

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ ALK Total Assay Kit

**Products: ALSU-TALK-A500, ALSU-TALK-A10K, ALSU-TALK-A50K,
ALSU-TALK-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize two distinct invariant epitopes on ALK. The protein detected by this kit corresponds to GenBank Accession NP_004295. ALK is also known as CD246; NBLST3; TFG/ALK. These antibodies recognize ALK of human origin. Other species should be tested on a case-by-case basis.

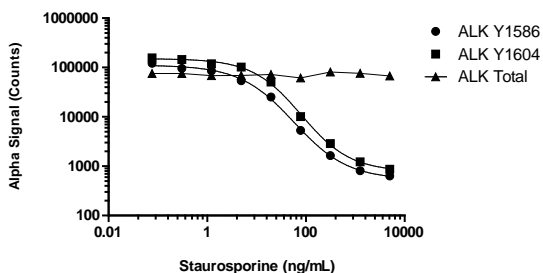
Control Lysate Information

Positive Control Lysate: Prepared from Karpas-299 cells cultured in T175 flasks to 10^6 cells/mL in medium containing 10% FBS, and lysed with 5X *SureFire Ultra* Lysis buffer at a final cell concentration of 0.3×10^6 cells/mL.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Karpas-299 cells were plated at 25K cells per well in a 96 well plate in HBSS + 0.1% BSA, then treated with staurosporine at the indicated concentrations for 2 hours at 37°C. Cells were then lysed with 5X Lysis buffer, and assayed for phospho (Tyr1586 or Tyr1604) or total ALK in separate wells of an assay plate with respective *SureFire Ultra* kits.

**Staurosporine inhibition of ALK phosphorylation in Karpas-299 cells
SureFire Ultra ALK assays**



For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Assay Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>