

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-PI3 Kinase p85 (Tyr458) (Mouse Specific) Assay Kit

ALSU-PP13K-A500, ALSU-PP13-A10K, ALSU-PP13K-A50K, ALSU-PP13K-A-HV

Kit-Specificity Information

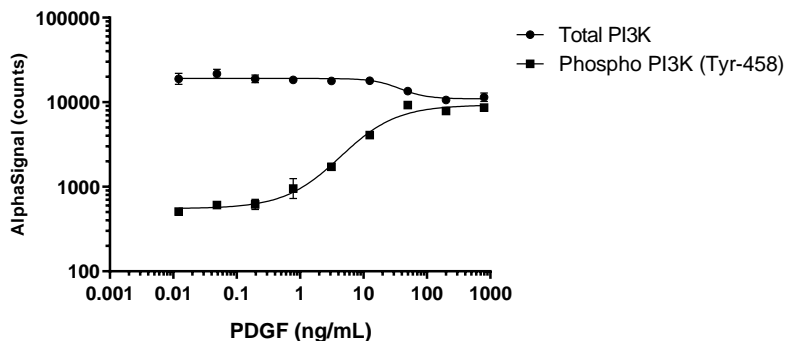
This assay kit contains antibodies which recognize the phospho-Tyr458 epitope, and a distal epitope, on PI3 Kinase p85. The protein detected by this kit corresponds to UniProt ID P27986. PI3 Kinase p85 is also known as Phosphoinositide 3-kinase. These antibodies recognize PI3 Kinase p85 of mouse origin. 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>.