

# MSD<sup>®</sup> sAPP Recombinant Protein Standards

## Recombinant Human sAPP $\beta$

**Contents:** 2.5  $\mu$ g recombinant sAPP $\beta$  protein

**Concentration:** 50  $\mu$ g/mL

**Volume:** 50  $\mu$ L

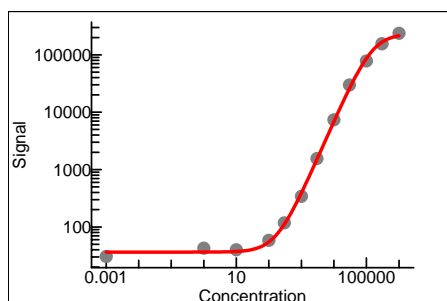
**Preparation:** Recombinant human sAPP $\beta$  protein was purified from overexpressing mammalian cells.

**Storage:** **Store at -80°C.** After initial thaw, protein should be divided into five single-use aliquots (9  $\mu$ L, allowing for 6  $\mu$ L recovery) to avoid additional freeze/thaw cycles.

**Quality Control:** Recombinant protein has been analyzed by SDS-PAGE and MSD MULTI-SPOT<sup>®</sup> Assays.

## MSD MULTI-SPOT<sup>®</sup> Assay Results

Typical titration curve for recombinant sAPP $\beta$  using the MSD MULTI-SPOT sAPP $\beta$  assay.



conc	Ave	StdDev	%CV	P/N
0	31	15	49	
1 pg/ml	43	15	34	1
10 pg/ml	40	8	20	1
100 pg/ml	59	13	23	2
300 pg/ml	118	12	10	4
1 ng/ml	342	16	5	11
3 ng/ml	1553	72	5	51
10 ng/ml	7350	541	7	240
30 ng/ml	29802	2262	8	972
100 ng/ml	78230	4371	6	2551
300 ng/ml	155928	6771	4	5085
1 ug/ml	237557	7991	3	7746

Detection limit (3 S.D. over background): 182 pg/ml

## SDS-PAGE

A 0.5 mg sample of each sAPP protein was run on a 4-12% Bis-Tris NuPAGE gel to demonstrate purity (>95%).

