# MSD®MULTI-ARRAY Assay System

### Phospho-Histone H3 (Ser10) Assay Base Kit

20-Plate Kit

K150EWA-3



www.mesoscale.com®

### MSD Phosphoprotein Assays

### Phospho-Histone H3 (Ser10) Assay Base Kit

This package insert must be read in its entirety before using this product.

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

MESO SCALE DISCOVERY<sup>®</sup> A division of Meso Scale Diagnostics, LLC. 1601 Research Boulevard Rockville, MD 20850-3173 USA www.mesoscale.com

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, MSD GOLD, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, U PLEX, S-PLEX, 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), U-PLEX (design), S-PLEX (design), V PLEX (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.

All other trademarks or service marks are the property of their respective owners.

# Table of Contents

Introduction	4
Principle of the Assay	4
Reagents Supplied	5
Required Material and Equipment – not supplied	5
Optional Material – not supplied	5
Safety	6
Reagent Preparation	6
Sample Preparation and Storage	8
Assay Protocol	9
Typical Data	11
Assay Components	11
Limitations of the Procedure	12
References	12
Appendix	13
Summary Protocol	14
Plate Diagrams	15

### **Contact Information**

**MSD** Customer Service

Phone: 1-240-314-2795 Fax: 1-301-990-2776 Email: CustomerService@mesoscale.com

### MSD Scientific Support

Phone:1-240-314-2798Fax:1-240-632-2219 attn: Scientific SupportEmail:ScientificSupport@mesoscale.com

### Introduction

**Histone H3**, along with H2a, H2b, H1/H5, and H4, is one of the five main classes of histones, which are found in the nucleus of eukaryotic cells, and are the main building blocks of chromatin. Histone H3 has an amino terminal tail and post-translational modification of this tail—an epigenetic modification—is involved in chromatin relaxation and transcriptional regulation.<sup>1</sup> Histones undergo many different types of post translational modifications, such as acetylation, methylation, phosphorylation, ubiquitylation, and sumoylation.<sup>2</sup> These modifications can change the chromatin structure and allow for binding of transcription factors to the DNA.<sup>3</sup>

Acetylation of lysine residues on histones is frequently associated with actively transcribed regions of the genome, whereas methylation of lysine is seen both at transcribed and transcriptionally repressed areas of the genome.<sup>4</sup> Epigenetic modifications, of which histone post-translational modification is one example, have been shown to play a role in cancer, embryonic stem cell function, and mammalian development.<sup>5</sup>

### Principle of the Assay

MSD phosphoprotein assays provide a rapid and convenient method for measuring the levels of protein targets within a single, small-volume sample. The assays are available in both singleplex and multiplex formats. In a singleplex assay, an antibody for a specific protein target is coated on one electrode (or "spot") per well. In a multiplex assay, an array of capture antibodies against different targets is patterned on distinct spots in the same well. The Phospho-Histone H3 (Ser10) Assay is a sandwich immunoassay (Figure 1). MSD provides a plate that has been pre-coated with capture antibody for total Histone H3. The user adds the sample and a solution containing the detection antibody—anti-phospho-Histone H3 (Ser10) conjugated with an electrochemiluminescent compound, MSD SULFO-TAG<sup>™</sup> label—over the course of one or more incubation periods. Analyte in the sample binds to the capture antibody immobilized on the working electrode surface; recruitment of the conjugated detection antibody by bound analyte completes the sandwich. The user adds an MSD read buffer that provides the appropriate chemical environment for electrochemiluminescence and loads the plate into an MSD instrument for analysis. Inside the MSD instrument, a voltage applied to the plate electrodes causes the labels bound to the electrode surface to emit light. The instrument measures intensity of emitted light to provide a quantitative measure of phosphorylated Histone H3 present in the sample.



**Figure 1**. Spot diagram showing placement of analyte capture antibody. The numbering convention for the different spots is maintained in the software visualization tools, on the plate packaging, and in the data files. A unique bar code label on each plate allows complete traceability back to MSD manufacturing records.

