxylase 2	Glutamate decarboxylase 2	GAD2/GAD65	Q05329
45	GDF-15	Growth/differentiation factor 15	Growth/differentiation factor 15	GDF15	Q99988
46	GlnRS/QARS	Glutamyl-tRNA synthetase	Glutamine--tRNA ligase	QARS1	P47897
47	GlyRS/EJ	Glycyl-tRNA synthetase	Glycine--tRNA ligase	GARS1	P41250
48	GP2	Pancreatic secretory granule membrane major glycoprotein GP2	Pancreatic secretory granule membrane major glycoprotein GP2	GP2	P55259
49	GP210	Nuclear pore membrane glycoprotein 210	Nuclear pore membrane glycoprotein 210	NUP210/gp210	Q8TEM1
50	Hemocyanin	Hemocyanin	Hemocyanin		
51	Heparan Sulphate	Heparan Sulphate	Basement membrane-specific heparan sulfate proteoglycan core protein	HSPG2	P98160

52	Histone H2A	Histone H2A	Histone H2A type 1	H2AC11	P0C0S8
	Histone H2B	Histone H2B	Histone H2B	Histone H2B type 1-C/E/F/G/I	P62807
53	Histone H3	Histone H3	Histone H3.1	H3C1	P68431
	Histone H4	Histone H4	Histone H4	H4C1	P62805
54	HMGR	HMG-CoA reductase	3-hydroxy-3-methylglutaryl-coenzyme A	HMGR	P04035
55	IA2	Insulinoma-associated protein (IA-2)	islet cell antigen 512 (ICA 512); Receptor-type tyrosine-protein phosphatase-like N	PTPRN	Q16849
56	IARS/OJ/IleRS	Isoleucyl-tRNA synthetase	Isoleucine--tRNA ligase, cytoplasmic	IARS1	P41252
57	IFN-A1	Interferon alpha-1/13	Interferon alpha-1/13	IFNA1	P01562
58	IFN-B	Interferon beta	Interferon beta	IFNB1	P01574
59	Intrinsic factor (IF)	Gastric intrinsic factor	Cobalamin binding intrinsic factor	CBLIF	P27352
60	Jo-1	Histidyl-tRNA Synthetase (Jo-1)	Histidine--tRNA ligase, cytoplasmic	HARS1	P12081
61	KU (P70/P80)	Human KU protein, composed of P70 and P80 subunits	XRCC5 and XRCC6	XRCC5	P12956
				XRCC6	P13010
62	La/SSB	Lupus La protein homolog	Lupus La protein homolog	SSB	P10881
63	Laminin	Laminin	Laminin subunit gamma-1	LAMC1	P11047
64	LC1 (liver cytosol)	LC1 (liver cytosol)	Formimidoyltransferase-cyclodeaminase	FTCD	Q95954
65	M2	Mitochondrial 2-oxo acid dehydrogenase multienzyme	PDC-E2, OGDC-E2, and BCOADC-E2	DLAT	P10515
66	MDA5	melanoma differentiation-associated protein 5 (MDA5; CADM 140)	Interferon-induced helicase C domain-containing protein 1	IFIH1	Q9BYX4
67	Mi-2	nucleosome remodeling deacetylase (NuRD) complex, subunit	Chromodomain-helicase-DNA-binding protein	CHD3	Q12873
				CHD4	Q14839
68	MJ/NXP-2/MORC3	MORC family CW-type zinc finger protein 3	Nuclear matrix protein 2	MORC3	Q14149
69	MPO	Myeloperoxidase	Myeloperoxidase	MPO	P05164
70	Muscarinic receptor	Muscarinic receptor	Muscarinic acetylcholine receptor M5	CHRM5	P08912
71	MuSK	Muscle-specific tyrosine-protein kinase receptor	Muscle, skeletal receptor tyrosine-protein kinase	MUSK	Q15146
72	Myelin basic protein	Myelin basic protein (MBP)	Myelin basic protein	MBP	P02686
73	Myosin	Myosin	Myosin Calcium activated from bovine muscle	MYH	
74	Myosin heavy & light chain	Myosin heavy and light chain	Myosin Heavy and Light Chain from human heart	MYH1	P12882
					P05976
					P08590
					P10916
75	Neuropilin-1	Neuropilin-1	Neuropilin-1	NRP1	Q14786
76	NT5C1A	Cytosolic 5'-nucleotidase 1A	Cytosolic 5'-nucleotidase 1A	NT5C1A	Q98X13
77	Nucleolin	Nucleolin	Nucleolin	NCL	P19338
78	Nucleosome antigen	Nucleosome antigen	Nucleosome (Native. Purified from bovine tissues)		
