

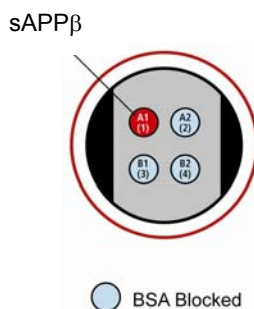
MSD[®] 96-Well MULTI-ARRAY[®] Mouse/Rat sAPP β Assay

This protocol has been optimized for detection of soluble amyloid precursor protein β in mouse and rat cerebrospinal fluid.

Storage

MSD Materials

| | |
|---|---------------|
| <input type="checkbox"/> Read Buffer T (with surfactant), 4X | RT |
| <input type="checkbox"/> Blocker A | RT |
| <input type="checkbox"/> 10% Sodium Dodecyl Sulfate (SDS) | RT |
| <input type="checkbox"/> Nonidet P-40 (NP-40) | RT |
| <input type="checkbox"/> MULTI-SPOT [®] 96-well 4 Spot sAPP β plate(s) | 2-8°C |
| <input type="checkbox"/> Tris Wash Buffer (10X) | 2-8°C |
| <input type="checkbox"/> SULFO-TAG [™] Anti-APP (22C11) Antibody (50X) | 2-8°C |
| <input type="checkbox"/> 1M Dithiothreitol (DTT) | -20°C |
| <input type="checkbox"/> GF1 Assay Diluent | -20°C |
| <input type="checkbox"/> sAPP β Calibrator (50 μ g/mL) | \leq -70 °C |



The SECTOR[®] Imager data file will identify spots according to their well location, not by the coated capture antibody name.

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



Notes:

Other Materials & Equipment (not supplied)

- Deionized water for diluting Wash Buffer and Read Buffer
- Automated plate washer, Multidrop[®], or other efficient multi-channel pipetting equipment for washing 96-well plates
- Appropriate liquid handling equipment for desired throughput that must accurately dispense 10 – 150 μ L into a 96-well micro plate
- Microcentrifuge tubes (e.g. VWR catalog# 20170-662)
- Adhesive Plate Seals

Protocol at a Glance

The following protocol is optimized for quantifying sAPP β in rodent cerebrospinal fluid. The protocol can be completed in approximately 3 to 3½ hours if each reagent is prepared during the preceding blocking or incubation step.

Read the entire detailed instructions before beginning work.

Step 1. Add Blocking solution, incubate 1 hour, wash.

Step 2. Add 25 μ L of Denaturing Assay Solution
Add 25 μ L of samples or Calibrator, incubate 1 hour, wash.

Step 3. Add 25 μ L of Detection Antibody, incubate 1 hour, wash

Step 4. Add 150 μ L of Read Buffer and analyze plate.

Detailed Instructions

Prepare a stock of 1X Tris Wash Buffer:

- a) 1X Tris Wash Buffer will be used throughout the assay to make other reagents as well as to wash plates. Approximately 250 mL per plate is required – more if using an automatic plate washer.
- b) In a 250 mL bottle combine:
 - 25 mL 10X Tris Wash Buffer
 - 225 mL deionized water

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

Prepare Blocker A Solution:

- a) Prepare 20 mL per plate.
- b) In a 50 mL tube combine:
 - 20 mL 1X Tris Wash Buffer
 - 600 mg Blocker A

Solutions containing Blocker A should be kept at 4°C and discarded after 14 days.



Notes:

Prepare Denaturing Assay Solution:

- a) Prepare 6 mL per plate.
- b) In a 15 mL tube combine:
 - 5730 μ L GF1 Assay Diluent
 - 120 μ L 10% SDS
 - 120 μ L NP-40
 - 30 μ L 1M DTT

Denaturing Assay Solution should be made immediately prior to use, kept at 4°C and discarded after use.

Prepare Antibody Dilution Buffer:

- a) Prepare 3 mL per plate.
- b) In a 15 mL tube combine:
 - 1 mL Blocker Solution A
 - 2 mL 1X Tris Wash Buffer

Prepare sAPP β Calibrators:

- a) Prepare highest Calibrator by adding 10 μ L of sAPP β (50 μ g/mL) to 490 μ L of Denaturing Assay Solution.
- b) Prepare the next Calibrator by transferring 100 μ L of the highest Calibrator to 200 μ L of Denaturing Assay Solution.
- c) Repeat 3-fold serial dilutions 5 additional times to generate 7 Calibrators.
- d) This yields the following Calibrator concentrations:

Calibrator dilutions should be kept at 4°C and discarded after use.

| <u>Calibrator</u> | <u>sAPPβ (ng/mL)</u> |
|-------------------|---------------------------------------|
| Cal 7 | 1,000 |
| Cal 6 | 333 |
| Cal 5 | 111 |
| Cal 4 | 37 |
| Cal 3 | 12 |
| Cal 2 | 4.1 |
| Cal 1 | 1.3 |
| Cal 0 | 0.0 |

- e) Use Denaturing Assay Solution for Cal 0. These Calibrators will be sufficient to run an 8-point calibration curve in triplicate for one plate. Do not store diluted Calibrators.



Prepare the 1X Detection Antibody Solution:

Notes:

- a) In a 15 mL tube combine:
 - 60 μ L of 50X SULFO-TAG Anti-APP (22C11) Antibody
 - 2940 μ L of Antibody Dilution Buffer
- b) This will yield 3 mL of diluted Detection Antibody Solution at the working concentration with sufficient volume for one plate.

Dilute Read Buffer:

- In a 50 mL tube, combine (per plate):
- 10 mL 4X Read Buffer T
 - 10 mL deionized water

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

STEP 1 ***Block Plate:***

- a) Add 150 μ L/well of Blocker A Solution.
- b) Incubate with shaking for 1 hour at room temperature.
- c) Wash plate three times with 1X Tris Wash Buffer.

STEP 2 ***Start the Incubation:***

- a) Add 25 μ L/well of Denaturing Assay Solution.
- b) Immediately add 25 μ L/well of Calibrator or sample.
- c) Incubate with shaking for 1 hour at room temperature.
- d) Wash plate three times with 1X Tris Wash Buffer.

STEP 3 ***Detection Antibody Addition:***

- a) Add 25 μ L/well of 1X Detection Antibody Solution.
- b) Incubate with shaking for 1 hour at room temperature.
- c) Wash plate three times with 1X Tris Wash Buffer.

STEP 4 ***Read Plate:***

- a) Add 150 μ L/well of 2X Read Buffer T.
- b) Read plate on SECTOR Imager immediately after Read Buffer addition and analyze data.

Avoid bubbles while adding the Read Buffer; it will interfere with accurate reading of the plate.

