

OmicsArray™ Alzheimer's Disease Autoantigen Array			Catalog No. PA023-V1
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	Ig control 1	IgG, IgM, IgA and IgE (human and mouse)	
22	Ig control 2		
23	Ig control 3		
24	Ig control 4		
25	anti-Ig control 1	anti-IgG, anti-IgM, anti-IgA and anti-IgE (human and mouse)	
26	anti-Ig control 2		
27	anti-Ig control 3		
28	anti-Ig control 4		