

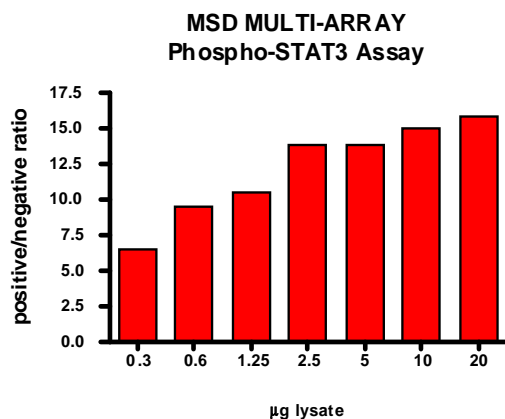
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-STAT3 (Tyr 705)

Catalog No:	C11DI-1
Contents:	2 x 100 µg MSDLY0019 pSTAT3 Negative Cell Lysate Cell lysate from growing HeLa cell monolayers 2 x 100 µg MSDLY0048 pSTAT3 Positive Cell Lysate Cell lysate from growing HeLa cell monolayers pretreated with 1 mM sodium vanadate for 4 hours followed by Oncostatin M (40 ng/mL) for 5 minutes to stimulate STAT3 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-ARRAY® Assays.

MSD MULTI-ARRAY Assay Results

The figure below illustrates a typical titration for MSDLY0019 (pSTAT3 negative) and MSDLY0048 (pSTAT3 positive) cell lysates with the MSD MULTI-ARRAY Phospho-STAT3 Whole Cell Lysate Kit. Results are presented as a ratio of the signals obtained with pSTAT3 positive and pSTAT3 negative lysates. In the phospho-STAT3 assay, the phospho-STAT3 signal ratio increases 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

MSDLY0019 and MSDLY0048 whole cell lysates (20 µg each) were analyzed by Western Blot with phospho specific (tyrosine 705) and total STAT3 antibodies.

