

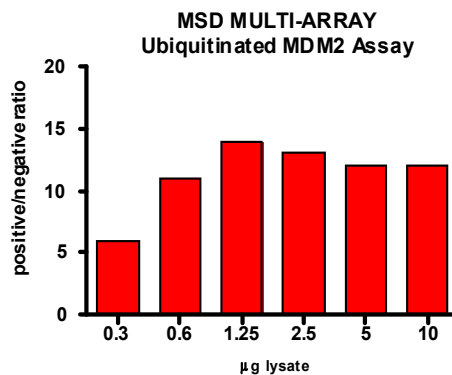
Meso Scale Discovery® Whole Cell Lysate Set

Ubiquitinated MDM2

Catalog No:	C12FJ-1
Contents:	2 x 100 µg MSDLY0058 Ubiquitinated MDM2 Negative Cell Lysate Cell lysate from growing HCT116 cells 2 x 100 µg MSDLY0059 Ubiquitinated MDM2 Positive Cell Lysate Cell lysate from HCT116 cells treated with 1 µM doxorubicin for 21 hours and 1 µM epoxomicin for 6 hours to accumulate ubiquitinated MDM2
Concentration:	2 mg/mL in MSD Complete Tris Lysis Buffer supplemented with 20 mM EDTA and 20 mM NEM
Volume:	2 vials (50 µL) negative lysate 2 vials (50 µL) positive lysate
Preparation:	Following cell treatment, HCT116 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.

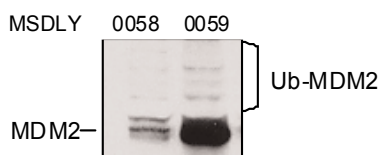
MSD MULTI-ARRAY Assay Results

The figure below illustrates typical lysate titrations for MSDLY0058 (ubiquitinated MDM2 negative) and MSDLY0059 (ubiquitinated MDM2 positive) cell lysates using the MSD MULTI-ARRAY Ubiquitinated MDM2 Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with ubiquitinated MDM2 positive and ubiquitinated MDM2 negative lysates. The ubiquitinated MDM2 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

MSDLY0058 and MSDLY0059 whole cell lysates (20 µg each) were analyzed by Western Blot with total MDM2 antibodies.



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

20212-v1-2008Mar