

Research use only. Not for use in diagnostic procedures.

AlphaLISA® SureFire® Ultra™

## Human and Mouse SQSTM1 p62 Total Detection Kit

Product number: ALSU-TSQSTM-A500, ALSU-TSQSTM-A10K,

ALSU-TSQSTM-A50K, ALSU-TSQSTM-A-HV



## Kit specificity:

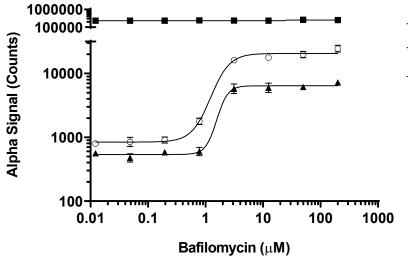
This assay kit contains antibodies which recognize distinct epitopes on SQSTM1 p62. The protein detected by this kit corresponds to UniProt ID Q13501. SQSTM1 p62 is also known as sequestosome 1 and EBIAP. These antibodies recognize SQSTM1 p62 of human and mouse origin. Other species should be tested on a case-by-case basis.

## **Control lysate information:**

Positive Control Lysate: Prepared from HeLa cells cultured to confluence in T175 flasks in 10% FBS containing medium, treated with 1  $\mu$ M Okadaic acid for 2 hours and lysed with 5 mL of Lysis Buffer.

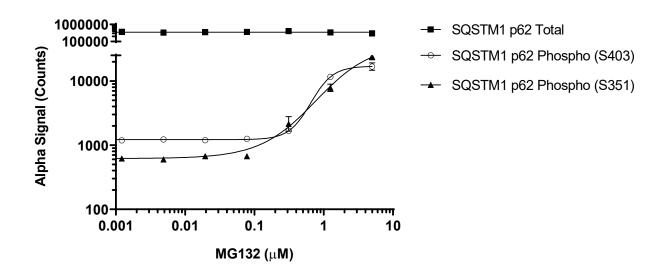
## Representative data

Data obtained with a 2-plate, 2-incubation protocol. A549 cells were seeded at 20K cells/well in a 96 well plate and incubated overnight. Cells were treated with Bafilomycin at the indicated concentrations for 24 hours. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser351 and Ser403) and Total SQSTM1 p62 using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



- SQSTM1 p62 Total
- ★ SQSTM1 p62 Phospho (Ser351)
- → SQSTM1 p62 Phospho (Ser403)

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with MG132 at the indicated concentrations for 16 hours. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser351 and Ser403) and Total SQSTM1 p62 using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF"



