

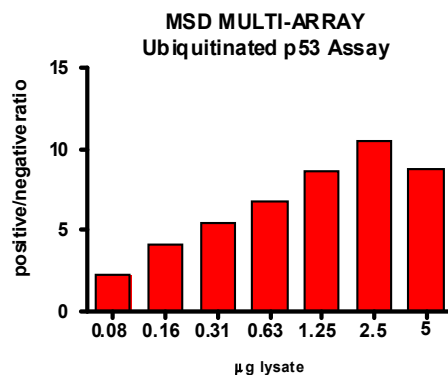
Meso Scale Discovery® Whole Cell Lysate Set

Ubiquitinated p53

Catalog No:	C11FK-1
Contents:	2 x 100 µg MSDLY0058 Ubiquitinated p53 Negative Cell Lysate Cell lysate from growing HCT116 cells 2 x 100 µg MSDLY0059 Ubiquitinated p53 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 p53
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. Ubiquitinated p53 lysates will lose more than 50% of activity after a single round of freeze/thaw.
Quality Control:	Lysates have been tested for performance in Western Blot and MSD MULTI-ARRAY® Assays.

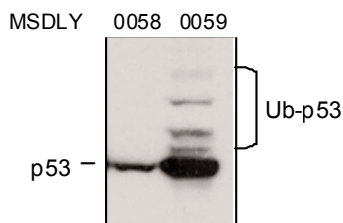
MSD MULTI-ARRAY Assay Results

The figure below illustrates typical lysate titrations for MSDLY0058 (ubiquitinated p53 negative) and MSDLY0059 (ubiquitinated p53 positive) cell lysates using the MSD MULTI-ARRAY Ubiquitinated p53 Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with ubiquitinated p53 positive and ubiquitinated p53 negative lysates. The ubiquitinated p53 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 p53 antibodies.



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

20213-v1-2008Mar