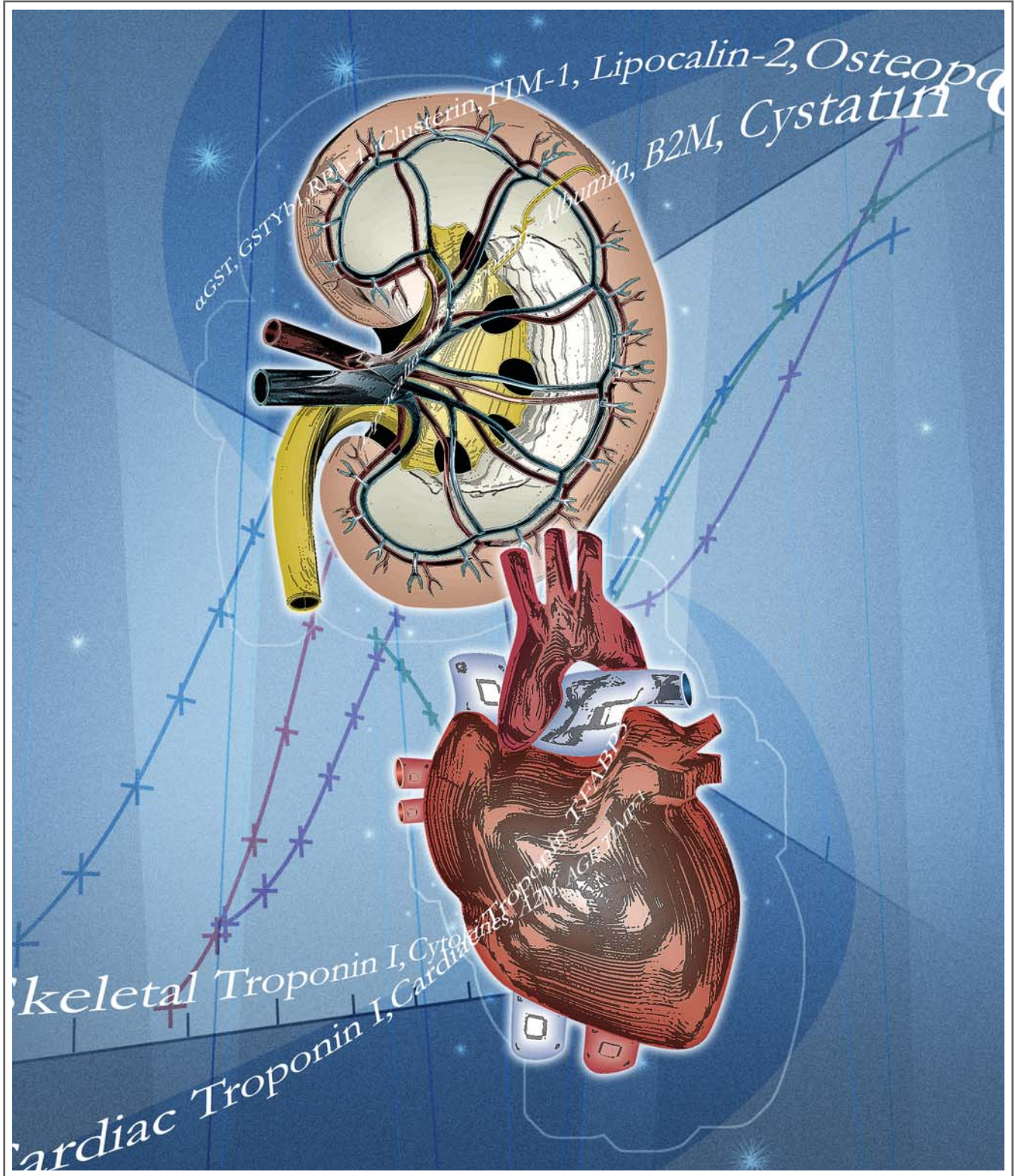
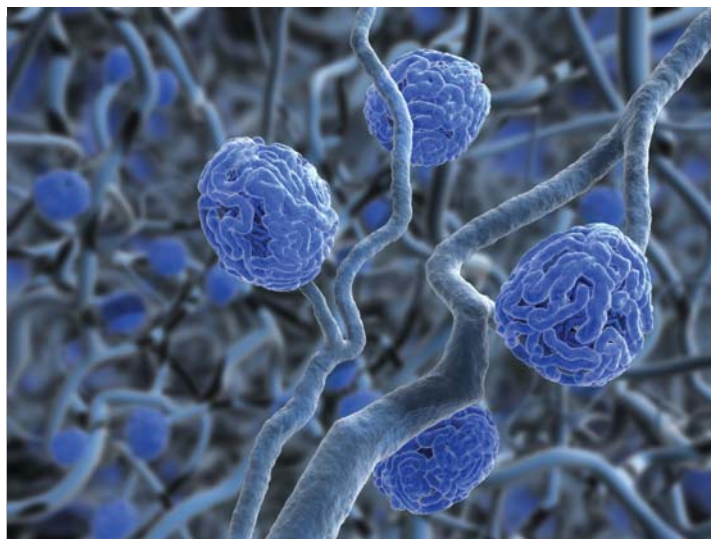


# TOXICOLOGY APPLICATIONS



# Meso Scale Discovery® Toxicology Applications

Preclinical toxicity determination for drug candidates is an area that increasingly requires measurement of several protein biomarkers. The best biomarkers are specific to particular organs or tissue types. Researchers require an assay system with consistent performance, high sensitivity and large dynamic range. MSD's® technology provides all of these characteristics plus multiplexing to save time and precious samples.



Glomeruli in the human kidney

■ The MSD Advantage	1
■ MSD Technology	2
■ Nephrotoxicity	3
■ Myotoxicity	8
■ Reproductive Toxicity	11
■ Inflammation and Growth Factors	12
■ Cytokines	15
■ Intracellular Signaling	17
■ Metabolic Disease	19
■ References	20
■ Select Catalog Numbers	21



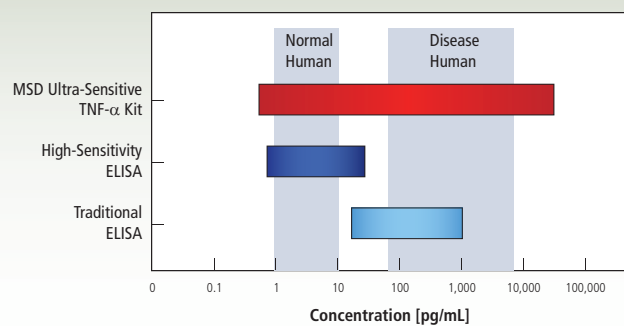
Meso Scale Discovery  
A division of Meso Scale Diagnostics, LLC.

9238 Gaither Road  
Gaithersburg, MD 20877 USA  
Phone: 1.301.947.2085  
Fax: 1.301.990.2776  
Email: [customerservice@mesoscale.com](mailto:customerservice@mesoscale.com)  
[www.mesoscale.com](http://www.mesoscale.com)

# The MSD Advantage

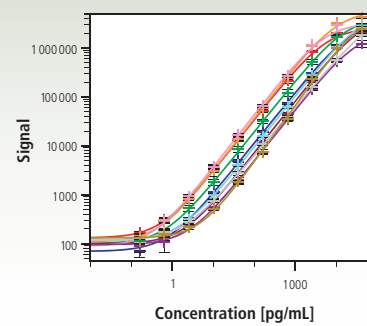
MESO SCALE DISCOVERY'S MULTI-ARRAY® Technology enables the detection of biomarkers in single and multiplex formats. Using the power of our patented electrochemiluminescence detection, MSD offers assays that have ultra-low detection limits, provide up to five logs of linear dynamic range, use minimal sample and handle difficult matrices easily. The extraordinary analytical performance of MSD assays translates into valuable advantages. Our sensitivity and dynamic range enable the development of assays that can measure native levels of biomarkers in normal and diseased samples without multiple dilutions, saving time, resources, sample and cost. Our multiplex panels allow measurements of multiple analytes from a single sample, without extra work or time. Our simple protocols and streamlined formats help validations go smoothly. These advantages—and many others—are why so many people use MSD assays. It's not only what MSD assays can do... it's what you can do with MSD assays.

## Rapid, Robust, Reproducible



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

## Multiplexing



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

## Assay Solutions and Services



- Over 300 single analyte kits
- Over 200 multiple analyte kits
- Customized multiplexed kits
- Qualified kits
- ELISA conversion packs
- QuickPlex® packs
- Prototype Printing Services
- On-site scientific support
- Contract assay development

## SECTOR® Instruments



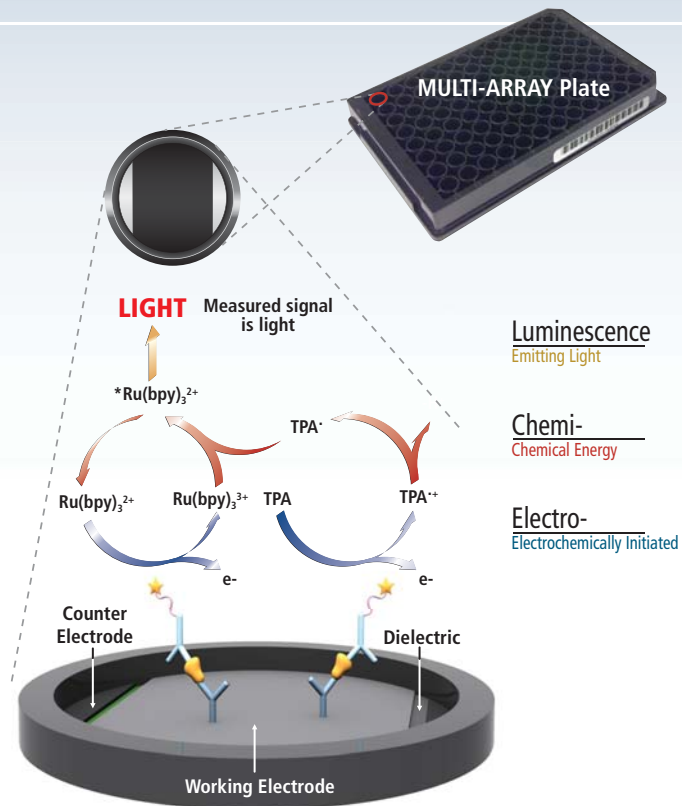
- Ultra-fast read time
- No fluidics
- No calibration required
- Reliable measurements
- Integrated data analysis tool
- Comprehensive validation packages
- 21 CFR 11 compliant software
- Comprehensive service plans
- On-site assistance

# MSD Technology

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

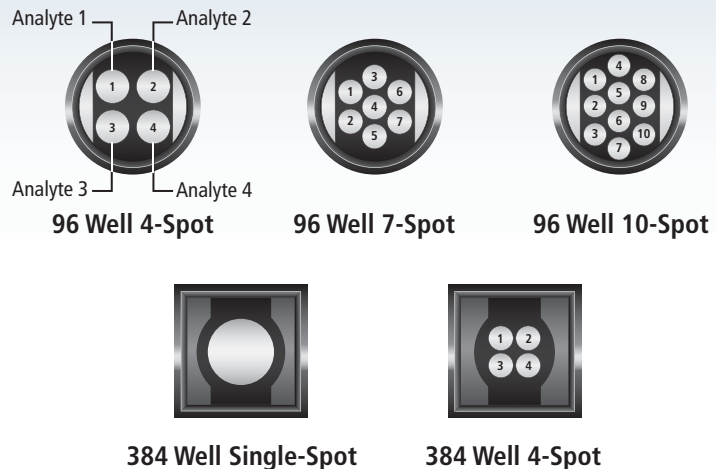
## Electrochemiluminescence Features:

- 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 are conveniently conjugated to biological molecules
- Emission at ~620 nm - eliminating problems with color quenching
- Signal amplification - multiple excitation cycles of each label enhance light levels and improve sensitivity
- Flexible surface coatings to suit most any biology
- 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® software provides customers with a powerful tool for data analysis



