

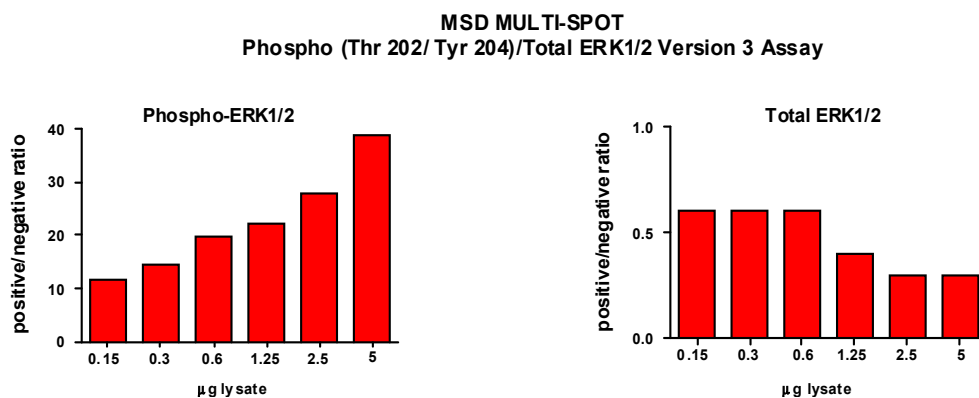
# Meso Scale Discovery® Whole Cell Lysate Set

## Phospho-ERK1/2

<b>Catalog No:</b>	C11CM-1
<b>Contents:</b>	2 x 100 µg MSDLY0004 pERK1/2 Negative Control Cell Lysate Cell lysate from growing Jurkat cells treated with LY294002 (50 µM) for 2.5 hours  2 x 100 µg MSDLY0003 pERK1/2 Positive Control Cell Lysate Cell lysate from growing Jurkat cells treated with PMA (200 nM) for 15 minutes
<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, Jurkat 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.

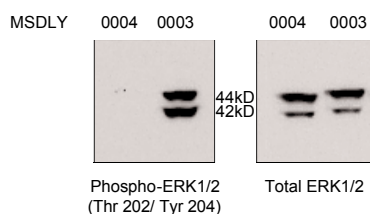
### MSD MULTI-SPOT Assay Results

The figure below illustrates a typical titration for MSDLY0004 (pERK1/2 negative) and MSDLY0003 (pERK1/2 positive) cell lysates with the MSD MULTI-SPOT Phospho/Total ERK1/2 (Version 3) Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with pERK1/2 positive and pERK1/2 negative lysates. The phospho-ERK1/2 signal ratios increase with the amount of lysates, and the signal ratios for total ERK1/2 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

MSDLY0004 and MSDLY0003 whole cell lysates (20 µg each) were analyzed by Western Blot with phospho-specific ERK1/2 and total ERK1/2 antibodies.



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

20039-v3-2008May