

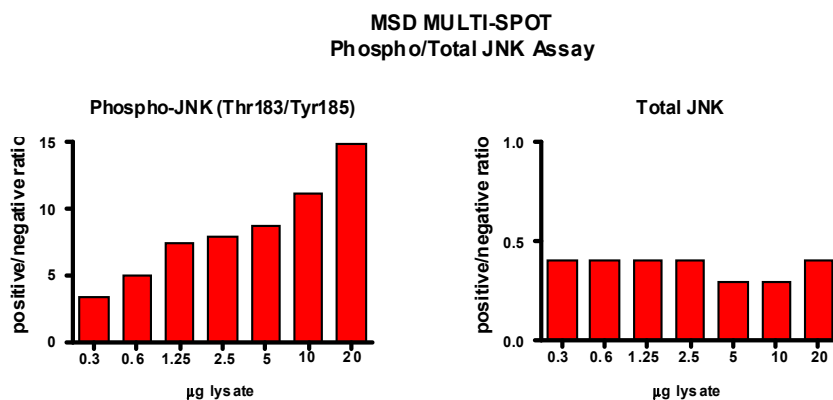
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-JNK

Catalog No:	C11CU-1
Contents:	2 x 100 µg MSDLY0009 pJNK Negative Control Cell Lysate Cell lysate from growing HEK293 cells treated with 1 µM rapamycin for 3 hours to inhibit JNK phosphorylation 2 x 100 µg MSDLY0007 pJNK Positive Control Cell Lysate Cell lysate from serum-deprived HEK293 cells harvested 30 minutes after UV irradiation (40 mJ/cm ²) to stimulate JNK 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.

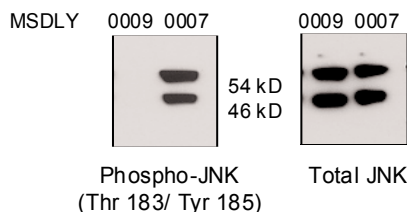
MSD MULTI-SPOT Assay Results

The figure below illustrates typical lysate titrations for MSDLY0009 (pJNK negative) and MSDLY0007 (pJNK positive) cell lysates using the MSD MULTI-SPOT Phospho/Total JNK Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with pJNK positive and pJNK negative lysates. The phospho-JNK signal ratios increase with the amount of lysate. The signal ratios for the total JNK remain at a constant level 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 MSDLY0007 whole cell lysates (20 µg each) were analyzed by Western Blot with Phospho-specific JNK and Total JNK antibodies.



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

20191-v2-2008May