

OmicsArray™ General Autoimmune Antigen Array, 82-plex				Catalog No. PA018
Antigen#	Antigene ID	Antigen Name	UniProt ID	Symbol
1	Actin Bovine muscle	Actin, alpha skeletal muscle	P68138	ACTA1
2	Actin bovine cardiac	Actin, alpha cardiac muscle 1	Q3ZC07	ACTC1
3	Alpha KGDH	Alpha-ketoglutarate dehydrogenase	Q02218	OGDH
4	BCOADC E2	Branched chain 2-oxo-acid dehydrogenase complex component E2	P11182	DBT
5	B2GPI	Beta-2-glycoprotein 1 (non recombinant Human protein)	P02749	APOH
6	Histone H4	Histone H4	P62805	H4C1
7	Histone H2A	Histone H2A type 2-A	Q6FI13	H2AC18; H2AC19
8	Histone H2b/F2b	Histone H2B type 2-F	Q5QNW6	H2BC18
9	Histone H3	Histone H3 Bovine non recombinant	P68432	H3.1
10	BPI	Bactericidal permeability-increasing protein	P17213	BPI
11	Myosin-6	Myosin-6	P13533	MYH6
12	Myosin Calcium activated from bovine muscle	Unconventional myosin-VI	E1BPK6	MYO6
13	Myosin bovine cardiac muscle	Myosin-6	F1N2G0	MYH6
14	CENP-B	Major centromere autoantigen B	P07199	CENPB
15	Collagen I	Collagen alpha-1(I) chain	P02452	COL1A1
16	Collagen III	Collagen alpha-1(III) chain	P02461	COL3A1
17	Collagen V	Collagen alpha-1(V) chain	P20908	COL5A1
18	Desmin	Desmin	P17661	DES
19	SCL70 (Topo I)	DNA topoisomerase 1	P11387	TOP1
20	Enolase	Alpha-enolase	P06733	ENO1
21	Alpha Enolase (citrullinated)	Alpha-enolase	P06733	ENO1
22	Glycyl tRNA Synthetase (EJ)	Glycine--tRNA ligase	P41250	GARS1
23	Histone H4	Histone H4 (1-103 aa)	P62805	H4C1
24	HSP27	Heat shock protein beta-1	P04792	HSPB1
25	HSP47	Serpin H1	P50454	SERPINH1
26	Insulin	Insulin	P01308	INS
27	SSB (La)	Lupus La protein	P05455	SSB
28	LC1	Formimidoyltransferase-cyclodeaminase	Q95954	FTCD
29	M2	Mitochondrial 2-oxo acid dehydrogenase multienzyme complexes	P10515 , P36957 , P11182	DLST, DLAT, DBT
30	MPO	Myeloperoxidase	P05164	MPO
31	Mucarinic Rceptor 3	Muscarinic acetylcholine receptor M3	P20309	CHRM3
32	Nucleolin	Nucleolin	P19338	NCL
33	Nucleosome	Native nucleosome antigen from bovine		N/A
34	PDC-E2	Dihydrolipoylysine-residue acetyltransferase component of pyruvate dehydrogenase complex, mitochondrial	P10515	DLAT
35	PDHA1	Pyruvate dehydrogenase E1 component subunit alpha	P08559	PDHA1
36	PM/ScI-75	Exosome complex component RRP45	Q06265	EXOSC9
37	SSA/Ro-60	RNA-binding protein RO60	P10155	RO60
38	Sm (Smith) antigen	Smith antigen (Sm Ag) nonrecombinant from Bovine	Q3ZC10	SNRPD1

39	SmD	Small nuclear ribonucleoprotein Sm D1	P62314	SNRPD1
40	SmD3	Small nuclear ribonucleoprotein Sm D3	P62318	SNRPD3
41	snRNP C	U1 small nuclear ribonucleoprotein C	P09234	SNRPC
42	Total histone	Histone from bovine (calf) thymus		N/A
43	CENP-A	Histone H3-like centromeric protein A	P49450	CENPA
44	Collagen type IV	Collagen alpha-1(IV) chain	P02462	COL4A1
45	dsDNA	double strand DNA		N/A
46	Jo-1	Histidine--tRNA ligase, cytoplasmic	B4E1C5	HARS1
47	M5 Muscarinic receptor	Muscarinic acetylcholine receptor M5	P08912	CHRM5
48	PM/ScI-100	Exosome complex component 10	Q01780	EXOSC10
49	PR3	Myeloblastin	P24158	PRTN3
50	SSA/Ro-52	E3 ubiquitin-protein ligase TRIM21	P19474	TRIM21
51	SmD1	Small nuclear ribonucleoprotein Sm D1	P62314	SNRPD1
52	SmD2	Small nuclear ribonucleoprotein Sm D2	P62316	SNRPD2
53	Ig control 1	Control		
54	Ig control 2	Control		
55	Ig control 3	Control		
56	Ig control 4	Control		
57	anti-Ig control 1	Control		
58	anti-Ig control 2	Control		
59	anti-Ig control 3	Control		
60	anti-Ig control 4	Control		