

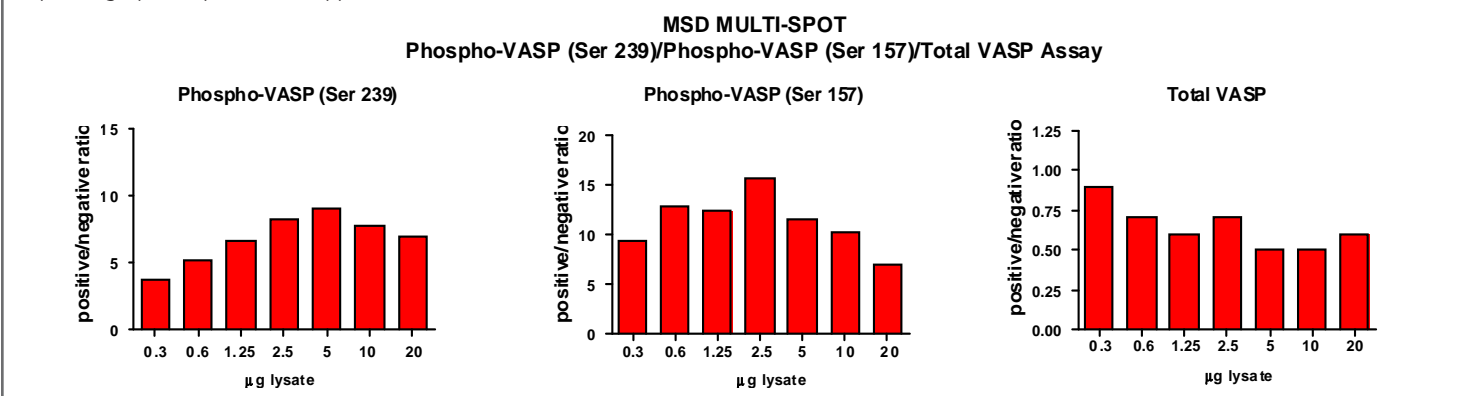
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-VASP (Ser 157/239)

Catalog No:	C11FG-1
Contents:	2 x 100 µg MSDLY0053 pVASP Negative Control Cell Lysate Cell lysate from serum deprived A431 cells 2 x 100 µg MSDLY0054 pVASP Positive Control Cell Lysate Cell lysate from serum deprived A431 cells treated with 100 µM forskolin and 100 nM calyculin A for approximately 1 hour to stimulate VASP 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, A431 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.

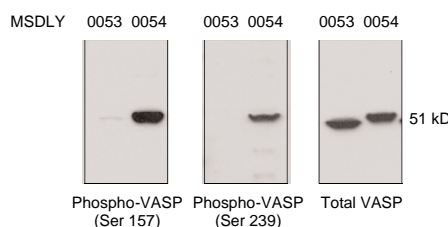
MSD MULTI-SPOT Assay Results

The figure below illustrates typical titrations for MSDLY0053 (pVASP negative) and MSDLY0054 (pVASP positive) cell lysates with the MSD MULTI-SPOT Phospho-VASP (Ser 239)/Phospho-VASP (Ser 157)/Total VASP Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with pVASP positive and pVASP negative lysates. The Phospho-VASP (Ser 239), Phospho-VASP (Ser 157), and Total VASP signal ratios remain close to 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

MSDLY0053 and MSDLY0054 whole cell lysates (20 µg each) were analyzed by Western Blot with phospho serine 239, phospho serine 157 and total VASP antibodies.



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

20052-v1-2010Feb