

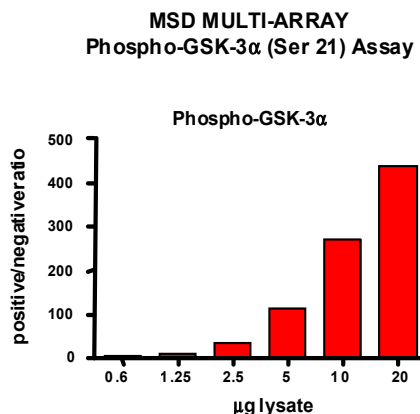
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-GSK-3 α

Catalog No:	C11CO-1
Contents:	2 x 100 μ g MSDLY0006 pGSK-3 α Negative Control Cell Lysate Cell lysate from growing Jurkat cells treated with 50 μ M LY294002 and 1 μ M staurosporine for 2.5 hours 2 x 100 μ g MSDLY0001 pGSK-3 α Positive Control Cell Lysate Cell lysate from growing Jurkat cells
Concentration:	2 mg/mL in MSD® Complete Tris Lysis Buffer
Volume:	2 vials (50 μ L) negative lysate 2 vials (50 μ L) positive lysate
Preparation:	Following cell treatment, Jurkat cell lysates were prepared on ice in MSD Complete Tris Lysis Buffer. Cell debris was cleared by centrifugation.
Storage:	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).
Quality Control:	Lysates have been tested for performance in Western Blot and MSD MULTI-ARRAY® Assays.

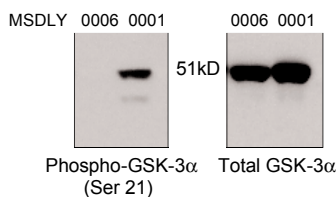
MSD MULTI-ARRAY Assay Results

The figure below illustrates typical lysate titrations for MSDLY0006 (pGSK-3 α negative) and MSDLY0001 (pGSK-3 α positive) cell lysates using the MSD MULTI-ARRAY Phospho-GSK-3 α Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with pGSK3 α positive and pGSK-3 α negative lysates. The phospho-GSK-3 α signal ratios increase with the amount of lysate. 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 μ g each) were analyzed by Western Blot with Phospho (Ser 21)-specific and Total GSK-3 α antibodies.



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

20195-v2-2008May