# MSD® TECHNOLOGY PLATFORM





# **MESO SCALE DISCOVERY® (MSD) Technology**

Г

MESO SCALE DISCOVERY'S MULTI-ARRAY® technology enables the detection of biomarkers in single and multiplex formats utilizing the next generation of electrochemiluminescence detection. The combination of electrochemiluminescence and patterned arrays bring speed and high density of information to research through miniaturization, organization, and parallel processing of biological assays. The MSD product line includes a diverse menu of single and multiplex assay kits for profiling biomarkers, cell signaling pathways, and other applications, as well as a suite of plates and reagents for assay development.



Antibodies on a carbon surface

| Company Overview                              | 1  |
|---|----|
| MSD Technology                                | 2  |
| MSD - Spot the Difference                     | 3  |
| Measurement of Biomarkers in Complex Samples  | 4  |
| Assessment of Toxic Effects                   | 5  |
| Rapid Quantification of Phosphoproteins       | 6  |
| Study of Compound Inhibition                  | 7  |
| Custom Assays                                 | 8  |
| Prototype Printing Services                   | 9  |
| Evaluation of Protein Therapeutics            | 10 |
| SECTOR Instruments                            | 11 |
| MSD Resources at Your Service                 | 12 |
| Publications using MSD MULTI-ARRAY Technology | 13 |



MESO SCALE DISCOVERY A division of Meso Scale Diagnostics, LLC.

1601 Research Blvd. Rockville, MD 20850 USA Phone: 1.240.314.2795 Fax: 1.301.990.2776 Email: customerservice@mesoscale.com www.mesoscale.com

Spot the Difference

# **Company Overview**

MESO SCALE DISCOVERY offers a unique, multiplexed immunoassay platform for the quantification of proteins in biological samples. With over 400 convenient assay kits and assay customization capabilities, MSD has enabled scientists to make accurate and precise determinations of levels of cytokines, phosphoproteins, and other biomarkers in different matrices. High quality data can be obtained in less time on the MSD platform with minimal effort and low cost. MSD's MULTI-ARRAY technology has been adopted by major pharmaceutical companies, clinical research organizations, biotech companies, personalized medicine companies, and academic and government institutions.

The graphic below illustrates the broad spectrum of applications of the MSD technology.



The benefits of MULTI-ARRAY technology are well recognized. Here are just a few peer-reviewed publications highlighting some of the advantages of the platform.

- high inter-laboratory reproducibility, low matrix effects and high reliability (Fichorova, R.N., et al. Anal Chem. 2008 80(12): 4741-51.)<sup>1</sup>
- reproducibility and cost-effectiveness over Western blot (Gowan, S.M., et al. Assay and Drug Dev Technol. 2007 5(3): 391-401.)<sup>2</sup>
- novel sandwich assay that provides sensitivity, accuracy, reproducibility and absolute quantification (Cao, L., et al. Cancer Research. 2008 68: 8039-8048.)<sup>3</sup>
- most suitable to detect low affinity anti-drug antibodies (Liang, M., et al. Assay and Drug Dev Technol. 2007 5(5): 1-8.)<sup>4</sup>

# **MSD** Technology

MSD's electrochemiluminescence detection technology uses SULFO-TAG<sup>™</sup> labels, which emit light upon electrochemical stimulation initiated at the electrode surfaces of MULTI-ARRAY and MULTI-SPOT® microplates.

#### **Electrochemiluminescence Features:** Minimal background signals and high signal to background ratios - the stimulation mechanism (electricity) is decoupled from the signal (light) Proximity - only labels bound near the electrode surface are detected, enabling non-washed assays Flexibility - labels are stable, non-radioactive, and LIGHT Measured signal is light conveniently conjugated to biological molecules \*Ru(bpy)<sub>3</sub><sup>2+</sup> Emission at ~620 nm - eliminates problems with color TPA quenching Signal amplification - multiple excitation cycles of each label Ru(bpy)<sub>3</sub><sup>3+</sup> TPA'+ TPA Ru(bpy)<sub>3</sub><sup>2+</sup> enhance light levels and improve sensitivity e-Flexible surface coatings to suit most any biology Counter

