

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

## Recombinant Human Swedish sAPP $\beta$

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

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

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

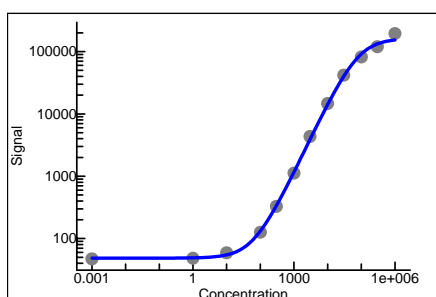
**Preparation:** Recombinant human sAPP $\beta$  protein containing the Swedish amino acid substitutions (670, 671) was purified from overexpressing mammalian cells.

**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 Swedish sAPP $\beta$  using the MSD MULTI-SPOT Swedish sAPP $\beta$  assay.



Conc	Ave	StDev	%CV	P/N
0	47	9	19	
1 pg/ml	48	11	24	1
10 pg/ml	59	22	37	1
100 pg/ml	126	18	14	3
300 pg/ml	329	15	4	7
1 ng/ml	1120	37	3	24
3 ng/ml	4345	176	4	92
10 ng/ml	14664	356	2	310
30 ng/ml	41785	125	0	883
100 ng/ml	81860	20236	25	1729
300 ng/ml	119720	3060	3	2529
1 ug/ml	194794	19973	10	4115

Detection limit (3 S.D. over background): 39 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%).

