

# AlphaLISA® SureFire® Ultra™

## Human p-SHP-1 (Tyr536) Detection Kit

**Product number:** 

ALSU-PSHP1-A500, ALSU-PSHP1-A10K,

ALSU-PSHP1-A50K, ALSU-PSHP1-A-HV



### Kit specificity:

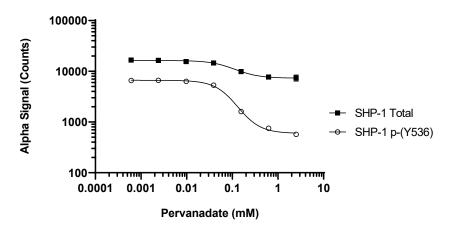
This assay kit contains antibodies which recognize the phospho-Tyr536 epitope and a distal epitope on SHP-1. The protein detected by this kit corresponds to UniProt ID P29350. SHP-1 is also known as Tyrosine-protein phosphatase non-receptor type 6 and Hematopoietic cell protein-tyrosine phosphatase. These antibodies recognize SHP-1 of human origin. Other species should be tested on a case-by-case basis.

#### **Control lysate information:**

Positive Control Lysate: Prepared from Jurkat cells, cultured in 10% FBS containing media. Cells were harvested and washed in HBSS + 0.1% BSA, spun down at 300 RCF for 5 minutes and lysed with Lysis Buffer at 3.2 x 10<sup>6</sup> cells/mL.

#### **Representative data**

Data obtained with a 2-plate, 2-incubation protocol. Jurkat cells were washed and resuspended in HBSS at 2 x 10<sup>6</sup> cells/mL. Cells were treated with Pervanadate at the indicated concentrations for 30 minutes. Cells were lysed with Lysis Buffer at 0.8 x 10<sup>6</sup> cells/mL and assayed separately for Phospho (Tyr536) and Total SHP-1 using respective *SureFire Ultra* kits. Equivalent to approximately 8,000 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"





Revvity, Inc. 940 Winter Street Waltham, MA 02451 USA (800) 762-4000 www.revvity.com For a complete listing of our global offices, visit <u>www.revvity.com</u> Copyright ©2023, Revvity, Inc. All rights reserved.

VersionTGRSU228.2