

Storage

## MSD<sup>®</sup> MULTI-SPOT Human (6E10) Abeta Triplex Assay

| Materials | Included |
|-----------|----------|
| materials | monaucu  |

| MULTI-SPOT <sup>®</sup> Abeta Peptide 3-Plex Plate        | 2–8°C  |
|---|--------|
| SULFO-TAG <sup>TM</sup> 6E10 Detection Antibody $(50X)^1$ | 2–8°C  |
| Aβ1-38 Peptide  | ≤-70°C |
| Aβ1-40 Peptide  | ≤-70°C |
| Aβ1-42 Peptide  | ≤-70°C |
| Tris Wash Buffer (10X)                                    | 2–8°C  |
| Blocker G (100X)  | 2–8°C  |
| Blocker A (dry powder)                                    | 2–8°C  |
| Read Buffer T (4X)  | RT     |



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

#### Safety

Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling kit components. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines. Additional safety information is available in the product safety data sheet, which can be obtained from MSD Customer Service.

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



<sup>&</sup>lt;sup>1</sup> SULFO-TAG conjugated detection antibodies should be stored in the dark.

#### Other Materials & Equipment (not supplied)

- Deionized water for diluting concentrated buffers
- $\square \quad 500 \text{ mL bottle}$
- $\square \quad 50 \text{ mL tubes}$
- $\square 15 \text{ mL tubes}$
- □ Adhesive plate seals
- □ Microtiter plate shaker
- Various microcentrifuge tubes for making serial dilutions of lysates (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 μL and 150 μL into a 96-well microplate

### Protocol at a Glance

- 1. Add Blocker A solution; incubate 1 hour; wash.
- 2. Add detection antibody.
- 3. Add samples or calibrators; incubate 2 hours; wash.
- 4. Add Read Buffer T and analyze plate.

This protocol can be completed in approximately 3 to 3 1/2 hours if each reagent is prepared during the preceding incubation. Alternatively, all reagents with the exception of diluted peptides can be prepared ahead of time. This lengthens the overall time required for the assay but frees up time during incubation steps.

#### **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 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
- $\square$  315 mL deionized water

Prepare 1% Blocker A Solution. You will need 50 mL per plate.

In a 50 mL tube, combine:

- □ 50 mL 1X Tris Wash Buffer
- □ 500 mg Blocker A



*Read the entire detailed instructions before beginning work.* 

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

Notes:

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



# **Prepare Detection Antibody Solution.** You will need 3.0 mL per plate at a 1X final concentration.

In a 15 mL tube, combine:

- **Ο** 60 μL 50X SULFO-TAG 6E10 Detection Antibody
- □ 30 µL 100X Blocker G
- **□** 2910 μL 1% Blocker A Solution

*NOTE:* For serum and plasma samples Blocker G should be omitted from the Detection Antibody Solution

**Prepare Read Buffer T.** You will need 20 mL per plate at a final 2X concentration.

In a 50 mL tube, combine:

- □ 10 mL 4X Read Buffer T
- □ 10 mL deionized water

| <b>Prepare Aβ1-38, Aβ1-40, and Aβ1-42 Peptide Standards.</b> MSD              |
|---|
| supplies individual vials of $A\beta$ peptide stock calibrators. For the lot- |
| specific concentrations of each calibrator, refer to the certificate of       |
| analysis (C of A) supplied with the kit. You can also find a copy of the C    |

of A at www.mesoscale.com.

To prepare 7 calibrator solutions plus a zero calibrator for up to 4 replicates:

- a) Thaw the stock calibrators at room temperature and mix well by vortexing.
- b) Prepare individual diluted stocks of the A $\beta$ 1-38 and A $\beta$ 1-40 peptides as follows:
  - 20 μL of the supplied Aβ1-38 or Aβ1-40 stock calibrator
    180 μL of 1% Blocker A Solution.

Mix well by vortexing.

- c) Combine the following to prepare the highest calibrator:
  - $\Box$  22.5 µL of pre-diluted A $\beta$ 1-38 Peptide
  - **\Box** 50 µL of pre-diluted A $\beta$ 1-40 Peptide
  - $\square \quad 15 \ \mu L \ of \ supplied \ A\beta 1-42 \ Peptide$

- d) Prepare the next calibrator by transferring 100  $\mu$ L of the highest calibrator to 200  $\mu$ L of 1% Blocker A Solution. Mix well by vortexing. Repeat 3-fold serial dilutions 5 additional times to generate 7 calibrators.
- e) Use 1% Blocker A Solution as the zero calibrator.



#### Notes:

*Excess diluted read buffer may be stored in a tightly sealed container at room temperature.* 

To avoid the possibility of aggregation and/or sticking of the peptides to the dilution tubes, the dilutions should be prepared immediately before use.

Thorough mixing of stock and diluted kit reagents is required.

Do not store diluted calibrators.

Both peptide standards and samples should be assayed in duplicate.



|        | Begin with a MULTI-SPOT 96-well 4-Spot Abeta Peptide 3-Plex Plate.<br>No pre-treatment is necessary.                  | Notes:   |
|--------|---|--|
| STEP 1 | Add 150 µL/well of 1% Blocker A Solution.   |  |
|        | <b>Incubate</b> with shaking at room temperature for 1 hour. During this time, prepare peptide standards and samples. |  |
| STEP 2 | <b>Wash</b> plate(s) four times with at least 150 $\mu$ L/well of 1X Tris Wash Buffer.                                | CSF samples should be assayed<br>'neat' or diluted no more than                                    |
|        | Add 25 $\mu$ L/well of detection antibody solution.   | 2-fold for optimal peptide<br>sensitivity  |
|        | Add 25 $\mu$ L/well of the samples prepared during Step 1 incubation.   | Shaking the plate accelerates analyte capture.   |
|        | <b>Incubate</b> with shaking for 2 hours at room temperature.   |  |
| STEP 3 | Wash plate(s) four times with at least 150 $\mu$ L/well of 1X Tris Wash Buffer  | Add read buffer carefully using<br>reverse pipetting technique.<br>Bubbles in the read buffer will |
|        | Add 150 $\mu$ L/well of diluted Read Buffer T.  | interfere with reliable imaging<br>of the plate if carried into the<br>wells.                      |
|        | Analyze with MSD instrument.  |  |
|        |   | Read plate(s) immediately after adding read buffer.  |

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, V-PLEX, STREPTAVIDIN GOLD, MESO, www.mesoscale.com, SMALL SPOT (design), 96 WELL 1, 4, 7, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC. The 6E10 antibody used in MSD Aβ assays is supplied by Covance Research Products, Inc ©2013 Meso Scale Diagnostics, LLC. All rights reserved.

