

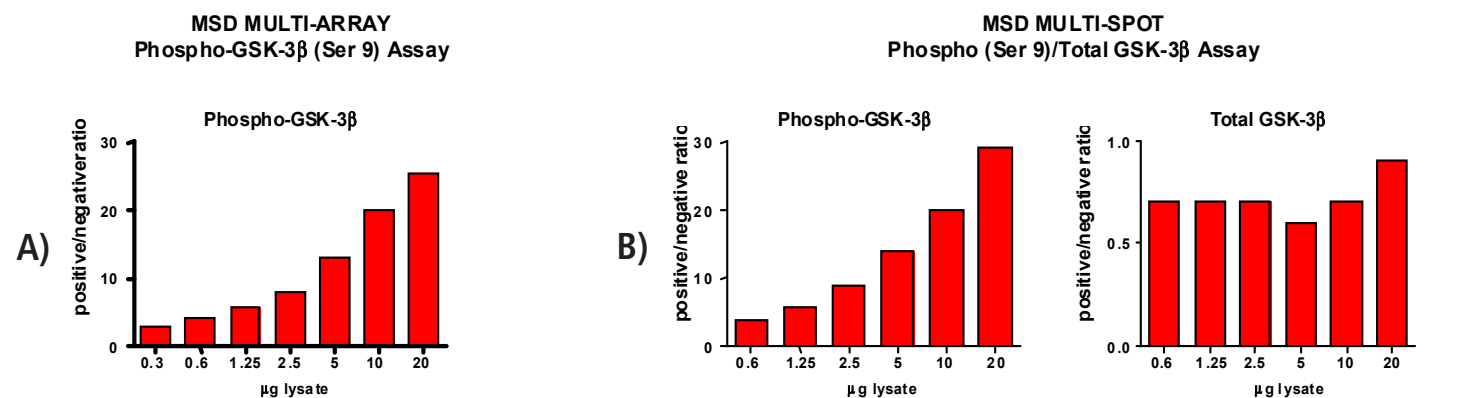
# Meso Scale Discovery® Whole Cell Lysate Set

## Phospho-GSK-3β

<b>Catalog No:</b>	C11CQ-1
<b>Contents:</b>	2 x 100 µg MSDLY0006 pGSK-3β Negative Control Cell Lysate Cell lysate from growing Jurkat cells treated with 50 µM LY294002 and 1 µM staurosporine for 2.5 hours to inhibit GSK-3β phosphorylation  2 x 100 µg MSDLY0001 pGSK-3β Positive Control Cell Lysate Cell lysate from growing Jurkat cells
<b>Concentration:</b>	2 mg/mL in MSD® Complete Tris Lysis Buffer
<b>Volume:</b>	2 vials (50 µL) negative lysate    2 vials (50 µL) positive lysate
<b>Preparation:</b>	Following cell treatment, Jurkat cell lysates were prepared on ice in MSD Complete Tris Lysis Buffer. Cell debris was cleared by centrifugation.
<b>Storage:</b>	Lysates should be stored at -80°C. Lysates will retain approximately 90% of activity after a single round of freeze thaw if handled properly (thawed on ice and immediately refrozen in smaller aliquots).
<b>Quality Control:</b>	Lysates have been tested for performance in Western Blot and MSD Assays.

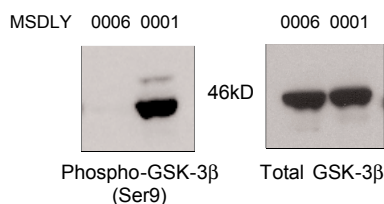
### MSD MULTI-SPOT Assay Results

Figures A and B (below) illustrate typical titrations for MSDLY0006 (pGSK-3β negative) and MSDLY0001 (pGSK-3β positive) cell lysates. Figure A depicts the MSD MULTI-ARRAY Phospho-GSK-3β (Ser 9) Whole Cell Lysate Kit; Figure B depicts the MSD MULTI-SPOT Phospho (Ser 9)/Total GSK-3β Whole Cell Lysate Kit. The results are presented as a ratio of pGSK-3β positive to pGSK-3β negative lysate signals. In both assays, the phospho-GSK-3β signal ratios increase with the titration of lysates. Total GSK-3β signal ratios in the Phospho/Total GSK-3β Assay remain at a constant level close to one throughout the titration. The representative results shown below are for demonstration purposes only and individual results may vary depending upon experimental application.



### Traditional Western Blot Results

MSDLY0006 and MSDLY0001 whole cell lysates (20 µg each) were analyzed by Western Blot with Phospho (Ser 9)-specific and total GSK-3β antibodies.



FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC OR THERAPEUTIC PROCEDURES.

20196-v1-2008May