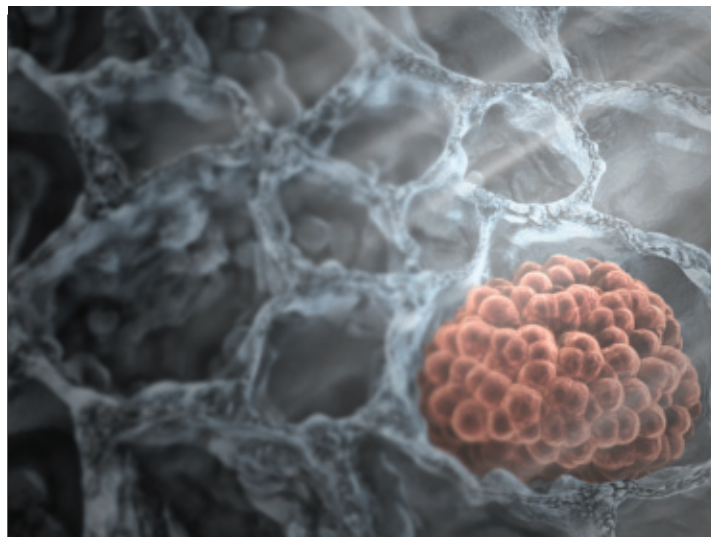


ONCOLOGY APPLICATIONS



MESO SCALE DISCOVERY® Oncology Applications

MESO SCALE DISCOVERY (MSD®) offers a diverse product line of biomarker assays including many oncology-focused kits. We feature phospho-protein and intracellular signaling assay kits designed for use with whole cell lysates and other complex matrices. Clinical biomarker assays for analysis of serum, plasma, and tissue culture samples are also available from MSD. Whether your focus of study is angiogenesis, apoptosis, or another aspect of oncology research, MSD can enable you with the tools and technology to enhance all stages of cancer therapeutics development.



Cancerous tumor in the lung



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
www.mesoscale.com



■ MSD Technology	1
■ MSD Multiplex Phosphoprotein Assays	2
■ The Challenge	3
■ The Transition from Discovery	4
■ Monitoring the Periphery	5
■ Recombinant Phosphoprotein Standards	6
■ Cell Signaling Markers	7
■ Inflammation and Cancer	8
■ Biomarker Profiling	9
■ Cytokines and Chemokines Markers	10
■ Prototype Printing Services	11
■ References	12
■ The MSD Advantage	13

MSD Technology

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

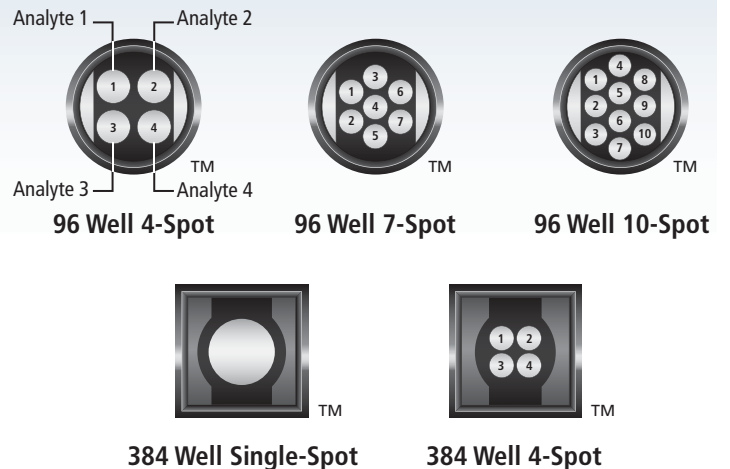
Electrochemiluminescence Features:

- 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 conveniently conjugated to biological molecules
- Emission at ~620 nm - eliminates 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:

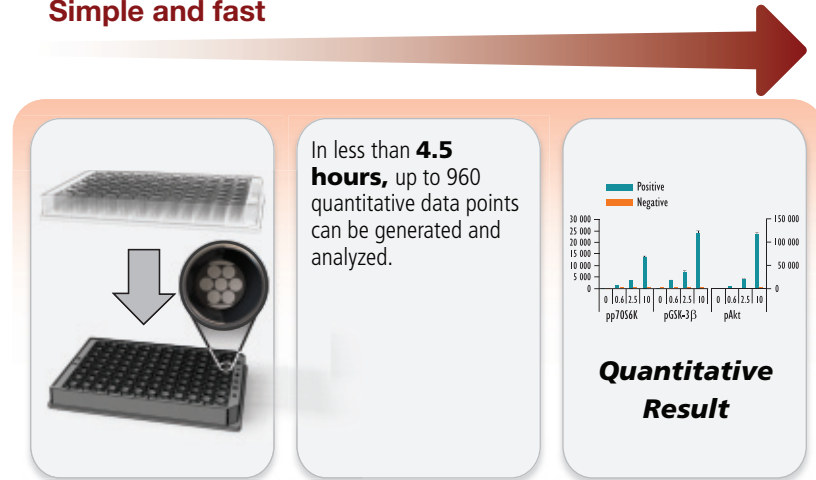
- Capability to simultaneously measure multiple analytes in the same well
- High density arrays for high throughput multiplexing of biomarkers
- The unique bar code label on each plate enables complete traceability back to MSD manufacturing records
- The MSD DISCOVERY WORKBENCH® software provides customers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



MSD Multiplex Phosphoprotein Assays: Speeding Discovery, Advancing Oncology

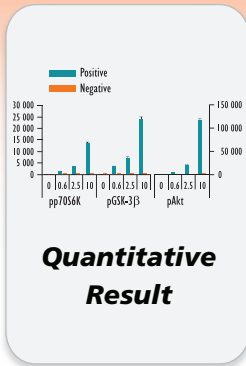
MSD

Simple and fast



Sample Preparation
Lysate preparation in 96-well plate and direct transfer to MSD plate

In less than **4.5 hours**, up to 960 quantitative data points can be generated and analyzed.

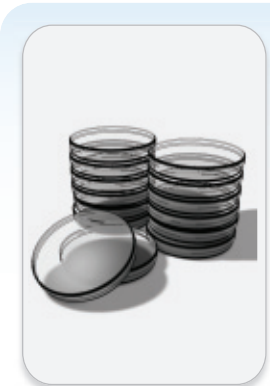


- reduce cost and time
- gain sensitivity
- reduce cell culture volume
- analyze wide range of matrices
- conserve precious samples
- gain with MSD multiplexing
- gain in throughput

- Multiple phosphoproteins as well as total pools can be accurately quantified in a single well from as few as 1000 cells
- % Phosphorylation of protein can also be calculated

Western Blot

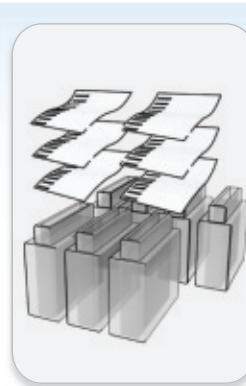
Laborious and time consuming



Sample Preparation

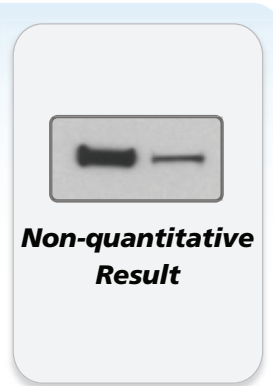


Gel Electrophoresis
6 gels ~ equivalent to one 96-well plate



Blotting and transfer to membrane

With Western blot, more than **10 hours**, and **up to 2 days** may be required to analyze 10-fold less data.