# Nephrotoxicity

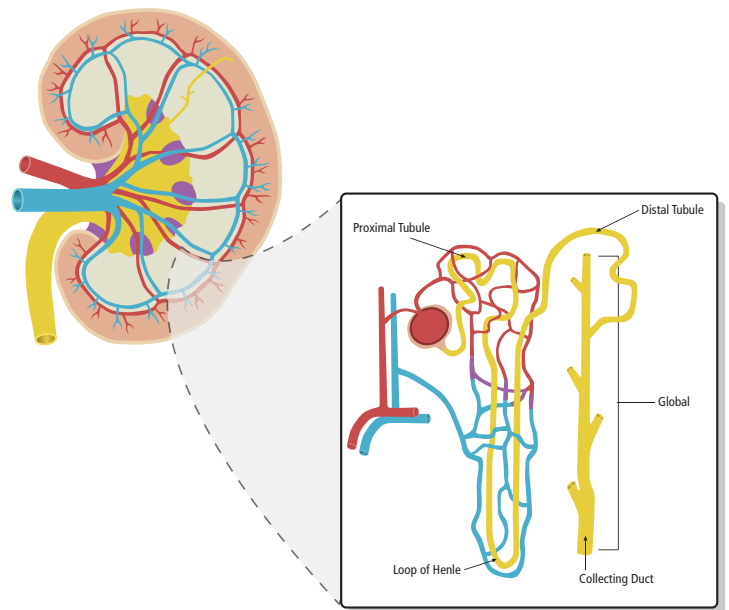
## Representative Nephrotoxicity Products for Toxicology Applications

Clusterin (rat)*	NGAL (human)†	B2M (rat)†
TIM-1/KIM-1/HAVCR (rat)*	VEGF (rat)	IL-18 (human)†
Argutus AKI Test® (rat)* (αGST, GSTYb1, RPA-1)	Kidney Injury Panel 1 (rat)* (Lipocalin-2, Osteopontin, Albumin, TIM-1)	Kidney Injury Panel 2 (rat)† (Lipocalin-2, Osteopontin, Albumin, TIM-1, αGST, Cystatin C, Clusterin)

\* Qualified Kit    † In Development    ‡ Prototype

For a complete list of products please check our website at [www.mesoscale.com](http://www.mesoscale.com)  
Please inquire about other human injury markers.

Traditional clinical markers for kidney injury, such as BUN and serum creatinine, are not sensitive enough to detect subtle kidney damage and often do not correlate to damage measured by histopathology. The Food and Drug Administration (FDA) and The European Agency for the Evaluation of Medicinal Products (EMA) have acknowledged the need for novel biomarkers to identify kidney injury earlier in the drug discovery process. MSD, working with members of the Critical Path Institute (C-Path) and Health and Environmental Sciences Institute (HESI) consortiums, has developed singleplex and multiplex panels to rapidly screen large numbers of compounds for drug induced kidney toxicity. Our Argutus AKI Test, which includes αGST, GSTYb1, and RPA-1, was developed in conjunction with Argutus Medical (formerly Biotrin International). HESI samples tested on this panel have been shown to correlate well to previous measurements. Our Kidney Injury Panel 1 (rat) includes assays for albumin, TIM-1 (a.k.a. KIM-1 or HAVCR), NGAL (a.k.a. Lipocalin-2) and osteopontin. All four analytes have been shown by C-Path Predictive Safety Testing Consortium (PSTC) members to be indicative of early detection of kidney damage. Many of these analytes have been submitted for approval to the FDA as biomarkers for pre-clinical nephrotoxicity measurements. MSD continues to be at the forefront of developing new panels and singleplex assays for toxicity applications.

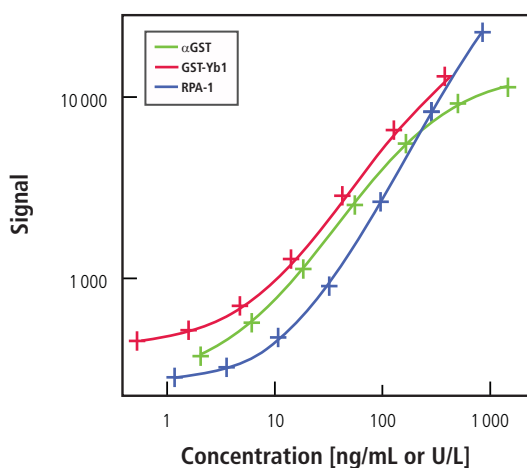


# Nephrotoxicity

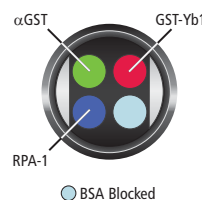
## Argutus AKI Test®

The Argutus AKI (Acute Kidney Injury) Test measures three biomarkers of drug-induced nephrotoxicity:  $\alpha$ GST, GST Yb1, and RPA-1. The HESI recently submitted these biomarkers to the FDA as predictive biomarkers. By combining measurements of  $\alpha$ GST (proximal tubular injury), GST Yb1 (distal tubular injury) and RPA-1 (collecting duct injury), the specific location of kidney injury can be inferred.<sup>1,2</sup> Collectively, these biomarkers offer a quantitative assessment of the location and extent of acute kidney injury. We qualified this panel according to the principles outlined in "Fit-for-Purpose Method Development and Validation for Successful Biomarker Measurement" by Lee, J.W., et. al<sup>3</sup>. The qualification procedure involved measurement of control samples on multiple days, establishment of limits of quantitation, spike recovery, dilutional linearity, and measurement of control and treated samples.

### $\alpha$ GST, GSTYb1, and RPA-1



	Analyte		
	$\alpha$ GST (ng/mL)	GSTYb1 (ng/mL)	RPA-1 (U/L)
LLOD	0.87	0.71	2.0
LLOQ	6.6	12.4	30
ULOQ	196	333	800



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

The LLOQ (Lower Limit of Quantitation) is determined as the lowest concentration where the %CV of the calculated concentration is less than 20% and the percent recovery of the standard is between 80% and 120%. For RPA-1, the percent recovery is between 75% and 125%.

The ULOQ (Upper Limit of Quantitation) is determined as the highest concentration where the %CV of the calculated concentration is less than 20% and the percent recovery of the standard is between 80% and 120%. For RPA-1, the percent recovery is between 75% and 125%.

### Catalog Numbers for Argutus AKI Test Qualified Kit

	Kit Size	Catalog Numbers
Rat	1 Plate	K15156C-1
	5 Plate	K15156C-2
	25 Plate	K15156C-4

### Precision: Multi-Day Study

Three levels of controls (high, mid, and low) were prepared for precision studies. Controls were made by spiking papilla medulla extract into rat urine.

	Control	Plates	Concentration	Average Intra-plate % CV	Inter-day % CV
$\alpha$ GST (ng/mL)	High	8	79.1	2.9	6.5
	Mid	8	61.7	3.3	6.6
GSTYb1 (ng/mL)	High	8	86.5	4.1	5.1
	Low	8	28.4	4.3	5.7
RPA-1 (U/L)	High	8	471	4.7	5.6
	Mid	8	89.6	5.5	6.3

### Samples

Rat urine samples from known injury inducing drugs were run on the Argutus AKI Test at 5-fold dilution. Measurements in italics were below the assay LLOQ at a 5-fold dilution. Measurements in bold were made at a 20-fold dilution of the urine samples.

MSD Assay Kits			$\alpha$ GST		GSTYb1		RPA-1	
Nephro-toxicant	Associated Biomarker	Sample ID	Concentration (ng/mL)	Concentration CV	Concentration (ng/mL)	Concentration CV	Concentration (U/L)	Concentration CV
Control	None	B671	11.7	2.9	<i>13.4</i>	35.6	714	1.4
		B672	26.8	16.0	<i>16.7</i>	20.5	1382	2.6
NPAA	RPA-1	B673	<i>7.1</i>	<i>3.1</i>	<i>29.0</i>	<i>13.1</i>	<b>8700</b>	<b>9.8</b>
		B674	<i>7.3</i>	<i>13.3</i>	<i>14.9</i>	<i>34.7</i>	<b>5482</b>	<b>5.5</b>
Tenidap	Clusterin	B675	20.3	7.6	<i>23.4</i>	7.2	4363	9.1
		B676	7.8	7.8	<i>19.5</i>	37.6	1657	5.2
Cisplatin	Alpha-GST	B680	69.2	3.9	<i>38.0</i>	36.6	757	6.3
		B681	176	2.6	<i>43.3</i>	30.2	704	4.4

### Example Protocol

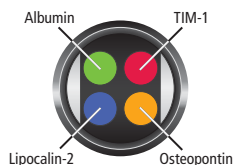
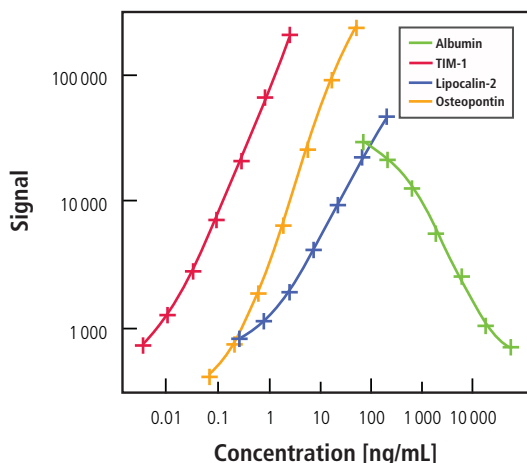
1. Add 150  $\mu$ L Blocking Solution, incubate 1 hour at RT.
2. Wash with PBS-T. Add 25  $\mu$ L Diluent, add 25  $\mu$ L Calibrator/Sample, incubate 2 hours at RT.
3. Wash with PBS-T. Add 25  $\mu$ L Detection Antibody, incubate 2 hours at RT.
4. Wash with PBS-T. Add 150  $\mu$ L Read Buffer T, read.

# Nephrotoxicity

## Kidney Injury Panel 1 (rat)

Our Kidney Injury Panel 1 (rat) measures albumin, TIM-1, lipocalin-2 (NGAL), and osteopontin (OPN) in urine samples. Albumin is very abundant (10's of  $\mu\text{g/mL}$ ); we measure it by a competitive immunoassay to enable multiplexing with lower-abundance biomarkers. The other analytes are measured by sandwich immunoassays. Albumin and TIM-1 are now biomarkers that the FDA and EMEA will consider during their drug review process.<sup>4</sup> Lipocalin-2 and osteopontin have also been implicated as important biomarkers of nephrotoxicity and acute kidney injury and have been studied extensively by CPATH members.<sup>5,6,7,8</sup> We qualified this panel according to the principles outlined in "Fit-for-Purpose Method Development and Validation for Successful Biomarker Measurement" by Lee, J.W., et. al<sup>3</sup>. The qualification procedure involved measurement of control samples on multiple days, establishment of limits of quantitation, spike recovery, dilutional linearity, and measurement of control and treated samples.

### Albumin, TIM-1, Lipocalin-2, and Osteopontin



	Analyte		
	TIM-1	Lipocalin-2	Osteopontin
LLOD (ng/mL)	0.0013	0.250	0.066
LLOQ (ng/mL)	0.02	1.56	0.39
ULOQ (ng/mL)	1.88	150	37.5

	Analyte
	Albumin
LLOD (ng/mL)	108
LLOQ (ng/mL)	781
Upper Limit (ng/mL)	28481

LLOD is defined as 80% of the maximum signal for albumin.

The upper limit of albumin was based on 20% of the signal over the highest calibrator. The LLOQ for albumin is the point on the standard curve where the %CV of the standard is less than 20% and the percent recovery of the standard was between 80% and 120%.