Spot the Difference

### **Reagents Supplied**

Product Description	Storage	Quantity per Kit K150EWA-3
MULTI-SPOT 96-Well 4-Spot Histone H3 Plate(s) N450EWA-1	2–8°C	20 plates
SULFO-TAG Anti-Phospho-Histone H3 (Ser10) Antibody <sup>1</sup> (50X)	2-8°C	4 vials (375 μL ea)
Read Buffer T (4X) R92TC-3 (50 mL), R92TC-2 (200 mL)	RT	1 bottle (200 mL)

### Additional Materials and Equipment

- Appropriately sized tubes and bottles for reagent preparation
- Polypropylene microcentrifuge tubes for preparing dilutions
- Liquid handling equipment suitable for dispensing 10 to 150 µL/well into a 96-well microtiter plate
- Plate washing equipment: automated plate washer or multichannel pipette
- Microtiter plate shaker (rotary) capable of shaking at 500–1,000 rpm
- Adhesive plate seals
- Deionized water
- Vortex mixer

### Optional Material — not supplied

Phosphoprotein Reagent Support Pack (K0000D-3)

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

## Safety

Use safe laboratory practices: wear gloves, safety glasses, and lab coats when handling assay components. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines.

Additional product-specific safety information is available in the safety data sheet (SDS), which can be obtained from MSD Customer Service or at <u>www.mesoscale.com</u>.

### **Reagent Preparation**

#### Note

The instructions below will prepare the reagents needed as described in the assay protocol. All supplemental reagents (inhibitors, buffers, and blocking reagents) are available for purchase in the MSD Phosphoprotein Reagent Support Pack, or alternatively can be purchased and prepared separately by the end user.

#### Prepare Tris Wash Buffer

Dilute 10X stock of Tris Wash Buffer to 1X as shown below. Tris Wash Buffer (1X) will be used throughout the assay to make additional reagents and wash plates. Approximately 350 mL per plate is required—more if using an automatic plate washer.

For one plate, combine:

- □ 35 mL of Tris Wash Buffer (10X)
- □ 315 mL deionized water

Excess Tris Wash Buffer may be stored at room temperature in a tightly sealed container for later use.

#### **Prepare Blocking Solution**

For one plate, combine:

- □ 600 mg Blocker A (dry powder)
- 20 mL 1X Tris Wash Buffer

#### **Prepare Antibody Dilution Buffer**

For one plate, combine:

- **Ο** 30 μL 10% Blocker D-R
- □ 1 mL blocking solution
- □ 1.97 mL 1X Tris Wash Buffer

Set aside on ice.



#### Prepare Complete Lysis Buffer

To 10 mL of Tris Lysis Buffer, add the following supplemental materials to prepare the complete lysis buffer (sufficient for 2-3 plates):

- □ 100 µL Protease Inhibitor Solution (100X stock)
- □ 100 µL Phosphatase Inhibitor Solution I (100X stock)
- □ 100 µL Phosphatase Inhibitor Solution II (100X stock)

The complete lysis buffer should be ice cold before use.

#### **Prepare Detection Antibody Solution**

For one plate, combine:

- □ 2.94 mL antibody dilution buffer
- □ 60 µL 50X SULFO-TAG Anti-Phospho-Histone H3 (Ser10) Antibody (1X final concentration)

#### **Prepare Read Buffer T**

For one plate, combine:

- □ 5.0 mL Read Buffer T (4X)
- □ 15 mL deionized water

Diluted read buffer (1X) may be stored at room temperature in a tightly sealed container for later use.

#### **Prepare MSD Plate**

This plate has been pre-coated with antibody for the analyte shown in Figure 1. The plate can be used as delivered; no additional preparation (e.g., pre-wetting) is required. The plate has also been exposed to a proprietary stabilizing treatment to ensure the integrity and stability of the immobilized antibodies.



### Sample Preparation and Storage

This cell lysis protocol is provided as a reference. Specific cell types or targets may benefit from alternative buffer components or techniques, depending upon the particular research application. Most lysis buffers are compatible with MSD MULTI-SPOT plates, although high concentrations of denaturing detergents (>0.1%) and reducing agents (DTT >1mM) should be avoided. Please contact MSD Scientific Support with any questions regarding lysate preparation options.

All manipulations should be performed on ice. The amount of complete lysis buffer required will vary depending on scale of preparation and type of cells. Larger cells (e.g. NIH3T3, HeLa) should be lysed at concentrations of  $1-5 \times 10^6$  cells per mL of lysis buffer. Smaller cells (e.g. Jurkat) should be lysed at concentrations of  $1-5 \times 10^7$  cells per mL of lysis buffer.