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

The Challenge:

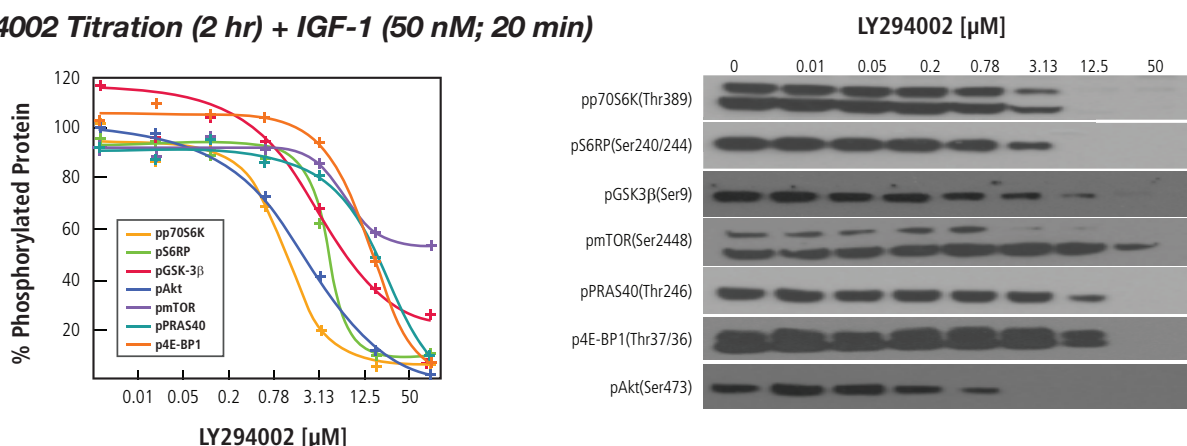
Rapid Discovery to Advance Promising Compounds in Oncology

Recent advances in understanding the molecular pathogenesis of cancer present challenges and opportunities in oncology research. For example, somatic mutations that result in up-regulation of the PI3 kinase (PI3K) signaling pathway are common in cancers and numerous discovery programs seek to identify specific inhibitors. Discovery teams demand a rapid screening approach to accurately assign IC_{50} and correctly rank order compounds with respect to their potency and in vivo efficacy. One research group replaced 6 weeks of western blot screening of 150 compounds with less than one week of testing using MSD multiplex assays. In just a few days, eight-point titrations of 150 compounds, in triplicate, against three phosphoproteins (10 800 data points) were completed.

“Compared to MSD assays, western blotting is time-consuming, with most antibodies requiring overnight incubations. Western blotting assays using two antibodies require 2 days, whereas the MSD assay can be completed and analyzed within 1 day.” (Gowan et al.)¹

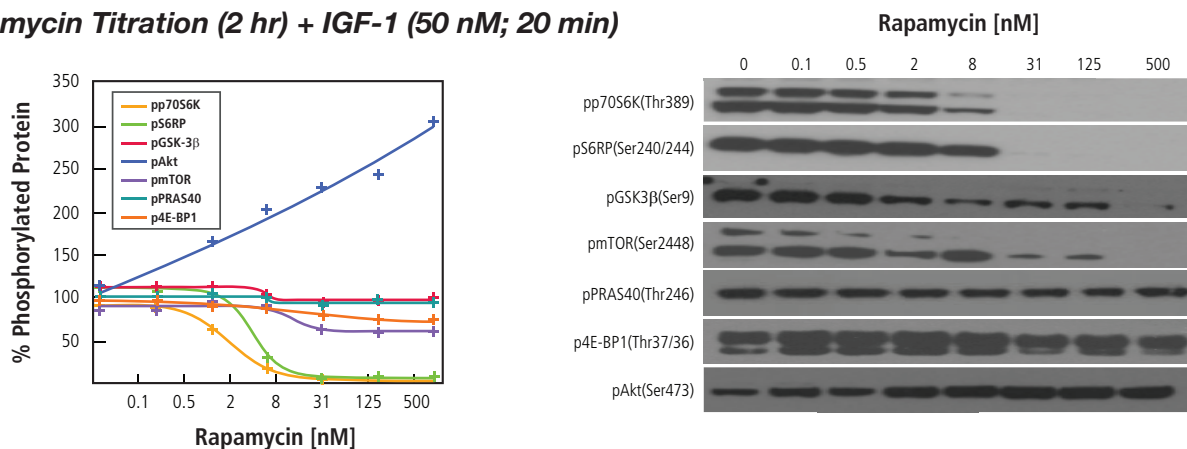
The figures below illustrate the dose-dependent effects of inhibitors of Akt signaling pathway using human MCF7 tumor cell lines stimulated with IGF-1. MCF7 cells were treated with the indicated amount of inhibitor followed by a 20 minute stimulation with IGF-1 (50 nM).

LY294002 Titration (2 hr) + IGF-1 (50 nM; 20 min)



- $\frac{1}{4}$ of whole cell lysate preparation from a well of a 96-well plate is used for each MSD data point. A 2-fold larger amount of lysate was used to complete traditional western blot analysis.

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



By using MSD's multiplexed assays containing targets of the PI3K pathway, Raynaud et al.² presented supportive data for the development of GDC-0941, a potent inhibitor of PI3K in glioblastoma cell line. GDC-0941 is now in clinical trials for solid breast tumor malignancies and non-Hodgkins lymphoma.

The Transition from Discovery: Multiplex Assays for Pharmacodynamics in Human Xenografts

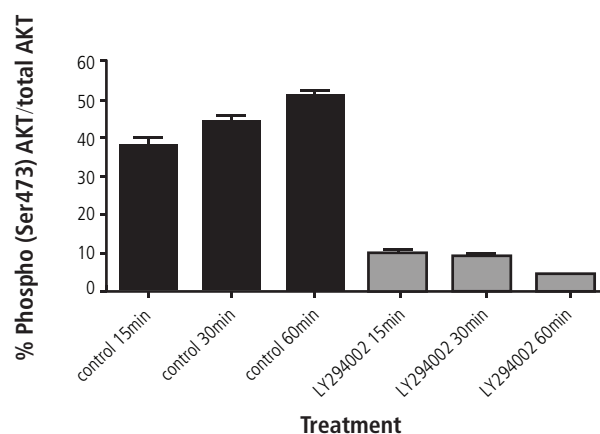
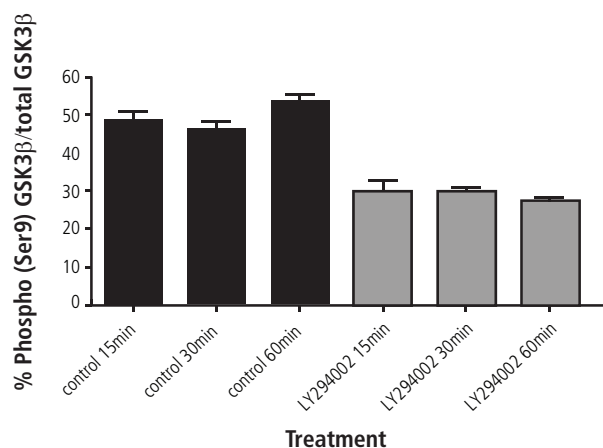
The same MSD assays that power compound characterization in tumor cell lines have proven sensitive and robust enough for the study of compound activity in xenografts. Gowan et al¹ selected MSD over all other available methods such as enzyme-linked immunosorbent assay (ELISA), western blotting, immunohistochemistry, reverse transcription-polymerase chain reaction, flow cytometry, magnetic resonance imaging/spectroscopy, and gene expression profiling. The MSD Phospho/Total Akt and Phospho/Total GSK-3 β multiplex assays were used to measure both in vitro and in vivo changes in response to LY294002 within two different tumor types.