### Catalog Numbers for Kidney Injury Panel 1 Qualified Kit

	Kit Size	Catalog Numbers
Rat	1 Plate	K15162C-1
	5 Plate	K15162C-2
	25 Plate	K15162C-4

### Precision: Multi-Day Study

A multi-day, multi-plate study over 13 plates was performed to show reproducibility. In addition to the standard curves, control samples (high, mid and low) were measured on each plate. Each sample was run in triplicate. The average intra-plate %CV and inter-plate %CV of the concentrations are shown below.

	Control	Plates	Concentration (ng/mL)	Average Intra-plate % CV	Inter-plate % CV
Albumin	High	13	12040	3.5	5.8
	Mid	13	4137	3.4	4.5
	Low	13	1765	3.1	7.2
TIM-1	High	13	0.960	3.5	4.4
	Mid	13	0.329	3.8	5.7
	Low	13	0.133	4.2	9.9
Lipocalin-2	High	13	194	8.4	10.4
	Mid	13	91.5	6.4	9.0
	Low	13	42.3	5.6	8.2
Osteopontin	High	13	15.6	5.0	6.4
	Mid	13	6.06	4.5	6.4

### Example Protocol

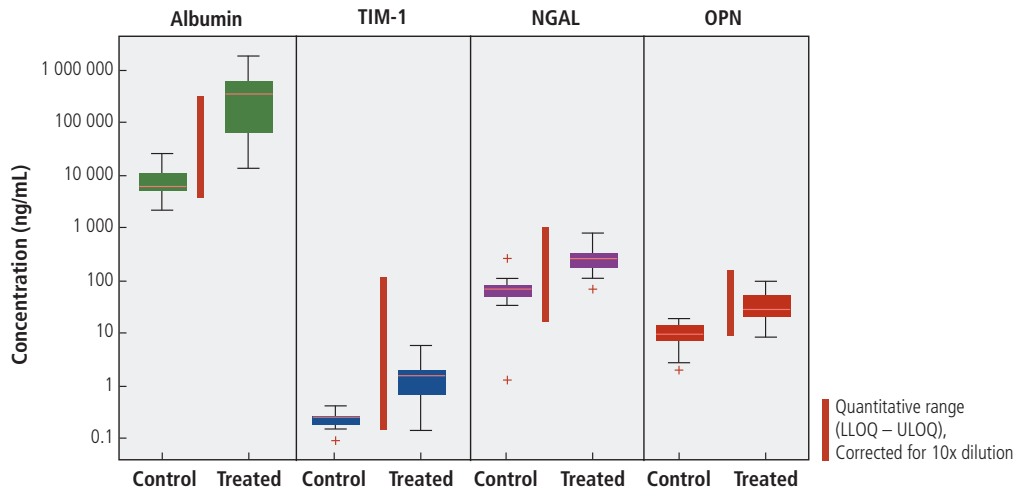
1. Add 150  $\mu\text{L}$  Blocking Solution, incubate 1 hour at RT.
2. Wash with PBS-T. Add 50  $\mu\text{L}$  sample or standard premixed with Albumin tracer, incubate 2 hours at RT.
3. Wash with PBS-T. Add 25  $\mu\text{L}$  Detection Antibody, incubate 2 hours at RT.
4. Wash with PBS-T. Add 150  $\mu\text{L}$  Read Buffer T, read.

# Nephrotoxicity

## Kidney Injury Panel 1 (rat)

### Samples

Rat urine samples were assayed at 10-fold dilution on the Kidney Injury Panel 1. The “treated” samples tested were from rats exposed to known nephrotoxicants prior to sample collection. Some treated samples that assayed above the ULOQ were assayed again at 40-fold dilution. For all of the analytes on the panel, significant correspondence between histo-pathology score and abundance was observed (data not shown).



### Catalog Numbers for Kidney Injury Panel 1 Qualified Kit

	Kit Size	Catalog Numbers
Rat	1 Plate	K15162C-1
	5 Plate	K15162C-2
	25 Plate	K15162C-4

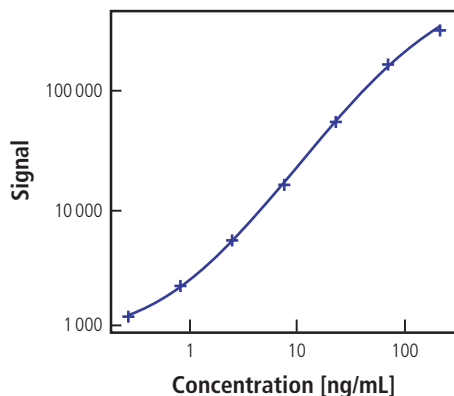
### Example Protocol

1. Add 150  $\mu$ L Blocking Solution, incubate 1 hour at RT.
2. Wash with PBS-T. Add 50  $\mu$ L sample or standard premixed with Albumin tracer, incubate 2 hours at RT.
3. Wash with PBS-T. Add 25  $\mu$ L Detection Antibody, incubate 2 hours at RT.
4. Wash with PBS-T. Add 150  $\mu$ L Read Buffer T, read.

## Rat Clusterin

Urinary clusterin has been accepted by the FDA and EMEA as a biomarker to be considered during drug review.<sup>4</sup> The MSD assay for clusterin in rat urine samples is currently available as a singleplex qualified kit that offers excellent sensitivity and dynamic range. We qualified this panel according to the principles outlined in “Fit-for-Purpose Method Development and Validation for Successful Biomarker Measurement” by Lee, J.W. et al<sup>1</sup>. The qualification procedure involved measurement of control samples on multiple days, establishment of limits of quantitation, spike recovery, dilutional linearity, and measurement of control and treated samples.

### Clusterin



	Clusterin (ng/mL)
LLOD	0.017
LLOQ	0.82
ULOQ	150

### Catalog Numbers for Clusterin Qualified Kit

	Kit Size	Catalog Numbers
Rat	1 Plate	K153HXC-1
	5 Plate	K153HXC-2
	25 Plate	K153HXC-4

### Example Protocol

1. Add 150  $\mu$ L Blocking Solution, incubate 1 hour at RT.
2. Wash with PBS-T. Add 25  $\mu$ L Assay Diluent, add 25  $\mu$ L Calibrator/ Sample, incubate 2 hours at RT.
3. Wash with PBS-T. Add 25  $\mu$ L Detection Antibody, incubate 2 hours at RT.
4. Wash with PBS-T. Add 150  $\mu$ L Read Buffer T, read.



# Nephrotoxicity

## Rat Clusterin

### Precision: Multi-Day Study

High, mid, and low controls were made by spiking recombinant protein into stock calibrator diluent. The controls were run in quadruplicate on each of 10 plates run across three days. The controls were run at a 10-fold dilution.

	Control	Plates	Concentration (ng/mL)	Average Intra-plate % CV	Inter-day % CV
Clusterin	High	10	139	7.27	10.4
	Mid	10	17.5	4.30	6.06
	Low	10	1.90	4.80	6.57

## Spike Recovery

Rat urine samples were spiked with calibrators at multiple values throughout the range of the assay. For the Argutus AKI Test panel and the Rat Clusterin assay, the Calibrators were spiked into 5-fold diluted rat urine. For the Kidney Injury Panel 1 panel the spiked samples were tested at a 10-fold dilution into diluent. The recombinant Osteopontin may be bound to proteins in the urine, making it under-recover.

$\% \text{ Recovery} = \text{measured} / \text{expected} \times 100$

Analyte	% Recovery in Urine Sample			Kit
	High Conc.	Mid Conc.	Low Conc.	
$\alpha$ GST	100	94	105	Argutus AKI Test
GST-Yb1	95	105	94	Argutus AKI Test
RPA-1	95	102	98	Argutus AKI Test
Albumin	102	101	98	Kidney Injury Panel 1 (rat)
TIM-1	95	101	104	Kidney Injury Panel 1 (rat)
Lipocalin-2	103	106	99	Kidney Injury Panel 1 (rat)
Osteopontin	63	53	38	Kidney Injury Panel 1 (rat)
Clusterin	94	93	97	Rat Clusterin

## Dilutional Linearity

Serial dilutions of rat urine samples were tested to assess linearity. Percent recovery is calculated as the measured concentration divided by the concentration for the previous dilution (expected).

$\% \text{ Recovery} = (\text{measured} \times \text{dilution factor}) / \text{expected} \times 100$

Analyte	% Dilutional Linearity in Urine Sample						Kit
	5	10	20	40	80	160	
$\alpha$ GST	102	101	-	-	-	-	Argutus AKI Test
GST-Yb1	105	84	86	-	-	-	Argutus AKI Test
RPA-1	86	80	89	96	100	94	Argutus AKI Test
Albumin	-	105	103	100	-	-	Kidney Injury Panel 1 (rat)
TIM-1	-	109	101	98	-	-	Kidney Injury Panel 1 (rat)
Lipocalin-2	-	145	112	103	-	-	Kidney Injury Panel 1 (rat)
Osteopontin	-	193	187	179	-	-	Kidney Injury Panel 1 (rat)
Clusterin	-	125	119	107	106	-	Rat Clusterin

# Myotoxicity

## Representative Myotoxicity Products for Toxicology Applications

Skeletal Troponin I (sTnI) (rat)\*

Cardiac Injury Panel 2 (rat)\*  
(cTnI, cTnT and FABP3)

Cardiac Injury Panel 3 (rat)\*  
(cTnI, cTnT, FABP3, Myl3)

Muscle Injury Panel 1 (rat)\*  
(cTnI, cTnT, sTnI, FABP3, Myl3)

Muscle Injury Panel 2 (rat)\*  
(Parvalbumin, TIMP-1, CK)

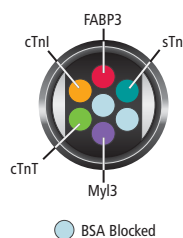
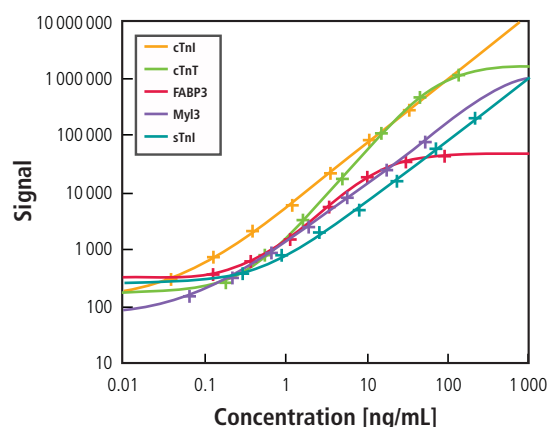
\* Qualified Kit † In Development

For a complete list of products please check our website at [www.mesoscale.com](http://www.mesoscale.com)

## Muscle Injury Panel 1 (rat)

Muscle Injury Panel 1 (rat) measures five important biomarkers of drug-induced myotoxicity.<sup>9,10</sup> The biomarkers on the panel were selected for their ability to collectively distinguish myocardial damage from damage to skeletal muscle or other tissues. Using these biomarkers, damage to skeletal muscle can be additionally stratified into slow and fast-twitch types. The selectivity of the assays available from MSD is demonstrated in the specificity section below. We qualified this panel according to the principles outlined in "Fit-for-Purpose Method Development and Validation for Successful Biomarker Measurement" by Lee, J.W., et al<sup>3</sup>. The qualification procedure involved measurement of control samples on multiple days, establishment of limits of quantitation, spike recovery, dilutional linearity, and measurement of control and treated samples.

### cTnI, cTnT, FABP3, Myl3, and sTnI



### Catalog Numbers for Muscle Injury Panel 1 Qualified Kit

	Kit Size	Catalog Numbers
Rat	1 Plate	K15181C-1
	5 Plate	K15181C-2
	25 Plate	K15181C-4

	Analyte				
	cTnI	cTnT	FABP3	Myl3	sTnI
LLOD (ng/mL)	0.010	0.079	0.069	0.023	0.148
LLOQ (ng/mL)	0.098	0.488	0.391	0.054	0.781
ULOQ (ng/mL)	20.0	100	15.0	44.0	160

### Samples

Serum, heparin plasma, and EDTA plasma samples collected from normal Sprague-Dawley rats were tested at 2-fold dilution on the Muscle Injury Panel 1 (rat). Shown here are the median and range of concentrations for each sample set. Skeletal Troponin I was below the quantitative range for all samples.