Analysis of proteins in their activated state (i.e. phosphorylated) usually requires stimulation prior to cell lysis. Verification of cell stimulation and sample preparation should be performed prior to using this kit. Phosphate Buffered Saline (PBS) should be ice-cold prior to use.

#### **Suspension Cells**

Pellet cells by centrifugation at 500 x g for 3 minutes at 2-8°C. Discard supernatant and wash the pellet once with cold PBS. Pellet the cells again, discard supernatant and resuspend in complete lysis buffer at  $1 - 5 \times 10^7$  cells per mL. Incubate on ice for 30 minutes. A shorter incubation time of 15 minutes may be adequate for many targets. Clear cellular debris from the lysate by centrifugation greater than or equal to  $10000 \times g$ , at 2-8°C for 10 minutes. Discard the pellet and determine protein concentration in the lysate using a detergent compatible protein assay such as BCA. Unused lysates should be aliquoted and quickly frozen in a dry ice-ethanol bath and stored at  $\leq$ -70°C.

#### **Adherent Cells**

All volumes are determined for cells plated in 15 cm dishes. Remove media from the plates and wash cells one time with 5 mL cold PBS. Add 2 mL PBS to the plates and scrape the cells from the surface of the dish and transfer into 15 mL conical tubes. Pellet the cells by centrifugation at 500 x g for 3 minutes at 2-8°C. Discard supernatant and resuspend cells in 0.5 - 2 mL of complete lysis buffer per dish. Incubate on ice for 30 minutes. A shorter incubation time of 15 minutes may be adequate for many targets. Clear cellular debris from the lysate by centrifugation greater than or equal to 10000 x g, at 2-8°C for 10 minutes. Discard the pellet and determine protein concentration in the lysate using a detergent compatible protein assay such as BCA. Unused lysates should be aliquoted and quickly frozen in a dry ice-ethanol bath and stored at  $\leq$ -70°C.

Refer to Appendix I for cell lysate preparation protocol modifications that accommodate the use of 96-well culture plates.

## Assay Protocol

The following protocol describes the most conservative approach to achieving optimal results with the MULTI-ARRAY<sup>®</sup> Phospho-Histone H3 (Ser10) Assay. The entire assay, including plate analysis on the MSD reader, can be completed in 3.5 hours. Once desired results are achieved, the protocol can be streamlined to eliminate multiple incubations and wash steps. Samples may be prepared for testing in the manner outlined in the Sample Preparation and Storage section.

#### 1. Block Plate and Prepare Samples:

- Add 150 μL of blocking solution into each well. Seal the plate with an adhesive plate seal, and incubate for 1 hour with vigorous shaking (300–1000 rpm) at room temperature.
- b. Prepare complete lysis buffer just prior to sample dilution.

**Note:** Samples, including cell lysates, etc., may be used neat or after dilution.

- MSD plates are compatible with most sample matrices. Avoid reagents that will denature the capture antibodies (e.g. high concentrations of reducing agents such as DTT should be avoided, and also SDS and other ionic detergents should be 0.1% or less in the sample applied to the well).
- Depending on the stability of the target in the matrix, additional protease and phosphatase inhibitors may be required in the matrix or diluent.
- If working with purified protein, only a few nanograms per well will generally provide a strong assay signal. Purified recombinant proteins may exhibit differences in both signal and background as compared to native proteins in cell lysates.
- > Keep diluted samples on ice until use
- c. Prepare positive and negative cell lysates:
  - (if purchased separately).
  - Thaw cell lysate samples on ice, and dilute them immediately before use. Keep on ice during all manipulations, and discard all remaining thawed, unused material.
  - Dilute cell lysate in complete lysis buffer to a final concentration of 0.8 µg/µL. This will deliver 20 µg/well lysate in 25 µL. A dilution series may also be prepared if desired.

#### Notes

Read entire protocol prior to beginning the assay.

Solutions containing MSD Blocker A should be stored at 2-8°C and discarded after 14 days.

Complete lysis buffer should be kept ice-cold during all experimental manipulations.

The sensitivity of MSD immunoassays rivals that of ELISAs and Western blots. The amount of sample required for a given assay will depend on the abundance of the analyte in the matrix and the affinities of the antibodies used.

