

Alpha SureFire® Ultra™ Multiplex

## Human p-ALK (Tyr1604) + Total ALK Detection Kit

**Product number:** MPSU-PTALK-K500, MPSU PTALK-K10K,  
MPSU-PTALK-K50K, MPSU-PTALK-K-HV



### Kit specificity:

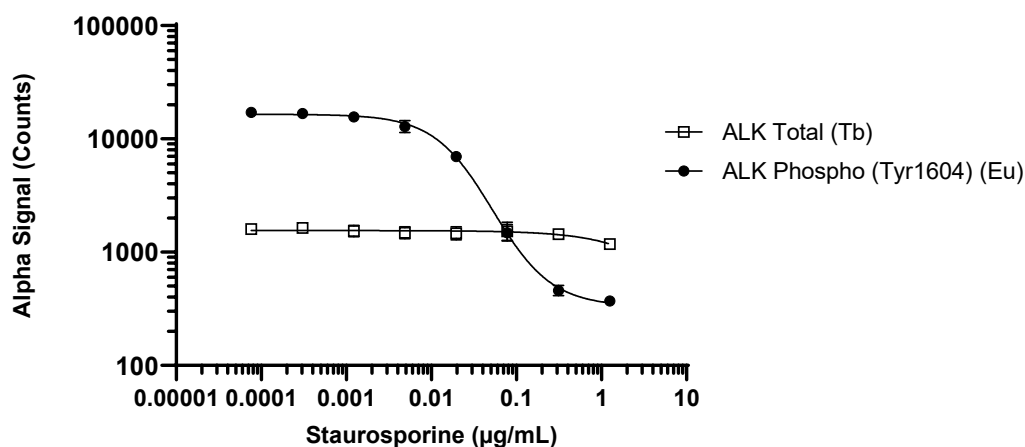
This assay kit contains antibodies which recognize the phospho-Tyr1604 epitope and a distal epitope on ALK plus antibodies to detect Total ALK. The protein detected by this kit corresponds to UniProt ID Q9UM73. These antibodies recognize ALK of human origin. Other species should be tested on a case-by-case basis.

### Control lysate information:

Positive Control Lysate: Prepared from KARPAS 299 cells cultured in 10% FBS containing medium, washed in HBSS + 0.1% BSA, spun down at 300 RCF for 5 minutes and lysed with Lysis Buffer at  $1.0 \times 10^6$  cells/mL.

### Representative data

Data obtained with a 2-plate, 2-incubation protocol. KARPAS 299 cells were seeded at 200K cells/well in a 96 well plate in HBSS + 0.1% BSA, then treated with Staurosporine at the indicated concentrations for 2 hours. Cells were lysed with 5X Lysis Buffer and assayed for Phospho (Tyr1604) and Total ALK in the same wells of an assay plate using the SureFire Ultra Multiplex kit. Equivalent to approximately 8,000 cells/data point.



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"