Sample	Statistic	cTnI	cTnT	FABP3	Myl3	sTnI
Serum	Median (ng/mL)	1.59	0.83	19.84	0.61	< 1.56
	Range (ng/mL)	0.44 - 3.53	< 0.976 - 2.04	5.28 - > 30	0.25 - 1.05	< 1.56
	N	10	10	10	10	10
Heparin Plasma	Median (ng/mL)	1.81	1.06	28.35	0.55	< 1.56
	Range (ng/mL)	0.37 - 4.10	< 0.976 - 2.79	4.12 - > 30	0.16 - 1.16	< 1.56
	N	10	10	10	10	10
EDTA Plasma	Median (ng/mL)	2.31	1.28	35.63	0.81	< 1.56
	Range (ng/mL)	1.45 - 4.05	< 0.976 - 2.80	19.6 - > 30	0.43 - 1.17	< 1.56
	N	10	10	10	10	10

### Example Protocol

1. Add 150  $\mu$ L Blocking Solution, incubate 1 hour at RT.
2. Wash with PBS-T. Add 25  $\mu$ L Antibody Diluent, add 25  $\mu$ L Calibrator/Sample, incubate 2 hours at RT.
3. Wash with PBS-T. Add 25  $\mu$ L Detection Antibody, incubate 2 hours at RT.
4. Wash with PBS-T. Add 150  $\mu$ L Read Buffer T, read.

# Myotoxicity

## Muscle Injury Panel 1 (rat)

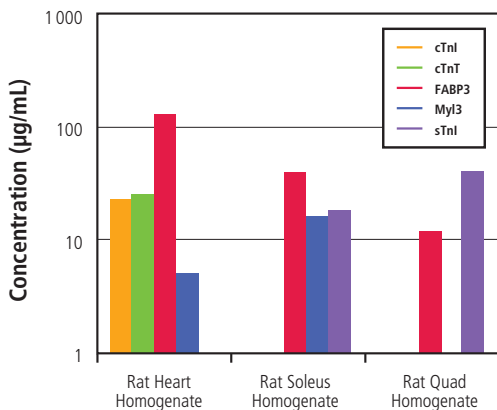
### Precision: Multi-Day Study

Control samples were measured on 9 plates across 3 days. The controls were run in triplicate or quadruplicate on each plate. Normal rat serum, rat soleus homogenate, and assay Calibrators are used to make control samples. The high control contains 25% normal rat serum and Calibrators. The mid control contains rat soleus homogenate and Calibrators. The low control contains only the assay Calibrators. The average intra-plate %CV and inter-plate %CV of the concentrations are shown below.

	Control	Plates	Concentration (ng/mL)	Average Intra-plate % CV	Inter-plate % CV
cTnI	High	9	9.36	3.8	5.9
	Mid	9	1.31	3.9	6.0
	Low	9	0.28	4.0	5.1
cTnT	High	9	35.4	3.9	5.0
	Mid	9	6.46	2.5	4.0
	Low	9	1.06	3.3	4.6
FABP3	High	9	13.9	5.9	7.5
	Mid	9	8.44	5.5	7.6
	Low	9	2.85	2.8	4.8
MyI3	High	9	30.3	5.9	7.0
	Mid	9	3.01	3.8	6.9
	Low	9	0.32	6.0	7.1
sTnI	High	9	108	4.3	5.8
	Mid	9	17.4	2.4	3.1
	Low	9	2.88	3.5	10.5

### Specificity

Tissue homogenates from heart, fast twitch, and slow twitch muscle were tested at 100X, 1000X, and 10000X sample dilution. The assays for cardiac troponins were positive for cardiac homogenates and negative for other muscle homogenates, demonstrating specificity for cardiac tissue. The assay for skeletal Troponin I was specific for fast and slow twitch skeletal muscle. The assay measured FABP3 in cardiac muscle and skeletal muscle. The slow twitch muscle was positive for MyI3, while approximately 200X less MyI3 was measured in fast twitch. Concentrations below are corrected for dilution.



Sample Group	cTnI		cTnT		FABP3		MyI3		sTnI	
	Sample Dilution	Conc. (µg/mL)	Sample Dilution	Conc. (µg/mL)	Sample Dilution	Conc. (µg/mL)	Sample Dilution	Conc. (µg/mL)	Sample Dilution	Conc. (µg/mL)
Rat Heart Homogenate	1000	22.6	1000	25.1	10000	125.2	1000	5.0	100	< LLOD
Rat Soleus Homogenate (slow twitch)	100	< LLOD	100	< LLOD	10000	38.8	1000	16.4	1000	18.1
Rat Quad Homogenate (fast twitch)	100	< LLOD	100	< LLOD	1000	12.2	100	0.08	1000	40.9

# Myotoxicity

## Spike Recovery

Rat serum, EDTA plasma and heparin plasma samples were spiked with calibrators at multiple values throughout the range of the assay. For the Rat sTnI assay, the Calibrators were spiked into 2-fold diluted samples. For the Muscle Injury Panel 2 assay, the Calibrators were spiked into 20-fold diluted plasma and 100-fold diluted serum samples.

$$\% \text{ Recovery} = \text{measured} / \text{expected} \times 100$$

Analyte	% Recovery									Kit
	Serum			EDTA Plasma			Heparin Plasma			
	High Conc.	Mid Conc.	Low Conc.	High Conc.	Mid Conc.	Low Conc.	High Conc.	Mid Conc.	Low Conc.	
cTnI	100	106	105	96	111	97	103	104	93	Cardiac Injury Panel 3 (rat)
	116	116	111	128	116	108	134	123	105	Muscle Injury Panel 1 (rat)
cTnT	88	89	89	83	90	99	113	112	106	Cardiac Injury Panel 3 (rat)
	77	86	86	89	78	71	100	104	95	Muscle Injury Panel 1 (rat)
FABP3	108	96	105	100	108	104	103	97	115	Cardiac Injury Panel 3 (rat)
	114	113	98	100	93	91	90	99	87	Muscle Injury Panel 1 (rat)
MyI3	110	106	108	120	114	104	112	105	96	Cardiac Injury Panel 3 (rat)
	115	120	105	136	131	107	138	128	107	Muscle Injury Panel 1 (rat)
sTnI	108	106	106	100	96	91	109	101	96	Rat Skeletal Troponin I (sTnI)
	101	109	100	107	108	99	110	107	100	Muscle Injury Panel 1 (rat)
Parvalbumin	119	123	120	122	114	112	136	121	117	Muscle Injury Panel 2 (rat)
TIMP-1	111	97	90	123	105	101	123	101	97	Muscle Injury Panel 2 (rat)
CK	93	107	123	96	108	107	92	99	105	Muscle Injury Panel 2 (rat)

## Dilutional Linearity

Serial dilutions of rat serum, EDTA plasma and heparin plasma samples were tested to assess linearity. For Cardiac Injury Panel 3 markers iso-proterenol treated rat serum was used. Percent recovery is calculated as the measured concentration divided by the concentration for the previous dilution (expected).

$$\% \text{ Recovery} = (\text{measured} \times \text{dilution factor}) / \text{expected} \times 100$$

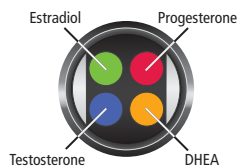
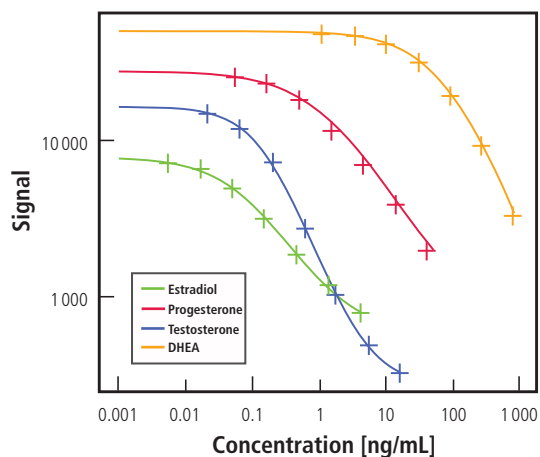
Analyte	% Dilutional Linearity									Kit
	Serum			EDTA Plasma			Heparin Plasma			
	2-fold	4-fold	8-fold	2-fold	4-fold	8-fold	2-fold	4-fold	8-fold	
cTnI	82	84	87	90	91	93	97	90	94	Cardiac Injury Panel 3 (rat)
	101	90	92	108	109	<LLOD	98	96	89	Muscle Injury Panel 1 (rat)
cTnT	99	83	77	104	88	<LLOD	92	93	<LLOQ	Cardiac Injury Panel 3 (rat)
	<LLOD	<LLOD	<LLOD	<LLOD	<LLOD	<LLOD	101	<LLOD	<LLOD	Muscle Injury Panel 1 (rat)
FABP3	>ULOQ	NA	107	>ULOQ	NA	81	>ULOQ	NA	85	Cardiac Injury Panel 3 (rat)
	NA	77	88	93	100	<LLOD	NA	93	104	Muscle Injury Panel 1 (rat)
MyI3	91	86	82	82	83	91	99	91	90	Cardiac Injury Panel 3 (rat)
	86	92	96	99	94	89	102	96	107	Muscle Injury Panel 1 (rat)
sTnI	79	87	92	104	91	92	88	95	97	Muscle Injury Panel 1 (rat)
	20-fold	40-fold	80-fold	20-fold	40-fold	80-fold	20-fold	40-fold	80-fold	
Parvalbumin	-	-	-	75	82	97	-	-	-	Muscle Injury Panel 2 (rat)
TIMP-1	-	-	-	102	87	<LLOQ	-	-	-	Muscle Injury Panel 2 (rat)
CK	-	-	-	82	110	<LLOQ	-	-	-	Muscle Injury Panel 2 (rat)

# Reproductive Toxicity

## Steroid Hormone Panel 1 (human, mouse, rat)

Toxicity to the endocrine system is a common mode of failure for drugs in the clinic, and steroid hormones are recognized as important biomarkers in this area. The Steroid Hormone Panel 1 is designed to measure critical steroid hormones in serum and plasma from human, mouse, and rat.

### Estradiol, Progesterone, Testosterone and DHEA



	Analyte			
	Estradiol (ng/mL)	Progesterone (ng/mL)	Testosterone (ng/mL)	DHEA (ng/mL)
LLOD	0.023	0.306	0.046	13.7
LLOQ	0.049	0.617	0.198	29.6
ULOQ	4.0	50.0	5.33	800

### Catalog Numbers for Steroid Hormone Panel 1

	Kit Size	Catalog Numbers
Human/	1 Plate	K15178C-1
Mouse/	5 Plate	K15178C-2
Rat	25 Plate	K15178C-4

### Example Protocol

1. Add 150  $\mu$ L Blocking Solution, incubate 1 hour at RT.
2. Wash, add 50  $\mu$ L standard/sample, incubate 2 hours at RT.
3. Add 15  $\mu$ L tracer, incubate 30 min at RT.
4. Wash, add 150  $\mu$ L Read Buffer T, read.

