

The antigen comparison of PA001-V2 and PA001-V3							
Ag#	ID	Protein Name	Present in V2	Present in V3	Alternative name /Synonyms	Gene symbol	UniProt ID
1	Aggrecan	Aggrecan	Y	Y	cartilage-specific proteoglycan core protein (CSPCP) or chondroitin sulfate proteoglycan 1	ACAN	<a href="#">P16112</a>
2	AGTR	Type-1 angiotensin II receptor	Y	Y	Type-1 angiotensin II receptor	AGTR1	<a href="#">P30556</a>
3	Alpha Fodrin	Alpha Fodrin	Y	Y	Spectrin alpha chain, non-erythrocytic 1	SPTAN1	<a href="#">Q13813</a>
4	Alpha-actinine	Alpha-actinine	Y	Y	Alpha-actinin-1; F-actin cross-linking protein	ACTN1	<a href="#">P12814</a>
5	Amyloid	Amyloid	Y	Y	Amyloid-beta precursor; proteinAlzheimer disease amyloid protein; Protease nexin-II	APP	<a href="#">P05067</a>
6	AQP4	Aquaporin-4	Y	Y	Aquaporin-4	AQP4	<a href="#">P55087</a>
7	B2 glycoprotein 1	B2-glycoprotein 1 (B2GPI)	Y	Y	Apolipoprotein H (APO-H)	APOH	<a href="#">P02749</a>
8	B2-microglobulin	B2-microglobulin	Y	Y	Beta-2-microglobulin	B2M	<a href="#">P61769</a>
9	B-actin	B-actin	Y	Y	Actin, cytoplasmic 1	ACTB	<a href="#">P60709</a>
10	BPI	Bactericidal permeability-increasing protein	Y	Y	Bactericidal permeability-increasing protein	BPI	<a href="#">P17213</a>
11	Calprotectin	Calprotectin	Y	Y	Protein S100, calgranulin, leucocyte L1 complex, migration inhibitory factor-related protein, MRP8/14, S100 calcium-binding protein	S100A8	<a href="#">P05109</a>
						S100A9	<a href="#">P06702</a>
12	Cardiolipin	Cardiolipin	Y	Y	Diphosphatidylglycerol	Cardiolipin sodium salt from bovine	<a href="#">c0563</a>
13	CD4	T-cell surface glycoprotein CD4	Y	Y	T-cell surface glycoprotein CD4	CD4	<a href="#">P01730</a>
14	CD8	T-cell surface glycoprotein CD8 beta chain	Y	Y	T-cell surface glycoprotein CD8 beta chain	CD8B	<a href="#">P10966</a>
15	CENP-A	Centromere protein A	Y	Y	Histone H3-like centromeric protein A, Centromere protein A	CENPA	<a href="#">P49450</a>
16	CENP-B	Centromere protein B	Y	Y	Centromere protein B	CENPB	<a href="#">P07199</a>
17	Chondroitin Sulfate C	Chondroitin Sulfate C	Y	Y	Chondroitin Sulfate Proteoglycan (CSPG), sulfated glycosaminoglycan polysaccharide	Chondroitin sulfate sodium salt from	<a href="#">Q6UVK1</a>
18	COL17A1	Collagen alpha-1(XVII) chain	Y	Y	COL17A1; collagen, type XVII, alpha 1; BPAG2	COL17A1	<a href="#">Q9UMD9</a>
19	Collagen I	Collagen type I	Y	Y	Collagen alpha-1(I) chain	COL1A1	<a href="#">P02452</a>
20	Collagen II	Collagen Type II	Y	Y	Collagen alpha-1(II) chain	COL2A1	<a href="#">P02458</a>
21	Collagen III	Collagen type III	Y	Y	Collagen alpha-1(III) chain/COL3A1	COL3A1	<a href="#">P02461</a>
22	Collagen IV	Collagen type IV	Y	Y	Collagen alpha-1(IV) chain	COL4A1	<a href="#">P02462</a>
23	Collagen V	Collagen type V	Y	Y	Collagen alpha-1(V) chain	COL5A1	<a href="#">P20908</a>
