

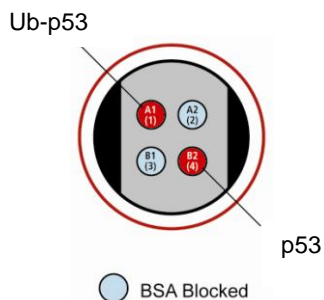
MSD[®] 96-Well MULTI-SPOT[®] Ubiquitinated, Total p53 Assay

The following assay protocol has been optimized for quantifying ubiquitinated and total p53 in whole cell lysate.

Storage

Materials Included

<input type="checkbox"/> Read Buffer T (with surfactant), 4X	RT
<input type="checkbox"/> Blocker A	RT
<input type="checkbox"/> Blocker B	RT
<input type="checkbox"/> EDTA (0.5M, pH 8.0)	RT
<input type="checkbox"/> MULTI-SPOT 96-well 4 Spot Ub-, Total p53 Plate(s)	2-8°C
<input type="checkbox"/> SULFO-TAG [™] Anti-Total p53 (50X)	2-8°C
<input type="checkbox"/> Tris Wash Buffer (10X)	2-8°C
<input type="checkbox"/> Tris Lysis Buffer (1X)	2-8°C
<input type="checkbox"/> Phosphatase Inhibitor I (100X)	2-8°C
<input type="checkbox"/> Phosphatase Inhibitor II (100X)	2-8°C
<input type="checkbox"/> Blocker D-R (10%)	≤-10°C ¹
<input type="checkbox"/> Protease Inhibitor Solution(50X)	≤-10°C



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

Meso Scale Discovery, Meso Scale Diagnostics, www.mesoscale.com, MSD, MSD (design), Discovery Workbench, Quickplex, Multi-Array, Multi-Spot, Sulfo-Tag and Sector are trademarks of Meso Scale Diagnostics, LLC.
© 2009 Meso Scale Discovery a division of Meso Scale Diagnostics, LLC. All rights reserved.

¹ Blocker D-R can tolerate up to 5 freeze-thaw cycles. Alternatively, an aliquot of blocker D-R can be stored at 2-8°C for up to 1 month.

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
- One 250 mL bottle
- Two 50 mL tube
- One 15 mL tube
- 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 micro plate

Read the entire detailed instructions before beginning work.

Protocol at a Glance

The following protocol describes the most conservative approach toward achieving highly sensitive results using MSD technology to quantify phosphoproteins. The protocol can be completed in approximately 3 to 3 1/2 hours if each reagent is prepared during the preceding incubation. All reagents with the exception of diluted lysates can also be prepared ahead of time. This lengthens the overall time required for the assay but frees up time during incubation steps.

Once desired results are achieved, the protocol can be streamlined to eliminate multiple incubation and wash steps to increase throughput.

1. Add Blocking Solution, incubate 1 hour, wash.
2. Add lysate, incubate 1 hour at room temperature, wash.
3. Add Detection Antibody, incubate 1 hour, wash.
4. Add 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 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 Tris Wash Buffer may be prepared at once and stored at room temperature for later use.



Notes:

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

Prepare Blocking Solution-A:

- a) Prepare 20 mL per plate.
- b) In a 50 ml tube combine:
 - 20 mL 1X Tris Wash Buffer
 - Blocker A 600mg (3%)

Prepare Antibody Dilution Buffer:

- a) Prepare 3 mL per plate.
- b) In a 15 mL tube combine:
 - 1 mL Blocking Solution-A
 - 1.97 mL 1X Tris Wash Buffer
 - 30 µL 10% Blocker D-R

Begin with a MULTI-SPOT 96-well 4 Spot Ub-, Total p53 Plate.
No pre-treatment is necessary.

STEP 1

Add 150 µL/well of Blocking Solution.

Plates may also be blocked overnight at 4°C.

Incubate with shaking at room temperature for 1 hour. Prepare Complete Tris Lysis Buffer, and dilute cell lysates during this time.

Prepare Complete Tris Lysis Buffer:

- a) To 10 mL of Tris Lysis Buffer, add the following:
 - 100 µL Phosphatase Inhibitor I (100X stock)
 - 100 µL Phosphatase Inhibitor II (100X stock)
 - 200 µL Protease Inhibitor Solution (50X stock)
 - 400 µL 0.5M EDTA, pH 8.0
- b) Keep Complete Tris Lysis Buffer on ice until use.

Complete Tris Lysis Buffer should be made each day of experimentation.

Prepare 5% Blocking Solution-B:

- a) Prepare 1 mL 5% Blocking Solution-B. Combine the following:
 - 50 mg Blocker B
 - 1 mL Complete Lysis Buffer
- b) Mix well to completely dissolve Blocker B.
- c) Keep 5% Blocking Solution-B on ice until use.

Prepare samples or positive and negative cell lysates:

(Note: Recommendations for cell lysate handling are provided; however, the suggested concentrations listed below may need to be adjusted depending upon specific samples tested.)



Notes:

- a) Thaw cell lysate samples on ice and dilute immediately before use. Keep on ice during all manipulations and discard all remaining thawed unused material.
- b) Dilute lysates and 5% Blocking Solution-B together in Complete Lysis Buffer to prepare a sample containing 0.1 µg/µL of lysate (2.5 µg/well in 25 µL) and 0.5% Blocking Solution-B.
- c) *For example*, to prepare 200 µL of 0.1 µg/µL lysate-0.5% Blocking Solution-B, combine the following: 20 µL 5% Blocking Solution-B, 10 µL 2 µg/µL lysate and 170 µL Complete Lysis Buffer.
- d) Additional dilutions may be prepared from the 0.1 µg/µL lysate sample if desired. Use a stock of Complete Lysis Buffer-0.5% Blocking Solution-B (dilute 5% Blocking Solution-B 1:10 in Complete Lysis Buffer) for any further lysate dilutions.

Wash plates three times with Wash Buffer.

STEP 2

Dispense 25 µL/well of diluted lysates.

Incubate with shaking for 1 hour at room temperature. Prepare SULFO-TAG Anti-Total p53 Detection Antibody Solution during this time.

Shaking a 96-well MSD MULTI-ARRAY[®] or MULTI-SPOT plate accelerates capture at the working electrode.

Prepare Detection Antibody Solution:

- a) Prepare 3.0 mL per plate.
- b) In a 15 mL tube combine:
 - 2.94 mL cold Antibody Dilution Buffer
 - 60 µL 50X Anti-Total p53 Antibody (Final concentration: 1X)

Wash plates three times with Wash Buffer.

STEP 3

Add 25 µL/well of Detection Antibody Solution.

Incubate with shaking at room temperature for 1 hour. Prepare Read Buffer during this time.

Dilute Read Buffer:

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

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

Wash plates three times with Wash Buffer.

Bubbles in the Read Buffer will interfere with reliable imaging of the plate if carried into the wells.

STEP 4

Add 150 µL/well of diluted Read Buffer T (with surfactant).

Analyze with SECTOR Imager plate reader.

Plates can be imaged immediately following the addition of Read Buffer.

