

OmicsArray™ Autoimmune Myocarditis and Myositis (AMM) Antigen Array, 120-plex antigen list - Cat.#PA016

Antigen Number	Antigen ID	Antigen Name	UniProt ID
1	20S P	Human 20S Proteasome Protein	
2	ACO2	Mitochondrial aconitase	Q99798
3	ACTA1	Alpha actin (Skeletal)	P68133
4	ACTB	Beta actin	P60709
5	ACTC1	Alpha actin (cardiac), Actin Alpha 1, Cardiac Muscle	P68032
6	ACTN1	Alpha-actinin	P12814
7	ADRA1A	Alpha-1A adrenergic receptor	P35348
8	ADRA2A	Alpha-2A adrenergic receptor	P08913
9	ADRB1	Beta-1 adrenergic receptor	P08588
10	ADRB2	Beta-2 adrenergic receptor	P07550
11	ANT2/SLC25A5	Adenine Nucleotide Translocator 2	P05141
12	ARS	Aminoacyl-tRNA Synthetase	P55000
13	AsnRS/KS	Asparaginyl-tRNA Synthetase (KS)	O43776
14	AT1R	Angiotensin II type 1-receptor	P30556
15	ATP1A1	ATPase, Na+/K+ transporting, alpha 1	P05023
16	ATP1B1	ATPase Na+/K+ Transporting Subunit Beta 1	P05026
17	ATP1B4	ATPase Na+/K+ Transporting Subunit Beta 4	Q9UN42
18	ATP2C1	ATPase secretory pathway Ca2+ transporting 1	P98194
19	B2GPI	Beta 2 Glycoprotein I	P02749
20	B2M	Beta-2 microglobulin	P61769
21	BCKDHA	Branched chain alpha-ketoacid dehydrogenase dihydrolipoyl transacylase	P12694
22	BCOADC-E2	branched chain 2-oxo-acid dehydrogenase complex	P11182
23	BHM	Bovine Heart Mitochondria	
24	CALR	Calreticulin	P27797
25	Cardiac Myosin	Myosin (bovine cardiac muscle)	
26	Cardiolipin	Cardiolipin (CL)	
27	CD4	CD4	P01730
28	CD8	CD8	P01732 P10966
29	CDH1	E-Cadherin/Cadherin-1	P12830
30	CDH2	Cadherin 2/N-cadherin	P19022
31	CDH8	Cadherin 8 (CDH8)	P55286
32	CDH9	Cadherin 9 (CDH9)	Q9ULB4
33	CHRM2	Muscarinic acetylcholine receptor M2	P08172
34	CHRM3	Cholinergic Receptor Muscarinic 3 (CHRM3)	P20309
35	cMyBP-C	Cardiac Myosin Binding Protein-C	Q14896
36	cTnI	Cardiac Troponin I	P19429
37	CTNNA1	alpha catenin	P35221
38	CTNNB1	beta catenin	P35222
39	CXADR/CAR	coxsackie and adenovirus receptor	P78310
40	CXCR3	CXC Motif Chemokine Receptor 3	P49682
41	DES	Desmin	P17661

42	DLG4	Disks Large Homolog 4/DLG4/PSD95	P78352
43	DSC1	Desmocollin-1	Q08554
44	DSC2	Desmocollin-2	Q02487
45	DSC3	Desmocollin-3	Q14574
46	dsDNA	dsDNA	
47	DSG1	Desmoglein 1 (DSG1)	Q02413
48	DSG2	Desmoglein 2 (DSG2)	Q14126
49	DSG3	Desmoglein 3 (DSG3)	P32926
50	DSG4	Desmoglein 4 (DSG4)	Q86SJ6
51	EDNRA	Endothelin receptor type A	P25101
52	EDNRB	Endothelin receptor type B	P24530
53	FARS2	Phenylalanyl-tRNA Synthetase 2	O95363
54	GAPDH	Glyceraldehyde-3-phosphate dehydrogenase	P00357
55	GJA1	Connexin 43/GJA1	P17302
56	GlnRS/QARS	Glutaminyl-tRNA synthetase (QARS)	P47897
57	GlyRS/EJ	Glycyl-tRNA Synthetase (EJ)	P41250
58	HMGCR	3-hydroxy-3-methylglutaryl-coenzyme A reductase	P04035
59	HSP60/HSPD1	Heat shock protein-60/chaperonins (Cpn)/HSPD1	P10809
60	HSP70/HSP1A1	Heat shock protein-70/Hu HSP70/HSPA1A Protein	P09446
61	HTR4	5-hydroxytryptamine receptor 4	Q13639
62	IARS/OJ/IleRS	Isoleucyl tRNA Synthetase (IARS/OJ)	Q9NSE4
63	IFN-a	Interferon-alpha	P01563
64	IFN-b	Interferon-beta	P01574
65	ITGB1	Integrin Subunit Beta 1	P05556
66	ITGB4	Integrin, beta 4 (ITGB4) /CD104	P16144
67	Jo-1	Histidyl-tRNA synthetase (Jo-1)	P12081
68	JUP	Junction Plakoglobin	P14923
69	KCNIP2	Kv channel-interacting protein2.6	Q9NS61
70	KCNJ5/Kir3.4/GIRK4	potassium channel subunit 3.4 (Kir3.4)	P48544
71	KCNK2/TREK1/TWIK1	Potassium channel subfamily K member 2	O95069
72	KCNQ1	Potassium voltage-gated channel subfamily KQT member 1	P51787
73	KRT24	Keratin 24	Q2M2I5
74	Ku (p70/p80)	Ku (p70/p80)	P13010
75	La/SSB	La	P05455
76	Laminin	Laminin subunit beta-1	P07942
77	M1MR/CHRM	M1 Muscarinic receptor (human)	P11229
78	M2	M2 (Mitochondrial 2-oxo acid dehydrogenase multienzyme complexes): PDC-E2, OGDC-E2, and BCOADC-E2	P10515
			P36957
			P11182
79	M3MR	M3 Muscarinic receptor (human) membrane preparation	P20309
80	M4MR	M4 Muscarinic receptor (human) membrane preparation	P08173
81	M5MR	M5 Muscarinic receptor, Human	P08912
82	MuSK	Muscle-specific tyrosine-protein kinase receptor (MuSK)	Q15146
83	MDA-5/IFIH1	melanoma differentiation-associated protein 5	Q9BYX4
84	Mi-2	Nucleosome remodeling-deacetylase (NuRD) complex	Q14839