- Carbon electrode plate surface has 10X greater binding capacity than polystyrene
- Custom surface coatings and patterns



## **MULTI-ARRAY and MULTI-SPOT Features:**

- Capability to simultaneously measure multiple analytes in the same well
- High density arrays for high throughput multiplexing of biomarkers
- The unique bar code label on each plate enables complete traceability back to MSD manufacturing records
- The MSD DISCOVERY WORKBENCH<sup>®</sup> software provides customers with a powerful tool for data analysis







96 Well 10-Spot





384 Well Single-Spot

384 Well 4-Spot



# **MSD - SPOT THE DIFFERENCE®**

MESO SCALE DISCOVERY's unique spot patterns are the hallmark of its MULTI-ARRAY technology for the detection of biomarkers in single and multiplex formats. MSD offers an innovative platform for immunoassays that have ultra-low detection limits, provide up to five logs of linear dynamic range, use minimal sample, and handle difficult matrices easily. Combined, these advantages enable the measurement of native levels of biomarkers in normal and diseased samples without multiple dilutions. MSD's simple protocols and streamlined assay formats from MSD reduce workflow without compromising quality. The result is an increase in productivity while eliminating unnecessary costs.



- Large dynamic range
- High sensitivity
- High precision
- Low background
- Conserves sample volume
- Simple protocols
- Reduces matrix effects
- Eliminates multiple dilutions

## **Assay Solutions and Services**

- Over 300 single analyte kits
- Over 200 multiple analyte kits
- Customized multiplexed kits
- ELISA conversion packs
- QUICKPLEX<sup>®</sup> packs
- Prototype Printing Services
- Complimentary on-site scientific support
- Contract assay development

### Multiplexing



- Multiple analytes in one well
- No compromise in performance or speed
- Catalog assay panels for rapid delivery
- Custom panels available

## **SECTOR®** Instruments

- Ultra-fast read time
- No fluidics
- No user calibration required
- Reliable measurements
- Integrated data analysis tool
- Comprehensive validation packages
- Software support for 21 CFR Part 11 compliance
- Comprehensive service plans



# **Measurement of Biomarkers in Complex Samples**

Clinical samples push the limits of traditional methods for immunoassays (e.g. ELISA, RIA). Complex matrices (e.g. sputum, vaginal fluids, etc.), widely ranging concentations of analytes, and limited sample volume can make assays intractable. MSD's assays improve sensitivity, expand the dynamic range, enable measurement of multiple analytes from a single sample (i.e. multiplexing), and work well in difficult sample types. The MSD platform has also proven to be well-suited for use in regulated work environments with available packages for IQ/OQ/PQ validation and software designed to support 21 CFR Part 11 and GLP compliance. Visit www.mesoscale.com for a complete list of assays and reagents.

Catalog Numbers for Human ProInflammatory 9-Plex Ultra-Sensitive Kit

|       | Kit Size | Catalog Numbers |
|-------|----------|-----------------|
|       | 1 Plate  | K15007C-1       |
| Human | 5 Plate  | K15007C-2       |
|       | 25 Plate | K15007C-4       |

#### Representative Data from Clinical Samples using MSD Human ProInflammatory 9-Plex



- Wide dynamic range enables biomarker measurement in controls and diseased samples with minimal dilution
- Highly adaptable assays allow quantification of analytes in complex sample matrices
- High sensitivity and multiplexing capability facilitate analysis of several biomarkers in a single sample

|              | Analyte |       |          |       |        |       |      |       |       |
|--------------|---------|-------|----------|-------|--------|-------|------|-------|-------|
| Analyte      | IL-2    | IL-8  | IL-12p70 | IL-1β | GM-CSF | IFN-γ | IL-6 | IL-10 | TNF-α |
| LLOD (pg/mL) | 0.35    | 0.090 | 1.4      | 0.36  | 0.34   | 0.53  | 0.27 | 0.21  | 0.50  |

LLOD (Lower Limit of Detection) is defined as 2.5 SD above the background.





In the above study<sup>\*</sup>, the MSD Human ProInflammatory 9-plex Ultra-Sensitive Assay was used to analyze a total of 127 human sera samples, which included diseased pools and controls. The upper end of the calibrator curve for this panel was 10000 pg/mL for all cytokines and the lower limit of detection (LLOD) was determined as 2.5 standard deviations above the background.

