

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ PI3 Kinase p85 Total Assay Kit

ALSU-TP13K-A500, ALSU-TP13K -A10K, ALSU-TP13K-A50K, ALSU-TP13K-A-HV

Kit-Specificity Information

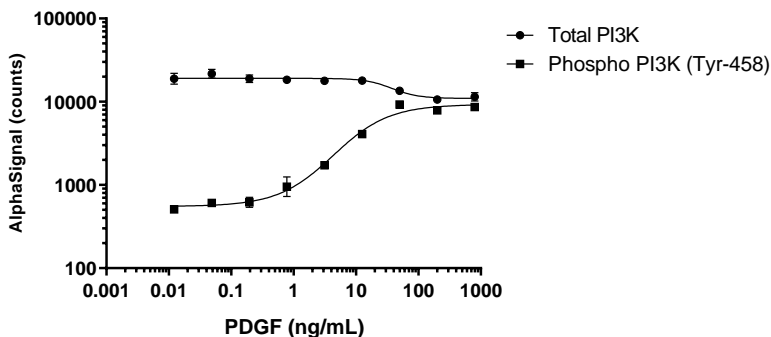
This assay kit contains antibodies which recognize invariant epitopes on PI3 Kinase p85. The protein detected by this kit corresponds to UniProt ID P27986. PI3 Kinase p85 is also known as Phosphoinositide 3-kinase. These antibodies recognize PI3 Kinase p85 of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from NIH/3T3 cells, cultured to 80-90% confluence in T175 flasks in 10% FBS containing medium, medium removed and then treated with 50 ng/mL PDGF-BB for 5 minutes and lysed in 5 mL of 1X *SureFire Ultra* Lysis buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. NIH3T3 cells seeded at 40k/well DMEM + 10% FBS in a 96 well plate and incubated overnight. Cells were treated for 10 mins with varying concentrations of PDGF in DMEM + 10% FCS. Cells were lysed in *SureFire Ultra* Lysis Buffer and assayed for Phospho (Tyr-458) and Total PI3 Kinase p85 using respective *SureFire Ultra* assay kits. Equivalent to approximately 4,000 cells/datapoint.



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>.