Major Obstacles with Conventional Methods

- Large sample volume
- High matrix interference
- Several days of procedure and analysis
- Qualitative versus quantitative

MSD Solutions and Customer's Comments

- **Reduced sample volume**
"...possibility of detailed and accurate analyses of valuable patient material as several proteins can be screened from a small sample within 1 day."¹
- **Rapid and robust**
"The assays are robust (coefficient of variation for phospho-Akt 13.4%) and offer significant advantages (in terms of speed and quantitation) over western blots."¹
- **Low background**
"The assay background signal...detected with MSD technology was very low...and the peak signal was very high... resulting in a superior signal-to-noise ratio..."³
- **Matrix compatibility**
"MSD protein profiling system offers a sensitive, adaptable platform for sandwich immunoassays in microplates with one to 10 carbon electrode arrays per well."¹
- **Absolute quantification**
"...novel sandwich assay (ECL) as it provides sensitivity...accuracy, reproducibility and absolute quantification..."⁴
- **One technology and one platform for different stages of drug discovery and development**
"This optimized procedure can be used for both in vitro and in vivo analysis, unlike an established fixed-cell ELISA with a time resolved fluorescent end point."¹



Inhibition of phosphorylation on Akt and GSK3 β phosphoprotein in human tumor xenograft treated with LY294002. Approximately 3×10^6 cells were injected bilaterally subcutaneously in the flanks of female NCr athymic mice. The effect of LY294002 at 100 mg/kg (two doses, 30 min apart) was compared to solvent controls at 15 min, 30 min, and 1 h after administration. Tumors were excised and immediately snapfrozen in liquid nitrogen. Samples were stored at -80°C until processed. Tumors were ground and lysed in the optimal lysis buffer (buffer B). Protein concentrations were measured, and samples (10 μg) were analyzed using MSD phosphoprotein assays.¹

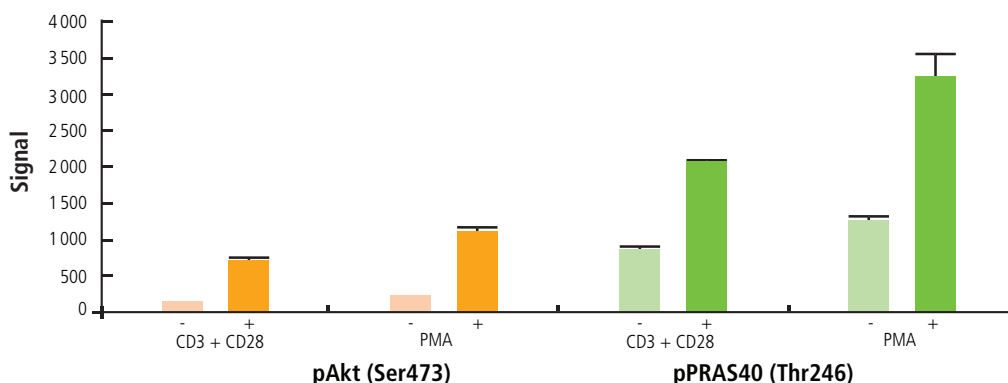
Monitoring the Periphery: Peripheral Blood Mononuclear Cells (PBMC) as Readout for Drug Efficacy

Chemical entities that inhibit kinases have been the subject of intense drug discovery and development owing to the documented role of these proteins in tumorigenesis. Pharmacodynamic studies ideally use the substrates of these kinases, phosphoproteins, to report drug activity. In many cases, the tumors that are the target of these therapies are unavailable or inaccessible. However, a number of the pathways that are the objective of therapeutic intervention are constitutively active in PBMCs. Tan et al.⁵ have reviewed the concept of using peripheral cells to monitor drug activity.

The data below illustrate the utility of MSD phosphoprotein assays for monitoring drug activity against the PI3 kinase pathway using PBMCs.

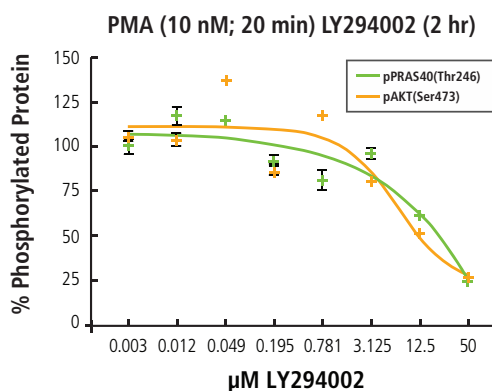
Dose-dependent inhibition of Akt and PRAS40 in PBMCs with a potent PI3 Kinase inhibitor

Human PBMCs were purified from whole blood. Whole cell lysates were prepared and examined using the MSD MULTI-ARRAY Phospho-Akt (Ser473) and Phospho-PRAS40 (Thr246) Whole Cell Lysate Kits.

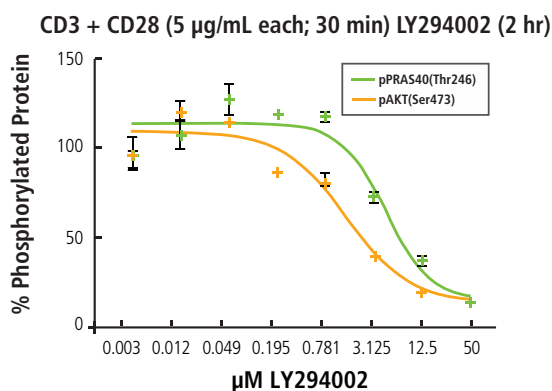


Human PBMCs were stimulated with either PMA (10 nM; 20 min) or CD3 + CD28 antibodies (5 µg/mL each; 30 min).

Human PBMCs were treated with serial dilutions of the PI3 Kinase inhibitor LY294002. The cells were then stimulated with either PMA (10 nM; 20 min) or CD3 + CD28 antibodies (5 µg/mL each; 30 min).



IC₅₀ curves for LY294002 inhibition of activated Akt and PRAS40 proteins after stimulation with PMA.



IC₅₀ curves for LY294002 inhibition of activated Akt and PRAS40 proteins after stimulation with a mixture of CD3 and CD28 antibodies.

"This technology...is cost-effective compared with western blots." (Gowan et al.)¹

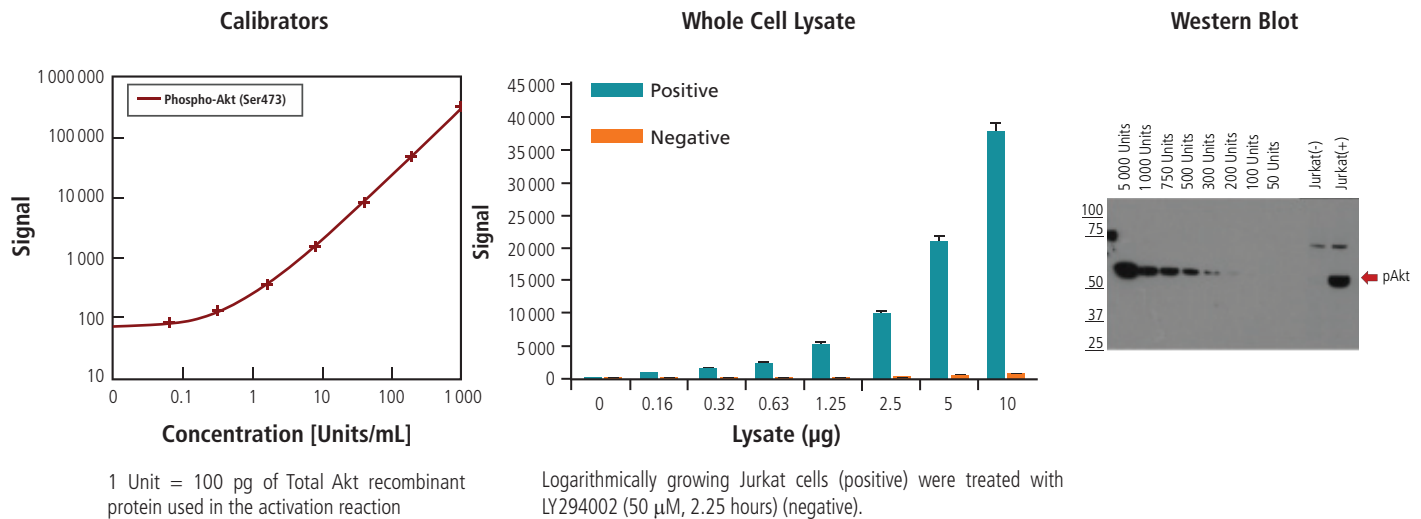
MSD multiplex panels have also been used by Ghosh, et al.⁶ to profile cytokine response in PBMCs treated with toll-like receptor agonists individually or in combination. The cell-free supernatants were collected and tested for IFN- α , IFN- β , IFN- γ , IL-12p70, and TNF- α using MULTI-SPOT plates on the MESO SCALE DISCOVERY platform.

