

AlphaLISA® SureFire® Ultra™

Human p-IRAK4 (Thr345/Ser346) Detection Kit

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



Kit specificity:

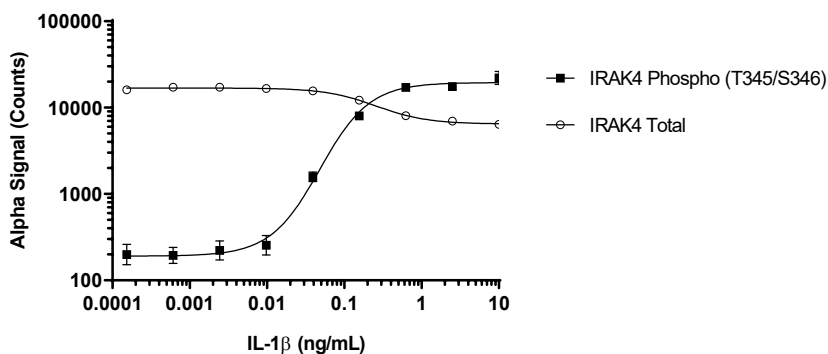
This assay kit contains antibodies which recognize the phospho-Thr345/Ser346 epitope and a distal epitope on IRAK4. The protein detected by this kit corresponds to UniProt ID Q9NWZ3. IRAK4 is also known as Interleukin-1 receptor associated kinase 4. These antibodies recognize IRAK4 of human origin. Other 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 2×10^6 cells/mL and then treated with 10 ng/mL IL-1 β for 10 minutes. Cells were spun down at 300 RCF for 5 minutes and lysed in Lysis Buffer at 1×10^6 cells/mL.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Karpas299 cells were seeded at 200K cells/well in 10% FBS containing medium and grown for 24 hours. Cells were harvested, washed in HBSS + 0.1% BSA and plated at 200K cells/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 separately for Phospho (Thr345/Ser346) and Total IRAK4 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"