79	Nup 62	Nucleoporin Nuclear pore glycoprotein p62	Nucleoporin Nuclear pore glycoprotein p62	Nup62	P37198
80	PCNA	Proliferating cell nuclear antigen	Proliferating cell nuclear antigen	PCNA	P12004
81	Phosphatidyl-serine	1,2-Diacyl-sn-glycero-3-phospho-L-	1,2-Diacyl-sn-glycero-3-phospho-L-serine	DGPS	
82	PL-12	Alanyl-tRNA Synthetase (PL-12)	Alanyl-tRNA Synthetase (PL-12)	AARS1	P49588
83	PL-7	Threonyl-tRNA Synthetase (PL-7)	Threonyl-tRNA Synthetase (PL-7)	TARS1	P26639
84	PLP1	Myelin proteolipid protein	Lipophilin	PLP1	P60201
85	PM/Scl 100	Exosome component 10	Exosome component 10	EXOSC10	Q01780
86	PM/Scl-75	Exosome complex component	Exosome complex component RRP45	EXOSC9	Q06265
87	PR3	Proteinase 3/Myeloblastin	Proteinase 3/Myeloblastin	PRTN3	P24158
88	Proteoglycan	Proteoglycan 4	Proteoglycan 4	PRG4	F1MDS0
89	Prothrombin protein	Prothrombin	Prothrombin	F2	P00734
90	Ribo Phosphoprotein P0	Ribo Phosphoprotein P0	Ribosomal Phosphoprotein P0	RPLP0	P05388
91	Ribo Phosphoprotein P1	Ribo Phosphoprotein P1	Ribosomal Phosphoprotein P1	RPLP1	P05386
92	Ribo Phosphoprotein P2	Ribo Phosphoprotein P2	Ribosomal Phosphoprotein P2	RPLP2	P05387
93	Ro/SSA (52 Kda)	E3 ubiquitin-protein ligase TRIM21	E3 ubiquitin-protein ligase TRIM21	TRIM21	P19474
94	Ro/SSA (60 Kda)	RNA-binding protein RO60	RNA-binding protein RO60	RO60	P10155
95	S100	S100 Calcium Binding Protein B	S100 Calcium Binding Protein B	S100B	P04271
96	SAE1	SUMO-activating enzyme subunit 1	Ubiquitin-like 1-activating enzyme E1A	SAE1	Q9UBF0
97	SAE2/UBA2	SUMO-activating enzyme subunit 2	Ubiquitin-like modifier-activating enzyme 2	UBA2	Q9UBT2
98	Scl-70/Topoisomerase I	DNA Topoisomerase I (Scl-70; full	DNA Topoisomerase I (Scl-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	SRP14	Signal recognition particle 14 kDa protein	18 kDa Alu RNA-binding protein	SRP14	P37108
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	TGF-B1	Transforming growth factor beta-1 proprotein,	Transforming growth factor beta-1 proprotein,	TGFB1; TGFB	P01137
107	Thyroglobulin	Native Thyroglobulin Isolated from human thyroid glands	Native Thyroglobulin Isolated from human thyroid glands	TG	P01266
108	TIF1 gama	E3 ubiquitin-protein ligase TRIM33	E3 ubiquitin-protein ligase TRIM33	TRIM33; KIAA1113, RFG7, TIF1G	Q9UPN9
109	TNF-a	Tumor necrosis factor	Tumor necrosis factor	TNF TNFA, TNFSF2	P01375
110	TPO	Thyroid peroxidase; TPO	Thyroid peroxidase; TPO	TPO	P07202
111	Troponin-I	Troponin I, cardiac muscle	Troponin I, cardiac muscle	TNNI3 TNNC1	P19429
112	TrpRS/WARS	Tryptophanyl-tRNA	Tryptophan--tRNA ligase, mitochondrial	WARS2	Q9UGM6
113	TTG	Tissue transglutaminase (Ttg)/Protein-glutamine gamma-glutamyltransferase 2 (TGase-2)	Tissue transglutaminase (Ttg)/Protein-glutamine gamma-glutamyltransferase 2 (TGase-2)	TGM2	P21980
114	TyrRS/YARS/HA	Tyrosyl-tRNA synthetase (TyrRS)	Tyrosine--tRNA ligase, mitochondrial	YARS2	Q9Y274
115	U1-snRNP 68/70	U1 small nuclear ribonucleoprotein 70 kD, U1 snRNP 70 kDa, U1-70K,	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, Sm protein B/B', Sm-B/B', SmB/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	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		