24	Collagen VI	Collagen type VI	Y	Y	Collagen alpha-1(VI) chain	COL6A1	<a href="#">P12109</a>
25	complement C1q	complement C1q	Y	Y	Complement C1q subcomponent subunit A	C1QA	<a href="#">P02745</a>
26	complement C3/C3a	complement C3/C3a	Y	Y	Complement C3	C3	<a href="#">P01024</a>
27	Core Histone	Core Histone	Y	Y	H2A, H2B, H3 and H4		
28	CRP	C-reactive protein	Y	Y	C-reactive protein	CRP	<a href="#">P02741</a>
29	dsDNA	double strand DNA	Y	Y	double strand DNA		
30	EBNA1	Epstein-Barr nuclear antigen 1	Y	N	Epstein-Barr nuclear antigen 1	EBNA1	<a href="#">Q3KSS4</a>
31	Elastin	Elastin	Y	Y	Tropoelastin	ELN	<a href="#">P15502</a>
32	Factor B	Complement factor B	Y	Y	Complement factor B	CFB	<a href="#">P00751</a>
33	Factor D	Complement factor D	Y	Y	Complement factor D	CFD	<a href="#">P00746</a>
34	Factor H	Complement factor H	Y	Y	Complement factor H	CFH	<a href="#">P08603</a>
35	Factor I	Complement factor I	Y	Y	Complement factor I	CFI	<a href="#">P05156</a>
36	Factor P	Complement factor P	Y	Y	Complement factor P/Properdin	CFP	<a href="#">P27918</a>

The antigen comparison of PA001-V2 and PA001-V3							
Ag#	ID	Protein Name	Present in V2	Present in V3	Alternative name /Synonyms	Gene symbol	UniProt ID
37	Fibrinogen IV	Fibrinogen IV	Y	Y	Fibrinogen alpha chain/Fibrinogen beta chain	FGA	<a href="#">P02672</a>
38	Fibrinogen S	Fibrinogen Type I-S	Y	Y	Fibrinogen gamma-B chain	FGB	<a href="#">P02676</a>
							<a href="#">P12799</a>
39	Fibronectin	Fibronectin	Y	Y	Fibronectin	FN1	<a href="#">P02751</a>
40	GAD65	Glutamate decarboxylase 2	Y	Y	Glutamate decarboxylase 2	GAD2/GAD65	<a href="#">Q05329</a>
41	GBM	Glomerular Basement Membrane	Y	N	Glomerular Basement Membrane		
42	GDF-15	Growth/differentiation factor 15	Y	Y	Growth/differentiation factor 15	GDF15	<a href="#">Q99988</a>
43	GM-CSF	Granulocyte-macrophage colony-stimulating factor	Y	N	Granulocyte-macrophage colony-stimulating factor	CSF2/GM-CSF	<a href="#">P04141</a>
44	GP2	Pancreatic secretory granule membrane major glycoprotein GP2	Y	Y	Pancreatic secretory granule membrane major glycoprotein GP2	GP2	<a href="#">P55259</a>
45	GP210	Nuclear pore membrane glycoprotein 210	Y	Y	Nuclear pore membrane glycoprotein 210	NUP210/gp210	<a href="#">Q8TEM1</a>
46	Hemocyanin	Hemocyanin	Y	Y	Hemocyanin		
47	Heparan Sulphate	Heparan Sulphate	Y	Y	Basement membrane-specific heparan sulfate proteoglycan core protein	HSPG2	<a href="#">P98160</a>
48	Heparin	Heparin	Y	N	Heparin		
49	Histone H1	Histone H1	Y	N	Histone H1 human	H1-0	<a href="#">P07305</a>
50	Histone H2A	Histone H2A	Y	Y	Histone H2A type 1	H2AC11	<a href="#">P0C058</a>
51	Histone H2B	Histone H2B	Y	Y	Histone H2B	Histone H2B type 1-C/E/F/G/I	<a href="#">P62807</a>
52	Histone H3	Histone H3	Y	Y	Histone H3.1	H3C1	<a href="#">P68431</a>
