

Assay Kits from MSD[®]

MSD offers many kits for measurement of cytokines and chemokines, cell signaling targets, bioprocess contaminants, 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

Download protocols for individual assays at 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.
- 384-well format is available in 50- plate kits.
- Customer must supply diluents and calibrators.

Material	Storage	Ultrasensitive Kit	Tissue Culture Kit	Base Kit
MULTI-ARRAY [®] or MULTI-SPOT [®] plate	2-8°C	X	X	X
SULFO-TAG [™] Detection Antibody (individual or blend)	2-8°C	X	X	X
Cytokine Calibrator (individual or blend)	≤-70°C	X	X	
Diluent 1	2-8°C		X	
Diluent 100	2-8°C		X	X
Blockers (assay specific)	2-8°C	X	X	
Read Buffer T (4X)	RT	X	X	X
Serum Cytokine Assay Diluent	≤-10°C	X		
Serum Cytokine Antibody Diluent	≤-10°C	X		

Cell Signaling Pathways (Phosphoproteins, Intracellular Markers, Ubiquitylation)

View a complete list of available markers at:

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

Download protocols for individual assays at 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

- Contents: Plates, Detection Antibody, Read Buffer T (4X).
- Available in 20- plate kits.
- 384-well format is available in 50- plate kits. Please call for availability.
- 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

Material	Storage	Whole Cell Lysate Kit	Base Kit
MULTI-SPOT Plate	2-8°C	X	X
SULFO-TAG Detection Antibody	2-8°C	X	X
Blocker A	2-8°C	X	
Assay Specific Blockers	2-8°C or RT	X	
Tris Wash Buffer (10X)	2-8°C	X	
Tris Lysis Buffer (1X)	2-8°C	X	
Phosphatase Inhibitor I (100X)	2-8°C	X	
Phosphatase Inhibitor II (100X)	2-8°C	X	
Protease Inhibitor Solution (100X)	2-8°C	X	
Read Buffer T (4X)	RT	X	X

Clinical Biomarker Assays (Alzheimer's, Bone, Cardiac, Kidney, 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.

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 diluents, blockers, inhibitors, and additives.

Bioprocess Contaminant Assays

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.

Complete Kit

- 96-well format is available in 1-, 5-, and 20- plate kits.
- Reagents included and protocol are formatted for Bio-Production samples.

Base Kit

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

Assay Development Packs

View technical and application notes for assay development 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

Cell Based Assays

For developing cell based assays

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	Cell Based Assays
MULTI-ARRAY 96-well High Bind Plates	10			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		
MULTI-ARRAY 96-well High Bind Avidin Gold Plates			5	5	
MULTI-ARRAY 96-well Streptavidin Gold Plates			5	10	
MULTI-ARRAY 96-well High Bind Streptavidin Plates			5		
SULFO-TAG NHS Ester, 150 nMoles	2		2	4	
Spin Columns	4		4	8	
SULFO-TAG Labeled Streptavidin	1	1		1	1
SULFO-TAG Labeled Anti Mouse Antibody	1	1	1		1
SULFO-TAG Labeled Anti Rabbit Antibody	1	1	1		1
SULFO-TAG Labeled Anti Human Antibody				1	
Blocker A Kit	1	1	1	1	1
Blocker B	1	1	1	1	1
Read Buffer T (4X)	1	1	1	1	2

QUICKPLEX® Multiplexing Assay Kits

Kit Contents

Materials Included	QUICKPLEX ABCD High Bind Kit	QUICKPLEX ABCD Kit	QUICKPLEX ABCsA High Bind Kit	QUICKPLEX ABCsA Kit
96-well 4-Spot High Bind QUICKPLEX ABCD Plates	5			
96-well 4-Spot QUICKPLEX ABCD Plates		5		
96-well 4-Spot High Bind QUICKPLEX ABCsA Plates			5	
96-well 4-Spot QUICKPLEX ABCsA Plates				5
Linker A NHS Ester	1	1	1	1
Linker B NHS Ester	1	1	1	1
Linker C NHS Ester	1	1	1	1
Linker D NHS Ester	1	1		
Biotin NHS Ester			1	1
SULFO-TAG NHS Ester	1	1	1	1
Spin Columns	8	8	8	8
Blocker A Kit	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

Complete Blocker Kit

- Contents:
 - Blockers: A, B, C, D-B, D-M, D-R, D-G

Contact Information

Telephone: 1.301.947.2085

Fax: 1.301.990.2776

Email: customerservice@mesoscale.com

Web: www.mesoscale.com

