

OmicsArray™ Alzheimer's Disease Autoantigen Array			Catalog No. PA023-V2
Antigen serial No.	Antigen ID	Antigen Name	UniProtID
1	Amyloid B	Amyloid β Protein Fragment 1-40	
2	Tau	Microtubule-associated protein Tau	P10636
3	pTau 181	Phospho-Tau (Thr181)	
4	ATCAY	Caytaxin	Q86WG3
5	NME7	Nucleoside diphosphate kinase homolog 7	Q9Y5B8
6	ITPR1	Inositol 1,4,5-triphosphate receptor type 1 (ITPR1)	Q14643
7	WDR1	WD repeat- containing protein 1 (WDR1)	O75083
8	ASRGL1	Isoaspartyl peptidase/L-asparaginase	Q7L266
9	CDKN1A	Cyclin-dependent kinase inhibitor 1	P38936
10	BACE1	Beta-secretase 1	P56817
11	BIM	Bcl-2-like protein 11	O43521
12	OMGP	Oligodendrocyte- myelin glycoprotein (OMG)	P23515
13	MOG	Myelin-oligodendrocyte glycoprotein (MOG)	Q16653
14	S100b	Protein S100-B	P04271
15	GFAP	Glial fibrillary acidic protein	P14136
16	YWHAG (14-3-3 γ)	14-3-3 protein gamma	P61981
17	NPTX2	Neuronal pentraxin-2	P97738
18	HIST1H3F	Histone H3.1	P68431
19	AGER	Advanced glycosylation end product-specific receptor	Q15109
20	AGTR1	Type-1 angiotensin II receptor	P30556
21	KDM4D	Lysine-specific demethylase 4D	Q6B016
22	PLD4	Phospholipase D4 (PLD4)	Q96BZ4
23	SRPX	Sushi repeat- containing protein (SRPX)	P78539
24	ATP2A2	Sarco/endoplasmic reticulum calcium ATPase 2 (ATP2A2)	P16615
25	PAIP2	Polyadenylate-binding protein-interacting protein 2	Q9BPZ3
26	DNAJC8	DnaJ homolog subfamily C member 8	O75937
27	hnRNP H	Heterogeneous nuclear ribonucleoprotein H (hnRNP H)	P31943
28	CBLN3	Cerebellin-3 (CBLN3)	Q6UW01
29	ADD2	Beta-adducin	P35612
30	SERF1A	Small EDRK-rich factor 1	O75920
31	CTSF	Cathepsin F (CTSF)	Q9UBX1
32	SERPINE2	Glia-derived nexin	P07093
33	FMOD	Fibromodulin (FMOD)	Q06828
34	CDH13	Cadherin-13 (CDH13)	P55290
35	NQO1	NAD(P)H dehydrogenase [quinone] 1 (NQO1)	P15559
36	SLC3A2	4F2 cell-surface antigen heavy chain (SLC3A2)	P08195
37	TIMP-2	Metalloproteinase inhibitor 2 (TIMP-2)	P16035
38	IVD	Isovaleryl-CoA dehydrogenase	P26440
39	ASXL1	Polycomb group protein ASXL1	Q8IXJ9
40	CYFIP1	Cytoplasmic FMR1-interacting protein 1	Q7L576
41	Ig control 1	IgG, IgM, IgA and IgE (human and mouse)	
42	Ig control 2		
43	Ig control 3		
44	Ig control 4		

45	anti-Ig control 1	anti-IgG, anti-IgM, anti-IgA and anti-IgE (human and mouse)	
46	anti-Ig control 2		
47	anti-Ig control 3		
48	anti-Ig control 4		