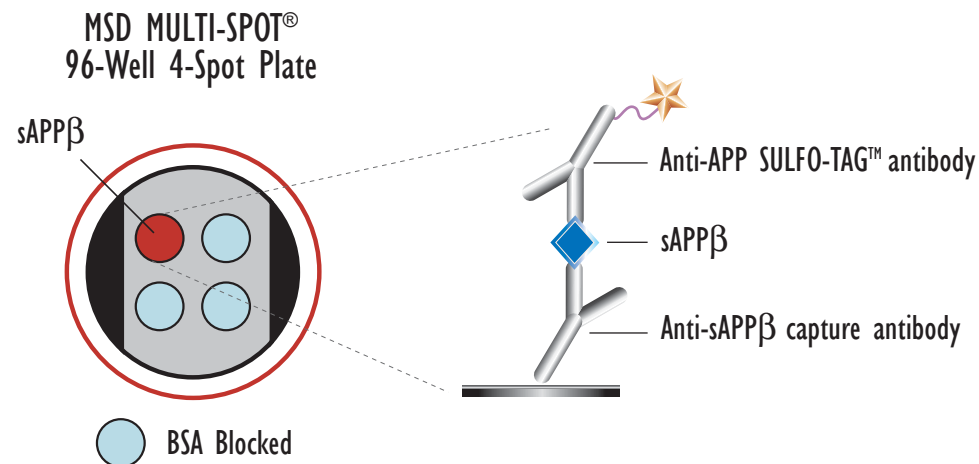
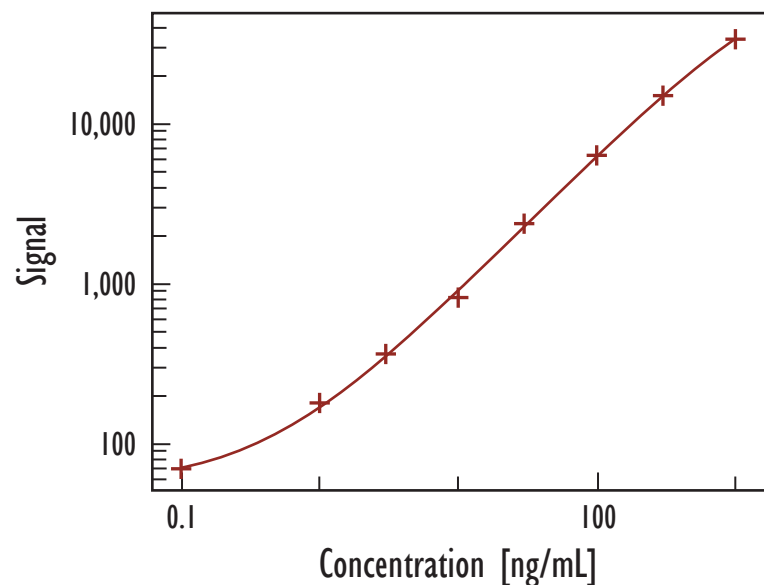


MULTI-ARRAY[®] Mouse/Rat sAPP β Assay

Detection of Secreted Fragment of Amyloid Precursor Protein β in Mouse and Rat Cerebrospinal Fluid



Recombinant human sAPP β was purified from mammalian cells (>95% pure) and diluted in denaturing assay diluent. Samples were added to MSD MULTI-SPOT 4 spot plates coated with anti-sAPP β antibody on one of four spatially distinct electrodes per well. sAPP β protein was detected with anti-APP antibody (22C11), labeled with MSD SULFO-TAG reagent.

Peptide (ng/mL)	sAPP β		
	Average	StdDev	%CV
0	71	7	10
1.4	181	12	7
4.1	369	27	7
12.3	824	36	4
37	2,387	37	2
111	6,349	1,047	16
333	14,923	321	2
1,000	33,336	3,291	10

sAPPβ LLOD	0.4 ng/mL
------------------------------------	------------------

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

Kit Size	SI2400	SI6000
1 plate	K150BTE-1	K110BTE-1
5 plates	K150BTE-2	K110BTE-2
20 plates	K150BTE-3	K110BTE-3
20 plates (Base)	K150BTA-3	K110BTA-3

MULTI-ARRAY[®] Mouse/Rat sAPP β Assay

Detection of Secreted Fragment of Amyloid Precursor Protein β in Mouse and Rat Cerebrospinal Fluid

Spike Recovery

- Measured analyte from spike recovery in 3 Rodent CSF samples

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

Spike (ng/mL)	Average % Recovery
500	110
250	99
125	90
62.5	94

Dilutional Linearity

- MSD recommends that Rodent CSF samples be run neat
- Measured analyte in 3 Rodent CSF samples followed by subsequent dilution

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

Dilution	Average % Recovery
1/2	78
1/4	72