

Meso Scale Discovery® Whole Cell Lysate Set

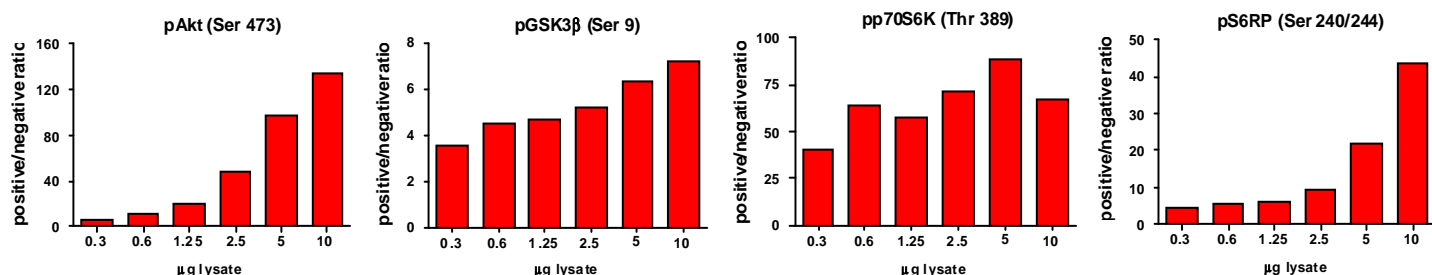
Phospho-Akt Signaling Panel 2

Catalog No:	C1177-1
Contents:	2 x 100 µg MSDLY0066 pAkt/ pGSK-3β/ pp70S6K / pS6RP Negative Cell Lysate Cell lysate from MCF-7 cells treated with 50 µM LY294002 for 2.5 hours 2 x 100 µg MSDLY0067 pAkt/ pGSK-3β/ pp70S6K / pS6RP Positive Cell Lysate Cell lysate from MCF-7 cells treated with 100 nM IGF-1 for 20 minutes to stimulate phosphorylation
Concentration:	2 mg/mL in MSD® Complete Tris Lysis Buffer
Volume:	2 vials (50 µL) negative lysate 2 vials (50 µL) 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-SPOT® Assays.

MSD MULTI-SPOT Assay Results

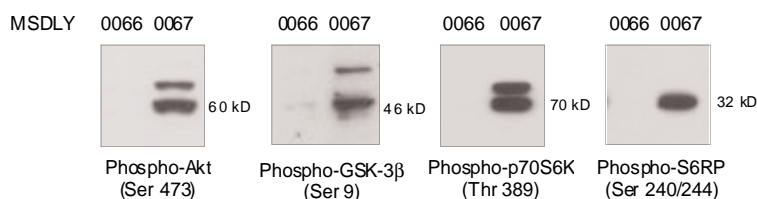
The figure below illustrates typical lysate titrations for MSDLY0066 (negative) and MSDLY0067 (positive) using the MSD MULTI-SPOT Akt Signaling Panel 2 Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with phospho-Akt/phospho-GSK-3β/phospho-p70S6K/phospho-S6RP positive and negative lysates. In all cases, the signal ratios increase with the amount of lysates. The representative results shown below are for demonstration purposes only and individual results may vary depending upon experimental application.

MSD MULTI-SPOT Assays
Akt Signaling Panel 2



Traditional Western Blot Results

MSDLY0066 and MSDLY0067 whole cell lysates (20 µg each) were analyzed by Western Blot with phospho-specific Akt (Ser 473), GSK-3β (Ser 9), p70S6K (Thr 389), pS6RP (Ser 240/244) antibodies.



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

20216-v1-2008Aug