

AlphaLISA® SureFire® Ultra™

Human p-IKKβ (Ser177/181) Detection Kit

Product number: ALSU-PIKKB-A500, ALSU PIKKB-A10K,
ALSU-PIKKB-A50K, ALSU-PIKKB-A-HV



Kit specificity:

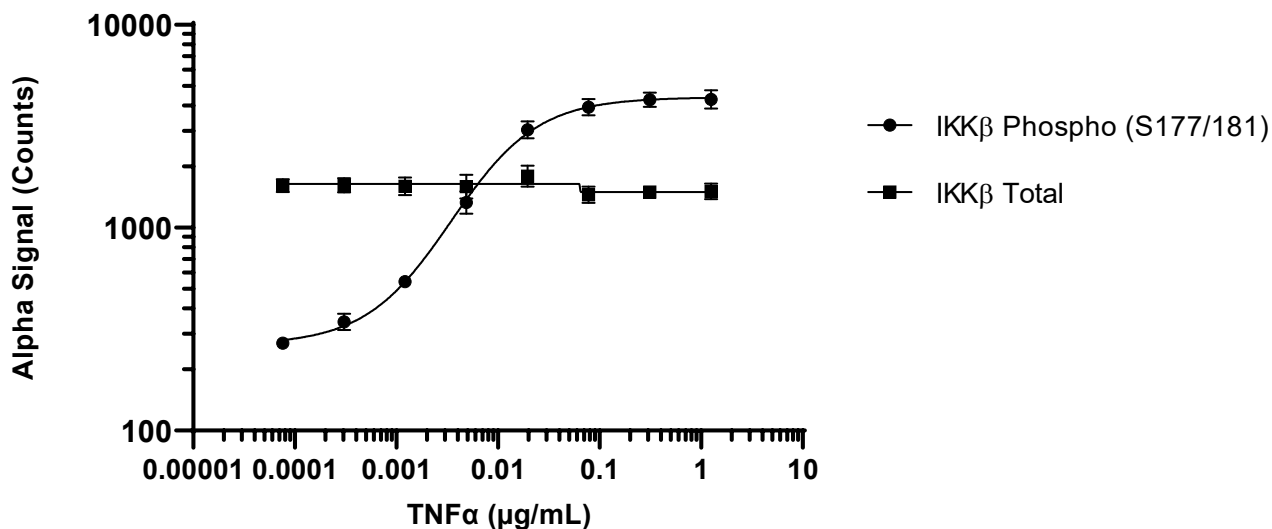
This assay kit contains antibodies which recognize the Ser177/181 epitope and a distal epitope on IKKβ. The protein detected by this kit corresponds to UniProt ID O14920. IKKβ is also known as Inhibitor of nuclear factor kappa-B kinase subunit beta. These antibodies recognize IKKβ of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

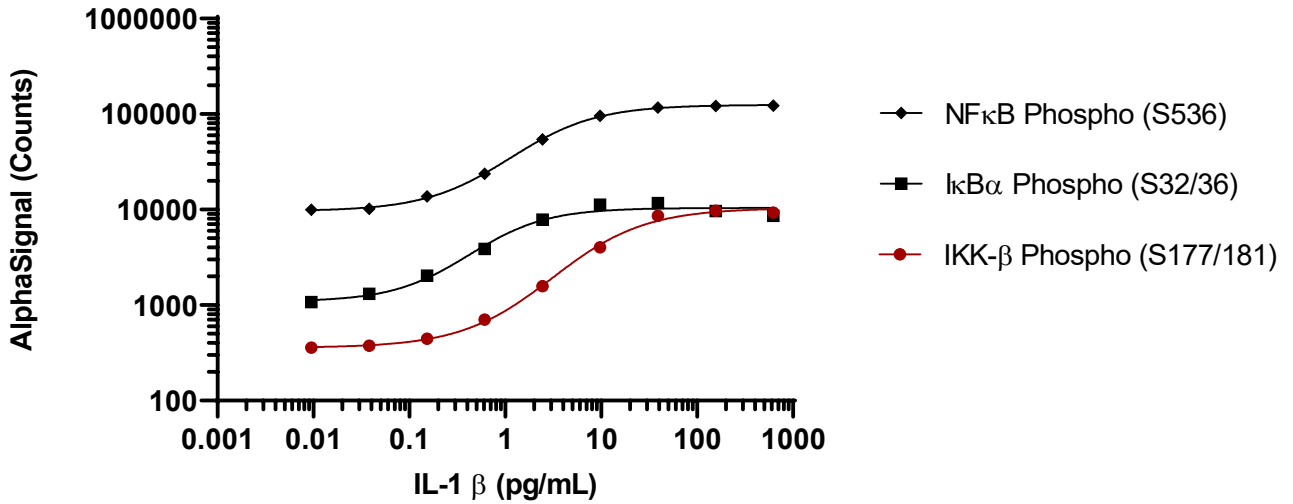
Positive Control Lysate: Prepared from KARPAS 299 cells cultured in 10% FBS containing media. Cells were harvested and washed with HBSS + 0.1% BSA, adjusted to 10 x10⁶ cells/mL and then treated with 2 ng/mL rhIL1-B for 5 minutes. Cells were spun down at 300 RCF for 5 minutes and lysed in Lysis Buffer at 8 x 10⁶ cells/mL.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. A549 cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with TNFα at the indicated concentrations for 10 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser177/181) and Total IKKβ, using respective SureFire Ultra kits. Equivalent to approximately 4,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. Karpas 299 cells were washed with HBSS + 0.1% BSA and seeded at 300K cells per well in a 96 well plate. Cells were treated with IL-1 β at the indicated concentrations for 10 minutes. Cells were lysed with 5x Lysis Buffer and assayed for various targets using respective SureFire Ultra kits. Equivalent to approximately 12,000 cells/datapoint (2,400 cells for NF κ B).



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"

