

MSD[®] 96-Well MULTI-SPOT sAPP α /sAPP β Assay

Base Catalog No: K15120E

I. Materials Included

Reagent	Storage	Catalog No.	Size	Quantity Supplied		
				1-Plate Kit	5-Plate Kit	20-Plate Kit
MULTI-SPOT [®] 96-well 4-spot sAPP α /sAPP β Plate	2–8 °C	N45120B-1	4-spot	1	5	20
APP Antibody (50X) (SULFO-TAG [™] Detection Antibody)	2–8 °C	D21EA-2	1 vial	1	1	4
sAPP α Calibrator (50 μ g/mL)	\leq -70 °C	C01BS-2	1 vial	1	5	20
sAPP β Calibrator (50 μ g/mL)	\leq -70 °C	C01BT-2	1 vial	1	5	20
Blocker A (dry powder)	RT	R93BA-4	15 g	1	1	1
Read Buffer T (4X)	RT	R92TC-3	50 mL	1	1	-
		R92TC-2	200 mL	-	-	1
Tris Wash Buffer (10X)	2–8 °C	R61TX-2	200 mL	1	1	-
		R61TX-1	1000 mL	-	-	1

RT = room temperature
Dash (-) = not applicable

II. Other Materials & Equipment (not supplied)

- Deionized water for diluting concentrated buffers
- 500 mL bottle
- 50 mL tubes
- 15 mL tubes
- Adhesive plate seals
- Microtiter plate shaker
- Various microcentrifuge tubes for making serial dilutions of supernatants (if desired)
- Automated plate washer or other efficient multi-channel pipetting equipment for washing 96-well plates
- Appropriate liquid handling equipment for desired throughput that must accurately dispense 25 μ L and 150 μ L into a 96-well micro plate
- Vortex mixer

Note: A spot map identifying the location of each assay can be found on the plate packaging. This information will be needed for data analysis.

FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.

III. Protocol at a Glance

1. Add blocking solution, incubate 1 hour, wash.
2. Add calibrator or samples, incubate 1 hour, wash.
3. Add detection antibody, incubate 1 hour, wash.
4. Add Read Buffer T and analyze the plate.

The following protocol is optimized for quantifying sAPP α and sAPP β . The protocol takes approximately 3 to 3½ hours to complete. All reagents can be prepared ahead of time. This lengthens the overall time required for the assay but frees up time during incubation steps.

Notes:

Read the entire detailed instructions before beginning work.

IV. Detailed Instructions

Prepare a stock of 1X Tris Wash Buffer. 1X Tris Wash Buffer is used throughout the assay to dilute other reagents and wash the plates. Approximately 350 mL per plate is required—more if using an automatic plate washer.

In a 500 mL bottle, combine:

- 35 mL 10X Tris Wash Buffer
- 315 mL deionized water

A larger amount of Tris Wash Buffer may be prepared at once and stored at room temperature for later use.

Prepare 3% Blocker A Solution. You will need 20 mL per plate.

In a 50 mL tube, combine:

- 20 mL 1X Tris Wash Buffer
- 600 mg Blocker A (3% w/v)

Solutions containing Blocker A should be kept at 2–8 °C and discarded after 14 days.

Prepare Antibody Dilution Buffer. You will need 3 mL per plate.

In a 15 mL tube, combine:

- 2 mL 1X Tris Wash Buffer
- 1 mL of 3% Blocker A solution

Prepare Detection Antibody Solution. You will need 3 mL per plate.

In a 15 mL tube, combine:

- 60 μ L 50X SULFO-TAG Anti-APP Detection Antibody
- 2.94 mL cold Antibody Dilution Buffer

Detection antibody solution should be stored in the dark at 2–8 °C.

Prepare Read Buffer T. MSD provides Read Buffer T as a 4X stock solution. The working solution is 1X. You will need 20 mL per plate at a 1X concentration.

In a 50 mL tube, combine:

- 15 mL deionized water
- 5 mL 4X Read Buffer T

Diluted read buffer may be kept in a tightly sealed container at room temperature for later use.

Prepare Standards

1000 ng/mL:	6 µL of 50 µg/mL sAPP α Calibrator plus 6 µL of 50 µg/mL sAPP β Calibrator solution plus 288 µL diluent
316 ng/mL:	100 µL of the 1000 ng/mL solution plus 216 µL diluent
100 ng/mL:	100 µL of the 316 ng/mL solution plus 216 µL diluent
32 ng/mL:	100 µL of the 100 ng/mL solution plus 216 µL diluent
10 ng/mL:	100 µL of the 32 ng/mL solution plus 216 µL diluent
3.2 ng/mL:	100 µL of the 10 ng/mL solution plus 216 µL diluent
1 ng/mL:	100 µL of the 3.2 ng/mL solution plus 216 µL diluent
0.32 ng/mL:	100 µL of the 1 ng/mL solution plus 216 µL diluent
0.10 ng/mL:	100 µL of the 0.32 pg/mL solution plus 216 µL diluent
0.032 ng/mL:	100 µL of the 100 pg/mL solution plus 216 µL diluent
0.010 ng/mL:	100 µL of the 0.032 pg/mL solution plus 216 µL diluent
0 ng/mL:	diluent alone

Begin with a MULTI-SPOT sAPP α /sAPP β plate. No pre-treatment is necessary.

STEP 1 Add Blocker A Solution

- Add 150 µL/well of 3% Blocker A solution to the plate(s).
- Incubate the plate(s) at room temperature with shaking for 1 hour.
- Wash the plate(s) three times with 300 µL/well of 1X Tris Wash Buffer.

