

**OmicsArray™ Cardiac Autoimmune Diseases (CAD) antigen array, 54-plex Antigen List -
Catalog No. PA009**

Antigen#	ID	Antigen Name
1	20S P	20S proteasome complex
2	ACO2	Mitochondrial aconitase
3	ACTA1	Alpha actin (Skeletal)
4	ACTB	Beta Actin
5	ACTC1	Actin Alpha 1, α -cardiac actin, Cardiac Muscle
6	ACTN1	α -Actinin
7	ATP1A1	ATPase, Na ⁺ /K ⁺ transporting, alpha 1
8	ATP1B1	ATPase Na ⁺ /K ⁺ Transporting Subunit Beta 1
9	ATP1B4	ATPase Na ⁺ /K ⁺ Transporting Subunit Beta 4
10	ATP2C1	ATPase secretory pathway Ca ²⁺ transporting 1
11	BCKDHA	branched chain alpha-ketoacid dehydrogenase
12	CALR	Calreticulin
13	CDH1	E-Cadherin/Cadherin-1
14	CDH2	Cadherin 2
15	CDH8	Cadherin 8 (CDH8)
16	CDH9	Cadherin 9 (CDH9)
17	CHRM2	Muscarinic acetylcholine receptor M2
18	CHRM3	Cholinergic Receptor Muscarinic 3 (CHRM3)
19	cMyBP-C	Cardiac Myosin Binding Protein-C
20	cTnl	Cardiac Troponin I
21	DLG4	Disks Large Homolog 4/DLG4/PSD95
22	DSC1	Desmocollin-1
23	DSC2	Desmocollin-2
24	DSC3	Desmocollin-3
25	DSG1	Desmoglein 1 (DSG1)
26	DSG2	Desmoglein 2 (DSG2)
27	DSG3	Desmoglein-3 (DSG-3)
28	DSG4	Desmoglein-4 (DSG-4)
29	GAPDH	Glyceraldehyde-3-phosphate dehydrogenase
30	GJA1	Connexin 43/GJA1
31	HSP60/HSPD1	heat shock protein-60/chaperonins
32	HSP70/HSP1A1	heat shock protein-70/Hu HSP70/HSPA1A Protei
33	ITGB1	Integrin Subunit Beta 1
34	ITGB4	Integrin, beta 4 (ITGB4) /CD104
35	JUP	Junction Plakoglobin
36	KCNIP2	Kv channel-interacting protein2.6
37	KRT24	Keratin 24
38	M1MR	M1 Muscarinic receptor (human)
39	M2	M2 (Mitochondrial 2-oxo acid dehydrogenase
40	M3MR	M3 Muscarinic receptor (human) membrane
41	M4MR	M4 Muscarinic receptor (human) membrane
42	M5MR	M5 Muscarinic receptor, Human

43	M5MR	M5 Muscarinic receptor (human) membrane
44	MYH1	Myosin Heavy Chain (from human heart)
45	MYH9	Human Myosin-9 (MYH9), aa2-241
46	MYL1	Myosin Light Chain from human heart
47	Myosin	Myosin Calcium activated from bovine muscle
48	Myosin Cardiac	Myosin (bovine cardiac muscle)
49	PYGB	Brain glycogen phosphorylase
50	SDHA	flavoprotein subunit of succinate
51	SPTBN4	Spectrin, beta, non-erythrocytic 4
52	Tn-ITC	Troponin I-T-C complex
53	TPM1	Tropomyosin 1
54	T-T complex	Tropomyosin/Troponin Complex, Bovine
55	Ig control 1	Ig control 1
56	Ig control 2	Ig control 2
57	Ig control 3	Ig control 3
58	Ig control 4	Ig control 4
59	anti-Ig control 1	anti-Ig control 1
60	anti-Ig control 2	anti-Ig control 2
61	anti-Ig control 3	anti-Ig control 3
62	anti-Ig control 4	anti-Ig control 4