

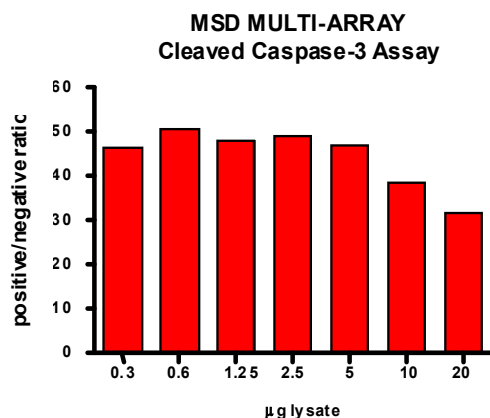
# Meso Scale Discovery® Whole Cell Lysate Set

## Cleaved Caspase-3

|                         |   |
|-------------------------|---|
| <b>Catalog No:</b>      | C11CF-1   |
| <b>Contents:</b>        | 2 x 100 µg MSDLY0001<br>Cleaved Caspase-3 Negative Cell Lysate<br>Cell lysate from growing Jurkat cells<br><br>2 x 100 µg MSDLY0002<br>Cleaved Caspase-3 Positive Cell Lysate<br>Cell lysate from growing Jurkat cells treated with 1 µM staurosporine for 4 hours to induce Caspase-3 cleavage |
| <b>Concentration:</b>   | 2 mg/mL in MSD® Complete Tris Lysis Buffer  |
| <b>Volume:</b>          | 2 vials (50 µL) negative lysate    2 vials (50 µL) positive lysate  |
| <b>Preparation:</b>     | Following cell treatment, Jurkat cell lysates were prepared on ice in MSD Complete Tris Lysis Buffer. Cell debris was cleared by centrifugation.  |
| <b>Storage:</b>         | 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).  |
| <b>Quality Control:</b> | Lysates have been tested for performance in Western Blot and MSD MULTI-ARRAY® Assays.   |

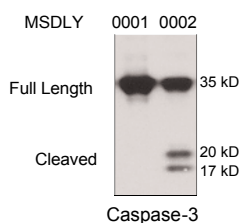
### MSD MULTI-ARRAY Assay Results

The figure below illustrates a typical titration for MSDLY0001 (cleaved Caspase-3 negative) and MSDLY0002 (cleaved Caspase-3 positive) cell lysates with the MSD MULTI-ARRAY Caspase-3 (Cleaved) Whole Cell Lysate Kit. Results are presented as a ratio of the signals obtained with cleaved Caspase-3 positive and cleaved Caspase-3 negative lysates. In the cleaved Caspase-3 singleplex assay, the cleaved Caspase-3 signal ratio remains constant with increasing amounts 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

MSDLY0001 and MSDLY0002 whole cell lysates (20 µg each) were analyzed by Western Blot with an antibody recognizing both full length and cleaved Caspase-3.



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

20046-v1-2008Mar