\* The Biomarker Reference Set for Cancers in Women (BRSCW) was provided by the National Cancer Institute on behalf of the Early Detection Research Network (EDRN).



# **Assessment of Toxic Effects**

In toxicity studies, compound or disease-induced changes are typically evaluated using a combination of histochemical endpoints and a number of potential biomarkers, each of which can indicate toxic change in tissues and organs. The best biomarkers are specific to particular organs or tissue types. Toxicologists require an assay system with consistent performance, high sensitivity, and large dynamic range. MSD technology provides all of these characteristics plus multiplexing to save time and precious samples. Featured here is the MSD Muscle Injury Panel I (rat) which measures cardiac Troponin I, skeletal Troponin I, cardiac Troponin T, FABP3, and Myl3.

For more detail on these and other relevant markers please see our Toxicology Brochure online at www.mesoscale.com.

| Catalog | <b>Numbers</b>    | for        |
|---------|-------------------|------------|
| Muscle  | <b>Injury Pan</b> | el 1 (rat) |

|     | Kit Size                       | Catalog Number                      |
|-----|--------------------------------|-------------------------------------|
| Rat | 1 Plate<br>5 Plate<br>25 Plate | K15181C-1<br>K15181C-2<br>K15181C-4 |

## **Muscle Injury Panel 1 (rat)**





LLOD (Lower Limit of Detection) is defined as 2.5 SD above the background.

LLOD (ng/mL) 0.0040 0.017

**MSD** has developed assays to rapidly screen for biomarkers of drug-induced toxicity in conjunction with the Critical Path Initiative (C-Path) and Health and Environmental Sciences (HESI) consortia.

## Specificity



 MSD assays for cardiac markers are positive for cardiac homogenates and negative for others, whereas skeletal marker assays are positive for skeletal muscle only

sTnl

0.16 0.015 0.047

**MSD** offers robust assays for markers of kidney injury, cardiotoxicity, and acute phase inflammatory response. Please visit www.mesoscale.com for a comprehensive list of toxicology assay kits.

# **Rapid Quantification of Phosphoproteins**



- Need expensive imaging system for semi-quantitative analysis
- Low throughput: stripping and reprobing are an inaccurate and undesirable approach for multiplexing

# **Study of Compound Inhibition**

The discovery of potent and highly selective small molecule compounds has proven to be useful in targeted therapy of cancers, cardiovascular diseases, and neurodegenerative disorders. They serve as valuable tools for deciphering the functions of many cell signaling pathways. MSD assays can be used to rapidly evaluate the potency of an inhibitor against cell signaling targets such as Akt. The following example demonstrates the use of the MSD platform for dose response studies. LY294002 is a highly selective inhibitor of phosphatidylinositol 3-kinase and blocks Akt phosphorylation. This inhibitory effect of LY294002 has been demonstrated below. Rapamycin inhibits mammalian target of rapamycin (mTOR), which is downstream of Akt in the cascade. The following plot illustrates the inhibition of mTOR signaling by rapamycin, which in turn leads to an increase of Akt phosphorylation by a negative feedback inhibition process.<sup>22</sup> MCF7 cells were used for the study.

#### Catalog Numbers for Akt Signaling Panel II

|        | Kit Size | Catalog Numbers |
|--------|----------|-----------------|
| Human/ | 1 Plate  | K15177D-1       |
| Mouse/ | 5 Plate  | K15177D-2       |
| Rat    | 20 Plate | K15177D-3       |



#### Akt Signaling: Percent Activated Protein Normalized to Total Protein



#### Rapamycin Titration (2 hr) + IGF-1 (50 nM; 20 min)





nM Rapamycin (+ 50 nM IGF-1)

#### p4c-br ((11157/30)

- Faster and more quantitative measurements than Western blot
- High throughput analysis enables a large number of samples to be tested in 96- and 384-well formats

pAkt(Ser473)

# **Custom Assays**

MSD supports you in every step of the research and drug development process. If you are studying a special combination of analytes, we can provide a multiplex panel to meet your needs. We will work with you to prepare a custom kit according to your preferences and provide you with a protocol for the assay. If you need help, our field application scientists and scientific support team are available to support you.

