

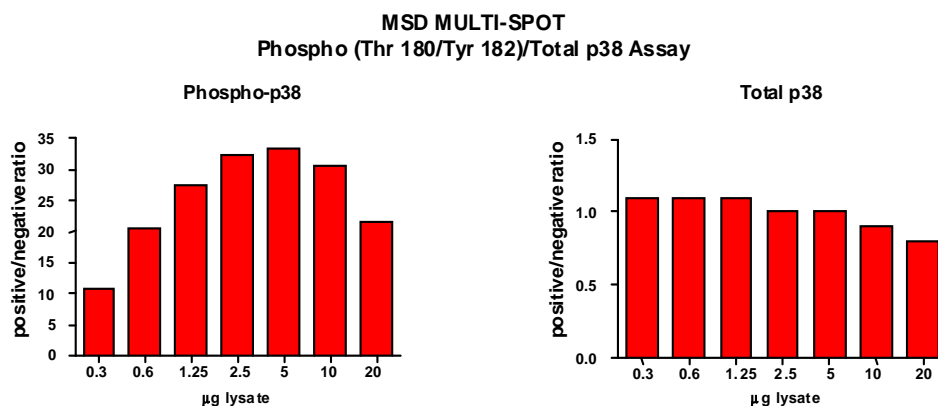
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-p38

Catalog No:	C11CY-1
Contents:	2 x 100 µg MSDLY0009 pp38 Negative Control Cell Lysate Cell lysate from growing HEK293 cells treated with 1 µM rapamycin for 3 hours to inhibit p38 phosphorylation 2 x 100 µg MSDLY0040 pp38 Positive Control Cell Lysate Cell lysate from growing HEK293 cells treated with 50 nM calyculin A for 30 minutes to stimulate p38 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.

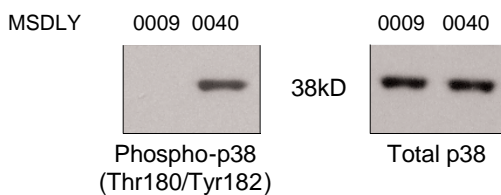
MSD MULTI-SPOT Assay Results

The figure below illustrates a typical titration for MSDLY0009 (pp38 negative) and MSDLY0040 (pp38 positive) cell lysates with the MSD MULTI-SPOT Phospho/Total p38 Whole Cell Lysate Kit. Results are presented as a ratio of the signals obtained with pp38 positive and pp38 negative lysates. The phospho-p38 signal ratios increase with the amount of lysates, and the signal ratios for total p38 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

MSDLY0009 and MSDLY0040 whole cell lysates (20 µg each) were analyzed by Western Blot with phospho-specific (Thr 180/ Tyr 182) and total p38 antibodies.



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

20051-v2-2008May