

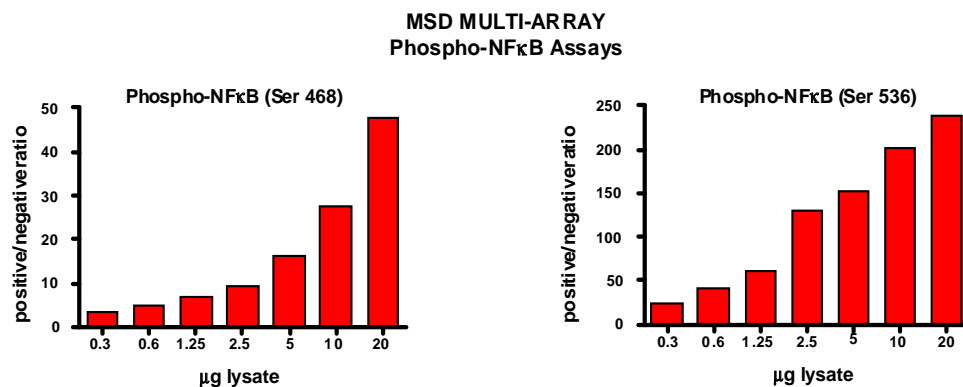
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-NFκB (Ser 468,536)

Catalog No:	C11ED-1
Contents:	2 x 100 µg MSDLY0001 pNFκB Kinase Negative Cell Lysate Cell lysate from growing Jurkat cells 2 x 100 µg MSDLY0051 pNFκB Kinase Positive Cell Lysate Cell lysate from Jurkat cells treated with 25 ng/mL TNF-α and 50 nM calyculin A for 15 minutes to stimulate NFκB 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, Jurkat 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.

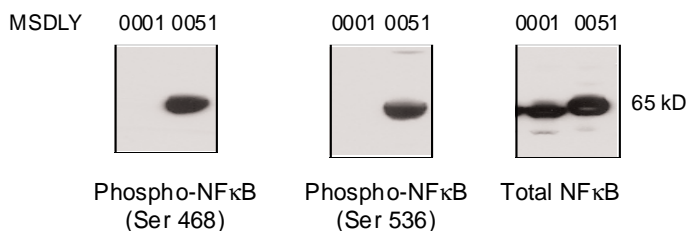
MSD MULTI-ARRAY Assay Results

The figure below illustrates typical lysate titrations for MSDLY0001 (pNFκB negative) and MSDLY0051 (pNFκB positive) cell lysates using the MSD MULTI-ARRAY Phospho-NFκB(Ser468) and Phospho-NFκB(Ser536) Whole Cell Lysate Kits. The results are presented as a ratio of the signals obtained with pNFκB positive and pNFκB negative lysates. The phospho-NFκB 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

MSDLY0001 and MSDLY0051 whole cell lysates (20 µg each) were analyzed by Western Blot with phospho-specific and total NFκB antibodies.



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

20207-v1-2010Feb