

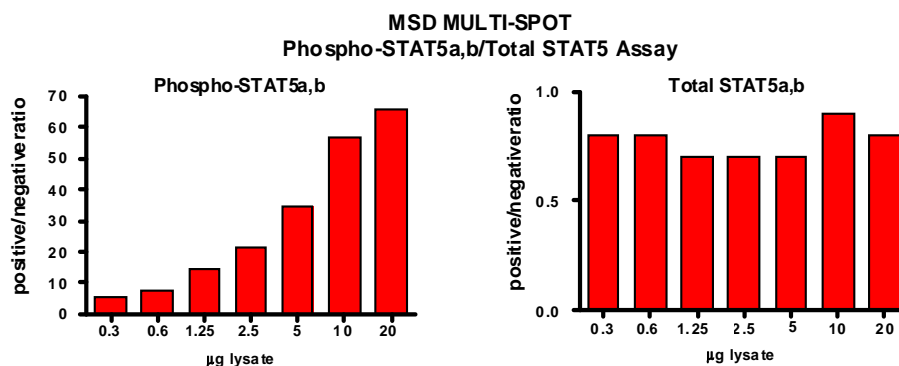
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-STAT5a,b (Tyr 694)

Catalog No:	C10IG-1
Contents:	2 x 100 µg MSDLY0019 pSTAT5a,b Negative Cell Lysate Cell lysate from growing HeLa cell monolayers 2 x 100 µg MSDLY0048 pSTAT5a,b Positive Cell Lysate Cell lysate from HeLa cell monolayers pre- treated with 1 mM Sodium Vanadate for 4 hours followed by Oncostatin M stimulation (40 ng/mL) for 5 minutes to stimulate STAT5 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 (pSTAT5a,b negative) and MSDLY0048 (pSTAT5a,b positive) cell lysates using the MSD MULTI-SPOT Phospho/Total STAT5a,b Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with pSTAT5a,b positive and pSTAT5a,b negative lysates. The phospho-STAT5a,b signal ratios increase with the amount of lysate. The signal ratios for total STAT5a,b 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 MSDLY0048 whole cell lysates (20 µg each) were analyzed by Western Blot with phospho-specific STAT5a,b and total STAT5a,b antibodies.