Recombinant Phosphoprotein Standards: Quantitative Assignment of Abundance

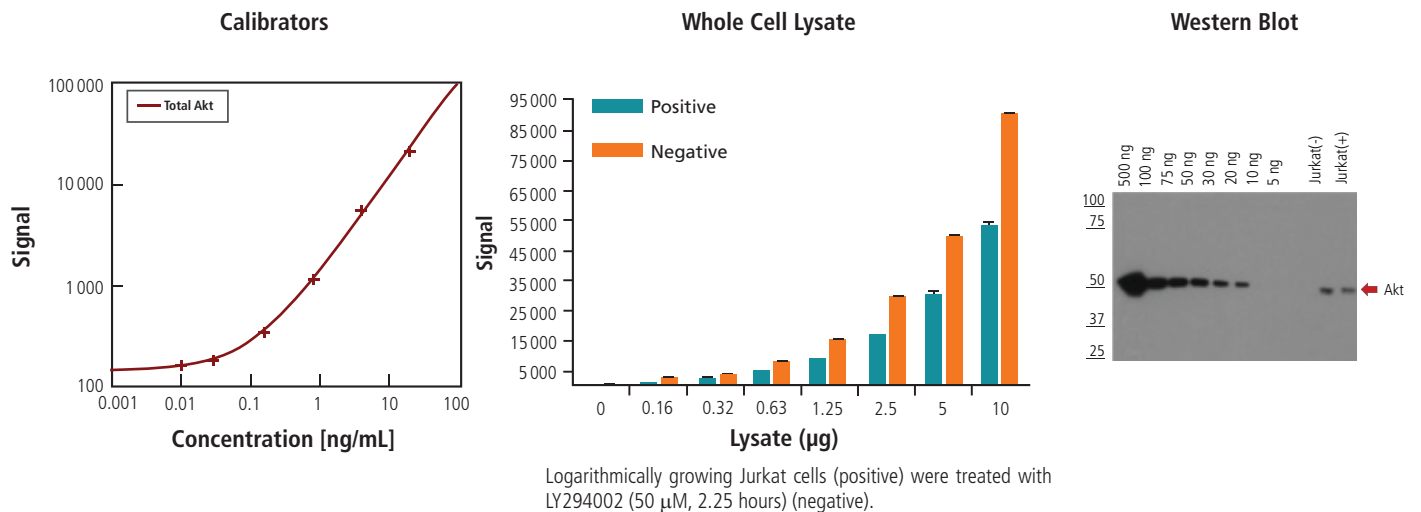
The absolute quantification of phosphoproteins in clinical samples will greatly add to their utility as prognostic, diagnostic, and pharmacodynamic markers. Such precision will require the use of purified, recombinant protein standards as calibrators, as is routinely practiced in concentration assignments for soluble proteins in bodily fluids. The production and handling of recombinant protein calibrators must assure that the epitopes of the capture and detection antibodies are readily recognized and that the post-translational phosphorylation is not lost. In the example below, MSD site-specific phosphorylated and total Akt recombinant standards are compared to MSD whole cell lysates and non-quantitative western blot analysis.

"...novel sandwich assay (ECL) as it provides sensitivity... accuracy, reproducibility and absolute quantification..." (Cao et al.)⁴

Phospho-Akt (Ser473) Assay



Total Akt Assay



Representative Cell Signaling Markers

Singleplex Kits

4E-BP1	pT37/46 [§] Total [‡]	GAPDH [‡]		Met	pY1349 Total	Smad1	pS463/465 [‡]
Akt	pS473 pT308 Total	GSK-3α	pS21 [‡] Total [‡]	mTOR	pS2448	SRC3	Total [‡]
Aurora A	pT288	GSK-3β	pS9 Total	NFκB	pS536 pS468 Total [‡]	STAT3	pY705
BAD	pS112 Total	HIF-1α	Total	p21	Total [‡]	STAT4	pY693 Total
c-Jun	pS63	Histone H3	pS10	p38	pT180/pY182 Total	STAT5a,b	pY694 Total
c-Kit	pY721 Total	HSP27	pS15 pS78 pS82 Total	p53	pS15 Total	Tau	pT231 pS202 pS262 [‡] pS396 [‡]
Caspase-3	p20/p17 (cleaved) Total	HSP70	Total	p70S6K	pT421/pS424 pT389 Total		Total pT181
EGFR	pY1173 pY1068 Total	IGF-1R	pY Total	PARP	D214 (cleaved)	Tuberin	pT1462 [‡]
eIF2α	pS51 [‡] Total [‡]	IR	pY Total	PDGFR-β	pY751	VASP	pS157 pS239 Total
EphB4	Total [‡]	IRS-1	pS312 Total	PDH E1α	pS293 [‡]	VEGFR2	pY1054 pY1175 Total
ErbB2	pY1248 Total	JAK2	pY1007/1008 [‡] Total [‡]	Pleckstrin	pS [‡]		
ErbB3 (HER3)	pY [‡]	JNK	pT183/pY185 Total	PRAS40	pT246 Total [§]		
ERK1/2	(pT202/pY204), (pT185/pY187) Total	MAPKAPK-2	pT334	Rb	pS608 pS780 Total		
ERK5	pT218/pY220	MEK1/2	pS217/221 Total	RON	pY1238/1239 [‡]		
FOXO3a	pT32 [§]	MEK2	Total	S6RP	pS235/236 pS240/244 Total		
FRS2	pY196 pY436						

[‡] Available through our Prototype Printing Services [§] Coming Soon

Multiplex Kits

Activated/Total Panels

Akt (pS473), Total Akt	NFκB (pS468), Total NFκB [‡]
BAD (pS112), Total BAD	NFκB (pS536), Total NFκB [‡]
Caspase-3 (Cl. p20/p17), Total Caspase-3	p38 (pT180/pY182), Total p38
EGFR (pY1173), Total EGFR	p53 (pS15), Total p53
eIF4E (pS209), Total [‡]	p70S6K (pT421/pS424), Total p70S6K
ErbB2 (pY1248), Total ErbB2	Rb (pS608), Total Rb
ERK1/2 (pT202/pY204)/(pT185/pY187), Total ERK1/2	Rb (pS780), Total Rb
GSK-3α (pS21), Total GSK-3α [‡]	S6RP (pS240/244), Total S6RP
GSK-3β (pS9), Total GSK-3β	STAT5a,b (pY694), Total STAT5a,b
HSP27 (pS15), Total HSP27	Tau (pT231), Total Tau
HSP27 (pS78), Total HSP27	VASP (pS157), VASP (pS239), Total VASP
HSP27 (pS82), Total HSP27	
JNK (pT183/pY185), Total JNK	
c-Kit (pY721), Total c-Kit	
MEK1/2 (pS217/221), Total MEK1/2	
Met (pY1349), Total Met	
mTOR (pS2448), Total mTOR	

