

OmicsArray™ Kidney Autoimmune Antigen Array, 51-plex				Catalog No. PA020
Antigen Number	Antigen ID	Antigen Name	Gene symbol	UniProt ID
1	ACO2	Aconitate hydratase, mitochondrial	ACO2	Q99798
2	Vcl	Vinculin	VCL	P18206
3	CALD1	Caldesmon	CALD1/CAD/CDM	Q05682
4	LMNA	Prelamin-A/C	LMNA/LMN1	P02545
5	SRSF9	Serine/arginine-rich splicing factor 9	SRSF9/SFRS9, SRP30C	Q13242
6	PPID	Peptidyl-prolyl cis-trans isomerase D	PPID/CYP40, CYPD	Q08752
7	CAPZB	F-actin-capping protein subunit beta	CAPZB	P47756
8	PSMA1	Proteasome subunit alpha type-1	PSMA1/HC2/NU/PROS30/PSC2	P25786
9	PRDX1	Peroxiredoxin-1	PRDX1/PAGA/ PAGB/ TDPX2	Q06830
10	PLA2R	Secretory phospholipase A2 receptor	PLA2R1/CLEC13C	Q13018
11	THSD7A	Thrombospondin type-1 domain-containing protein 7A	THSD7A/KIAA0960	Q9UPZ6
12	ANXA1	Annexin A1	ANXA1/ANX1/LPC1	P04083
13	ANXA2	Annexin A2	ANXA2/ANX2/ANX2L4/ CAL1H/LPC2D	P07355
14	NEPH1 (KIRREL1)	Kin of IRRE-like protein 1	KIRREL1/KIRREL, NEPH1	Q96J84
15	PTPRO	Receptor-type tyrosine-protein phosphatase O	PTPRO/GLEPP1/ PTPU2	Q16827
16	PODXL	Podocalyxin	PODXL/PCLP, PCLP1	O00592
17	TRPC6	Transient receptor protein 6 (TRP-6)	TRPC6/TRP6	Q9Y210
18	Nephrin	Renal glomerulus-specific cell adhesion receptor	NPHS1/NPHN	O60500
19	ITGA3	Integrin alpha 3 beta 1/VLA-3 Protein	ITGA3/MSK18	P26006
20	DAG1	Dystroglycan	DAG1	Q14118
21	NEPH2 (Kirrel3)	Kin of IRRE-like protein 3	KIRREL3/KIAA1867/NEPH2	Q8IZU9
22	AGTR1	Type-1 angiotensin II receptor	AGTR1/AGTR1A/AGTR1B/AT2R1/ AT2R2	P30556
23	Synpo	Synaptopodin	SYNPO/KIAA1029	Q8N3V7
24	NPHS2	Podocin	NPHS2	Q9NP85
25	CNTN1	Contactin 1	CNTN1	Q12860
26	GBM	Glomerular basement membrane protein		
27	NRP2	Neuropilin-2	NRP2/VEGF165R2	O60462
28	NRP1	Neuropilin-1	NRP1/NRP/VEGF165R	Q14786
29	Tubulin	Tubulin protein from Porcine brain	TUBA1A	P02550
30	Tubulin alpha	alpha Tubulin	TUBA1A/TUBA3	Q71U36
31	MPO	Myeloperoxidase (pANCA)	MPO	P05164
32	PR3	Leukocyte proteinase 3	PRTN3/MBN	P24158
33	dsDNA	double strand DNA		
34	ssDNA	single strand DNA		
35	Sm	U-small nuclear riboprotein complexes		
36	Sm/RNP	Bovine Sm/RNP ribonuclear protein complex		
37	RO52	E3 ubiquitin-protein ligase TRIM21	TRIM21/RNF81/ RO52/SSA1	P19474
38	RO60	RNA-binding protein RO60	RO60/SSA2/TROVE2	P10155
39	La/SSB	Lupus La protein	SSB	P05455
40	SSA/SSB	SSA/SSB antigen		
41	Scl70	Scl-70 (Topoisomerase I)	TOP1	P11387
42	PM/SCL75	Human PM/Scl 75 protein component of nuclear exosome		

43	PM/SCL100	Human PM/Scl 100 protein component of nuclear exosome		
44	Jo-1	Histidyl-tRNA synthetase (HRS/Jo-1)	HARS1/HARS/HRS	P12081
45	CENP-B	Centromere protein B (CENP-B)	CENPB	P07199
46	PCNA	Proliferating cell nuclear antigen (PCNA)	PCNA	P12004
47	M2	Mitochondrial 2-oxo acid dehydrogenase multienzyme complexes		
48	CRP	C-reactive protein (CRP)	CRP/PTX1	P02741
49	Ribo P	Ribosomal P proteins (P0, P1, P2)		
50	Histone	Total Histone		
51	Nucleosome	Nucleosome (non-recombinant/bovine)		
52	Buffer			
53	Ig control 1	Ig control		
54	Ig control 2	Ig control		
55	Ig control 3	Ig control		
56	Ig control 4	Ig control		
57	anti-Ig control 1	anti-Ig control		
58	anti-Ig control 2	anti-Ig control		
59	anti-Ig control 3	anti-Ig control		
60	anti-Ig control 4	anti-Ig control		