

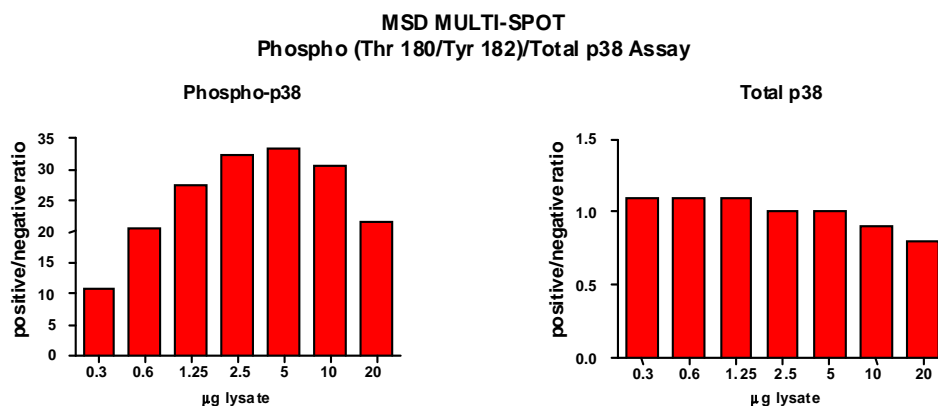
# Meso Scale Discovery® Whole Cell Lysate Set

## Phospho-p38

<b>Catalog No:</b>	C11CY-1
<b>Contents:</b>	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
<b>Concentration:</b>	2 mg/mL in MSD® Complete Tris Lysis Buffer
<b>Volume:</b>	2 vials (50 µL) negative lysate    2 vials (50 µL) positive lysate
<b>Preparation:</b>	Following cell treatment, HEK293 cell lysates were prepared on ice in MSD Complete Tris Lysis Buffer. Cell debris was cleared by centrifugation.
<b>Storage:</b>	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).
<b>Quality Control:</b>	Lysates have been tested for performance in Western Blot and MSD MULTI-SPOT® Assays.

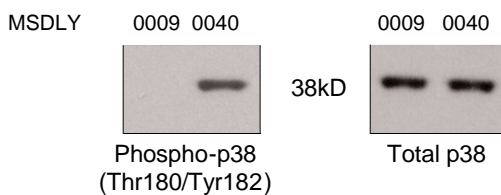
### MSD MULTI-SPOT Assay Results

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