## Design your multiplex in 3 easy steps



If you cannot find your assay in the MSD assay list, then tell us the antibody pairs you use in your ELISA, and we'll coat them for you through Prototype Printing Services.

## Up to 10 assays in 96-well plates

#### Mouse Metabolkine Multiplex Panel





|              | Analyte |        |         |       |        |          |       |
|--------------|---------|--------|---------|-------|--------|----------|-------|
|              | IL-6    | GM-CSF | Insulin | MCP-1 | Leptin | Resistin | TNF-α |
| LLOD (pg/mL) | 4.3     | 1.1    | 125     | 1.3   | 83     | 10       | 6.1   |

LLOD (Lower Limit of Detection) is defined as 2.5 SD over the background signal.

# Up to 4 assays in 384-well plates

#### Human Cytokine Multiplex





# **Prototype Printing Services**

MSD offers Prototype Printing Services to facilitate assay development by customers. Prototype Printing Services provide the customer with a rapid and convenient way to get MSD plates coated with materials of their choice.

## Types of Capture Materials coated on MSD plates

#### 1. Antibodies

Antibodies are readily coated on MSD plates. Customers routinely see significant improvements in assay performance when ELISAs are converted to the MSD format.



#### 2. Proteins and Peptides

The following example demonstrates the use of an MSD 10-Spot plate for antibody screening. Titration of hybridoma supernatants identifies the high binding clones, which are subsequently tested for cross-reactivity as well as specificity.



By screening antibodies against multiple antigens, cross-reactivity and specificity can be determined in the same well.

How to Order

STEP 2: Place your order Reference the quote

**STEP 1:** Obtain a quote from MSD for your analyte of interest by contacting MSD

Customer Service.

number on your order.

STEP 3: Receive your prototype assay.

#### 3. Carbohydrates and Polysaccharides

Carbohydrates have been successfully coated on MSD plates, which have enabled the testing of different Pneumococcal vaccines by Marchese, et al.<sup>5</sup> MSD plates coated with lipopolysaccharides have been used by Thompson, et al. for serodiagnosis of Brucellosis in ruminants.<sup>6</sup>

In addition to the above mentioned materials, MSD plates are highly amenable to coating with viral proteins, cell lysates, etc.

| Request        | a Quote   |
|----------------|---|
| Contact        | customer service to request a quote.            |
| tel:<br>Email: | 1.240.314.2795<br>customerservice@mesoscale.com |

# **Evaluation of Protein Therapeutics**

Immunogenicity testing is a crucial part of biopharmaceutical development. The EMA and FDA have mandated immunogenicity testing for biotherapeutics. More stringent recommendations regarding immunogenicity assay performance necessitates the development of more robust and tolerant assays. MSD has worked with leaders in the biotherapeutics field to fine-tune its technology to meet the needs of the immunogenicity community. MSD offers a suite of assay development materials and kits that provide superior solutions for each stage of drug development process. The typical steps in testing immunogenicity of therapeutic antibodies have been shown below. The development of cell-based neutralization assays on the MSD platform has also been featured below. Visit www.mesoscale.com for more information on immunogenicity assay development and a complete listing of materials and reagents.

#### Catalog Numbers for Immunogenicity **Development Pack**

Instrument SECTOR Imager SECTOR PR K13A04-1

**Catalog Numbers** K15A04-1



- MSD assays allow for higher free drug tolerance
- Fewer washes permits the detection of low affinity antibodies
- Flexible assay formats enable testing of many drug types including antibodies, proteins and peptides
- Rapid assay development makes the assay cost- and time-efficient

#### **Cell-based Neutralization Assay**





pot the Difference

# **SECTOR Instruments**

The SECTOR Imager 6000 and the SECTOR Imager 2400 offered by MSD are ideal for users seeking high information content and high throughput. Both instruments use ultra-low noise CCD cameras for ultimate sensitivity, wide dynamic range, and rapid read times.