Multiplex Kits

MAP Kinase Panel (Phosphoprotein) (pERK1/2, pJNK, pp38)
MAP Kinase Panel (Total Protein) (ERK1/2, JNK, p38)
EGFR Family Panel (pEGFR, pErbB2, pIGF-1R)
EGFR, ErbB2, ErbB3 (pY detection) [‡]
Akt Signaling Panel (Phosphoprotein) (pAkt, pp70S6K, pGSK-3β)
Akt Signaling Panel (Total Protein) (Akt, p70S6K, GSK-3β)
Akt Signaling Panel II (Phosphoprotein) (pAkt, pp70S6K, pGSK-3β, pS6RP)
Akt Signaling Panel II (Total Protein) (Akt, p70S6K, GSK-3β, S6RP) [‡]
Apoptosis Panel (pp53, p53, Cl. Caspase-3, Cl. PARP)
ERK-STAT3 Cascade Panel (pERK1/2, pMEK1/2, pSTAT3)
Insulin Signaling Panel (Phosphoprotein) (pIR, pIGF-1R, pIRS-1)
Insulin Signaling Panel (Total Protein) (IR, IGF-1R, IRS-1)
STAT Panel (pSTAT3, pSTAT4, pSTAT5a,b)

Ubiquitin Pathway Kits

MDM2 - p53 complex
Total MDM2
Ubiquitinated MDM2
Total p53
Ubiquitinated p53
Self-Ubiquitinated MDM2 (Activity Assay)
Multiplex Panels
MDM2 (Ub), Total MDM2
p53 (Ub), Total p53

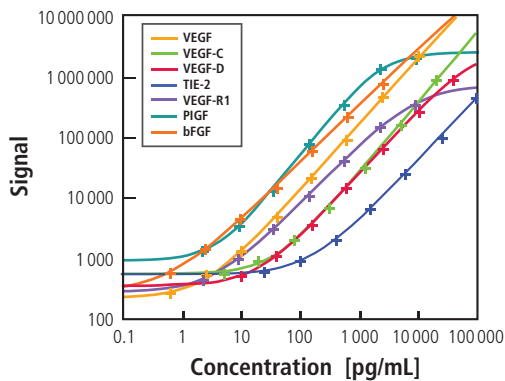
Inflammation and Cancer:

Evaluation of Angiogenic Biomarkers

Angiogenesis, the growth of new blood vessels, has emerged as a fundamental step in tumor growth, metastatic progression, and results in poor patient outcomes in several disease states including cancer. Understanding the critical regulators and pathways involved in the formation of new blood vessels is a complex process that involves not only neoplastic cells, but also endothelial cells, the basal membranes of neighboring capillaries and the stroma of the growing tumor mass. Key angiogenic factors, most notably VEGF, may serve as potential prognostic markers for disease activity or tumor progression and are often the target of inhibitors, receptor blockades or other chemotherapeutics.

MSD offers a variety of assays for pro-angiogenic and related proteins that provide a comprehensive assessment and understanding of the endothelial, vascular, and tumor cell microenvironment.

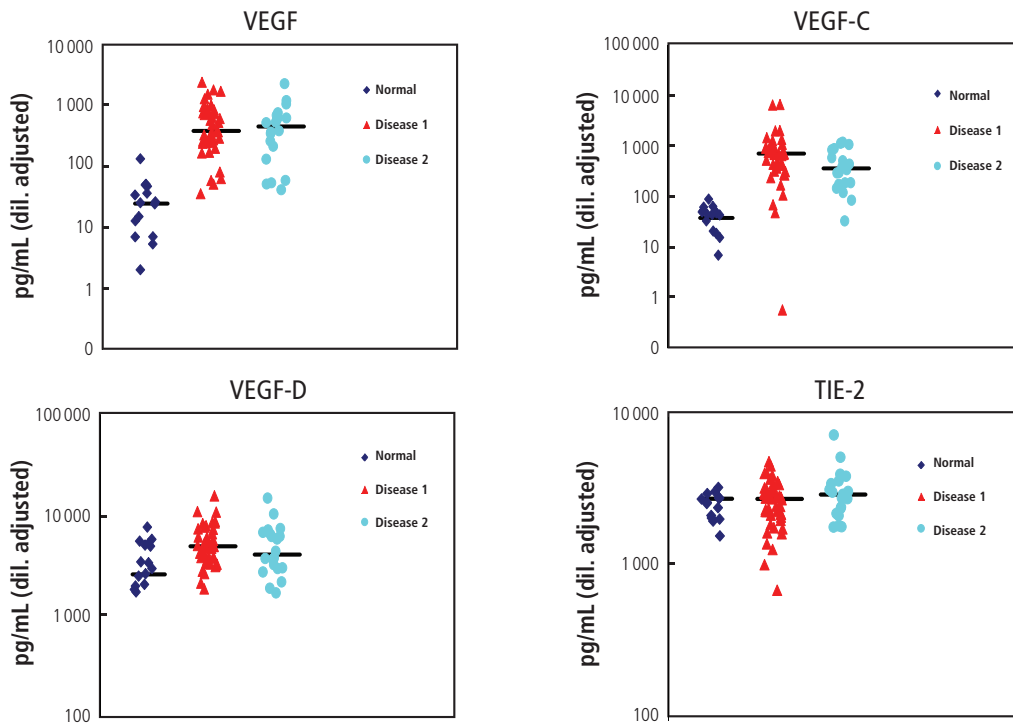
Human Angiogenesis Panel 1



LLOD (pg/mL)	Analyte						
	VEGF	VEGF-C	VEGF-D	TIE-2	VEGF-R1	PlGF	bFGF
	0.64	5.4	2.4	26	0.42	0.35	0.10

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

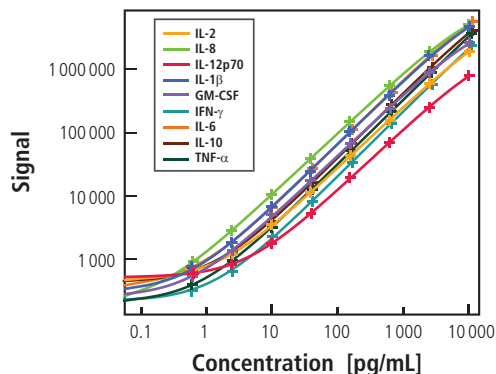
Illustrated below are biomarker levels in sera from normal and diseased populations tested at 2-fold dilution on the MSD Angiogenesis Panel 1. Some biomarkers may be elevated or suppressed in a given disease state whereas other biomarkers are not effected.



