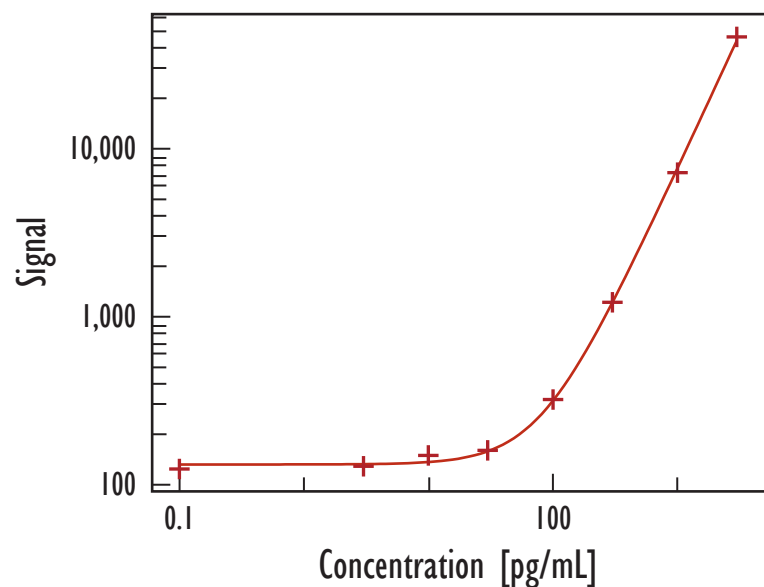


# MULTI-ARRAY<sup>®</sup> Human A $\beta$ 42 Ultra-Sensitive Assay

## Detection of Human Beta Amyloid 42 in Cerebral Spinal Fluid Samples

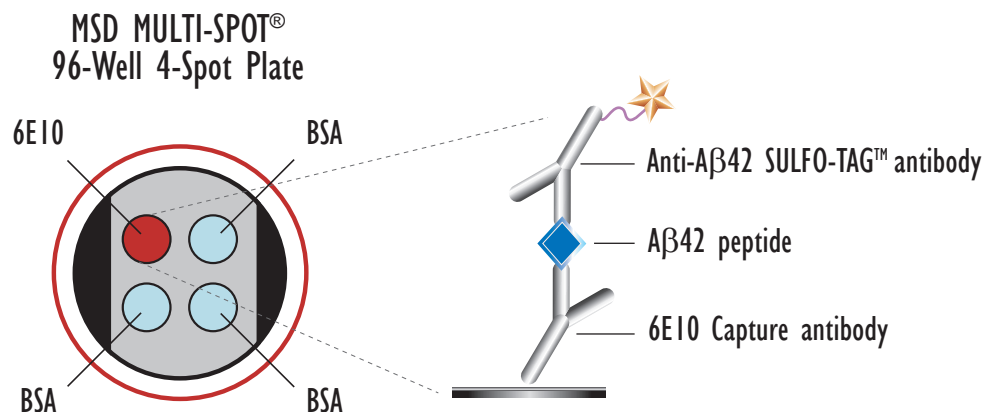
A $\beta$ 42



Peptide (pg/mL)	A $\beta$ 42		
	Average	StdDev	%CV
0	122	6	5
4.1	126	29	23
12.3	146	16	11
37	156	13	8
111	316	23	7
333	1,212	29	2
1,000	7,238	300	4
3,000	46,750	2,016	4

**A $\beta$ 42 Lower Limit of Detection 23 pg/mL**

LLOD (Lower Limit of Detection) is defined as 2.5x stdev above the background.



Synthetic A $\beta$ 42 peptides were diluted in 10% BSA/TBST and added to MSD MULTI-SPOT, 96-well 4 Spot plates coated with 6E10 antibody. A $\beta$ 42 peptides were detected with anti-A $\beta$ 42 labeled with MSD SULFO-TAG reagent. A representative titration curve is shown.

Kit Size	Catalog Number
1 plate	K151FUE-1
5 plates	K151FUE-2
20 plates	K151FUE-3
20 plates (Base)	K151FUA-3

# MULTI-ARRAY<sup>®</sup> Human A $\beta$ 42 Ultra-Sensitive Assay

## Detection of Human Beta Amyloid 42 in Cerebral Spinal Fluid Samples

### Spike Recovery

- Measured analyte from spike recovery in 3 pooled human CSF

$$\% \text{ recovery} = \frac{(\text{measured value} * 100)}{\text{expected value}}$$

Spike (pg/mL)	CSF Pool 1	CSF Pool 2	CSF Pool 3
500	104	105	116
250	112	107	112
125	110	105	112

### Dilution Linearity

- Measured analyte levels in 3 pooled human CSF samples followed by subsequent dilution

$$\% \text{ recovery} = \frac{(\text{measured value} * \text{dilution factor} * 100)}{\text{predicted value}}$$

Dilution Factor	CSF Pool 1	CSF Pool 2	CSF Pool 3
1/2	100	107	103
1/4	94	106	99