Samples and standards cannot be serially diluted in the MSD plate. Use microcentrifuge tubes or a separate 96-well polypropylene plate to prepare dilutions.  Wash and Add Samples: Wash the plate 3 times with 300 μL/well of Tris Wash Buffer. Add 25 μL of samples per well. Seal the plate with an adhesive plate seal, and incubate for 1 hour with vigorous shaking (300– 1000 rpm) at room temperature.

Prepare detection antibody solution during this time.

 Wash and Add Detection Antibody: Wash the plate 3 times with 300 μL/well of Tris Wash Buffer. Add 25 μL of detection antibody solution to each well of the MSD plate. Seal the plate with an adhesive plate seal, and incubate for 1 hour with vigorous shaking (300–1000 rpm) at room temperature.

Prepare 1X Read Buffer T during this time.

- 4. Wash and Read: Wash the plate 3 times with 300  $\mu$ L/well of Tris Wash Buffer. Add 150  $\mu$ L of 1X Read Buffer T to each well of the MSD plate. Analyze the plate on the MSD instrument:
  - a. Double click on DISCOVERY WORKBENCH<sup>®</sup> icon on computer desktop (if not already open).
  - b. Click the imager icon in upper left corner of screen (if not already open to plate reading screen).
  - c. From the pull down menu select "Read From Barcode."
  - d. If only reading one plate check "Return Plate to Input Stack." Then check "Read Plate(s)" checkbox and enter 1.
  - e. If reading multiple plates, check the "Read Plate(s)" checkbox and enter number of plates to be read in the text field. For example, if five plates need to be read, type in "5."
  - f. Click the "Run" button. The "Run Options" window will be displayed
  - g. If the data from each microplate is to be exported as individual files, select "Separate Files" in the "Export" area of the "Run Options" window. Select "Appended File" if all data from the entire stack run is to be exported to one file. Select "Default" in the "Export Format" area. Check the box to export default data file.
  - h. If desired, make selections to export a custom data file.
  - i. Browse and select the location to export data files.
  - j. Click OK to initiate the run.
  - k. Data will be automatically saved in the software database. Text versions of the requested data files will be exported to the designated folder.

#### Notes

Shaking a 96-well MSD MULTI-ARRAY or MULTI-SPOT plate during an incubation step will typically accelerate capture at the working electrode.

The lysate sample incubation time provided is optimized for the use of MSD cell lysates. Samples from other sources may require a longer incubation.

Excess diluted read buffer may be kept in a tightly sealed container at room temperature for later use.

Bubbles introduced during the read buffer addition will interfere with imaging of the plate and produce unreliable data.

Plate should be imaged within 5 minutes following the addition of read buffer. Due to the varying nature of each research application, assay stability should be investigated prior to allowing plates to sit with read buffer for extended periods.

An all-inclusive indelible copy of the data and associated instrument information will be saved on the internal database, regardless of data file export selection. Additional copies of the data can be exported in any layout at a later time using this database. Consult the instrument user manual for more information.



# Typical Data

Representative results for the MULTI-ARRAY Phospho-Histone H3 (Ser10) Assay are illustrated below. The signal and ratio values provided below are example data; individual results may vary depending upon the samples tested. Growing HeLa cells (negative) were treated with nocodazole (1 µg/mL; 19 hours) and calyculin A (50 nM; final 30 minutes) (positive). Whole cell lysates were added to MSD MULTI-SPOT 4-Spot plates coated with an anti-total Histone H3 antibody on one of the four spatially distinct electrodes per well. Phosphorylated Histone H3 was detected with anti-phospho-Histone H3 (Ser10) antibody conjugated with MSD SULFO-TAG reagent.



*Figure 2:* Sample data generated with the MULTI-ARRAY Phospho-Histone H3 (Ser10) Assay. Increased signal is observed with the titration of pHistone H3 positive cell lysate. Signal for negative lysate remains low throughout the titration.

### Assay Components

The capture and detection antibodies used in this assay are listed below. They cross-react with human, mouse, and rat whole cell lysates.

	Source Species		
Analyte	MSD Capture Antibody	MSD Detection Antibody	
Phospho-Histone H3	Rabbit Polyclonal	Rabbit Monoclonal	



### Limitations of the Procedure

The following points should be noted with the MULTI-ARRAY Phospho-Histone H3 (Ser10) Assay to maximize assay sensitivity and performance.

- > A no-wash assay format may be employed, however lower sensitivity may be observed.
- > All buffers containing phosphate should be avoided when detecting phosphoproteins.
- > Due to the unstable nature of phosphoproteins, cell lysates should be thawed immediately prior to use, and any remaining thawed material should be subsequently discarded.

