

Research use only. Not for use in diagnostic procedures.

AlphaLISA® SureFire® Ultra™

Human IKKα Total Detection Kit

Product number: ALSU-TIKKA-A500, ALSU-TIKKA-A10K,

ALSU-TIKKA-A50K, ALSU-TIKKA-A-HV



Kit specificity:

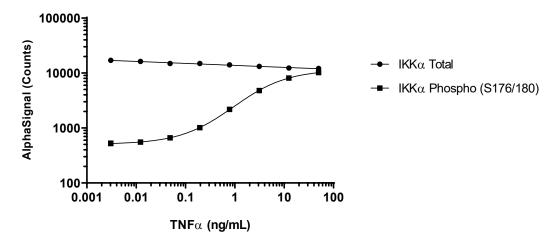
This assay kit contains antibodies which recognize distinct epitopes on IKK α . The protein detected by this kit corresponds to UniProt ID O15111. IKK α is also known as Inhibitor of nuclear factor kappa-B kinase subunit alpha. These antibodies recognize IKK α of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control lysate: Prepared from recombinant p-IKK α protein (Abcam, ab102103) at a concentration of 250 ng/mL.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells seeded at 20K cells/well in 96 well plate and incubated for 2 days. Cells were treated with 40 nM Calyculin A and TNF α at the indicated concentrations for 10 minutes. Cells were lysed with Lysis buffer and assayed for Phospho (Ser176/180) and Total IKK α 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"