85	MJ/NXP-2/MORC3	nuclear matrix protein 2 (NXP-2), MORC3 (MORC family CW-type zinc finger 3)	Q14149
86	MPO	Myeloperoxidase (MPO)	P05164
87	MYH1	Myosin Heavy Chain (from human heart)	P12882
88	MYH6	Myosin-6	P13533
89	MYH9	Human Myosin-9 (MYH9), aa2-241	P35579
90	MYL1	Myosin Light Chain from human heart	P05976
91	Myosin	Myosin Calcium activated from bovine muscle (Myosin heavy chain and light chain)	
92	NT5C1A	Cytosolic 5'-nucleotidase 1A (NT5C1A)	Q9BXI3
93	OGDC-E2	2-oxoglutarate dehydrogenase complex	P36957
94	PAR1	Proteinase-activated receptor 1	P25116
95	PAR2	Proteinase-activated receptor 2	P55085
96	PDC-E2	Pyruvate dehydrogenase (PDH) complex, Mitochondria	P10515
97	PKP2	Plakophilin 2	Q99959
98	PL-12	Alanyl-tRNA Synthetase (PL-12)	Q5JTZ9
99	PL-7	Threonyl-tRNA Synthetase (PL-7)	Q9BW92
100	PM/Scl100	Polymyositis/scleroderma (PM/Scl) 100	Q01780
101	PM/Scl75	Polymyositis/scleroderma (PM/Scl) 75	Q06265
102	PYGB	Brain glycogen phosphorylase	P11216
103	RNP/SM	U-small nuclear riboprotein complexes (U-snRNPs)	
104	Ro-52	Ro/SSA 52Kd	P19474
105	Ro-60	Ro/SSA 60Kd	P10155
106	SAE1	Small ubiquitin-like modifier-activating enzyme A subunit (SAE1)	M0QX65
107	SAE2/UBA2	SUMO-1 activating enzyme B subunit (SAE2)/Ubiquitin-like modifier activating enzyme 2 (UBA2)	Q9UBT2
108	TOP1 (Scl-70)	DNA topoisomerase I (Scl-70)	P11387
109	SDHA	Flavoprotein subunit of succinate dehydrogenase	P31040
110	SPTBN4	Spectrin, beta, non-erythrocytic 4	Q9H254
111	SRP14	Signal recognition partical 14 (component of 7SL RNA complex)	P37108
112	SRP54	Signal recognition partical 54 (component of 7SL RNA complex)	P61011
113	ADAM17/TACE	Disintegrin and metalloproteinase domain-containing protein 17	P78536
114	TIF-1R	Transcriptional intermediary factor-1 Gamma (TIF-1γ) / TRIM33	Q9UPN9
115	Tn-ITC	Troponin I-T-C complex	
116	TNNC1	Troponin C	P02585
117	TPM1	Tropomyosin 1	P09493
118	TrpRS/WARS	Tryptophanyl-tRNA synthetase	P23381
119	T-T complex	Tropomyosin/Troponin Complex, Bovine	
120	TyrRS/YARS/HA	Tyrosyl-tRNA synthetase (YARS)	Q9Y274
121	Ig control 1	Ig control 1	
122	Ig control 2	Ig control 2	
123	Ig control 3	Ig control 3	
124	Ig control 4	Ig control 4	
125	anti-Ig control 1	anti-Ig control 1	
126	anti-Ig control 2	anti-Ig control 2	
127	anti-Ig control 3	anti-Ig control 3	

128	anti-Ig control 4	anti-Ig control 4	
-----	-------------------	-------------------	--