

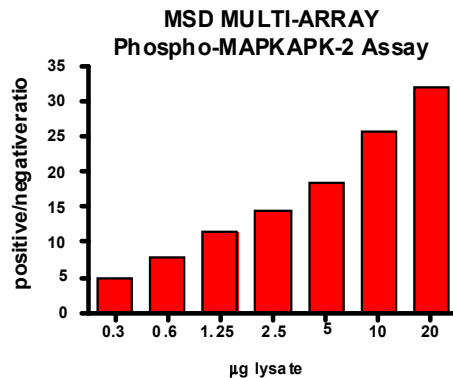
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-MAPKAPK-2

Catalog No:	C10FV-1
Contents:	2 x 100 µg MSDLY0009 pMAPKAPK-2 Negative Cell Lysate Cell lysate from HEK293 cells treated with 1 µM rapamycin for 3 hours to inhibit phosphorylation 2 x 100 µg MSDLY0040 pMAPKAPK-2 Positive Cell Lysate Cell lysate from HEK293 cells treated with 50 nM calyculin A for 30 minutes to stimulate MAPKAPK-2 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-ARRAY® Assays.

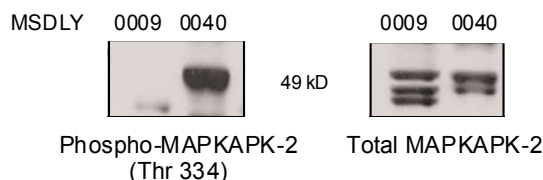
MSD MULTI-ARRAY Assay Results

The figure below illustrates typical lysate titrations for MSDLY0009 (pMAPKAPK-2 negative) and MSDLY0040 (pMAPKAPK-2 positive) cell lysates using the MSD MULTI-ARRAY Phospho-MAPKAPK-2 Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with pMAPKAPK-2 positive and pMAPKAPK-2 negative lysates. The phospho-MAPKAPK-2 signal ratios 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-specific MAPKAPK-2.



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

20205-v1-2008May