

AlphaLISA® SureFire® Ultra™

Human p-ALK (Tyr1604) Detection Kit

Product number: ALSU-PALK-C500, ALSU PALK-C10K,
ALSU-PALK-C50K, ALSU-PALK-C-HV



Kit specificity:

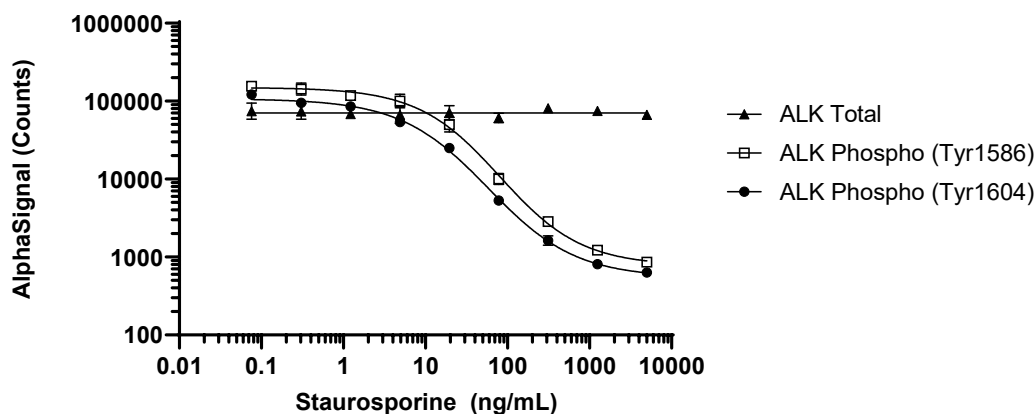
This assay kit contains antibodies which recognize the phospho-Tyr1604 epitope and a distal epitope on ALK. The protein detected by this kit corresponds to UniProt ID Q9UM73. ALK is also known as Anaplastic lymphoma kinase. These antibodies recognize ALK of human origin. Other species should be tested 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 washed in HBSS + 0.1% BSA, spun down at 300 RCF for 5 minutes and then lysed with Lysis Buffer at 1.0×10^6 cells/mL.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. KARPAS 299 cells were harvested, washed in HBSS + 0.1% BSA and seeded at 25K cells/mL in a 96 well plate. Cells were treated with Staurosporine at the indicated concentrations for 2 hours. Cells were lysed with 5x Lysis Buffer and assayed separately for Phospho and Total ALK using respective SureFire Ultra kits. Equivalent to approximately 1,000 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"