# Inflammation and Growth Factors

## Representative Inflammation and Growth Factor Markers

E-Cadherin (human)	HGF (human)	E-Selectin (human)
P-Cadherin (human)	sICAM-1 (human)	P-Selectin (human)
c-Kit (human)	sICAM-3 (human)	TGF-β1 (human)
CRP (human)	KDR (sVEGFR2) (human)	Thrombomodulin (human)
bFGF (human)	PIGF (human)	sVCAM-1 (human)
sFlt-1 (sVEGFR1) (human)	SAA (human)	VEGF (human, mouse/rat)

## Multiplex Panels

Vascular Injury Panel I (human) (sICAM-3, E-Selectin, P-Selectin, Thrombomodulin)	Inflammation Panel 1 (rat)* (Lipocalin-2, TSP-1, TIMP-1, MCP-1)	Growth Factor Panel II (human) (c-Kit, KDR)
Vascular Injury Panel II (human) (CRP, sICAM-1, sVCAM-1, SAA)	Growth Factor Panel I (human) (bFGF, VEGF, sFlt-1, PIGF)	Acute Phase Protein Panel 1 (rat)* (A2M, AGP)
c-Kit (pY721), Total c-Kit		

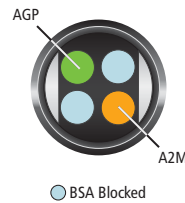
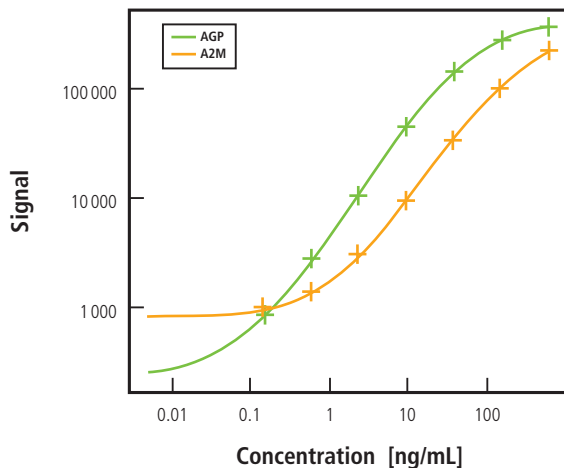
\* Qualified Kit

For a complete list of products please check our website at [www.mesoscale.com](http://www.mesoscale.com)

Circulating levels of cytokines and other inflammation markers have been studied in patients with chronic renal failure<sup>11</sup> and persistent inflammation has been the suspect of other pathophysiologic alterations in chronic kidney disease.<sup>12</sup> Biomarkers of inflammation have also emerged as potential preclinical indicators of heart failure.<sup>13</sup>

## Acute Phase Protein Panel 1 (rat)

### AGP and A2M



	Analyte	
	AGP (ng/mL)	A2M (ng/mL)
LLOD	0.012	0.036
LLOQ	0.146	1.00
ULOQ	150	400

### Catalog Numbers for Acute Phase Protein Panel 1 (rat) Qualified Kit

	Kit Size	Catalog Numbers
Rat	1 Plate	K15175C-1
	5 Plate	K15175C-2
	25 Plate	K15175C-4

### Precision: Multi-Day Study

High, mid, and low controls were defined that are rat serum samples diluted 10000-fold into assay diluent. Samples were chosen that span a large proportion of the dynamic range for each assay. These controls were run in triplicate on each of 9 plates run across 3 days of testing. The concentration values shown below are corrected for the 10000-fold dilution.

	Control	Plates	Concentration (ng/mL)	Average Intra-plate % CV	Inter-plate % CV
AGP	High	9	42706	5.3	9.9
	Mid	9	8427	4.2	10.1
	Low	9	408	3.4	7.1
A2M	High	9	215203	6.2	8.6
	Mid	9	20468	2.4	6.3
	Low	9	1870	3.1	4.0

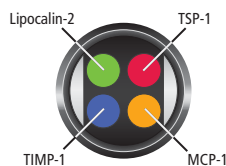
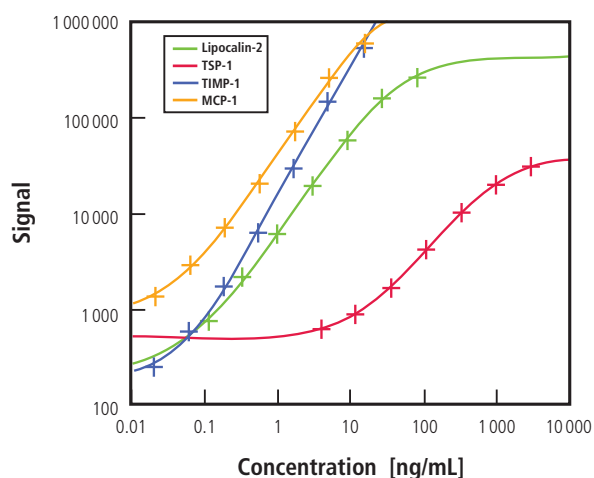
### Example Protocol

1. Add 25 µL Assay Diluent, incubate 30 min at RT.
2. Add 25 µL of Calibrators/Sample, incubate 2 hours at RT.
3. Wash with PBS-T. Add 25 µL of Detection Antibody, incubate 2 hours at RT.
4. Wash with PBS-T. Add 150 µL of Read Buffer, read.

# Inflammation and Growth Factors

## Rat Inflammation Panel 1

### Lipocalin-2, TSP-1, TIMP-1, and MCP-1



	Analyte			
	Lipocalin-2 (ng/mL)	TSP-1 (ng/mL)	TIMP-1 (ng/mL)	MCP-1 (ng/mL)
LLOD	0.0050	3.7	0.0046	0.0028
LLOQ	0.050	35	0.10	0.050
ULOQ	25	1500	12	12

### Example Protocol

1. Add 25  $\mu$ L Diluent, incubate 30 minutes at RT.
2. Add 25  $\mu$ L Calibrator/Sample, incubate 2 hours at RT.
3. Wash with PBS-T. Add 25  $\mu$ L Detection Antibody, incubate 2 hours at RT.
4. Wash with PBS-T. Add 150  $\mu$ L Read Buffer T, read.

### Precision: Multi-Day Study

Control samples of high, mid, and low levels of each analyte were measured on each plate. The controls were run in quadruplicate on each of 21 plates run across multiple days ( $n > 3$ ).

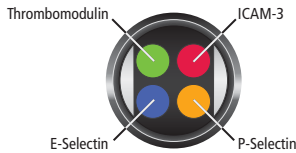
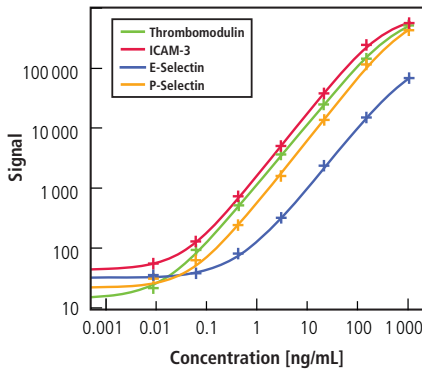
	Control	Plates	Concentration (ng/mL)	Average Intra-plate % CV	Inter-plate % CV
Lipocalin-2	High	21	12.8	3.7	6.0
	Mid	21	1.58	3.5	6.9
	Low	21	0.364	3.4	7.5
TSP-1	High	21	1083	9.8	12.5
	Mid	21	658	6.4	10.6
	Low	21	73.8	3.3	7.0
TIMP-1	High	21	6.55	6.8	9.0
	Mid	21	0.914	3.1	6.5
	Low	21	0.404	4.0	6.5
MCP-1	High	21	9.48	4.1	6.8
	Mid	21	0.836	5.8	8.7
	Low	21	0.0945	6.2	10.7

# Inflammation and Growth Factors

## Vascular Injury Panel I (human)

In pre-clinical safety studies, drug-induced vascular injury is of great concern. Endothelial cell dysfunction is considered one of the early indicators of vascular disorders. Drugs may trigger inflammatory signaling cascades and the upregulation of soluble adhesion molecules by activated endothelial cells. Since vascular injury involves multiple mediators and cell types, a multiplex panel of validated biomarkers may serve as a useful predictor of activation of inflammatory cascades in the development of vascular injury.

### Thrombomodulin, ICAM-3, E-Selectin, and P-Selectin



LLOD (ng/mL)	Analyte	
	Thrombomodulin	ICAM-3
	0.017	0.066

LLOD (ng/mL)	Analyte	
	E-Selectin	P-Selectin
	0.34	0.097

### Spike Recovery

	% Recovery			
	Thrombomodulin	ICAM-3	E-Selectin	P-Selectin
Serum	87	77	97	89
EDTA Plasma	69	71	69	85
Heparin Plasma	83	76	91	91

Pooled plasma and serum samples were spiked with known quantities of Thrombomodulin, ICAM-3, E-selectin, P-selectin at different levels and measured using MSD Vascular Injury Panel I.

### Catalog Numbers for Human Vascular Injury I

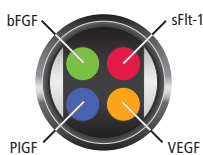
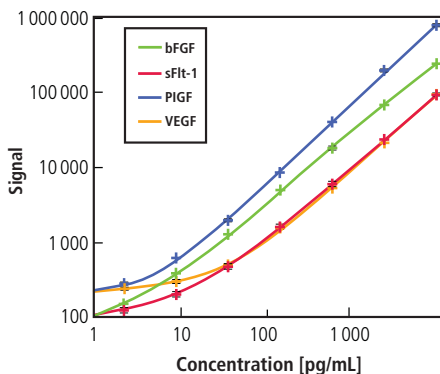
	Kit Size	Catalog Numbers
Human	1 Plate	K15135C-1
	5 Plate	K15135C-2
	20 Plate	K15135C-3

### Example Protocol

1. Add 150  $\mu$ L Blocking Solution, incubate 1 hour at RT.
2. Wash with PBS-T. Add 40  $\mu$ L Assay Diluent, add 10  $\mu$ L Calibrators/Sample, incubate 2 hours at RT.
3. Wash with PBS-T. Add 25  $\mu$ L of Detection Antibody, incubate 1 hour at RT.
4. Wash with PBS-T. Add 150  $\mu$ L of Read Buffer T, read.

## Growth Factor Panel I (human)

### bFGF, sFlt-1, PlGF, and VEGF



LLOD (pg/mL)	Analyte			
	bFGF	sFlt-1	PlGF	VEGF
	2.2	3.7	0.96	6.4

### Spike Recovery

	% Recovery			
	bFGF	sFlt-1	PlGF	VEGF
Serum	99	105	111	114
EDTA Plasma	77	94	98	95

Pooled plasma and serum samples were spiked with known quantities of bFGF, sFlt-1, PlGF, VEGF at different levels and measured using MSD Growth Factor Panel I.

### Catalog Numbers for Human Growth Factor I

	Kit Size	Catalog Numbers
Human	1 Plate	K15029C-1
	5 Plate	K15029C-2
	20 Plate	K15029C-3

### Example Protocol

1. Add 150  $\mu$ L Blocking Solution, incubate 1 hour at RT.
2. Wash with PBS-T. Add 25  $\mu$ L Assay Diluent, add 25  $\mu$ L Calibrators/Sample, incubate 2 hours at RT.
3. Wash with PBS-T. Add 25  $\mu$ L of Detection Antibody, incubate 2 hours at RT.
4. Wash with PBS-T. Add 150  $\mu$ L of Read Buffer T, read.