STEP 2 Add Sample or Calibrator

- Add 25 µL/well of samples or calibrator to the plate(s).
- Incubate the plate(s) at room temperature with shaking for 1 hour.
- Wash the plate(s) three times with 300 µL/well of 1X Tris Wash Buffer.

STEP 3 Add Detection Antibody

- Add 25 µL/well of detection antibody solution to the plate(s).
- Incubate the plate(s) at room temperature with shaking for 1 hour.
- Wash the plate(s) three times with 300 µL/well of 1X Tris Wash Buffer.

STEP 4 Read Plate

- Add 150 µL/well of 1X Read Buffer T to the plate(s).
- Incubate the plate(s) at room temperature (NO SHAKING) for 10 minutes.
- Analyze the plate(s) with a SECTOR® Imager instrument.

Notes:

The sAPP calibrators can be diluted in a solution of 1% Blocker A in 1X Tris Wash Buffer. If the calibration curve will be used for quantification of proteins in a complex matrix (culture supernatant, serum, CSF, etc.) a different diluent may be desired.

The pH changes that occur in a culture medium are detrimental to this assay, and it is recommended that culture medium samples be supplemented with HEPES buffer, pH 7.3 at a final concentration of 50 mM. Other matrices should be examined for pH effects also, supplemented with HEPES buffer as needed.

It is recommended that calibrators and samples be assayed in duplicate.

The sAPP α /sAPP β assay is sensitive to the use of denaturing reagents and to the heat generated during sonification or homogenization.

Shaking the plate accelerates analyte capture.

Bubbles in the read buffer will interfere with reliable imaging of the plate if carried into the wells.

The incubation in read buffer is essential for this assay.

The necessity of the incubation in read buffer may vary for different matrices.

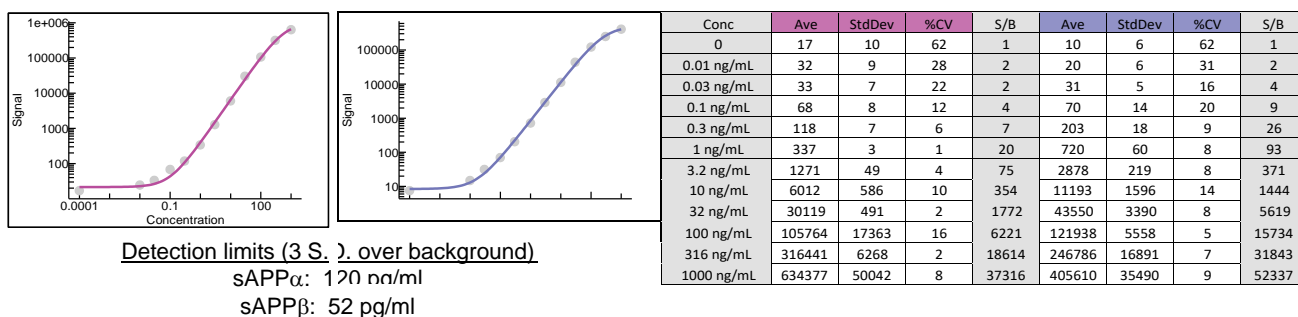
sAPP Calibrators

Recombinant Human sAPP α and sAPP β

Contents:	750 ng recombinant sAPP α and sAPP β proteins
Concentration:	50 μ g/mL
Volume:	15 μ L
Preparation:	Recombinant human sAPP proteins were purified from overexpressing mammalian cells.
Storage:	Store at ≤ -70 °C.
Quality Control:	Recombinant proteins have been analyzed by SDS-PAGE and MSD MULTI-SPOT Assays.

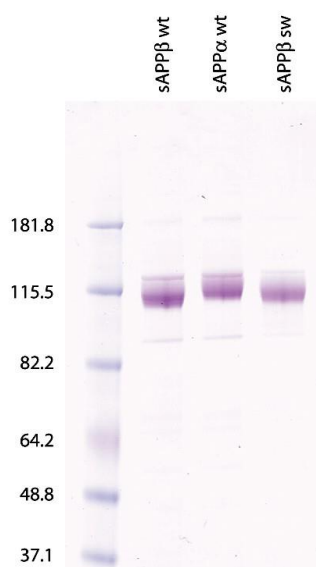
MSD MULTI-SPOT Assay Results

Typical titration curve for recombinant sAPP proteins using the MSD MULTI-SPOT sAPP α / β duplex assay.



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%).



MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, mesoscale.com, www.mesoscale.com, methodicalmind.com, www.methodicalmind.com, DISCOVERY WORKBENCH, InstrumentLink, MESO, MesoSphere, Methodical Mind, MSD GOLD, MULTI-ARRAY, MULTI-SPOT, QuickPlex, ProductLink, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, TeamLink, TrueSensitivity, TURBO-BOOST, TURBO-TAG, N-PLEX, R-PLEX, S-PLEX, T-PLEX, U-PLEX, V-PLEX, MSD (design), MSD (luminous design), Methodical Mind (design), 96 WELL SMALL-SPOT (design), 96 WELL 1-, 4-, 7-, 9-, & 10-SPOT (designs), 384 WELL 1- & 4-SPOT (designs), N-PLEX (design), R-PLEX (design), S-PLEX (design), T-PLEX (design), U-PLEX (design), V-PLEX (design), It's All About U, SPOT THE DIFFERENCE, The Biomarker Company, and The Methodical Mind Experience are trademarks and/or service marks owned by or licensed to Meso Scale Diagnostics, LLC. All other trademarks and service marks are the property of their respective owners..

©2012, 2020 Meso Scale Diagnostics, LLC. All rights reserved.