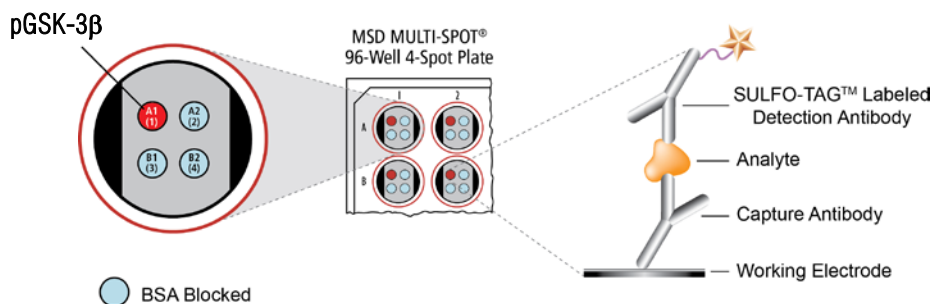


# MSD® Phospho-GSK-3β (Ser9) Assay Whole Cell Lysate Kit

For quantitative determination in human, mouse, and rat whole cell lysate samples

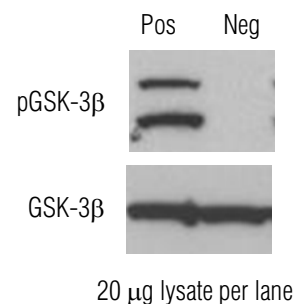
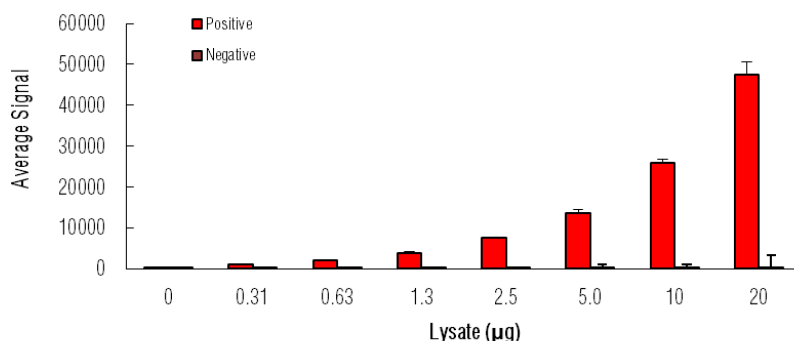


**Glycogen synthase kinase-3 (GSK-3)** is a serine/threonine protein kinase that is found in two cellular isoforms  $-\alpha$  and  $-\beta$ . GSK-3 has diverse cellular effects including involvement in metabolism, embryonic development, and cell survival. The two isoforms are regulated through phosphorylation, with inhibition as a result of growth factor and insulin-mediated phosphorylation by Akt on Ser 21 (GSK-3 $\alpha$ ) and Ser 9 (GSK-3 $\beta$ ). The inhibition of GSK-3 $\alpha$ /GSK-3 $\beta$  results in the dephosphorylation and activation of substrates such as glycogen synthase, eIF-2B, and C/EBP $\alpha$  causing increased protein and glycogen synthesis. Tyrosine (216) phosphorylation of GSK-3 $\beta$  results in its activation and the subsequent phosphorylation of various cellular proteins including cyclin D-1 and  $\beta$ -catenin. An important member of the Wnt signaling pathway, GSK-3 plays a role in cell fate in early embryonic development. GSK-3 $\beta$  has also been implicated in the progression of Alzheimer's disease through the phosphorylation of the microtubule-associated protein tau.

The MSD Phospho-GSK-3 $\beta$  (Ser9) Assay is available on 96-well 4-Spot plates. This datasheet outlines the performance of the assay.

## Typical Data

Representative results for the Phospho-GSK-3 $\beta$  (Ser9) Assay are illustrated below. The signal and ratio values provided below are example data; individual results may vary depending upon the samples tested. Western blot analyses of each lysate type were performed with Phospho-GSK-3 $\beta$  (Ser9) and total GSK-3 $\beta$  antibodies and are shown below for comparison. Logarithmically growing Jurkat cells (positive) were treated with LY294002 (50  $\mu$ M; 2.5 hours) and staurosporine (1  $\mu$ M; 2.5 hours) (negative). Whole cell lysates were added to MSD MULTI-SPOT® 4-Spot plates coated with anti-phospho-GSK-3 $\beta$  antibody on one of the four spatially distinct electrodes within a well. Phosphorylated GSK-3 $\beta$  was detected with anti-total GSK-3 $\beta$  antibody conjugated with MSD SULFO-TAG™ reagent.



