

AlphaLISA® SureFire® Ultra™

Human p-RIPK1 (Ser320) Detection Kit

Product number: ALSU-PRIPK1-B500, ALSU-PRIPK1-B10K,
ALSU-PRIPK1-B50K, ALSU-PRIPK1-B-HV



Kit specificity:

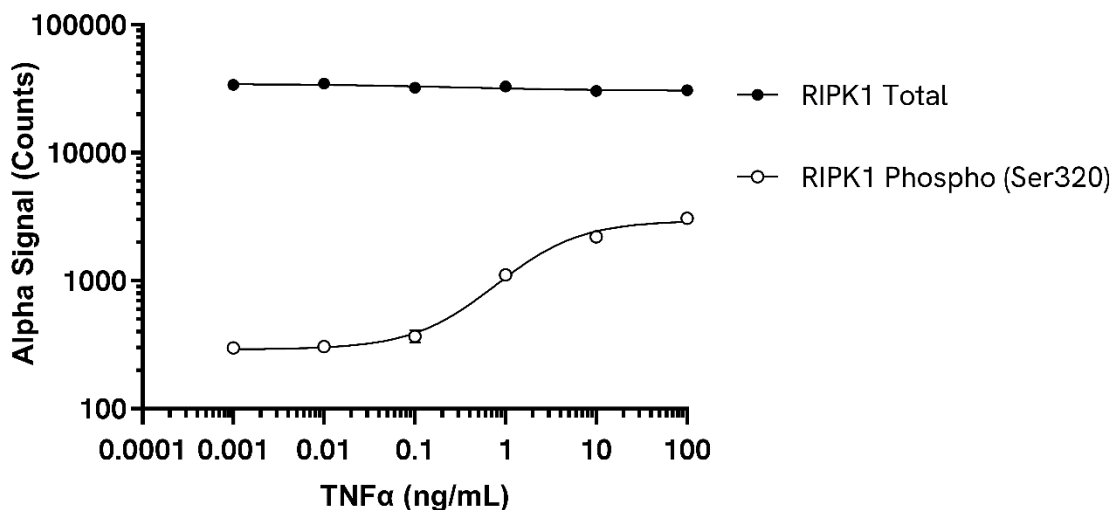
This assay kit contains antibodies which recognize phospho-Ser320 epitope and distal epitope on RIPK1. The protein detected by this kit corresponds to UniProt ID Q13546. RIPK1 is also known as Receptor-interacting serine/threonine-protein kinase 1 and Cell death protein RIP. These antibodies recognize RIPK1 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

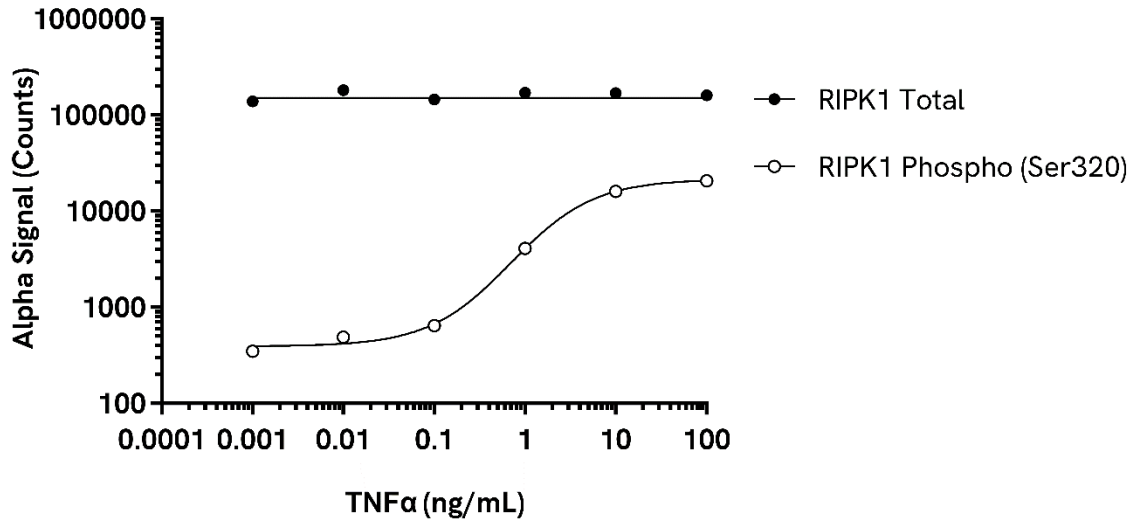
Positive Control Lysate: Prepared from RT4 cells, cultured to confluence in 10% FBS containing medium, then treated with 100 ng/mL of TNF α for 15 minutes and lysed with 5 mL of Lysis Buffer.

Representative data:

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with TNF α at the indicated concentrations for 15 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser320) and Total RIPK1 using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. RT4 cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with TNF α at the indicated concentrations for 15 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser320) and Total RIPK1 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"