MSD SECTOR PR<sup>®</sup> readers offer users a choice of multiplex and single readout capabilities in our popular and compact benchtop platform. These affordable readers provide a combination of speed, simplicity, and performance that makes them perfect for target validation, assay development, immunogenicity testing, and basic research applications. Assays developed on the SECTOR PR readers are fully portable to the SECTOR Imagers.

All the instruments use MSD DISCOVERY WORKBENCH software, which offers one-click assays, enhanced data export tools, partial-plate reading features, and supports 21 CFR Part 11 compliance.

## **SECTOR Imager Features**

- Single and multiplex assay formats
- Highly sensitive imaging systems
- No complicated fluidics
- Rapid read times (~1 minute/plate for SI6000, and ~3 minutes/plate for SI2400)
- Six logs dynamic range
- Non-washed assay formats
- Simple operation
- Workstation or automated operation
- Simultaneous bar code label reading on short and long sides of microplates

## **SECTOR PR Features**

- Photodiode array for fast and efficient detection
- No complicated fluidics
- Integrated barcode readers
- Multiplex and single array readouts

- Simple operation
- Five logs dynamic range
- Non-washed assay formats



**Catalog Numbers for** 

**SECTOR Instruments** 





| Model                   | Detection<br>Technology | Plate Read<br>Time | Multiplex<br>Capability | Automation<br>Integration |
|-------------------------|-------------------------|--------------------|-------------------------|---------------------------|
| SECTOR<br>Imager 6000   | CCD<br>Camera           | 70 seconds         | Yes                     | Yes                       |
| SECTOR<br>Imager 2400   | CCD<br>Camera           | 3.5 minutes        | Yes                     | Yes                       |
| SECTOR PR<br>400 Reader | Photodiode<br>array     | 2-5 minutes        | Yes                     | Optional                  |
| SECTOR PR<br>100 Reader | Photodiode<br>array     | 2 minutes          | No                      | Optional                  |

# **MSD** Resources at Your Service

## **Customer Support**

Phone: 1-240-314-2795 Fax: 1-301-990-2776 Email: CustomerService@mesoscale.com Web: www.mesoscale.com/support Hours of Operation: 5:00 AM to 8:00 PM, Monday – Friday, U.S. Eastern Time

## Scientific Support

Phone: 1-240-314-2798 Email: ScientificSupport@mesoscale.com Web: www.mesoscale.com/support Hours of Operation: 8:30 AM to 5:30 PM, Monday – Friday, U.S. Eastern Time

- Contract Assay Development Services
- Custom Assays and Prototype Printing Services
- On-Site Assistance

## **Field Service Engineers**

Phone: 1-301-947-2057 Email: InstrumentService@mesoscale.com Hours of Operation: 8:30 AM to 5:30 PM, Monday – Friday, U.S. Eastern Time After Hours: 1-301-767-5682

#### Literature

Browse our online library of product literature, technical application notes, and FAQs as well as an extensive list of research studies citing MSD technology.