### References

- 1. Strahl B, Allis CD. The language of covalent histone modifications. Nature. 2000 Jan 6;403(6765):41-5.
- 2. Peterson CL, Laniel MA. Histones and histone modifications. Curr Biol. 2004 Jul 27;14(14):R546-51.
- 3. Taverna SD, Li H, Ruthenburg AJ, Allis CD, Patel DJ. How chromatin-binding modules interpret histone modifications: lessons from professional pocket pickers. Nat Struct Mol Biol. 2007 Nov;14(11):1025-40. Epub 2007 Nov 5.
- 4. Verdone L, Caserta M, Mauro E. Role of histone acetylation in the control of gene expression. Biochem Cell Biol. 2005 Jun;83(3):344-53.
- 5. Spivakov M, Fisher AG. Fisher Epigenetic signatures of stem-cell identity. Nat Rev Genet. 2007 Apr;8(4):263-71.



## Appendix

#### 96-well Culture Plate Modifications

Successful adaptation to a 96-well culture format is cell type and target-dependent. The number of cells to be plated per well should be determined for each cell type. General recommended plating concentrations for adherent cells range from  $1 \times 10^4 - 5 \times 10^4$  cells per well and approximately  $2 \times 10^6$  cells per mL (50 - 75 µL per well) for suspension cells. These numbers are provided as a guide, and the optimal concentrations will vary depending upon cell line used.

#### **Suspension Cells**

For flat bottom plates, experiments should be designed such that the final volume per well is  $50 - 75 \mu$ L. Perform cell lysis using a 4X complete lysis buffer concentrate, supplemented with protease and phosphatase inhibitors at 4X concentrations. Add 4X complete lysis buffer directly to cells in the growth medium for a final 1X concentration in the well.

Note: With some effort, a 10X complete lysis buffer can also be prepared.

(For conical microwell plates, perform lysis by pelleting the cells, removing most of the growth medium, and adding a constant amount of 1X complete lysis buffer).

#### **Suspension Cells**

Plate cells on biologically treated tissue culture ware (such as BD BioCoat Cellware (Becton, Dickinson and Company, Franklin Lakes, NJ) to reduce variability due to cells lost as growth medium is removed. Treat cells as desired. Gently aspirate growth medium from microwell plate. A PBS wash step is not required and can introduce variability. Add 50-100 µL 1X complete lysis buffer per well.

Cell lysis time should be determined by the end user. Some targets are immediately available for detection. Other targets may require an incubation step at room temperature, 45°C, or on ice with gentle agitation.

Carefully pipet cell lysate onto prepared capture plate, and proceed with assay protocol.

It is important to transfer a constant volume and avoid pipetting too vigorously, as the introduction of air bubbles may result. (Targets can be captured from a volume greater than  $25 \ \mu$ L).

### **Summary Protocol**

#### MSD 96-well MULTI-ARRAY Phospho-Histone H3 (Ser10) Assay Kit

MSD provides this summary protocol for your convenience. Please read the entire detailed protocol prior to performing the MULTI-ARRAY Phospho-Histone H3 (Ser10) Assay.

#### Step 1 : Block Plate and Prepare Samples

Add 150 µL/well of blocking solution. Incubate at room temperature with vigorous shaking (300-1000 rpm) for 1 hour. Prepare complete lysis buffer just prior to sample dilution. Prepare positive and negative cell lysates and keep on ice until use.

#### Step 2 : Wash and Add Sample

Wash the plate 3 times with 300 µL/well of 1X Tris Wash Buffer. Dispense 25 µL/well samples. Incubate at room temperature with vigorous shaking (300-1000 rpm) for 1 hour.

#### Step 3 : Wash and Add Detection Antibody Solution

Wash the plate 3 times with 300  $\mu$ L/well of 1X Tris Wash Buffer. Dispense 25  $\mu$ L/well 1X detection antibody solution. Incubate at room temperature with vigorous shaking (300-1000 rpm) for 1 hour.

#### Step 4 : Wash and Read Plate

Wash the plate 3 times with 300  $\mu$ L/well of 1X Tris Wash Buffer. Dispense 150  $\mu$ L/well 1X Read Buffer T. Analyze plate on an MSD instrument within 5 minutes of read buffer addition.



### Plate Diagrams

