

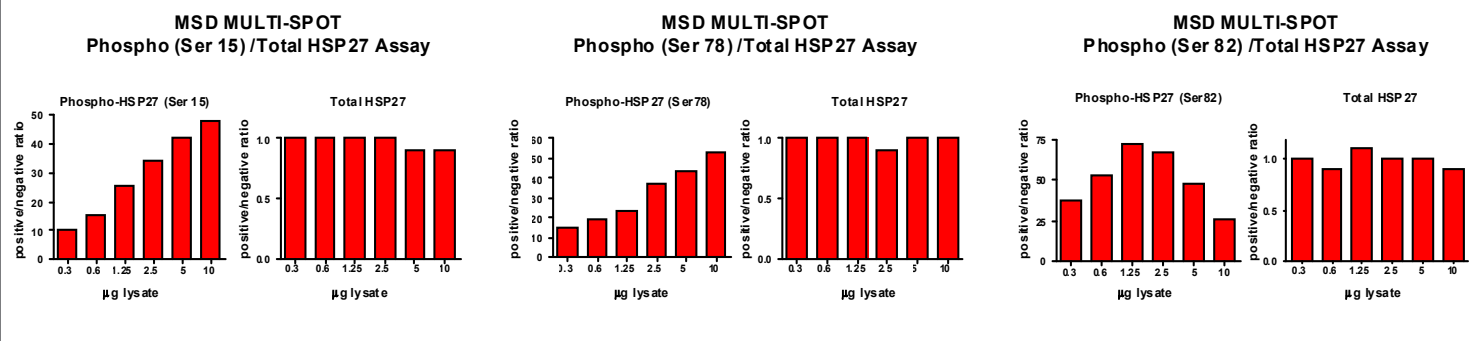
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-HSP27

Catalog No:	C11CS-1
Contents:	2 x 100 µg MSDLY0019 pHSP27 Negative Cell Lysate Cell lysate from growing HeLa cell monolayers 2 x 100 µg MSDLY0020 pHSP27 Positive Cell Lysate Cell lysate from growing HeLa cell monolayers treated with 0.4 M sorbitol for 30 minutes to stimulate HSP27 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, HeLa 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.

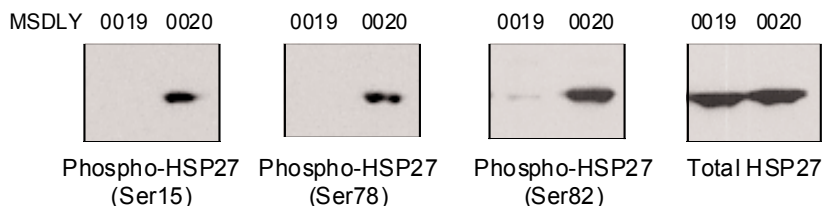
MSD MULTI-SPOT Assay Results

The figure below illustrates typical lysate titrations for MSDLY0019 (pHSP27 negative) and MSDLY0020 (pHSP27 positive) cell lysates using the MSD MULTI-SPOT Phospho (Ser15)/total HSP27, Phospho (Ser78)/total HSP27, Phospho (Ser82)/total HSP27 Whole Cell Lysate Kits. The results are presented as a ratio of the signals obtained with pHSP27 positive and pHSP27 negative lysates. The phospho HSP27 signal ratios increase with the amount of lysate. The signal ratios for total HSP27 remain at a constant level close to one 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

MSDLY0019 and MSDLY0020 whole cell lysates (20 µg each) were analyzed by Western Blot with Phospho-specific HSP27 and Total HSP27 antibodies.



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

20190-v2-2008May