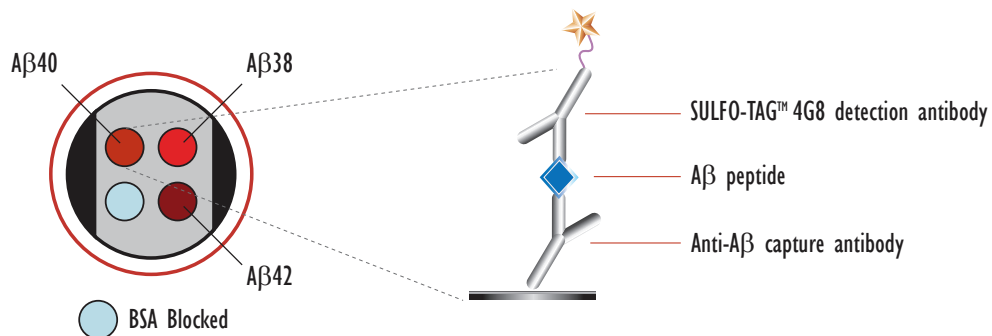
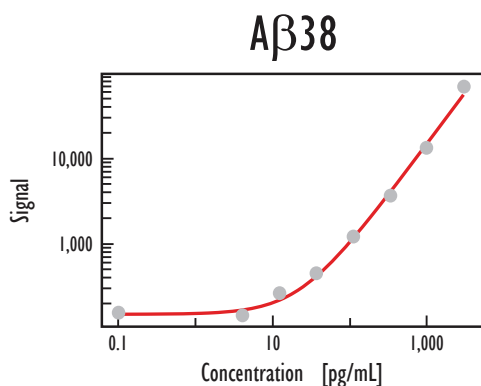


MULTI-SPOT[®] Human/Rodent (4G8) Abeta Triplex Ultra-Sensitive Assay

Detection of A β 38, A β 40, and A β 42

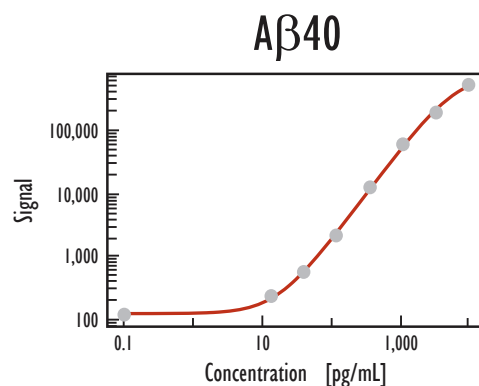


Kit Size	Catalog Number
1 plate	K15141E-1
5 plates	K15141E-2
20 plates	K15141E-3
20 plates (Base)	K15141A-3



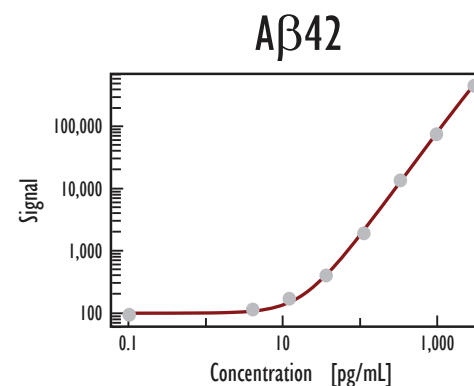
Peptide (pg/mL)	A β 38		
	Average	StdDev	%CV
0	154	17	11
4.1	142	32	23
12.3	256	16	6
37	443	25	6
111	1,187	17	1
333	3,693	65	2
1,000	13,498	1,128	8
3,000	69,717	5,664	8

A β 38 Detection Limit
(2.5 StdDev over background) **9 pg/mL**



Peptide (pg/mL)	A β 40		
	Average	StdDev	%CV
0	121	4	4
13.7	240	6	2
41	562	23	4
123	2,210	121	5
370	12,830	912	7
1,111	61,961	4,259	7
3,333	205,990	16,488	8
10,000	527,001	30,722	6

A β 40 Detection Limit
(2.5 StdDev over background) **5 pg/mL**



Peptide (pg/mL)	A β 42		
	Average	StdDev	%CV
0	92	14	15
4.1	116	6	6
12.3	165	3	2
37	402	30	7
111	1,939	139	7
333	13,557	1,355	10
1,000	86,511	10,396	12
3,000	437,890	24,665	6

A β 42 Detection Limit
(2.5 StdDev over background) **8 pg/mL**

Synthetic A β peptides are diluted in 1% BSA/TBST and incubated with SULFO-TAG 4G8 detection antibody on an MSD MULTI-SPOT, 96-well 4 Spot plate coated with A β peptide-specific antibodies. This kit and protocol can be adapted for use with cell culture medium, plasma and brain homogenates. Representative titrations of A β peptides using Kit Version 2 are shown.