

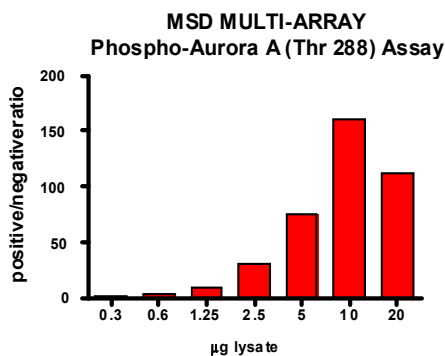
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-Aurora A (Thr 288)

Catalog No:	C10JC-1
Contents:	2 x 100 µg MSDLY0019 pAurora A Negative Cell Lysate Cell lysate from growing HeLa cell monolayers 2 x 100 µg MSDLY0060 pAurora A Positive Cell Lysate Cell lysate from HeLa cell monolayers treated with 1 µg/mL nocodazole for 19 hours followed by a 50 nM calyculin A treatment for 30 minutes to stimulate Aurora A 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.

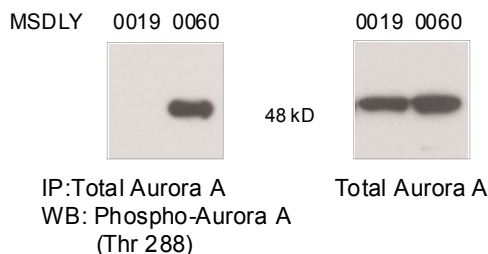
MSD MULTI-ARRAY Assay Results

The figure below illustrates typical lysate titrations for MSDLY0019 (pAurora A negative) and MSDLY0060 (pAurora A positive) cell lysates using the MSD MULTI-ARRAY Phospho-Aurora A (Thr288) Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with pAurora A positive and pAurora A negative lysates. The phospho-Aurora A signal ratios increase with the amount of lysate. 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 MSDLY0060 whole cell lysates (200 µg each) were analyzed by immunoprecipitation using a total antibody for IP and phospho-specific antibody for western blot. Whole cell lysates (20 µg) were analyzed using Total Aurora A antibody.



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

20200-v1-2008May