Biomarker Profiling: Analysis of Clinical Samples

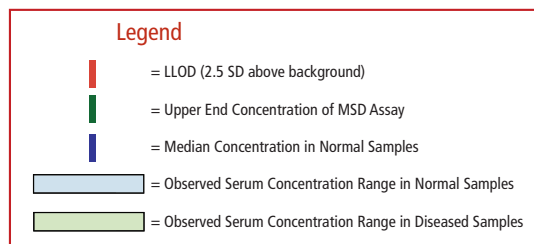
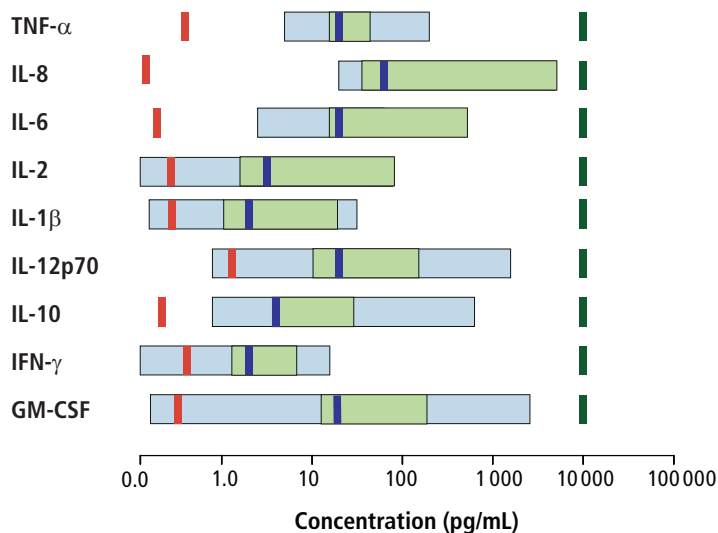
Clinical samples push the limits of traditional methods for immunoassays (e.g. ELISA, RIA). Complex matrices (e.g. sputum, vaginal fluids, etc.), widely ranging concentrations of analytes, and limited sample volume can make assays intractable. MSD's assays improve sensitivity, expand the dynamic range,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

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



LLOD (pg/mL)	Analyte								
	IL-2	IL-8	IL-12p70	IL-1β	GM-CSF	IFN-γ	IL-6	IL-10	TNF-α
	0.35	0.09	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*, a total of 127 human serum samples were tested, which included diseased pools and controls. The upper end of the calibrator curve for this panel is 10 000 pg/mL for all cytokines and the lower limit of detection (LLOD) was determined as 2.5 standard deviations above the background.

The MSD MULTI-ARRAY platform has been used by Bafadhel, et al.⁸ to study a panel of 11 cytokines and chemokines in a complex matrix such as sputum samples from subjects with chronic obstructive pulmonary disorder (COPD). The rapid assays enabled the measurement of multiple cytokines from several samples with minimal dilution and indicated that IL-5 in sputum is a repeatable and reproducible measurement of COPD.

"Reproducible interlaboratory measurement of all cytokines was found only for the MSD assays." (Fichorova et al.)⁹

* 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).

Representative Cytokines and Other Oncology Markers

Human				Mouse		Rat		Canine
BDNF	IL-6	MCP-1	RANTES	GM-CSF	MIP-1 α [‡]	CINC-2 [‡]	TIMP-1 [‡]	GM-CSF [‡]
CTACK [‡]	IL-6R	MCP-2 [‡]	SDF-1 α [‡]	IFN- γ	MIP-1 β [‡]	CINC-3 [‡]	TNF- α	IFN- γ [‡]
Eotaxin	IL-7	MCP-3 [‡]	TARC	IL-1 β	MIP-2 [‡]	CNTF-1 [‡]		IL-2
Eotaxin-3	IL-8	MCP-4	TGF- β 1	IL-2	MIP-3 α [‡]	GM-CSF		IL-6
G-CSF	IL-10	MDC	TGF- β 2 [‡]	IL-4	MMP-9 [‡]	IFN- γ		IL-8
GM-CSF	IL-12	MIG	TGF- β 3 [‡]	IL-5	RANTES	IL-1 α		IL-10
ENA-78 [‡]	IL-12p40	MIP-1 α	TIMP-1 [*]	IL-6	TNF- α	IL-1 β		MCP-1 [‡]
GRO- α	IL-12p70	MIP-1 β	TIMP-2 [‡]	IL-10	TNF-RI	IL-2 [‡]		TNF- α
IFN- α 2a	IL-13	MIP-3 α	TNF- α	IL-12	TNF-RII	IL-4		
IFN- β	IL-15	MIP-3 β [‡]	TNF-RI	IL-12p40		IL-5		
IFN- γ	IL-16	MMP-1	TNF-RII	IL-12p70		IL-6		
IL-1 α	IL-17	MMP-2	TSLP [‡]	IL-13		IL-10 [‡]		
IL-1 β	IL-18 [‡]	MMP-3	TWEAK [‡]	IL-17		IL-13		
IL-1ra [‡]	IL-23 [‡]	MMP-8 [‡]		KC/GRO/CINC (CXCL1)		KC/GRO/CINC (CXCL1)		
IL-2	IP-10	MMP-9		MCP-1		MCP-1		
IL-4	I-TAC	MMP-10		M-CSF [‡]		MIP-3 α		
IL-5	M-CSF	NGAL [‡]						

[‡] Available through our Prototype Printing Services ^{*} Qualified Kit

Multiplex Panels		Panel	Analytes
Human TH1/TH2 7-Plex	IFN- γ , IL-2, IL-4, IL-5, IL-10, IL-12p70, IL-13	Human Chemokine 9-Plex	Eotaxin, Eotaxin-3, IL-8, IP-10, MCP-1, MCP-4, MDC, MIP-1 β , TARC
Human TH1/TH2 10-Plex	IFN- γ , IL-1 β , IL-2, IL-4, IL-5, IL-8, IL-10, IL-12p70, IL-13, TNF- α	Human Demonstration 4-Plex	IL-1 β , IL-2, IL-6, TNF- α
Human Proinflammatory 4-Plex I	IFN- γ , IL-1 β , IL-6, TNF- α	Human Demonstration 7-Plex	GM-CSF, IL-1 β , IL-2, IL-5, IL-6, IL-8, TNF- α
Human Proinflammatory 4-Plex II	IL-1 β , IL-6, IL-8, TNF- α	Human Demonstration 10-Plex	GM-CSF, IL-1 β , IL-2, IL-4, IL-5, IL-6, IL-8, IL-10, IL-12p70, TNF- α
Human Proinflammatory 7-Plex	IFN- γ , IL-1 β , IL-6, IL-8, IL-10, IL-12p70, TNF- α	Mouse TH1/TH2 9-Plex	IFN- γ , IL-1 β , IL-2, IL-4, IL-5, KC/GRO/CINC, IL-10, IL-12 total, TNF- α
Human Proinflammatory 9-Plex	GM-CSF, IFN- γ , IL-1 β , IL-2, IL-6, IL-8, IL-10, IL-12p70, TNF- α	Mouse Proinflammatory 7-Plex	IFN- γ , IL-1 β , IL-6, IL-10, IL-12p70, KC/GRO/CINC, TNF- α
Human MMP 2-Plex	MMP-2, MMP-10	Rat Demonstration 7-Plex	IFN- γ , IL-1 β , IL-4, IL-5, IL-13, KC/GRO/CINC, TNF- α
Human MMP 3-Plex	MMP-1, MMP-3, MMP-9	Canine Proinflammatory Panel 3	IL-2, IL-6, IL-8, TNF- α
Human Chemokine 7-Plex	Eotaxin, IL-8, IP-10, MCP-1, MCP-4, MIP-1 β , TARC		