53	Histone H4	Histone H4	Y	Y	Histone H4	H4C1	<a href="#">P62805</a>
54	IA2	Insulinoma-associated protein (IA-2)	Y	Y	islet cell antigen 512 (ICA 512); Receptor-type tyrosine-protein phosphatase-like N	PTPRN	<a href="#">Q16849</a>
55	IFN-A1	Interferon alpha-1/13	Y	Y	Interferon alpha-1/13	IFNA1	<a href="#">P01562</a>
56	IFN-B	Interferon beta	Y	Y	Interferon beta	IFNB1	<a href="#">P01574</a>
57	IFN-E	Interferon epsilon	Y	N	Interferon epsilon	IFNE	<a href="#">Q86WN2</a>
58	IFN-L2	Interferon lambda-2	Y	N	Interferon lambda-2	IFNL2	<a href="#">Q8IZJ0</a>
59	IFN-r	Interferon gamma	Y	N	Interferon gamma	IFNG	<a href="#">P01579</a>
60	IL-12	Interleukin-12 subunit alpha/beta	Y	N	Interleukin-12 subunit alpha/beta	IL12A	<a href="#">P29459</a>
						IL12B	<a href="#">P29460</a>
61	IL-15	Interleukin-15	Y	N	Interleukin-15	IL15	<a href="#">P40933</a>
62	Insulin	Insulin	Y	N	Insulin	INS	<a href="#">P01308</a>
63	Jo-1	Histidyl-tRNA Synthetase (Jo-1)	Y	Y	Histidine--tRNA ligase, cytoplasmic	HARS1	<a href="#">P12081</a>
64	KU (P70/P80)	Human KU protein, composed of P70 and P80 subunits	Y	Y	XRCC5 and XRCC6	XRCC6	<a href="#">P12956</a>
						XRCC5	<a href="#">P13010</a>
65	La/SSB	Lupus La protein homolog	Y	Y	Lupus La protein homolog	SSB	<a href="#">P10881</a>
66	Laminin	Laminin	Y	Y	Laminin subunit gamma-1	LAMC1	<a href="#">P11047</a>
67	LC1 (liver cytosol )	LC1 (liver cytosol )	Y	Y	Formimidoyltransferase-cyclodeaminase	FTCD	<a href="#">Q95954</a>
68	LPS	Lipopolysaccharide	Y	N	Lipopolysaccharide		
69	M2	Mitochondrial 2-oxo acid dehydrogenase multienzyme complexes	Y	Y	PDC-E2, OGDC-E2, and BCOADC-E2	DLAT	<a href="#">P10515</a>
70	MDA5	melanoma differentiation-associated protein 5 (MDA5; CADM 140)	Y	Y	Interferon-induced helicase C domain-containing protein 1	IFIH1	<a href="#">Q9BYX4</a>

The antigen comparison of PA001-V2 and PA001-V3							
Ag#	ID	Protein Name	Present in V2	Present in V3	Alternative name /Synonyms	Gene symbol	UniProt ID
71	Mi-2	nucleosome remodeling deacetylase (NuRD) complex, subunit	Y	Y	Chromodomain-helicase-DNA-binding protein	CHD3	<a href="#">Q12873</a>
						CHD4	<a href="#">Q14839</a>
72	Mitochondrial antigen	Mitochondrial antigen	Y	N	Pyruvate dehydrogenase complex, mitochondrial	PDHA	<a href="#">Q9N1X8</a>
						PDHB	<a href="#">P11966</a>
73	MPO	Myeloperoxidase	Y	Y	Myeloperoxidase	MPO	<a href="#">P05164</a>
74	Muscarinic receptor	Muscarinic receptor	Y	Y	Muscarinic acetylcholine receptor M5	CHRM5	<a href="#">P08912</a>
75	Myelin basic protein (MBP)	Myelin basic protein (MBP)	Y	Y	Myelin basic protein	MBP	<a href="#">P02686</a>
76	Myosin	Myosin	Y	Y	Myosin Calcium activated from bovine muscle	MYH	
77	Myosin heavy & light chain	Myosin heavy and light chain	Y	Y	Myosin Heavy and Light Chain from human heart	MYH1	<a href="#">P12882</a>
							<a href="#">P05976</a>
