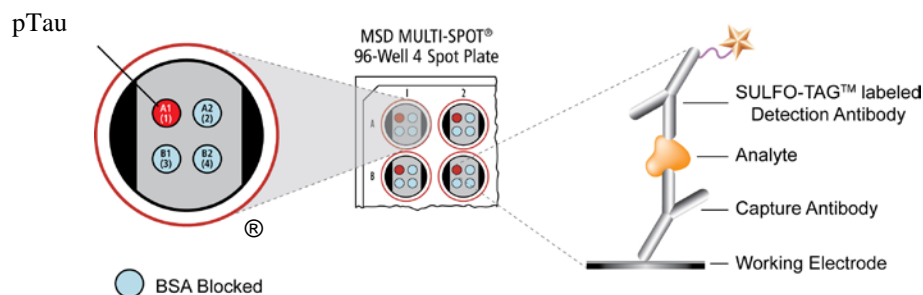


MSD[®] MULTI-SPOT Phospho-Tau (Thr 231) Assay

Storage

Materials Included

❑ MULTI-SPOT [®] 96-Well 4-Spot Phospho-Tau (Thr231) Plate(s)	2–8°C
❑ SULFO-TAG [™] Anti-Total Tau Antibody ¹	2–8°C
❑ Tau441 Calibrator (phosphorylated)	≤-70°C
❑ Tris Wash Buffer (10X)	2–8°C
❑ Blocker D–B (10%) ²	≤-10°C
❑ Blocker A (dry powder)	RT
❑ 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.

¹ SULFO-TAG conjugated detection antibodies should be stored in the dark.

² Blocker D–B can tolerate at least 5 freeze–thaw cycles. Alternatively, aliquots of the blocker can be stored at 2–8°C for up to 1 month.

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



Other Materials & Equipment (not supplied)

Notes:

- 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 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 μ L and 150 μ L into a 96-well microplate

Protocol at a Glance

Read the entire detailed instructions before beginning work.

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

The 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 lysates 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.

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

In a 500 mL bottle, combine:

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

Prepare 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

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



Prepare antibody dilution buffer. You will need 3.0 mL per plate.

In a 15 mL tube, combine:

- ❑ 1 mL Blocker A Solution
- ❑ 1.97 mL 1X Tris Wash Buffer
- ❑ 30 μ L Blocker D–B (10%)

Prepare detection antibody solution. You will need 3.0 mL per plate at a 10 nM final concentration.

In a 15 mL tube, combine:

- ❑ 91 μ L Anti-Total Tau Antibody
- ❑ 2909 μ L cold antibody dilution buffer

Prepare Read Buffer T. You will need 20 mL per plate at a final 1X concentration.

In a 50 mL tube, combine:

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

Prepare dilutions of Tau441 Calibrator (phosphorylated). To prepare 7 calibrator solutions plus a zero calibrator for up to 4 replicates:

- a) Thaw the stock calibrator and mix well by vortexing.
- b) Prepare the highest calibrator by diluting the supplied calibrator with your selected assay diluent.
 - a. Add 10 μ L of Tau441 Calibrator to 490 μ L of assay diluent.
 - b. Mix well by vortexing.
- c) Prepare the next calibrator by transferring 100 μ L of the highest calibrator to 300 μ L of assay diluent. Mix well by vortexing. Repeat 4-fold dilutions 5 additional times to generate 7 calibrators.
- d) Use assay diluent 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 to the dilution tubes, the dilutions should be prepared immediately before use.

The calibrator and all diluted samples should be prepared in a diluent that mimics the sample matrix as closely as possible (e.g., cell culture medium, lysis buffer, immunodepleted CSF, etc...).

The diluent used must contain sufficient protein to prevent non-specific sticking of tau to the assay well. In the absence of an optimized diluent, 10% MSD Blocker A in 1X Tris Wash Buffer is recommended.

Samples derived from biological fluids and/or tissues may require independent manipulations not described here.



This yields the following calibrator concentrations:

Calibrator	Phospho-Tau (Thr231) (units/well)	Dilution Factor
Calibrator-01	50	
Calibrator-02	13	4
Calibrator-03	3.1	4
Calibrator-04	0.78	4
Calibrator-05	0.20	4
Calibrator-06	0.049	4
Calibrator-07	0.012	4
Calibrator-08	0	n/a

Note: The source vial of calibrator specifies the units of phosphorylated tau (Thr231) per μg of total tau.

Begin with a MULTI-SPOT 96-Well 4-Spot Phospho-Tau (Thr231) Plate. No pre-treatment is necessary.

Notes:

STEP 1 Add 150 μL /well of Blocker A Solution.

Incubate with shaking at room temperature for 1 hour. During this time, prepare calibration curve and samples.

Shaking the plate accelerates analyte capture.

STEP 2 Wash plate(s) four times with at least 150 μL /well of 1X Tris Wash Buffer.

Add 25 μL /well of calibrator or samples prepared during Step 1 incubation.

Incubate with shaking for 1 hour at room temperature.

STEP 3 Wash plate(s) four times with at least 150 μL /well of 1X Tris Wash Buffer

Add 25 μL /well of detection antibody solution.

Incubate with shaking for 1 hour at room temperature.

STEP 4 Wash plate(s) four times with at least 150 μL /well of 1X Tris Wash Buffer.

Add 150 μL /well of diluted Read Buffer T.

Analyze with MSD instrument.

Add read buffer carefully using reverse pipetting technique. Bubbles in the read buffer will interfere with reliable imaging of the plate if carried into the wells.

Read plate(s) immediately after adding read buffer.