Research use only. Not for use in diagnostic procedures.

AlphaLISA® SureFire® Ultra™ Human p-VAV1 (Tyr174) Detection Kit

Product number:

ALSU-PVAV1-A500, ALSU-PVAV1-A10K,

ALSU-PVAV1-A50K, ALSU-PVAV1-A-HV



Kit specificity:

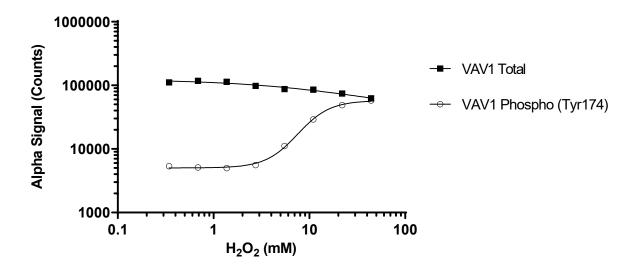
This assay kit contains antibodies which recognize the phospho-Tyr174 epitope and a distal epitope on VAV1. The protein detected by this kit corresponds to UniProt ID P15498. These antibodies recognize VAV1 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

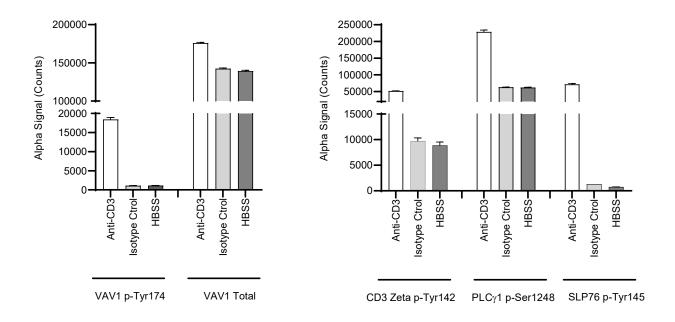
Positive Control Lysate: Prepared from THP1 cells, seeded at 200 K cells/mL and grown for two days in T175 flasks in 10% FBS containing medium. Cells were harvested, washed in HBSS + 0.1% BSA and adjusted to 2x10⁶ cells/mL. Cells were treated with 2 mM pervanadate for 10 minutes and lysed with the addition of 5X Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Jurkat cells were plated at 200K cells/well in a 96 well plate in HBSS + 0.1% BSA, then treated with H_2O_2 at the indicated concentrations for 2 minutes. Cells were lysed with 5X Lysis Buffer and assayed for Phospho (Tyr174) and Total VAV1 using respective *SureFire Ultra* kits. Equivalent to approximately 8,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. Jurkat cells were seeded at 5x10⁶ cells/well in a 12 well plate in HBSS + 0.1% BSA. Cells were incubated for 1 hour at 37°C then treated with 2.5 µg/mL of anti-CD3 Ab (BD cat# 555329), a mouse IgG Isotype control or HBSS for 2 minutes. Cells were lysed with 5X Lysis Buffer and assayed for Phospho (Tyr174) and Total VAV1 as well as other related targets 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"

ſe√

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.