							<a href="#">P08590</a>
							<a href="#">P10916</a>
78	Neuropilin-1	Neuropilin-1	Y	Y	Neuropilin-1	NRP1	<a href="#">Q14786</a>
79	Nucleolin	Nucleolin	Y	Y	Nucleolin	NCL	<a href="#">P19338</a>
80	Nucleosome antigen	Nucleosome antigen	Y	Y	Nucleosome (Native. Purified from bovine tissues)		
81	Nup 62	Nucleoporin Nuclear pore glycoprotein p62	Y	Y	Nucleoporin Nuclear pore glycoprotein p62	Nup62	<a href="#">P37198</a>
82	PCNA	Proliferating cell nuclear antigen	Y	Y	Proliferating cell nuclear antigen	PCNA	<a href="#">P12004</a>
83	Phosphatidyl-serine	1,2-Diacyl-sn-glycero-3-phospho-L-serine	Y	Y	1,2-Diacyl-sn-glycero-3-phospho-L-serine	DGPS	
84	PL-12	Alanyl-tRNA Synthetase (PL-12)	Y	Y	Alanyl-tRNA Synthetase (PL-12)	AARS1	<a href="#">P49588</a>
85	PL-7	Threonyl-tRNA Synthetase (PL-7)	Y	Y	Threonyl-tRNA Synthetase (PL-7)	TARS1	<a href="#">P26639</a>
86	PM/Sc1 100	Exosome component 10	Y	Y	Exosome component 10	EXOSC10	<a href="#">Q01780</a>
87	PM/Sc1-75	Exosome complex component RRP45	Y	Y	Exosome complex component RRP45	EXOSC9	<a href="#">Q06265</a>
88	PR3	Proteinase 3/Myeloblastin	Y	Y	Proteinase 3/Myeloblastin	PRTN3	<a href="#">P24158</a>
89	Proteoglycan	Proteoglycan 4	Y	Y	Proteoglycan 4	PRG4	<a href="#">F1MDS0</a>
90	Prothrombin protein	Prothrombin	Y	Y	Prothrombin	F2	<a href="#">P00734</a>
91	Rheumatoid factor (RF)	Rheumatoid factor (RF)	Y	N	Human IgG Fc fragment		
92	Ribo Phosphoprotein P0	Ribo Phosphoprotein P0	Y	Y	Ribosomal Phosphoprotein P0	RPLP0	<a href="#">P05388</a>
93	Ribo Phosphoprotein P1	Ribo Phosphoprotein P1	Y	Y	Ribosomal Phosphoprotein P1	RPLP1	<a href="#">P05386</a>
94	Ribo Phosphoprotein P2	Ribo Phosphoprotein P2	Y	Y	Ribosomal Phosphoprotein P2	RPLP2	<a href="#">P05387</a>
95	Ro/SSA (52 Kda)	E3 ubiquitin-protein ligase TRIM21	Y	Y	E3 ubiquitin-protein ligase TRIM21	TRIM21	<a href="#">P19474</a>
96	Ro/SSA (60 Kda)	RNA-binding protein RO60	Y	Y	RNA-binding protein RO60	RO60	<a href="#">P10155</a>
97	S100	S100 Calcium Binding Protein B	Y	Y	S100 Calcium Binding Protein B	S100B	<a href="#">P04271</a>
98	Sc1-70/Topoisomerase I	DNA Topoisomerase I (Sc1-70; full length)	Y	Y	DNA Topoisomerase I (Sc1-70; full length)	TOP1	<a href="#">P11387</a>
99	Sm	Small nuclear ribonucleoprotein Sm from Bovine tissue	Y	Y	Small nuclear ribonucleoprotein Sm from Bovine tissue	Sm	
100	Sm/RNP	RNP/Sm ribonucleoprotein complex native protein from bovine tissue	Y	Y	RNP/Sm ribonucleoprotein complex native protein from bovine tissue	RNP/Sm	
101	SmD	Small nuclear ribonucleoprotein Sm, recombinant D1,D2,D3 mixture	Y	Y	Small nuclear ribonucleoprotein Sm, recombinant D1,D2,D3 mixture	SNRPD1	<a href="#">P62314</a>
102	SP100	Nuclear dot-associated Sp100 protein; Speckled 100 kDa	Y	Y	Nuclear dot-associated Sp100 protein; Speckled 100 kDa	SP100	<a href="#">P23497</a>

