

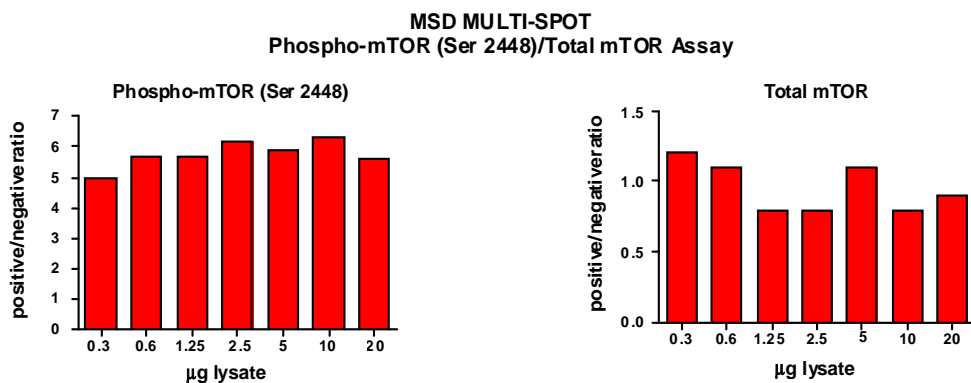
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-mTOR (Ser 2448)

Catalog No:	C10JE-1
Contents:	2 x 100 µg MSDLY0068 pmTOR Negative Cell Lysate Cell lysate from HEK293 cells treated with 100nM wortmannin for 3 hours to reduce basal mTOR phosphorylation 2 x 100 µg MSDLY0069 pmTOR Positive Cell Lysate Cell lysate from HEK293 cells treated with 1 µM PMA for 30 minutes to stimulate mTOR 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, HEK293 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.

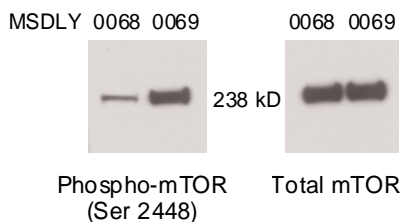
MSD MULTI-SPOT Assay Results

The figure below illustrates typical lysate titrations for MSDLY0068 (pmTOR negative) and MSDLY0069 (pmTOR positive) cell lysates using the MSD MULTI-SPOT Phospho (Ser2448)/Total mTOR Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with pmTOR positive and pmTOR negative lysates. The phospho-mTOR signal ratios remain constant with increasing amounts of lysate. The total mTOR remains 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

MSDLY0068 and MSDLY0069 whole cell lysates (20 µg each) were analyzed by Western Blot with phospho-specific and total mTOR antibodies.



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

20206-v1-2010Feb