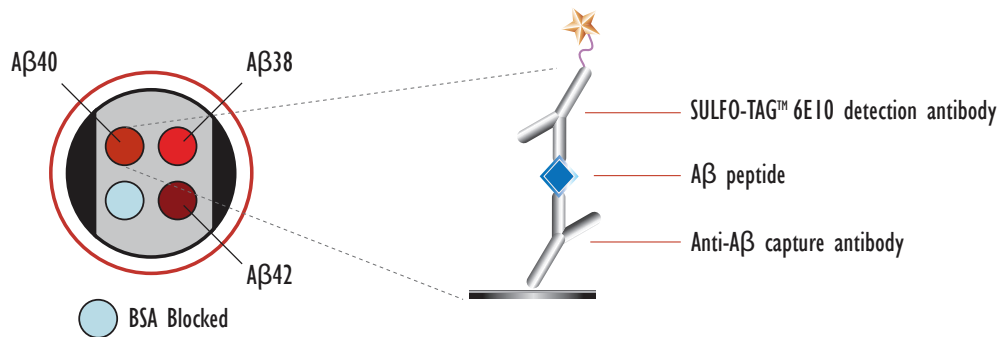
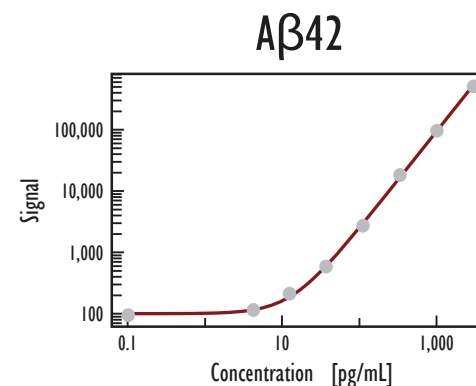
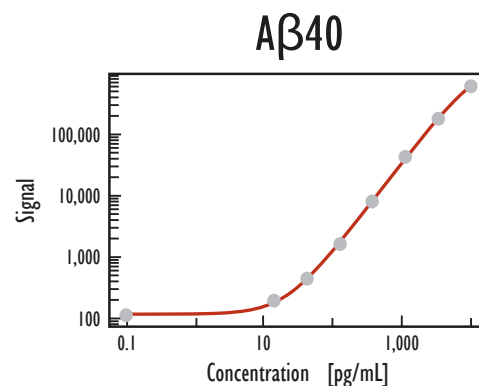
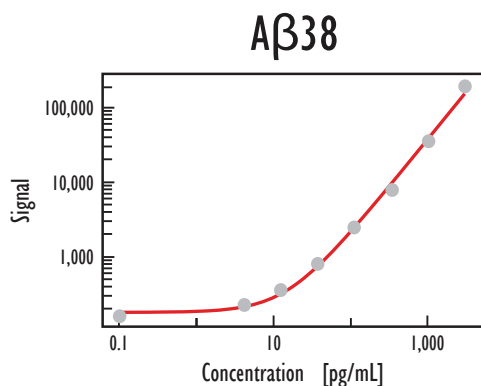


MULTI-SPOT[®] Human (6E10) Abeta Triplex Assay

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



Kit Size	Catalog Number
1 plate	K15148E-1
5 plates	K15148E-2
20 plates	K15148E-3
20 plates (Base)	K15148A-3



Peptide (pg/mL)	A β 38		
	Average	StdDev	%CV
0	158	18	11
4.1	221	20	9
12.3	348	28	8
37	782	49	6
111	2,541	383	15
333	8,028	1,454	18
1,000	35,809	4,218	12
3,000	207,371	23,013	11

A β 38 Detection Limit
(2.5 StdDev over background) **4 pg/mL**

Peptide (pg/mL)	A β 40		
	Average	StdDev	%CV
0	110	10	9
13.7	194	5	2
41	436	17	4
123	1,610	106	7
370	8,056	814	10
1,111	44,103	3,141	7
3,333	180,916	10,668	6
10,000	636,315	18,173	3

A β 40 Detection Limit
(2.5 StdDev over background) **6 pg/mL**

Peptide (pg/mL)	A β 42		
	Average	StdDev	%CV
0	93	8	9
4.1	119	4	4
12.3	217	12	6
37	576	96	17
111	2,753	297	11
333	18,099	1,889	10
1,000	95,493	25,992	27
3,000	536,356	47,048	9

A β 42 Detection Limit
(2.5 StdDev over background) **5 pg/mL**

Synthetic A β peptides are diluted in 1% BSA/TBST and incubated with SULFO-TAG 6E10 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.