

# Assay Kits from MSD

MSD offers many kits for measurement of cytokines and chemokines, cell signaling targets, and clinical biomarkers.

## Cytokines and Chemokines

View a complete list of available markers at:

[http://www.mesoscale.com/CatalogSystemWeb/WebRoot/products/assays/cytokines\\_singleplex.aspx](http://www.mesoscale.com/CatalogSystemWeb/WebRoot/products/assays/cytokines_singleplex.aspx)

Download protocols for individual assays at [www.mesoscale.com](http://www.mesoscale.com).

Kits are packaged according to sample matrix.

### Tissue Culture Kit

- 96-well format is available in 1-, 5-, and 25- plate kits.
- 384-well format is available in 5- and 50- plate kits. Please call for availability.
- Accommodates simple matrices.

### Ultra-Sensitive Kit

- 96-well format available in 1-, 5-, 25- plate kits.
- 384-well format available as a custom order.
- Accommodates simple and complex matrices.

### Base Kit

- 96-well format is available in 25- and 250- plate kits.
- Customer must supply diluents and calibrator.

Material	Storage	Ultrasensitive Kit	Tissue Culture Kit	Base Kit
MULTI-ARRAY® or MULTI-SPOT® plate	4°C	X	X	X
SULFO-TAG™ Detection Antibody (individual or blend)	4°C	X	X	X
Cytokine Calibrator (individual or blend)	-80°C	X	X	
Calibrator Diluent	4°C	X	X	
Antibody Diluent	4°C	X	X	X
Blockers (assay specific)	4°C	X	X	
Read Buffer T (4X)	RT	X	X	X
Serum Cytokine Assay Diluent	-20°C	X		
Serum Cytokine Antibody Diluent	-20°C	X		

# Cell Signaling Pathways (Phosphoproteins, Intracellular Markers, Ubiquitinylation)

View a complete list of available markers at:

[http://www.mesoscale.com/CatalogSystemWeb/WebRoot/products/assays/phosphoprotein\\_singleplex.aspx](http://www.mesoscale.com/CatalogSystemWeb/WebRoot/products/assays/phosphoprotein_singleplex.aspx)

Download protocols for individual assays at [www.mesoscale.com](http://www.mesoscale.com).

## Whole Cell Lysate Kit

- 96-well format is available in 1-, 5-, and 20- plate kits
- 384-well format is available in 5- and 50- plate kits. Please call for availability.
- Accommodates whole cell lysate samples. For alternative matrices, contact your scientific support representative.
- Customer must supply controls or separately purchase MSD positive/negative lysate sets.

## Base Kit

- Available in 20- plate kits.
- Customer must supply buffers, blockers, inhibitors, additives and controls. Information on preparation can be accessed at: [http://www.mesoscale.com/CatalogSystemWeb/WebRoot/literature/notes/pdf/Phosphoprotein\\_materials\\_list.pdf](http://www.mesoscale.com/CatalogSystemWeb/WebRoot/literature/notes/pdf/Phosphoprotein_materials_list.pdf)

Material	Storage	Whole Cell Lysate Kit	Base Kit
MULTI-SPOT Plate	4°C	X	X
SULFO-TAG Detection Antibody	4°C	X	X
Blocker A	4°C	X	
Assay Specific Blockers	4°C or RT	X	
Tris Wash Buffer (10X)	4°C	X	
Tris Lysis Buffer (1X)	4°C	X	
Phosphatase Inhibitor I (100X)	4°C	X	
Phosphatase Inhibitor II (100X)	4°C	X	
Protease Inhibitor Solution (50X)	-20°C	X	
Read Buffer T (4X)	RT	X	X

# Clinical Biomarker Assays (Alzheimer's, Metabolic, Hypoxia, Vascular/Growth)

View a complete list of available markers at:

<http://www.mesoscale.com/CatalogSystemWeb/WebRoot/products/assays.aspx>

Download protocols for individual assays at [www.mesoscale.com](http://www.mesoscale.com).

## Matrix-Specific Kit

- 96-well format is available in 1-, 5-, and 20- plate kits.
- Diluents included and protocol is matrix specific.

## Base Kit

- Contents: Plates, Detection Antibody, Read Buffer T (4X).
- Available in 20- plate kits.
- Customer supplies their own buffers, blockers, inhibitors, and additives.

# Assay Development Packs

View application notes for each assay development pack at:

<http://www.mesoscale.com/CatalogSystemWeb/WebRoot/literature/notes.aspx>

## ELISA Conversion Pack I (uncoated plates)

*For developing immunoassays by spotting antibody directly on the electrode surface*

## ELISA Conversion Pack II (anti-species plates)

*For developing immunoassays utilizing mouse or rabbit antibodies*

## ELISA Conversion Pack III (avidin and SA plates)

*For developing immunoassays utilizing biotinylated antibodies*

## Immunogenicity Development Pack

*For developing immunoassays for immunogenicity and bioproduction applications*

### Kit Contents

Materials Included	ELISA Conversion Pack I (uncoated plates)	ELISA Conversion Pack II (anti-species plates)	ELISA Conversion Pack III (biotin capture)	Immunogenicity Development Pack
MULTI-ARRAY 96-well High Bind Plates	10			10
MULTI-ARRAY 96-well Plates	10			
MULTI-ARRAY 96-well Anti-Rabbit (Goat) Plates		5		
MULTI-ARRAY 96-well Anti-Mouse (Goat) Plates		5		
MULTI-ARRAY 96-well Avidin Plates			5	5
MULTI-ARRAY 96-well High Bind Avidin Plates			5	
MULTI-ARRAY 96-well Streptavidin Plates			5	5
MULTI-ARRAY 96-well High Bind Streptavidin Plates			5	5
SULFO-TAG NHS Ester, 150 nMoles	2		2	2
Spin Columns	4		4	4
SULFO-TAG Labeled Streptavidin	1	1		1
SULFO-TAG Labeled Anti Mouse Antibody	1	1	1	
SULFO-TAG Labeled Anti Rabbit Antibody	1	1	1	
SULFO-TAG Labeled Anti Human Antibody				1
Blocker A Kit	1	1	1	1
Blocker B	1	1	1	1
Read Buffer T (4X)	1	1	1	1

# Supplemental Reagents

Designed for use with 20 - 25 96-well plates or 10 384-well plates.

View additional buffers, diluents and blockers available for sale at:

<http://www.mesoscale.com/CatalogSystemWeb/WebRoot/products/reagents.aspx>

## Cytokine Reagent Support Packs

- Three varieties available for human, mouse and rat assays.
- Contents:
  - Diluents for tissue culture samples
  - Serum cytokine assay and antibody diluents
  - Blocker B (included with human pack only)
  - Blocker D-B (included with mouse pack only)

## Phosphoprotein Reagent Support Pack

- Contents:
  - Tris Lysis and Wash Buffers
  - Phosphatase and Protease Inhibitors
  - Blockers: A, B, D-B, D-M, D-R, D-G

## Contact Information

Telephone: 1.240.631.2522 x4685

Fax: 1.301.990.2776

Email: [customerservice@mesoscale.com](mailto:customerservice@mesoscale.com)

Web: [www.mesoscale.com](http://www.mesoscale.com)