Oncology & Vascular Markers				Multiplex Panels
AFP (human)	EPO (human, mouse/rat)	β -NGF (human) [‡]	TGF- β 3 (human) [‡]	Acute Phase Protein Panel 1 (rat) (AZM, AGP) [*]
Angiopoietin 1 (human) [‡]	sErbB2 (human)	Osteopontin (human)	TIMP-1 (human [*] , rat [‡])	c-Kit (pY721), Total c-Kit (human)
Angiopoietin 2 (human) [‡]	sFas (human) [‡]	p21 (human) [‡]	Thrombomodulin (human)	EPO, VEGF (human)
E-Cadherin (human)	FGF acidic (human) [‡]	uPA (human) [‡]	sVCAM-1 (human)	Growth Factor Panel I (human) (bFGF, VEGF, sFlt-1, PIGF)
P-Cadherin (human)	bFGF (human)	PEDF/Serpin F1 (human) [‡]	VEGF _(121/165) (human, mouse/rat)	Growth Factor Panel II (human) (c-Kit, KDR)
Cancer Antigen 125 (human)	sFlt-1 (sVEGFR1) (human)	PIGF (human)	VEGF-C (human) [‡]	Hypoxia Panel (mouse/rat) (EPO, VEGF)
Cancer Antigen 19.9 (human)	HGF (human)	PSA (human) [‡]	VEGF-D (human) [‡]	Hypoxia Panel (human) (EPO, IGFBP-1, VEGF)
Cancer Antigen 15.3 (human) [‡]	HIF-1 α (human)	SAA (human)		Vascular Injury Panel I (human) (sICAM-3, E-Selectin, P-Selectin, Thrombomodulin)
CEA (human)	IGFBP-1 (human)	E-Selectin (human)		Vascular Injury Panel II (human) (CRP, sICAM-1, sVCAM-1, SAA)
c-Kit (human)	sICAM-1 (human)	L-Selectin (rat) [‡]		
CRP (human)	sICAM-3 (human)	P-Selectin (human)		
EGF (human) [‡]	KDR (sVEGFR2) (human)	TGF- β 1 (human)		
sEGFR (human)	sMet (human) [‡]	TGF- β 2 (human) [‡]		
Endoglin (human) [‡]	MMP-9 (human)			

For a complete list of products please check our website at www.mesoscale.com

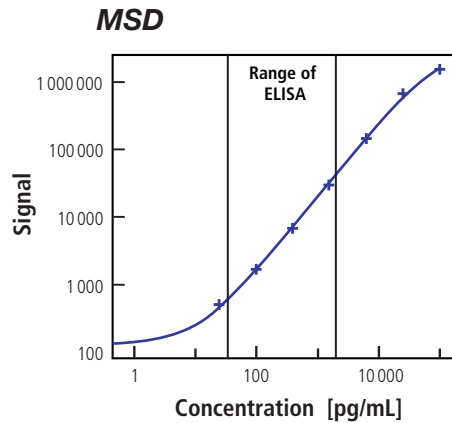
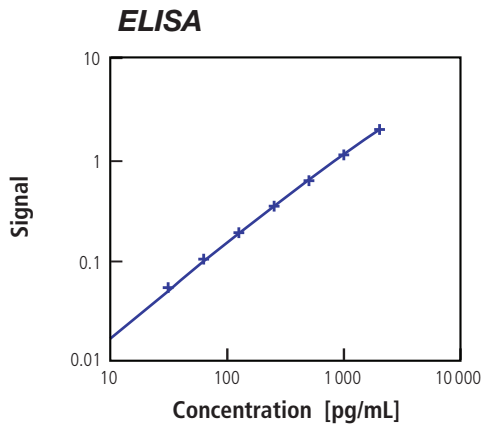
Prototype Printing Services

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. Coating antibodies expedites the conversion of ELISAs to the MSD platform with ease. An example of conversion of an ELISA to an MSD assay is demonstrated here. Customers routinely see significant improvements in assay performance when ELISAs are converted to the MSD format.

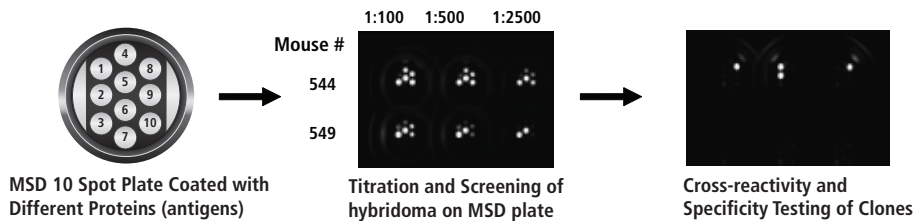


How to Order

- STEP 1:** Obtain a quote from MSD for your analyte of interest by contacting MSD Customer Service
- STEP 2:** Place your order Reference the quote number on your order.
- STEP 3:** Receive your Prototype assay

2. Proteins and Peptides

The MSD platform is very effective for screening antibodies. By testing antibodies against multiple antigens, customers can test for cross-reactivity and specificity in the same well. The example shown here 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.



3. Carbohydrates and Polysaccharides

Carbohydrates have been successfully coated on MSD plates, which has enabled the testing of different Pneumococcal vaccines by Marchese, et al.¹⁰ MSD plates coated with lipopolysaccharides have been used by Thompson, et al.¹¹ for serodiagnosis of Brucellosis in ruminants. Other than the above mentioned materials, MSD plates are highly amenable to coating with viral proteins, cell lysates, etc.

MSD's Prototype Printing Services Offer:

- 1. ELISA Conversion Packages:** MSD offers conversion packages for assay development.
- 2. Coating Optimization Package:** MSD can provide plates manufactured using several coating conditions on different plate surfaces as a means of determining the optimal conditions for a given assay.
- 3. On-site Scientific Support:** Support for assay development can be obtained from our Field Application Scientists and the Scientific Support team (ScientificSupport@mesoscale.com or 301-947-2025).
- 4. MSD Early Access Assays:** These are assays for which MSD has identified a working antibody pair and calibrator. MSD can provide suggestions for assay diluents and protocol.