# Cytokines

## Representative Cytokine Markers

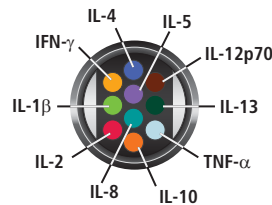
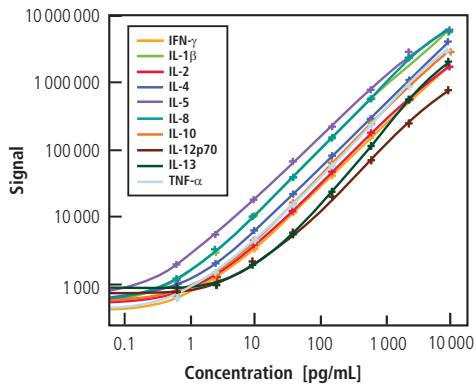
CINC-2 (rat) <sup>‡</sup>	IL-4 (human, mouse, rat)	IP-10 (human, mouse)	MMP-3 (human)
CINC-3 (rat) <sup>‡</sup>	IL-5 (human, mouse, rat)	I-TAC (human)	MMP-8 (human) <sup>‡</sup>
CTACK (human) <sup>‡</sup>	IL-6 (human, mouse, rat, canine)	KC/GRO/CINC (CXCL1) (mouse, rat)	MMP-9 (human, mouse) <sup>‡</sup>
ENA-78 (human) <sup>‡</sup>	IL-6R (human)	M-CSF (human, mouse) <sup>‡</sup>	MMP-10 (human)
Eotaxin (human)	IL-7 (human)	MCP-1 (human, mouse, rat, canine) <sup>‡</sup>	NGAL (human) <sup>‡</sup>
Eotaxin-3 (human)	IL-8 (human, canine)	MCP-2 (human) <sup>‡</sup>	RANTES (human, mouse)
G-CSF (human)	IL-10 (human, mouse, rat <sup>‡</sup> , canine <sup>‡</sup> )	MCP-4 (human)	SDF-1 $\alpha$ (human) <sup>‡</sup>
GM-CSF (human, mouse, rat, canine)	IL-12 (human, mouse)	MDC (human)	TARC (human)
GRO- $\alpha$ (human)	IL-12p40 (human, mouse)	MIG (human)	TGF- $\beta$ 1 (human)
IFN- $\alpha$ (human) <sup>‡</sup>	IL-12p70 (human, mouse)	MIP-1 $\alpha$ (human, mouse) <sup>‡</sup>	TIMP-1 (human, rat) <sup>‡</sup>
IFN- $\beta$ (human)	IL-13 (human, mouse, rat)	MIP-1 $\beta$ (human, mouse) <sup>‡</sup>	TIMP-2 (human) <sup>‡</sup>
IFN- $\gamma$ (human, mouse, rat, canine) <sup>‡</sup>	IL-15 (human)	MIP-2 (mouse) <sup>‡</sup>	TNF- $\alpha$ (human, mouse, rat, canine)
IL-1 $\alpha$ (human, rat)	IL-16 (human)	MIP-3 $\alpha$ (human, mouse, rat)	TNF-RI (human, mouse)
IL-1 $\beta$ (human, mouse, rat)	IL-17 (human, mouse)	MIP-3 $\beta$ (human) <sup>‡</sup>	TNF-RII (human, mouse)
IL-1ra (human) <sup>‡</sup>	IL-18 (human)	MMP-1 (human)	TSLP (human) <sup>‡</sup>
IL-2 (human, mouse, rat <sup>‡</sup> , canine)	IL-23 (human) <sup>‡</sup>	MMP-2 (human)	TWEAK (human) <sup>‡</sup>

<sup>‡</sup> Available through our Prototype Printing Services.

For a complete list of products please check our website at [www.mesoscale.com](http://www.mesoscale.com)

Multiplexed assays from MSD enable the detection of up to ten cytokines in a single well, in the same amount of time as a singleplex assay—using the same amount of sample!

## Human TH1/TH2 10-Plex Ultra-Sensitive



Analyte	Analyte									
	IFN- $\gamma$	IL-1 $\beta$	IL-2	IL-4	IL-5	IL-8	IL-10	IL-12p70	IL-13	TNF- $\alpha$
LLOD (pg/mL)	0.39	0.18	0.67	0.31	0.076	0.14	0.36	2.3	1.8	0.48

## Catalog Numbers for Human TH1 / TH2 10-Plex Ultra-Sensitive<sup>®</sup>

Human	Kit Size Catalog Numbers	
		1 Plate
	5 Plate	K15010C-2
	25 Plate	K15010C-4

<sup>®</sup> Visit [www.mesoscale.com](http://www.mesoscale.com) for Tissue Culture version of Human TH1 / TH2 10-Plex Kit.

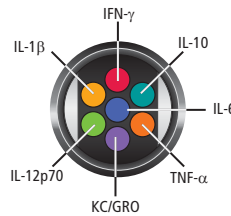
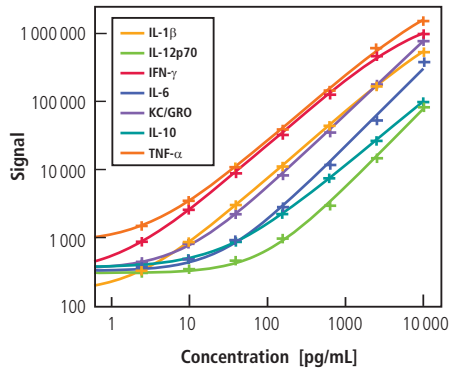
## Example Protocol

1. Add 25  $\mu$ L of Assay Diluent to plate, incubate 30 minutes.
2. Add 25  $\mu$ L of Calibrators/Sample, incubate 2 hours.
3. Wash with PBS-T. Add 25  $\mu$ L of Detection Antibody, incubate 1 hour.
4. Wash with PBS-T. Add 150  $\mu$ L Read Buffer T, Read.

# Cytokines

MSD offers several rodent cytokine assays in singleplex and multiplex formats. These assays enable the detection of up to 10 analytes in the same well. The wide dynamic range of the assays enables the measurement of native and disease state levels of cytokines in a single run. The panel on this page shows the sensitivity and range of one of our rodent cytokine assays. The panel was assayed with less than 25  $\mu\text{L}$  of sample.

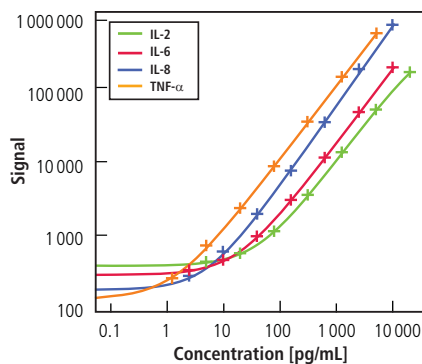
## Mouse ProInflammatory 7-Plex Ultra-Sensitive



LLOD (pg/mL)	Analyte						
	IL-1 $\beta$	IL-12p70	IFN- $\gamma$	IL-6	KC/GRO	IL-10	TNF- $\alpha$
	0.75	35	0.38	4.5	3.3	11	0.85

The exploration of the canine immune system has been driven by the dog's importance as an animal model for understanding the pathogenesis of several canine and human diseases. Dogs are used for drug toxicity trials, transplantation studies and in research of a variety of human diseases such as rheumatoid arthritis, cyclic neutropenia and systemic lupus erythematosus. The Canine ProInflammatory Panel 3 (IL-2, IL-6, IL-8 and TNF- $\alpha$ ) is a tool for researchers using canine models to study diseases, drug toxicity as well as vaccine development.

## Canine ProInflammatory Panel 3 Ultra-Sensitive



LLOD (pg/mL)	Analyte			
	IL-2	IL-6	IL-8	TNF- $\alpha$
	7.6	2.4	1.3	0.17

### Catalog Numbers for Mouse ProInflammatory 7-Plex Ultra-Sensitive <sup>®</sup>

Mouse	Kit Size	Catalog Numbers
	1 Plate	K15012C-1
5 Plate	K15012C-2	
25 Plate	K15012C-4	

<sup>®</sup> Visit [www.mesoscale.com](http://www.mesoscale.com) for Tissue Culture version of Mouse ProInflammatory 7-Plex Kit.

### Example Protocol

1. Add 25  $\mu\text{L}$  Assay Diluent, incubate 30 min at RT.
2. Add 10 to 25  $\mu\text{L}$  of Calibrator/ Sample incubate 2 hours at RT.
3. Wash with PBS-T. Add 25  $\mu\text{L}$  of Detection Antibody, incubate 1 to 2 hours at RT.
4. Wash with PBS-T. Add 150  $\mu\text{L}$  of Read Buffer, read.

### Catalog Numbers for Canine ProInflammatory Panel 3 Ultra-Sensitive

Canine	Kit Size	Catalog Numbers
	1 Plate	K15035C-1
5 Plate	K15035C-2	
25 Plate	K15035C-4	

### Example Protocol

1. Add 25  $\mu\text{L}$  Assay Diluent, incubate 30 min at RT.
2. Add 25  $\mu\text{L}$  of Calibrators/Sample, incubate 2 hours at RT.
3. Wash with PBS-T. Add 25  $\mu\text{L}$  of Detection Antibody, incubate 2 hours at RT.
4. Wash with PBS-T. Add 150  $\mu\text{L}$  of Read Buffer, read.

# Intracellular Signaling

## Representative Intracellular Signaling Markers

Akt (pS473) (human, mouse, rat)  
 Akt (pT308) (human, mouse)  
 Akt (Total) (human, mouse, rat)  
 GSK-3 $\alpha$  (pS21) (human)  
 GSK-3 $\beta$  (pS9) (human)

GSK-3 $\beta$  (Total) (human, mouse, rat)  
 IRS-1 (pS312) (human)  
 p70S6K (pT389) (human, mouse, rat)  
 p70S6K (pT421/pS424) (human, mouse, rat)  
 p70S6K (Total) (human, mouse, rat)

S6RP (pS235/236) (human, mouse, rat)  
 S6RP (pS240/244) (human, mouse, rat)  
 S6RP (Total) (human, mouse, rat)

## Multiplex Panels

Akt (pS473), Total Akt (human, mouse, rat)  
 GSK-3 $\beta$  (pS9), Total GSK-3 $\beta$  (human)  
 p70S6K (pT421/pS424), Total p70S6K (human, mouse, rat)

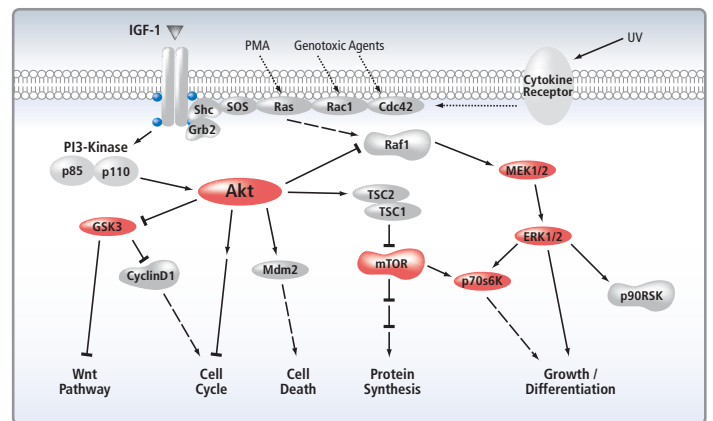
Akt Signaling Panel (Phospho Protein) (pAkt, pp70S6K, pGSK-3 $\beta$ ) (human, mouse, rat)  
 Akt Signaling Panel (Total Protein) (Akt, p70S6K, GSK-3 $\beta$ ) (human, mouse, rat)  
 Insulin Signaling Panel (Phospho Protein) (pIR, pIRS-1, pIGF-1R) (human)

Insulin Signaling Panel (Total Protein) (IR, IRS-1, IGF-1R) (human)

For a complete list of products please check our website at [www.mesoscale.com](http://www.mesoscale.com)

## Cell Signaling

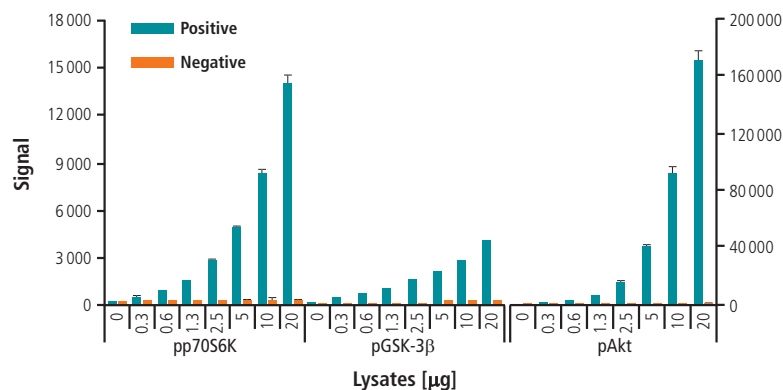
Insulin modulates blood glucose levels by driving glucose uptake and glycogen synthesis. Insulin signaling activates PI3 kinase which results in phosphorylation of Akt. This inhibits downstream substrates including GSK-3 $\beta$  thereby promoting protein synthesis. Akt also activates mTOR which stimulates protein synthesis through p70S6K. In addition, downstream targets of Akt coordinate glycogen synthesis, fatty acid oxidation and protein synthesis.



