

MSD[®] sAPP Recombinant Protein Standards

Recombinant Human sAPP α

Contents: 2.5 μ g recombinant sAPP α protein

Concentration: 50 μ g/mL

Volume: 50 μ L

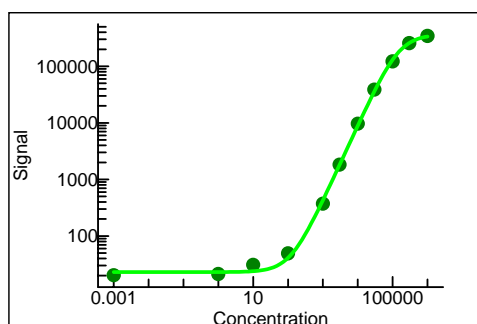
Preparation: Recombinant human sAPP α protein was purified from overexpressing mammalian cells.

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

Quality Control: Recombinant protein has been analyzed by SDS-PAGE and MSD MULTI-SPOT[®] Assays.

MSD MULTI-SPOT[®] Assay Results

Typical titration curve for recombinant sAPP α using the MSD MULTI-SPOT sAPP α Assay.



Conc	Ave	StdDev	%CV	S/B
0	20	22	109	
1 pg/ml	32	8	27	2
10 pg/ml	47	6	14	2
100 pg/ml	74	3	4	4
300 pg/ml	92	11	11	5
1 ng/ml	373	16	4	18
3 ng/ml	1827	181	10	90
10 ng/ml	9711	423	4	478
30 ng/ml	38780	971	3	1907
100 ng/ml	122396	1162	1	6019
300 ng/ml	257081	8794	3	12643
1 ug/ml	344233	13166	4	16930

Detection limit (3 S.D. over background): 254 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%).