**Fig. 1:** Sample data generated with the MULTI-ARRAY® Phospho-GSK-3 $\beta$  (Ser9) Assay. Increased signal is observed with the titration of pGSK-3 $\beta$  positive cell lysate. Signal for negative lysate remains low throughout the titration. The Phospho-GSK-3 $\beta$  (Ser9) Assay provides a quantitative measure of the data obtained with the traditional Western blot.

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

## Catalog Numbers

Phospho-GSK-3 $\beta$  (Ser9)  
Assay: Whole Cell Lysate Kit

### Kit size

1 plate	K150CQD-1
5 plates	K150CQD-2
20 plates	K150CQD-3

Phospho-GSK-3 $\beta$  Whole  
Cell Lysate Set

200 $\mu$ g	C11CQ-1
-------------	---------

## Ordering information

MSD Customer Service  
Phone: 1-301-947-2085  
Fax: 1-301-990-2776  
Email: CustomerService@mesoscale.com

## Company Address

MESO SCALE DISCOVERY®  
A division of  
Meso Scale Diagnostics, LLC.  
9238 Gaither Road  
Gaithersburg, MD 20877 USA

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

For Research Use Only.  
Not for use in diagnostic  
procedures.

# MSD Phosphoprotein Assays

## Lysate Titration

Data for pGSK-3 $\beta$  positive and negative Jurkat cell lysates using the MULTI-ARRAY Phospho-GSK-3 $\beta$  (Ser9) Assay are presented below.

Lysate ( $\mu$ g)	Positive			Negative			P/N
	Average Signal	StdDev	%CV	Average Signal	StdDev	%CV	
0	87	13	15.1	81	9	10.5	
0.31	1231	86	7.0	101	11	10.6	12
0.63	2215	116	5.3	117	6	5.1	19
1.3	3881	300	7.7	132	2	1.8	30
2.5	7579	213	2.8	161	3	1.6	47
5.0	13629	869	6.4	167	15	9.1	82
10	25872	871	3.4	191	14	7.2	135
20	47551	3109	6.5	211	12	5.6	226

## MSD Advantage

- **Multiplexing:** Multiple analytes can be measured in one well using typical sample amounts of 25  $\mu$ g/well or less without compromising speed or performance
- **Large dynamic range:** Linear range of up to five logs enables the measurement of native levels of biomarkers in normal and diseased samples without multiple dilutions
- **Minimal background:** The stimulation mechanism (electricity) is decoupled from the signal (light)
- **Simple protocols:** Only labels near the electrode surface are detected, enabling no-wash assays
- **Flexibility:** Labels are stable, non-radioactive, and conveniently conjugated to biological molecules
- **High sensitivity and precision:** Multiple excitation cycles of each label enhance light levels and improve sensitivity

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

## References using MSD's platform for the measurement of phosphoproteins

1. Beltran PJ, Chung YA, Moody G, Mitchell P, Cajulis E, Vonderfecht S, Kendall R, Radinsky R, Calzone FJ. Efficacy of Ganitumab (AMG 479), Alone and in Combination With Rapamycin, in Ewing's and Osteogenic Sarcoma Models. *J Pharmacol Exp Ther.* 2011 Jun;337(3):644-54.
2. Grimshaw KM, Hunter LJ, Yap TA, Heaton SP, Walton MI, Woodhead SJ, Fazal L, Reule M, Davies TG, Seavers LC, Lock V, Lyons JF, Thompson NT, Workman P, Garrett MD. AT7867 is a potent and oral inhibitor of AKT and p70 S6 kinase that induces pharmacodynamic changes and inhibits human tumor xenograft growth. *Mol Cancer Ther.* 2010 May;9(5):1100-10.
3. Guillard S, Clarke PA, Te Poele R, Mohri Z, Bjerke L, Valenti M, Raynaud F, Eccles SA, Workman P. Molecular pharmacology of phosphatidylinositol 3-kinase inhibition in human glioma. *Cell Cycle.* 2009 Feb 1;8(3):443-53.
4. Gowan SM, Hardcastle A, Hallsworth AE, Valenti MR, Hunter LJ, de Haven Brandon AK, Garrett MD, Raynaud F, Workman P, Aherne W, Eccles SA. Application of meso scale technology for the measurement of phosphoproteins in human tumor xenografts. *Assay Drug Dev Technol.* 2007 Jun;5(3):391-401.

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, 4-SPOT (DESIGN) and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.  
© 2011 Meso Scale Diagnostics, LLC. All rights reserved.

For Research Use Only. Not for use in diagnostic procedures.