Available AKT/mTOR pathway phosphoprotein targets

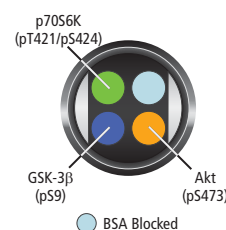
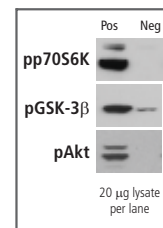
## Akt Signaling Panel

### Akt (pS473), GSK-3 $\beta$ (pS9) and p70S6K (pT421/pS424)



Logarithmically growing Jurkat cells were treated with PMA (200 nM, 15 min) (positive) or LY294002 (50  $\mu$ M) and staurosporine (1  $\mu$ M; 2.25 hr) (negative). Whole cell lysates were added to MSD MULTI-SPOT 96-Well 4 Spot plates coated with analyte-specific capture antibodies on spatially distinct electrodes in each well. Phosphorylated Akt, p70S6K, and GSK-3 $\beta$  were detected with multiplexed sandwich immunoassays using an anti-phospho-specific antibody as the capture or detection antibody in each sandwich.

## Traditional Western



## Catalog Numbers for Akt Signaling Panel

	Kit Size	Catalog Numbers
Human/	1 Plate	K15115D-1
Mouse/	5 Plate	K15115D-2
Rat	20 Plate	K15115D-3

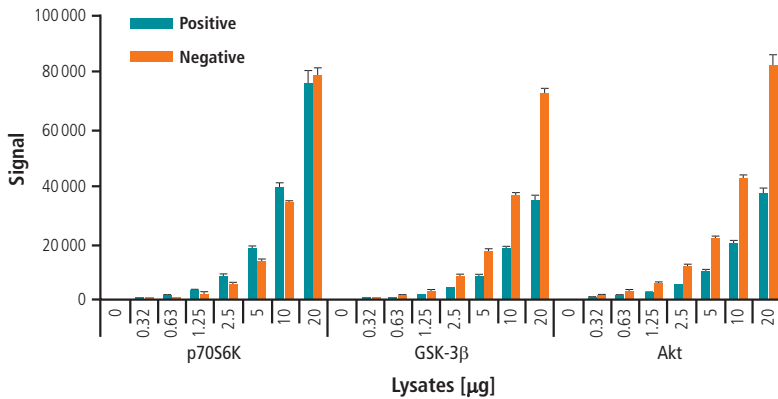
## Example Protocol

1. Add 150  $\mu$ L Blocking Solution, incubate 1 hour at RT.
2. Wash with TBS-T. Add 25  $\mu$ L lysates, incubate 1 hour at RT.
3. Wash with TBS-T. Add 25  $\mu$ L Detection Antibodies, incubate 1 hour at RT.
4. Wash with TBS-T. Add 150  $\mu$ L Read Buffer T, read.

# Intracellular Signaling

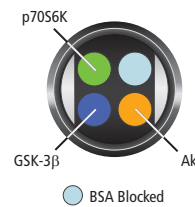
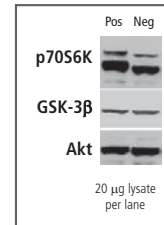
## Cell Signaling

### Akt Signaling (Total Protein) Panel Akt, GSK-3 $\beta$ and p70S6K



Logarithmically growing Jurkat cells were treated with PMA (200 nM, 15 min) (positive) or LY294002 (50  $\mu$ M) and staurosporine (1  $\mu$ M; 2.25 hr) (negative). Whole cell lysates were added to MSD MULTI-SPOT 96-Well 4 Spot plates coated with analyte-specific capture antibodies on spatially distinct electrodes in each well. Akt, p70S6K, and GSK-3 $\beta$  were detected in multiplex with anti-total-Akt, anti-total-p70S6K, and anti-total-GSK-3 $\beta$  antibodies. Detection antibodies for total proteins recognize epitopes distinct from the capture antibodies.

#### Traditional Western



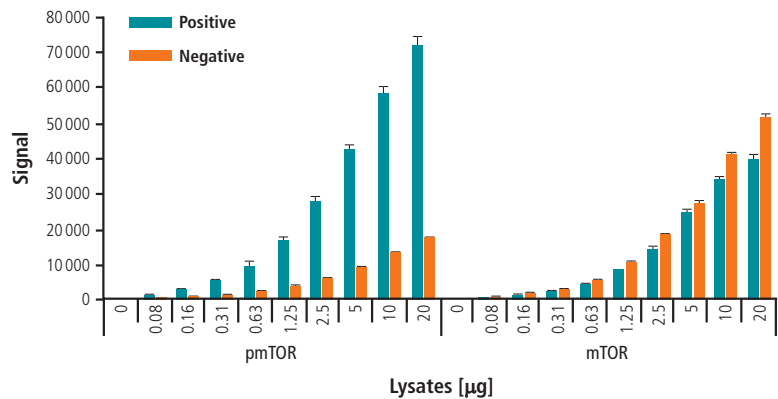
#### Catalog Numbers for Akt Signaling Panel (Total Protein)

	Kit Size	Catalog Numbers
Human/	1 Plate	K15133D-1
Mouse/	5 Plate	K15133D-2
Rat	20 Plate	K15133D-3

#### Example Protocol

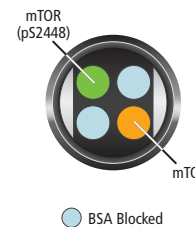
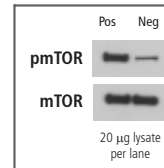
1. Add 150  $\mu$ L Blocking Solution, incubate 1 hour at RT.
2. Wash with TBS-T. Add 25  $\mu$ L lysates, incubate 1 hour at RT.
3. Wash with TBS-T. Add 25  $\mu$ L Detection Antibodies, incubate 1 hour at RT.
4. Wash with TBS-T. Add 150  $\mu$ L Read Buffer T, read.

### Phospho/Total mTOR



Growing HEK293 cells were treated with Wortmannin (100 nM, 3 h)(negative) or PMA (1  $\mu$ M, 30 min)(positive). Whole cell lysates were added to MSD MULTI-SPOT 4 Spot plates coated with anti-phospho-mTOR antibody and anti-total-mTOR antibody on two of the four spatially distinct electrodes per well. Phosphorylated and total mTOR were detected with anti-total-mTOR antibody labeled with MSD SULFO-TAG reagent.

#### Traditional Western



#### Catalog Numbers for Phospho/Total mTOR

	Kit Size	Catalog Numbers
Human/	1 Plate	K15170D-1
Mouse/	5 Plate	K15170D-2
Rat	20 Plate	K15170D-3

#### Example Protocol

1. Add 150  $\mu$ L Blocking Solution, incubate 1 hour at RT.
2. Wash with TBS-T. Add 25  $\mu$ L lysates, incubate 3 hours at RT.
3. Wash with TBS-T. Add 25  $\mu$ L Detection Antibody, incubate 1 hour at RT.
4. Wash with TBS-T. Add 150  $\mu$ L Read Buffer T, read.

# Metabolic Disease

## Representative Metabolic Markers

Adiponectin (human, mouse, rat)

Ghrelin (mouse/rat)

Active GLP-1 (7-36)amide (human, mouse/rat)

Active GLP-1 (human, mouse/rat)

GLP-1 (7-36)amide (human, mouse/rat)

Total GLP-1 (human, mouse/rat)

Glucagon (human, mouse/rat)

Insulin (human, mouse/rat)

Leptin (human, mouse, rat)

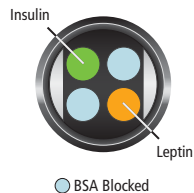
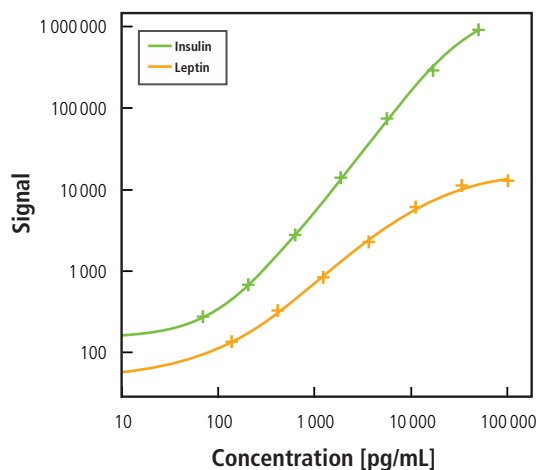
Resistin (human, mouse/rat)

For a complete list of products please check our website at [www.mesoscale.com](http://www.mesoscale.com)

Funicello, et al.<sup>14</sup> used the MSD multiplex Mouse Metabolic Panel to determine simultaneously plasma level of mouse insulin and leptin when studying the implications of proteases in adipogenesis.

## Mouse Metabolic Panel

### Mouse Insulin and Leptin



LLOD (pg/mL)	Analyte	
	Insulin	Leptin
	49	32

### Catalog Numbers for Insulin and Leptin

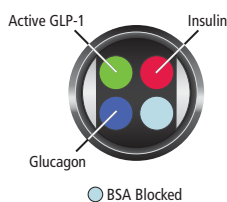
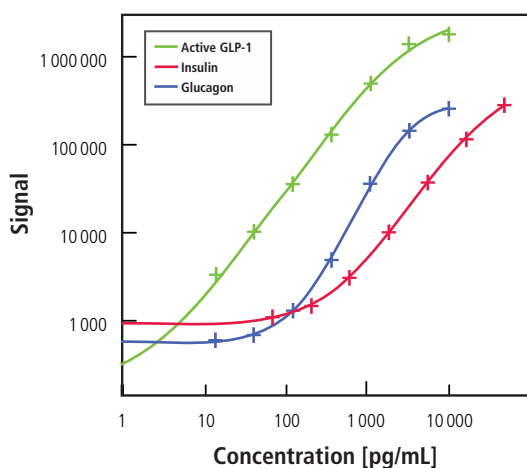
	Kit Size	Catalog Numbers
Human	1 Plate	K15164C-1
	5 Plate	K15164C-2
	20 Plate	K15164C-3
Mouse	1 Plate	K15124C-1
	5 Plate	K15124C-2
	20 Plate	K15124C-3
Rat	1 Plate	K15158C-1
	5 Plate	K15158C-2
	20 Plate	K15158C-3

### Example Protocol

1. Add 150  $\mu$ L Blocking Solution, incubate 1 hour at RT.
2. Wash with PBS-T. Add 40  $\mu$ L Detection Antibodies, add 10  $\mu$ L Calibrator/Sample, incubate 2 hours at RT.
3. Wash with PBS-T. Add 150  $\mu$ L Read Buffer T, read.

## Metabolic Multiplex

### Mouse/Rat Active GLP-1, Insulin, and Glucagon



LLOD (pg/mL)	Analyte		
	Active GLP-1	Insulin	Glucagon
	0.20	53	38

### Catalog Numbers for Active GLP-1, Insulin, and Glucagon

	Kit Size	Catalog Numbers
Mouse/Rat	1 Plate	K15184C-1
	5 Plate	K15184C-2
	20 Plate	K15184C-3






### Example Protocol

1. Add 150  $\mu$ L Blocking Solution, incubate 1 hour at RT.
2. Wash with PBS-T. Add 20  $\mu$ L Diluent, add 40  $\mu$ L Calibrator/Sample, incubate 2 hours at RT.
3. Wash with PBS-T. Add 25  $\mu$ L Detection Antibodies.
4. Wash with PBS-T. Add 150  $\mu$ L Read Buffer T, read.