References

- ¹ Gowan, S.M., Hardcastle, A., Hallsworth, A.E., Valenti, M.R., Hunter, L.J., de Haven Brandon, A.K., Garrett, M.D., Raynaud, F., Workman, P., Aherne, W., Eccles, S.A. **Application of Meso Scale Technology for the Measurement of Phosphoproteins in Human Tumor Xenografts.** *Assay and Drug Dev Technol.* 2007 Jun; Vol. 5 (3):391-401. 
- ² Raynaud, F.I., Eccles, S.A., Patel, S., Alix, S., Box, G., Chuckowree, I., Folkes, A., Gowan, S., De Haven Brandon, A., Di Stefano, F., Hayes, A., Henley, A.T., Lensun, L., Pergl-Wilson, G., Robson, A., Saghir, N., Zhyvoloup, A., McDonald, E., Sheldrake, P., Shuttleworth, S., Valenti, M., Wan, N.C., Clarke, P.A., Workman, P. **Biological properties of potent inhibitors of class I phosphatidylinositide 3-kinases: from PI-103 through PI-540, PI-620 to the oral agent GDC-0941.** *Mol Cancer Ther.* 2009 Jul; Vol. 8(7):1725-1738. 
- ³ Liang, M., Klakamp S.L., Funelas, C., Lu, H., Lam, B., Herl, C., Umble, A., Drake, A.W., Pak, M., Ageyeva, N., Pasumarthi, R., Roskos, L. **Detection of High- and Low-Affinity Antibodies Against a Human Monoclonal Antibody Using Various Technology Platforms.** *Assay and Drug Dev Technol.* 2007 Oct; Vol. 5(5): 1-8. 
- ⁴ Cao, L., Yu, Y., Darko, I., Currier, D., Mayeenuddin, L.H., Wan, X., Khanna, C., Helman, L.J. **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 Oct; Vol. 68:8039-8048. 
- ⁵ Tan, D.S.W., Thomas, G.V., Garrett, M.D., Banerji, U., De Bono, J.S., Kaye, S.B., Workman, P., **Biomarker-Driven Early Clinical Trials in Oncology A Paradigm Shift in Drug Development.** *The Cancer Journal.* 2009 Sep-Oct; Vol. 15, Number 5. 
- ⁶ Ghosh, T.K., Michelson, D.J., Solberg, J.C., Lipson, K.E., Inglefield, J.R., Alkan, S.S. **TLR-TLR cross talk in human PBMC resulting in synergistic and antagonistic regulation of type-1 and 2 interferons, IL-12 and TNF-alpha.** *Int Immunopharmacol.* 2007 Aug; Vol. 7(8):1111-21. 
- ⁷ Kelly, R.J., Rajan, A., Force, J., Keen, C., Cao, L., Yu, Y., Choyke, P.L., Turkbey, B., Raffeld, M., Xi, L., Steinberg, S.M., Wright, J.J., Lopez-Chavez, A., Kummar, S., Gutierrez, M., Giaccone, G. **Evaluation of KRAS mutations, angiogenic biomarkers and DCE-MRI in patients with advanced non-small cell lung cancer receiving sorafenib.** *Clin Cancer Res.* 2011 Mar 1;17(5):1190-9. 
- ⁸ Bafadhel, M., Saha, S., Siva, R., McCormick, M., Monteiro, W., Rugman, P., Dodson, P., Pavord, I.D., Newbold, P., Brightling, C.E. **Sputum IL-5 Concentration Is Associated with a Sputum Eosinophilia and Attenuated by Corticosteroid Therapy in COPD.** *Respiration.* 2009; Vol. 78:256-262. 
- ⁹ Fichorova R.N., Richardson-Harman, N., Alfano, M., Belec, L., Carbonneil, C., Chen, S., Cosentino, L., Curtis, K., Dezzutti, C.S., Donoval, B., Doncel, G.F., Donaghay, M., Grivel, J.C., Guzman, E., Hayes, M., Herold, B., Hillier, S., Lackman-Smith, C., Landay, A., Margolis, L., Mayer, K.H., Pasicznyk, J.M., Pallansch-Cokonis, M., Poli, G., Reichelderfer, P., Roberts, P., Rodriguez, I., Saidi, H., Sassi, R.R., Shattock, R., Cummins, J.E. Jr. **Biological and technical variables affecting immunoassay recovery of cytokines from human serum and simulated vaginal fluid: a multicenter study.** *Anal Chem.* 2008 Jun; Vol. 80(12): 4741-51. 
- ¹⁰ Marchese, R.D., Puchalski, D., Miller, P., Antonello, J., Hammond, O., Green, T., Rubinstein, L.J., Caulfield, M.J., Sikkema, D. **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 Mar; Vol. 16 (3):387-96. 
- ¹¹ Thompson, W.L., Karpus W.J., Van Eldik, L.J. **MCP-1-deficient mice show reduced neuroinflammatory responses and increased peripheral inflammatory responses to peripheral endotoxin insult.** *J Neuroinflam.* 2008 Aug; Vol. 5:35. 

 References using MSD MULTI-ARRAY technology



The MSD Advantage

Assay Solutions and Services



- Over 300 single analyte kits
- Over 200 multiple analyte kits
- Customized multiplexed 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 user calibration required
- Reliable measurements
- Integrated data analysis tool
- Comprehensive validation packages
- Software support for 21 CFR Part 11 compliance
- Comprehensive service plans



Ordering Information

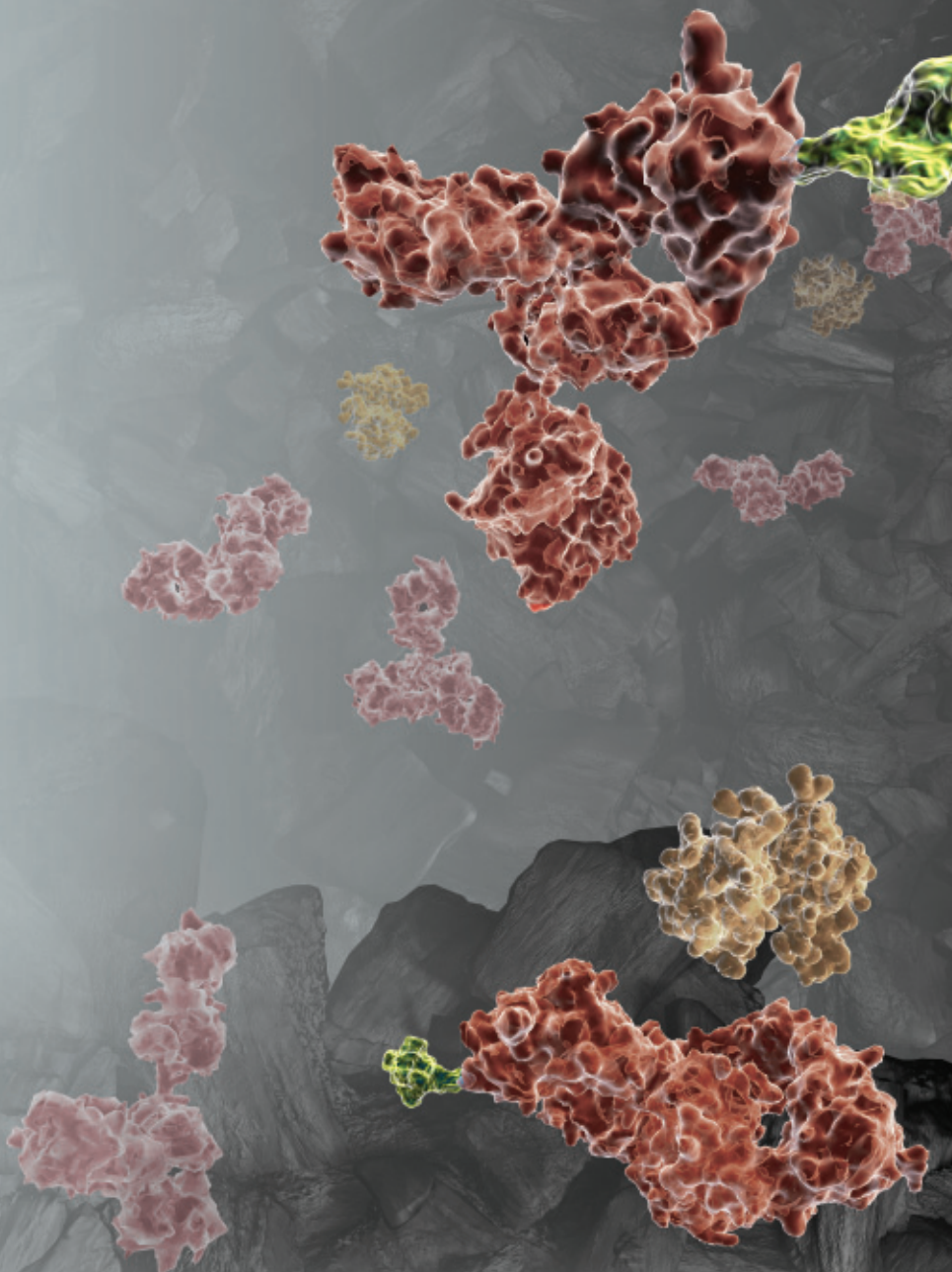
Purchase orders can be telephoned, faxed or emailed to the following:

TEL: 1.301.947.2085
FAX: 1.301.990.2776
EMAIL: customerservice@mesoscale.com

For a complete listing of our products,
visit our online catalog at

www.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

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, WWW.MESOSCALE.COM, MSD, MSD (DESIGN), DISCOVERY WORKBENCH, QUICKPLEX, MULTI-ARRAY, MULTI-SPOT, SULFO-TAG, SECTOR, SECTOR HTS, SECTOR PR, SMALL SPOT (DESIGN), 96 WELL 1, 4, 7, & 10-SPOT (DESIGNS), 384 WELL 1 & 4-SPOT (DESIGNS) and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.

© 2011 Meso Scale Diagnostics, LLC. All rights reserved.
19505-v4-2011.Jul

MESO SCALE DIAGNOSTICS®