- Brochures
- Customer Presentations
- Posters
- Product Inserts
- Publications





# Publications using MSD MULTI-ARRAY Technology

- <sup>1</sup> Fichorova, R.N., et al. Biological and technical variables affecting immunoassay recovery of cytokines from human serum and simulated vaginal fluid: a multicenter study. Anal Chem. 2008 80(12): 4741-51.
- <sup>2</sup> Gowan, S.M., et al. Application of Meso Scale Technology for the Measurement of Phosphoproteins in Human Tumor Xenografts. Assay and Drug Dev Technol. 2007 5(3): 391-401.
- <sup>3</sup> Cao, L., et al. Addiction to Elevated Insulin-like Growth Factor I Receptor and Initial Modulation of the AKT Pathway Define the Responsiveness of Rhabdomyosarcoma to the Targeting Antibody. Cancer Research. 2008 68: 8039-8048.
- <sup>4</sup> Liang, M., et al. Detection of High- and Low-Affinity Antibodies Against a Human Monoclonal Antibody Using Various Technology Platforms. Assay and Drug Dev Technol. 2007 5(5): 1-8.
- <sup>5</sup> Marchese, R.D., et al. Optimization and validation of a multiplex, electrochemiluminescence-based detection assay for the quantitation of immunoglobulin g serotype-specific antipneumococcal antibodies in human serum. Clin Vaccine Immunol. 2009 16(3): 387-96.
- <sup>6</sup> Thompson, I., et al. Competitive electrochemiluminescence (ECL) wash and no-washimmunoassays for the detection of serum antibodies to smooth Brucella strains. Clin Vaccine Immunol. 2009 16(5): 765-71.
- <sup>7</sup> Cludts, I., et al. Detection of neutralizing interleukin-17 antibodies in autoimmune polyendocrinopathy syndrome-1 (APS-1) patients using a novel non-cell based electrochemiluminescence assay. Cytokine. 2010 50(2): 129-37.
- <sup>8</sup> Lewczuk, P., et al. Soluble amyloid precursor proteins in the cerebrospinal fluid as novel potential biomarkers of Alzheimer's disease: a multicenter study. Molecular Psychiatry. 2010 15: 138–145.
- <sup>9</sup> Tai, E.S., *et al.* Insulin resistance is associated with a metabolic profile of altered protein metabolism in Chinese and Asian-Indian men. Diabetologia. 2010 53(4): 757-67.
- <sup>10</sup> Enewold, L., et al. Serum concentrations of cytokines and lung cancer survival in African Americans and Caucasians. Cancer Epidemiol Biomarkers Prev. 2009 18(1): 215-22.
- <sup>11</sup> Hansson, S.F., et al. Reduced levels of amyloid-beta-binding proteins in cerebrospinal fluid from Alzheimer's disease patients. J Alzheimers Dis. 2009 16(2): 389-97.
- <sup>12</sup> Lembo, A., et al. Use of serum biomarkers in a diagnostic test for irritable bowel syndrome. Aliment Pharmacol Ther. 2009 29(8): 834-42.
- <sup>13</sup> Loyet, K.M., et al. Technology comparisons for anti-therapeutic antibody and neutralizing antibody assays in the context of an anti-TNF pharmacokinetic study. J Immunol Methods. 2009 345(1-2): 17-289.
- <sup>14</sup> Mentor-Marcel, R.A., et al. Inflammation-associated serum and colon markers as indicators of dietary attenuation of colon carcinogenesis in ob/ob mice. Cancer Prev Res. 2009 2(1): 60-9.
- <sup>15</sup> Welge, V., et al. Combined CSF tau, p-tau181 and amyloid-beta 38/40/42 for diagnosing Alzheimer's disease. J Neural Transm. 2009 116(2): 203-12.
- <sup>16</sup> Leng, S.X., et al. ELISA and Multiplex Technologies for Cytokine Measurement in Inflammation and Aging Research. J Gerontol A Biol Sci Med Sci. 2008 63(8): 879-84.
- <sup>17</sup> Ryan, K.R., et al. Impaired dendritic cell maturation and cytokine production in patients with chronic mucocutanous candidiasis with or without APECED. Clin Exp Immunol. Vol. 2008 154(3): 406-14.
- <sup>18</sup> Lutterloh, E.C., et al. Inhibition of the RAGE products increases survival in experimental models of severe sepsis and systemic infection. Crit Care. 2007 11(6): 122.
- <sup>19</sup> Ghosh, T.K., et al. Toll-like receptor (TLR) 2-9 agonists-induced cytokines and chemokines: I. Comparison with T cell receptor-induced responses. Cell Immunol. 2006 243(1): 48-57.
- <sup>20</sup> Lu, Y., et al. A high throughput electrochemiluminescent cell-binding assay for therapeutic anti-CD20 antibody selection. J Immunological Methods. 2006 314(1-2): 74-9.
- <sup>21</sup> Moxness, M., et al. Immunogenicity Testing by Electrochemiluminescent Detection for Antibodies Directed against Therapeutic Human Monoclonal Antibodies. Clinical Chemistry. 2005 51: 1983-1985.

## **Other Cited References**

<sup>22</sup> Shi, Y., et al. Mammalian target of rapamycin inhibitors activate the AKT kinase in multiple myeloma cells by up-regulating the insulin-like growth factor receptor/insulin receptor substrate-1/phosphatidylinositol 3-kinase cascade. Mol Cancer Ther. 2005 4(10):1533-40.

