

Meso Scale Discovery® Whole Cell Lysate Set

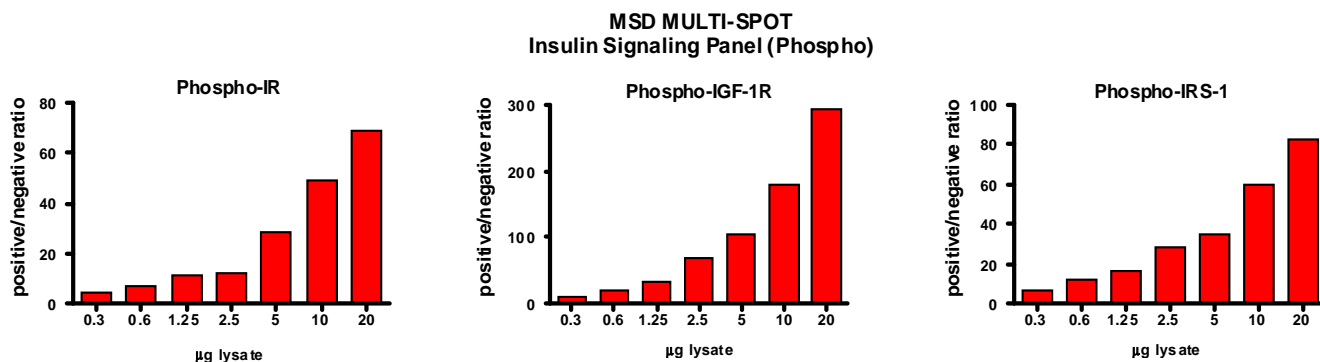
Insulin Signaling Panel

Catalog No:	C1151-1
Contents:	2 x 100 µg MSDLY0062 pIR/ pIGF-1R/ pIRS-1 Negative Cell Lysate Cell lysate from serum deprived MCF-7 cells 2 x 100 µg MSDLY0063 pIR/ pIGF-1R/ pIRS-1 Positive Cell Lysate Cell lysate from serum deprived MCF-7 cells treated with 100nM IGF-1 for 20 minutes to stimulate 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, MCF-7 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

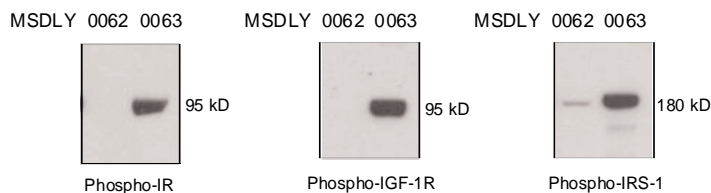
The figure below illustrates typical lysate titrations for MSDLY0062 (negative) and MSDLY0063 (positive) using the MSD MULTI-SPOT Insulin Signaling Panel (Phosphoprotein) Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with phospho-IR/phospho-IGF-1R/phospho-IRS-1 positive and negative lysates. In all cases, the 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.

Note: Lysate sets were also tested for performance on the MSD MULTI-SPOT Insulin Signaling Panel (Total protein) Whole Cell lysate kit (data not shown.)



Traditional Western Blot Results

MSDLY0062 and MSDLY0063 whole cell lysates (20 µg each) were analyzed by Western Blot with phospho (tyrosine)-specific IRS-1, IGF-1R and IR antibodies.



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

20215-v1-2008Aug