

OmicsArray™ Myasthenia Gravis (MG) Antigen Array Antigen List - Catalog No. PA017

Antigen serial Number	Antigen Name	Antigen Description	UniProtID
1	a-Actin	Alpha-cardiac actin	P68032
2	a-actinin	Alpha-actinin-1	P12814
3	AChE	Acetylcholinesterase	P22303
4	adiponectin	Adiponectin/Acrp30	Q15848
5	Agrin	Agrin	O00468
6	AITR/GITR/TNFRSF18	Glucocorticoid-induced TNFR-related protein	Q9Y5U5
7	b-actin	Beta-actin	P60709
8	CBG/Serpin A6	corticosteroid-binding globulin	P08185
9	CCL3/MIP1A	C-C motif chemokine 3, macrophage inflammatory protein 1α	P10147
10	CD40	TNF receptor superfamily member 5	P25942
11	CD40L	CD40 ligand	P29965
12	CHRM1/AChR M1	Muscarinic acetylcholine receptor M1	P11229
13	CHRM2/AChR M2	Muscarinic acetylcholine receptor M2	P08172
14	CHRM3/AChR M3	Muscarinic acetylcholine receptor M3	P20309
15	CHRM4	muscarinic acetylcholine receptor M4	P08173
16	CHRM5	muscarinic acetylcholine receptor M5	P08912
17	CHRNA1	Acetylcholine Receptor alpha 1/CHRNA1	P02708
18	CHRNA2	Acetylcholine Receptor alpha 2/CHRNA2	Q15822
19	CHRNA9	Neuronal acetylcholine receptor subunit alpha-9	Q9UGM1
20	CHRNA1 /AChR β	Acetylcholine receptor subunit beta/CHRNA1	P11230
21	CHRNA1 /AChR δ	Acetylcholine receptor subunit delta	Q07001
22	CHRNA1	Acetylcholine receptor subunit epsilon	Q04844
23	CHRNA1 /AChR γ	Acetylcholine Receptor Subunit Gamma (CHRNA1)	P07510
24	Collagen XIII	Collagen alpha-1(XIII) chain/COL13A1	Q5TAT6
25	Cortactin	Cortactin	Q14247
26	CXCL1	Growth-regulated alpha protein	P09341
27	Filamin A (FLNA)	Sapiens filamin A, alpha (actin binding protein 280) (FLNA)	P21333
28	HMGB1	High mobility group protein B1	P09429
29	HSP60/HSPD1	60 kDa heat shock protein, mitochondrial	P10809
30	HSP70 /HSPA1A	Heat shock 70 kDa protein 1A	P0DMV8
31	IFNA1	Interferon alpha-1	P01562
32	IFNA2	Interferon alpha-2	P01563
33	IFN-omega	Interferon omega	P05000
34	IL1a	Interleukin-1 alpha	P01583
35	IL1b	Interleukin-1 beta	P01584
36	IL2	Interleukin-2	P60568
37	IL4	Interleukin-4	P05112
38	IL6	Interleukin-6	P05231
39	IL12	Human IL-12 (p70)	P29460
40	IL17	Interleukin-17A	Q16552
41	Kappa light chain	Immunoglobulin kappa light chain	P0DOX7
42	KCNJ10	ATP-sensitive inward rectifier potassium channel 10	P78508
43	KCNQ1	Potassium voltage-gated channel subfamily KQT member 1	P51787

44	LRP4	Low density lipoprotein receptor-related protein 4	O75096
45	MMP2	Matrix metalloproteinase-2 (MMP-2), 72 kDa type IV collagenase	P08253
46	MMP3	Matrix metalloproteinase-3 (MMP-3), Stromelysin-1	P08254
47	MMP9	Matrix metalloproteinase-9	P14780
48	MuSk (ECD)	MuSK (extra-cellular domain)	O15146
49	MuSK (ICD)	MuSK (intra-cellular domain)	O15146
50	Myosin	Myosin-1	P12882
51	Myosin light chain	myosin light chain	P05976
52	PD-1	Programmed cell death protein 1	Q15116
53	RAGE	Advanced glycosylation end product-specific receptor	Q15109
54	Rapsyn	43 kDa receptor-associated protein of the synapse	Q13702
55	Resistin	Resistin	Q9HD89
56	RyR1	Ryanodine receptor 1 (749-1175aa)	P21817
57	RyR1	Ryanodine receptor 1 (2591-2939aa)	P21817
58	S100A12	Protein S100-A12	P80511
59	S100B	Protein S100-B	P04271
60	Titin (TTN)	Titin (14257-14543aa)	Q8WZ42
61	TNF alpha	Tumor necrosis factor	P01375
62	TNF beta	Lymphotoxin-alpha	P01374
63	Tropomyosin	Tropomyosin beta chain	P07951
64	UPAR	Urokinase plasminogen activator surface receptor	Q03405
65	Vimentin	Vimentin	P08670
66	Vinculin	Vinculin	P18206
67	Vitronectin	Vitronectin	P04004
68	ICOSL (CD275)/B7-H2	ICOS ligand, B7 homolog 2 (B7-H2)	O75144
69	Ig control 1	control	
70	Ig control 2	control	
71	Ig control 3	control	
72	Ig control 4	control	
73	anti-Ig control 1	control	
74	anti-Ig control 2	control	
75	anti-Ig control 3	control	
76	anti-Ig control 4	control	