

OmicsArray™ Rheumatoid Arthritis (RA) Antigen Array			Catalog No. PA022	
Antigen Number	Antigen ID	Antigen Name	Gene symbol	UniProt ID
1	CCP2	cyclic citrullinated peptide 2		
2	PAD2	Peptidylarginine deiminase II	PADI2	Q9Y2J8
3	PAD4	Peptidylarginine deiminase IV	PADI4	Q9UM07
4	Collagen type II	Collagen Alpha-1 type II	COL2A1	P02458
5	H2A	Histone H2A	H2AC7	P20671
6	H2B	Histone H2B	H2BC3	P33778
7	H3	Histone H3	H3Y1	P0DPK2
8	H4	Histone H4	H4C1; H4C2; H4C3; H4C4; H4C5; H4C6; H4C8; H4C9; H4C11; H4C12; H4C13; H4C14; H4C15; H4C16	P62805
9	Ro52 (SSA-52)	Ro/SSA 52KD, TRIM21	TRIM21	P19474
10	Ro60 (SSA-60)	Ro/SSA 60KD, TROVE2	RO60	P10155
11	Centromere B	Centromere protein B (CENP-B)	CENPB	P07199
12	IGFBP2	Insulin-like Growth Factor Binding Protein 2	IGFBP2	P18065
13	Fibrinogen	Fibrinogen	FGB	P02675
14	Cit-Fibrinogen	Human Fibrinogen (PAD2 Citrullinated)	FGB	P02676
15	α-Enolase	Alpha-enolase (ENO1)	ENO1	P06733
16	Cit-α-Enolase (CEP1)	Citrullinated α-Enolase (human)	ENO1	P06733
17	Cit-H2A	Citrullinated Histone H2A Type 1 (human)	H2AC7	P20671
18	Cit-H2B	Citrullinated Histone H2B	H2BC3	P33778
19	Cit-H3	Citrullinated Histone H3	H3Y1	P0DPK2
20	Cit-H4	Citrullinated Histone H4 (human)	H4C1; H4C2; H4C3; H4C4; H4C5; H4C6; H4C8; H4C9; H4C11; H4C12; H4C13; H4C14; H4C15; H4C16	P62805
21	Vimentin	Vimentin	VIM	P08670
22	Cit-Vimentin (Sa)	Citrullinated Vimentin	VIM	P08670
23	cit-LL37	Citrullinated LL37		
24	Cit-FLG	citrullinated Filaggrin	FLG	P20930
25	GNPTG	N-acetylglucosamine-1-phosphotransferase subunit gamma	GNPTG	Q9UJJ9
26	RF	Rheumatoid factor (RF)		
27	BRAF	Serine/threonine-protein kinase B-raf, v-Raf murine sarcoma viral oncogene homolog B1	BRAF	P15056
28	Acetylated ornithine (AcOrn)	Nα-Acetyl-L-ornithine		
29	N-acetylglucosamine -1-phosphatase transferase (OGT)	Human O-GlcNAcTransferase/OGT Protein	OGT	O15294
30	Malondialdehyde-acetaldehyde (MAA) modified protein	MAA modified albumin (MAA-Alb)		
31	Carbamylated Albumin	Carbamylated Albumin (Car-Alb)		
32	Carbamylated Fibrinogen	Carbamylated fibrinogen (Car-Fib)		
33	Ig control 1	Internal control		
34	Ig control 2	Internal control		
35	Ig control 3	Internal control		
36	Ig control 4	Internal control		
37	anti-Ig control 1	positive control		
38	anti-Ig control 2	positive control		
39	anti-Ig control 3	positive control		
40	anti-Ig control 4	positive control		