

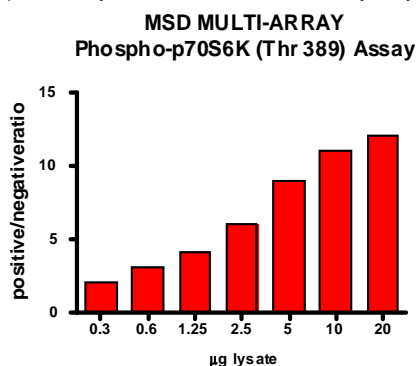
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-p70S6K (Thr 389)

Catalog No:	C11DN-1
Contents:	2 x 100 µg MSDLY0009 pp70S6K (Thr 389) Negative Cell Lysate Cell lysate from HEK293 cells treated with 1 µM rapamycin for 3 hours 2 x 100 µg MSDLY0040 pp70S6K (Thr 389) Positive Cell Lysate Cell lysate from HEK293 cells treated with 50 nM calyculin A for 30 minutes to stimulate p70S6K 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 Assays.

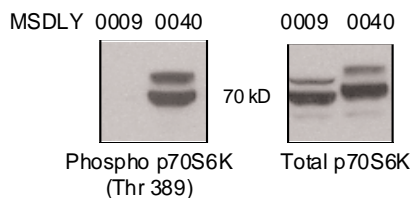
MSD MULTI-SPOT Assay Results

The figure below illustrates typical lysate titrations for MSDLY0009 (pp70S6K (Thr 389) negative) and MSDLY0040 (pp70S6K (Thr 389) positive) cell lysates using the MSD MULTI-ARRAY Phospho-p70S6K (Thr 389) Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with pp70S6K positive and pp70S6K negative lysates. The signal ratios for phospho-p70S6K (Thr 389) 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

MSDLY0009 and MSDLY0040 whole cell lysates (20 µg each) were analyzed by Western Blot with Phospho (Thr 389)-specific and Total p70S6K antibodies.



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

20218-v1-2008Aug