

# MSD<sup>®</sup> 384-Well MULTI-ARRAY<sup>®</sup> Total HIF-1 $\alpha$ Assay

Storage

## MSD Materials

<input type="checkbox"/> MULTI-SPOT <sup>®</sup> 384-well HIF-1 $\alpha$ Plate	2-8°C
<input type="checkbox"/> SULFO-TAG <sup>™</sup> Anti-Total HIF-1 $\alpha$ Antibody	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"/> Protease Inhibitor Solution (50X)	$\leq -10^{\circ}\text{C}$
<input type="checkbox"/> Blocker A	RT
<input type="checkbox"/> Read Buffer T , 4X (with surfactant)	RT

## Other Materials & Equipment (not supplied)

- ☐ Deionized water for diluting Wash Buffer and Read Buffer
- ☐ One 1 L bottle
- ☐ Two 50 mL tube
- ☐ One 15 mL tube
- ☐ Various microcentrifuge tubes for making serial dilutions of lysates (if desired)
- ☐ Adhesive plate seals
- ☐ Microtiter plate shaker
- ☐ Automated plate washer or other equipment for washing 384-well plates
- ☐ Appropriate liquid handling equipment for desired throughput that must accurately dispense 10  $\mu\text{L}$  and 35  $\mu\text{L}$  into a 384-well micro plate

Meso Scale Discovery, Meso Scale Diagnostics, [www.mesoscale.com](http://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.

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



## Protocol at a Glance

The following protocol describes the most conservative approach toward achieving highly sensitive results using MSD technology to quantify HIF-1 $\alpha$ . The protocol takes approximately 5 to 5 1/2 hours to complete 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 sample or lysate, incubate 2 hours, wash.
3. Add Detection Antibody, incubate 2 hours, 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 1 L per plate is required – more if using an automatic plate washer.
- b) In a 1 L bottle combine:
  - ☐ 100 mL 10X Tris Wash Buffer
  - ☐ 900 mL deionized water

### Prepare Blocking Solution-A:

- a) Prepare 20 mL per plate.
- b) In a 50 mL tube combine:
  - ☐ 20 mL 1X Tris Wash Buffer
  - ☐ 600 mg Blocker A (30 mg/mL or 3%)

### Prepare Antibody Dilution Buffer:

- a) Prepare 8 mL per plate.
- b) In a 15 mL tube combine:
  - ☐ 2.67 mL Blocking Solution A
  - ☐ 5.33 mL 1X Tris Wash Buffer

Begin with a MULTI-SPOT Custom Plate. No pre-treatment is necessary.

## Notes:

*Read the entire detailed instructions before beginning work.*

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

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



## STEP 1

Add 35  $\mu$ L/well of Blocking Solution-A.

Incubate with shaking at room temperature for 1 hour. Prepare Complete Lysis buffer, and prepare samples or dilute cell lysates during this time.

### ***Prepare Complete Lysis Buffer:***

- a) To 10 mL of Tris Lysis Buffer, add the following:
  - ☐ 200  $\mu$ L Protease Inhibitor Solution (50X stock)
- b) Keep Complete Lysis Buffer 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.)*

- 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 cell lysate in Complete Lysis Buffer to a final concentration of 0.5  $\mu$ g/ $\mu$ L. This will deliver 5  $\mu$ g/well in 10  $\mu$ L. A dilution series may also be prepared if desired.
- c) Keep diluted cell lysate on ice until use.

Wash plates four times with Tris Wash Buffer.

## STEP 2

Dispense 10  $\mu$ L/well of samples or diluted lysates.

Incubate with shaking at room temperature for 2 hours. Prepare Detection Antibody Cocktail during this time.

### ***Prepare Detection Antibody:***

- a) Dilute SULFO-TAG Anti-Total HIF-1 $\alpha$  Antibody to a final concentration of 10 nM.
- b) Use cold Antibody Dilution Buffer. Sufficient antibody is supplied to prepare 8 mL per plate.

Wash plates four times with Tris Wash Buffer.

## STEP 3

Add 10  $\mu$ L/well of Detection Antibody.

Incubate with shaking at room temperature for 2 hours. 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

Wash plates four times with Wash Buffer

## STEP 4

Add 35  $\mu$ L/well of diluted Read Buffer T (with surfactant).

Analyze with SECTOR<sup>®</sup> instrument.