The antigen comparison of PA001-V2 and PA001-V3							
Ag#	ID	Protein Name	Present in V2	Present in V3	Alternative name /Synonyms	Gene symbol	UniProt ID
103	Sphingomyelin	N-Acyl-D-sphingosine-1-phosphocholine, from egg yolk	Y	N	N-Acyl-D-sphingosine-1-phosphocholine, from egg yolk		
104	SRP54	Signal recognition particle 54 kDa protein, SRP54	Y	Y	Signal recognition particle 54 kDa protein, SRP54	SRP54	<a href="#">P61011</a>
105	ssDNA	Single strand DNA	Y	Y	Single strand DNA		
106	ssRNA	Single strand RNA	Y	N	Single strand RNA		
107	TGF-B1	Transforming growth factor beta-1 proprotein,	Y	Y	Transforming growth factor beta-1 proprotein,	TGFB1; TGFB	<a href="#">P01137</a>
108	Thyroglobulin	Native Thyroglobulin Isolated from human thyroid glands	Y	Y	Native Thyroglobulin Isolated from human thyroid glands	TG	<a href="#">P01266</a>
109	TIF1 gama	E3 ubiquitin-protein ligase TRIM33	Y	Y	E3 ubiquitin-protein ligase TRIM33	TRIM33; KIAA1113, RFG7, TIF1G	<a href="#">Q9UPN9</a>
110	TNF-a	Tumor necrosis factor	Y	Y	Tumor necrosis factor	TNF TNFA, TNFSF2	<a href="#">P01375</a>
111	TPO	Thyroid peroxidase; TPO	Y	Y	Thyroid peroxidase; TPO	TPO	<a href="#">P07202</a>
112	Troponin-I	Troponin I, cardiac muscle	Y	Y	Troponin I, cardiac muscle	TNNI3 TNNC1	<a href="#">P19429</a>
113	TTG	Tissue transglutaminase (Ttg)/Protein-glutamine gamma-glutamyltransferase	Y	Y	Tissue transglutaminase (Ttg)/Protein-glutamine gamma-glutamyltransferase 2 (TGase-2)	TGM2	<a href="#">P21980</a>
114	Tweek	Tumor necrosis factor ligand superfamily, member 12, TNFSF12, YNF-	Y	N	Tumor necrosis factor ligand superfamily, member 12, TNFSF12, YNF-related weak inducer of apoptosis, Apo-3	TNFSF12	<a href="#">Q43508</a>
115	U1-snRNP 68/70	U1 small nuclear ribonucleoprotein 70 kD, U1 snRNP 70 kDa, U1-70K, snRNP70	Y	Y	U1 small nuclear ribonucleoprotein 70 kD, U1 snRNP 70 kDa, U1-70K, snRNP70	SNRNP70 RNP1U1Z, RPU1, SNRP70,	<a href="#">P08621</a>
116	U1-snRNP A	U1 small nuclear ribonucleoprotein A, U1 snRNP A, U1-A, U1A	Y	Y	U1 small nuclear ribonucleoprotein A, U1 snRNP A, U1-A, U1A	SNRPA	<a href="#">P09012</a>
117	U1-snRNP B/B'	Small nuclear ribonucleoprotein-associated proteins B and B', snRNP-B,	Y	Y	Small nuclear ribonucleoprotein-associated proteins B and B', snRNP-B, Sm protein B/B', Sm-B/B', SmB/B'	SNRPB, COD, SNRPB1	<a href="#">P14678</a>
118	U1-snRNP C	U1 small nuclear ribonucleoprotein C, U1 snRNP C, U1-C, U1C	Y	Y	U1 small nuclear ribonucleoprotein C, U1 snRNP C, U1-C, U1C	SNRPC	<a href="#">P09234</a>
119	Vimentin	Vimentin	Y	Y	Vimentin	VIM	<a href="#">P08670</a>
120	Vitronectin	Vitronectin	Y	Y	Vitronectin, VN, S-protein, Serum-spreading factor, V75	VTN	<a href="#">P04004</a>
121	SRP14	Signal recognition particle 14 kDa protein	N	Y	18 kDa Alu RNA-binding protein	SRP14	<a href="#">P37108</a>
