

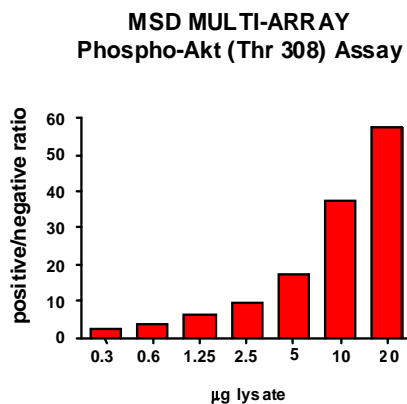
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-Akt (Thr 308)

Catalog No:	C11DY-1
Contents:	2 x 100 µg MSDLY0055 pAkt (Thr 308) Negative Control Cell Lysate Cell lysate from confluent NIH3T3 cells 2 x 100 µg MSDLY0056 pAkt (Thr 308) Positive Control Cell Lysate Cell lysate from confluent NIH3T3 cells treated with 100 ng/mL PDGF-BB for 5 minutes to stimulate Akt 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, NIH3T3 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.

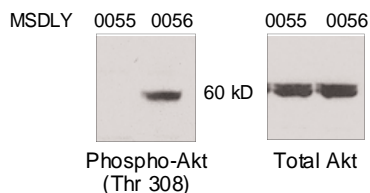
MSD MULTI-ARRAY Assay Results

The figure below illustrates a typical titration for MSDLY0055 (pAkt (Thr 308) negative) and MSDLY0056 (pAkt (Thr 308) positive) cell lysates with the MSD MULTI-ARRAY Phospho-Akt (Thr 308) Whole Cell Lysate Kit. Results are presented as a ratio of the signals obtained with pAkt (Thr 308) positive and pAkt (Thr 308) negative lysates. The phospho-Akt (Thr 308) signal ratios increase with the amount of lysates. The representative results shown below are for demonstration purposes only and individual results may vary depending upon experimental application.



Traditional Western Blot Results

MSDLY0055 and MSDLY0056 whole cell lysates (20 µg each) were analyzed by Western Blot with phospho-specific Akt threonine 308 and total Akt antibodies.



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

20049-v1-2008Aug