

Research use only. Not for use in diagnostic procedures.

AlphaLISA® SureFire® Ultra™ Human p-RIP3 (Ser227) Detection Kit

Product number:

ALSU-PRIP3-A500, ALSU-PRIP3-A10K,

ALSU-PRIP3-A50K, ALSU-PRIP3-A-HV



Kit specificity:

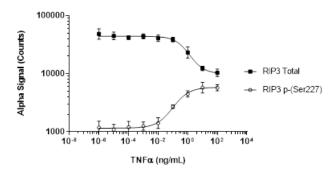
This assay kit contains antibodies which recognize the phospho-Ser227 epitope and a distal epitope on RIP3. The protein detected by this kit corresponds to UniProt ID Q9Y572. RIP3 is also known as RIP-like protein kinase 3 and Receptor-interacting protein 3. These antibodies recognize RIP3 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from U-937 cells cultured in 10% FBS containing medium. Cells were harvested and adjusted to 0.8 x 10⁶ cells/mL then treated with 20 μ M Z-VAD, 20 ng/mL rhTNF α and 200 nm SM-164 for 5 hours and 30 minutes. Cells washed with HBSS + 0.1% BSA, spun down at 300 RCF for 5 minutes and lysed in Lysis Buffer at 13.3 x 10⁶ cells/mL.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. U-937 cells were seeded at 400K cells/well in a 96 well plate then treated with TNF α at the indicated concentrations with 20 μ M Z-VAD and 200 nm SM-164 for 7 hours. Cells washed with HBSS, lysed with Lysis Buffer and assayed separately for Phospho (Ser227) and Total RIP3 using respective *SureFire Ultra* kits. Equivalent to approximately 20,000 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.