122	ARS	Aminoacyl tRNA synthetase	N	Y	Aminoacyl-tRNA Synthetase from Escherichia coli	alaS	<a href="#">P00957</a>
123	AsnRS/KS	Asparaginyl-tRNA Synthetase	N	Y	Human Asparaginyl-tRNA Synthetase (KS)	NARS1	<a href="#">Q43776</a>
124	Intrinsic factor (IF)	Gastric intrinsic factor	N	Y	Cobalamin binding intrinsic factor	CBIF	<a href="#">P27352</a>
125	PLP1	Myelin proteolipid protein	N	Y	Lipophilin	PLP1	<a href="#">P60201</a>
126	CCP	Cyclic citrullinated peptide	N	Y			
127	MuSK	Muscle-specific tyrosine-protein kinase receptor	N	Y	Muscle, skeletal receptor tyrosine-protein kinase	MUSK	<a href="#">Q15146</a>
128	DFS70	Dense fine speckles 70 kDa protein	N	Y	Transcriptional coactivator p75/p52	PSIP1	<a href="#">Q75457</a>
129	FARS2	Phenylalanine--tRNA ligase, mitochondrial	N	Y	Phenylalanyl-tRNA synthetase (PheRS)	FARS2	<a href="#">Q95363</a>
130	GlnRS/QARS	Glutaminyl-tRNA synthetase	N	Y	Glutamine--tRNA ligase	QARS1	<a href="#">P47897</a>
131	GlyRS/EJ	Glycyl-tRNA synthetase	N	Y	Glycine--tRNA ligase	GARS1	<a href="#">P41250</a>
132	IARS/OJ/IleRS	Isoleucyl-tRNA synthetase	N	Y	Isoleucine--tRNA ligase, cytoplasmic	IARS1	<a href="#">P41252</a>
133	MJ/NXP-2/MORC3	MORC family CW-type zinc finger protein 3	N	Y	Nuclear matrix protein 2	MORC3	<a href="#">Q14149</a>
134	NTSC1A	Cytosolic 5'-nucleotidase 1A	N	Y	Cytosolic 5'-nucleotidase IA	NTSC1A	<a href="#">Q98X13</a>
135	SAE1	SUMO-activating enzyme subunit 1	N	Y	Ubiquitin-like 1-activating enzyme E1A	SAE1	<a href="#">Q9UBE0</a>
136	SAE2/UBA2	SUMO-activating enzyme subunit 2	N	Y	Ubiquitin-like modifier-activating enzyme 2	UBA2	<a href="#">Q9UBT2</a>
137	HMGCR	HMG-CoA reductase	N	Y	3-hydroxy-3-methylglutaryl-coenzyme A reductase	HMGCR	<a href="#">P04035</a>
138	TrpRS/WARS	Tryptophanyl-tRNA synthetase (TrpRS)	N	Y	Tryptophan--tRNA ligase, mitochondrial	WARS2	<a href="#">Q9UGM6</a>
139	TyrRS/YARS/HA	Tyrosyl-tRNA synthetase (TyrRS)	N	Y	Tyrosine--tRNA ligase, mitochondrial	YARS2	<a href="#">Q9Y2Z4</a>

The antigen comparison of PA001-V2 and PA001-V3							
Ag#	ID	Protein Name	Present in V2	Present in V3	Alternative name /Synonyms	Gene symbol	UniProt ID
140	Ig control 1	Ig control 1	y	y	Human and mouse Ig control mix		
141	Ig control 2	Ig control 2	y	y	Human and mouse Ig control mix		
142	Ig control 3	Ig control 3	y	y	Human and mouse Ig control mix		
143	Ig control 4	Ig control 4	y	y	Human and mouse Ig control mix		
144	anti-Ig control 1	anti-Ig control 1	y	y	Anti-human and anti-mouse Ig control mix		
145	anti-Ig control 2	anti-Ig control 2	y	y	Anti-human and anti-mouse Ig control mix		
146	anti-Ig control 3	anti-Ig control 3	y	y	Anti-human and anti-mouse Ig control mix		
147	anti-Ig control 4	anti-Ig control 4	y	y	Anti-human and anti-mouse Ig control mix		