# References

- <sup>1</sup> Rozell, B., et al. **Glutathione transferases of classes alpha, mu and pi show selective expression in different regions of rat kidney.** Xenobiotica. 1993 23(8) 835-49.
- <sup>2</sup> Falkenberg, F.W., et al. **Urinary antigens as markers of papillary toxicity. I. Identification and characterization of rat kidney papillary antigens with monoclonal antibodies.** Arch. Toxicol. 1996 71(1-2) 80-92.
- <sup>3</sup> Lee, J.W., et al. **Fit-for-purpose method development and validation for successful biomarker measurement.** Pharm Res. 2006 Feb; 23(2):312-28.
- <sup>4</sup> <http://www.fda.gov/NewsEvents/Newsroom/PressAnnouncements/2008/ucm116911.htm>
- <sup>5</sup> Sieber, M., et al. **Comparative analysis of novel noninvasive renal biomarkers and metabonomic changes in a rat model of gentamicin nephrotoxicity.** Toxicol. Sci. 2009 109(2):336-49.
- <sup>6</sup> Waikar, S.S., and Bonventre, J.V. **Biomarkers for the diagnosis of acute kidney injury.** Nephron Clin. Pract. 2008 109(4):c192-7.
- <sup>7</sup> Denhardt, D.T., et al. **Role of osteopontin in cellular signaling and toxicant injury.** Annu. Rev. Pharmacol. Toxicol. 2001 41:723-49.
- <sup>8</sup> Lim, B.J., et al. **Osteopontin expression and microvascular injury in cyclosporine nephrotoxicity.** Pediatr. Nephrol. 2004 19(3):288-94.
- <sup>9</sup> Braunwald, E. **Biomarkers in Heart Failure.** N. Engl. J. Med. 2008 358:2148-2159.
- <sup>10</sup> O'Brien, P.J. **Cardiac troponin is the most effective translational safety biomarker for myocardial injury in cardiotoxicity.** Toxicology. 2008 245(3):206-18. 
- <sup>11</sup> Pecoits-Filho, R., et al. **Associations between circulating inflammatory markers and residual renal function in CRF patients.** Am. J. Kidney Dis. 2003 Jun;41(6):1212-8.
- <sup>12</sup> Carrero, J.J., and Stenvinkel, P. **Persistent inflammation as a catalyst for other risk factors in chronic kidney disease: a hypothesis proposal.** Clin. J. Am. Soc. Nephrol. 2009 Dec;4 Suppl 1:S49-55.
- <sup>13</sup> Lee, D.S., and Vasan, R.S. **Novel markers for heart failure diagnosis and prognosis.** Curr. Opin. Cardiol. 2005 May;20(3):201-10.
- <sup>14</sup> Funicello, M., et al. **Cathepsin K null mice show reduced adiposity during the rapid accumulation of fat stores.** PLoS ONE. 2007 Aug 1;2(1):e683. 

## Other references using MSD MULTI-ARRAY technology

- Sun, D., et al. **Electrochemiluminescent immunoassay for rat skeletal troponin I (Tnni2) in serum.** J. Pharmacol. Toxicol. Methods. 2010 61(1): 52-58. 
- Collings, F.B., and Vaidya, V.S. **Novel technologies for the discovery and quantitation of biomarkers of toxicity.** Toxicology. 2008 20;245(3):167-74. 
- Dieterle, F., and Marrer, E. **New technologies around biomarkers and their interplay with drug development.** Anal. Bioanal. Chem. 2008 390(1):141-54. 
- Kattla, J.J., et al. **Protein kinase B/Akt activity is involved in renal TGF- $\beta$ 1-driven epithelial-mesenchymal transition in vitro and in vivo.** Am. J. Physiol. Renal Physiol. 2008 295(1): F215-25. 
- O'Brien, P.J. **Cardiac troponin is the most effective translational safety biomarker for myocardial injury in cardiotoxicity.** Toxicology. 2008 245(3):206-18. 

# Select Catalog Numbers

## Toxicology Assays

For a complete listing of our products,  
visit our online catalog at  
[www.mesoscale.com](http://www.mesoscale.com)<sup>®</sup>

Assays	Description	Catalog Numbers
*TIM-1/KIM-1/HAVCR	Rat TIM-1/KIM-1/HAVCR (1 Plate) Rat TIM-1/KIM-1/HAVCR (5 Plates)	K153JHC-1 K153JHC-2
*Clusterin	Rat Clusterin Kit (1 Plate) Rat Clusterin Kit (5 Plates)	K153HXC-1 K153HXC-2
*Skeletal Troponin I (sTnI)	Rat Skeletal Troponin I Kit (1 Plate) Rat Skeletal Troponin I Kit (5 Plates)	K153IMC-1 K153IMC-2
*TIMP-1	Human TIMP-1 Kit (1 Plate) Human TIMP-1 Kit (5 Plates)	K151JFC-1 K151JFC-2
*Luteinizing Hormone (LH)	Human Luteinizing Hormone (LH) Kit (1 Plate) Human Luteinizing Hormone (LH) Kit (5 Plates)	K151ETC-1 K151ETC-2
*Argutus AKI Test <sup>®</sup> αGST, GSTYb1, RPA-1	Argutus AKI Test (rat) Kit (1 Plate) Argutus AKI Test (rat) Kit (5 Plates)	K15156C-1 K15156C-2
*Kidney Injury Panel 1 Albumin, TIM-1, Lipocalin-2, Osteopontin	Kidney Injury Panel 1 (rat) Kit (1 Plate) Kidney Injury Panel 1 (rat) Kit (5 Plates)	K15162C-1 K15162C-2
*Cardiac Injury Panel 2 cTnI, cTnT, FABP3	Cardiac Injury Panel 2 (rat) Kit (1 Plate) Cardiac Injury Panel 2 (rat) Kit (5 Plates)	K15155C-1 K15155C-2
*Cardiac Injury Panel 3 cTnI, cTnT, FABP3, Myl3	Cardiac Injury Panel 3 (rat) Kit (1 Plate) Cardiac Injury Panel 3 (rat) Kit (5 Plates)	K15161C-1 K15161C-2
*Muscle Injury Panel 1 cTnI, cTnT, FABP3, Myl3, sTnI	Muscle Injury Panel 1 (rat) Kit (1 Plate) Muscle Injury Panel 1 (rat) Kit (5 Plates)	K15181C-1 K15181C-2
*Acute Phase Protein Panel 1 A2M, AGP	Acute Phase Protein Panel 1 (rat) Kit (1 Plate) Acute Phase Protein Panel 1 (rat) Kit (5 Plates)	K15175C-1 K15175C-2
Steroid Hormone Panel 1 Estradiol, Progesterone, Testosterone, DHEA	Steroid Hormone Panel 1 (human, mouse, rat) Kit (1 Plate) Steroid Hormone Panel 1 (human, mouse, rat) Kit (5 Plates)	K15178C-1 K15178C-2

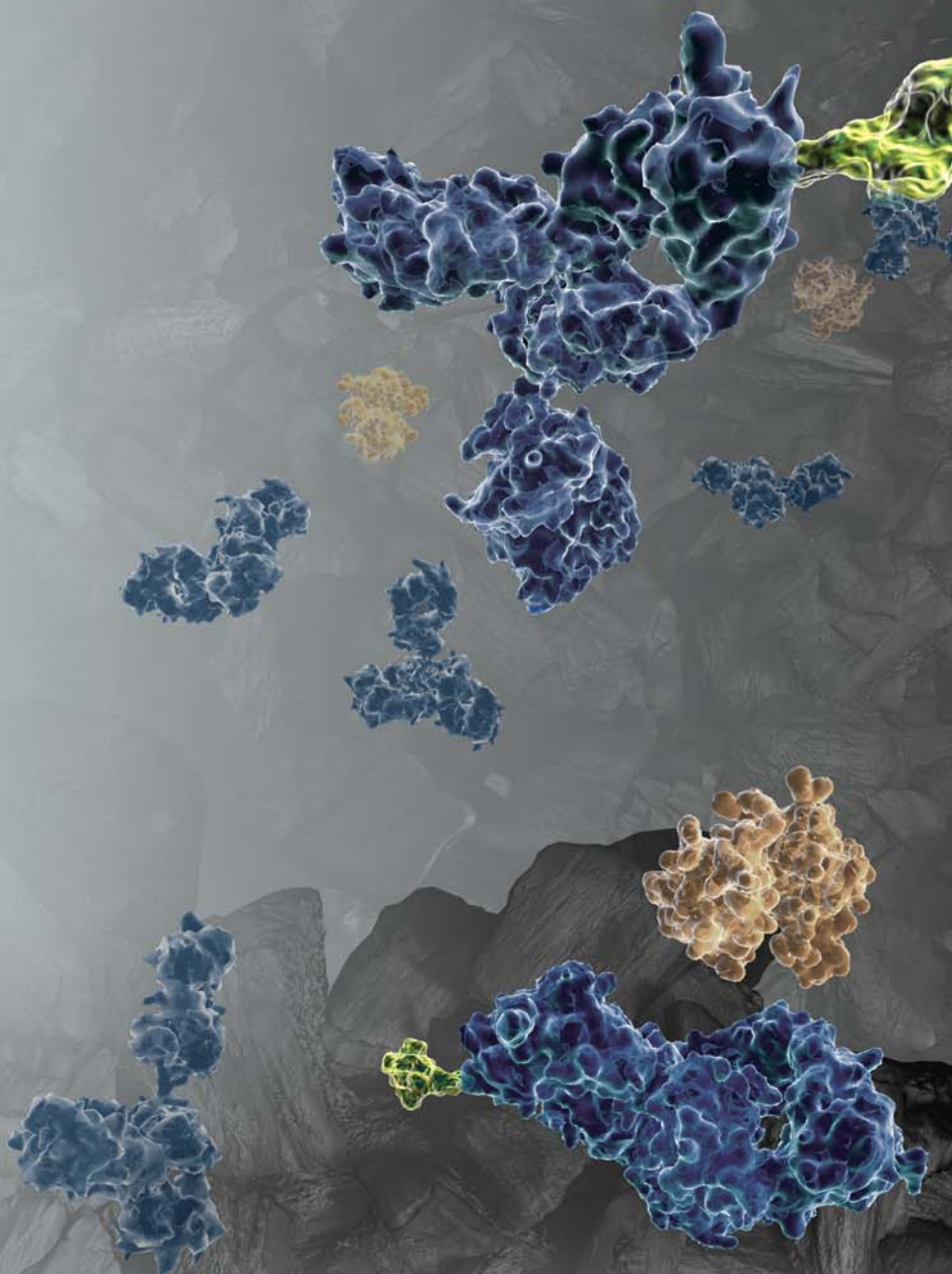
\* *Qualified Kit*

### Ordering Information

Purchase orders should be faxed or emailed to the following queries:

TEL: 1.301.947.2085  
FAX: 1.301.990.2776  
EMAIL: [customerservice@mesoscale.com](mailto:customerservice@mesoscale.com)

- Alzheimer's Disease
- BioProcess
- Cardiac
- Cell Signaling
- Clinical Immunology
- Cytokines
- Hypoxia
- Immunogenicity
- Inflammation
- Metabolic
- Oncology
- Toxicology
- Vascular



9238 Gaither Road Gaithersburg, MD 20877 USA Phone: 1.301.947.2085 Fax: 1.301.990.2776 Website: [www.mesoscale.com](http://www.mesoscale.com)

Meso Scale Discovery, Meso Scale Diagnostics, [www.mesoscale.com](http://www.mesoscale.com), MSD, MSD (design), MULTI-ARRAY, MULTI-SPOT and SULFO-TAG are trademarks and/or service marks of Meso Scale Diagnostics, LLC. AKI Test is a trademark of Argutus Medical Limited.

© 2011 Meso Scale Discovery, a division of Meso Scale Diagnostics, LLC. All rights reserved.  
19512-v3-2011Oct