

AlphaLISA® SureFire® Ultra™

# Human and Mouse PI3 Kinase p85 Total Detection Kit

**Product number:** ALSU-TP13K-A500, ALSU-TP13K-A10K,  
 ALSU-TP13K-A50K, ALSU-TP13K-A-HV



### Kit specificity:

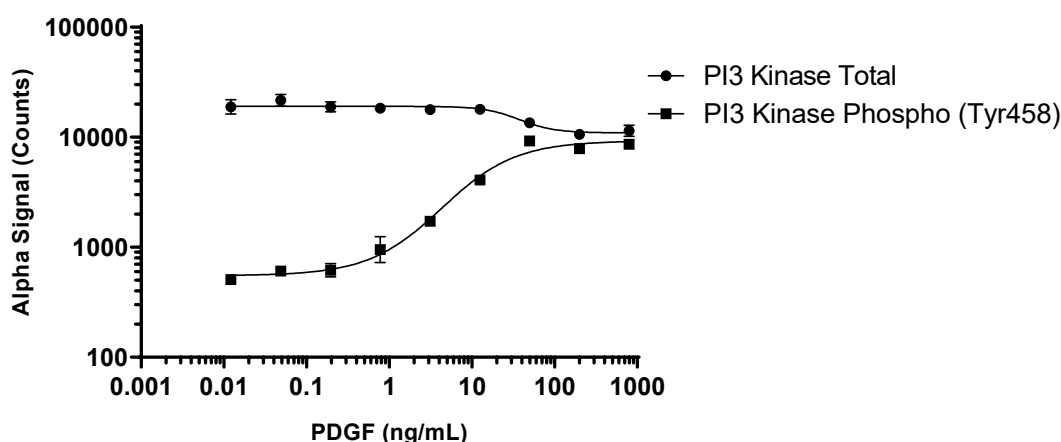
This assay kit contains antibodies which recognize distinct epitopes on PI3 Kinase p85. The protein detected by this kit corresponds to UniProt ID P27986. PI3 Kinase p85 is also known as Phosphoinositide 3-kinase. These antibodies recognize PI3 Kinase p85 of human and mouse origin. Other species should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

### Control lysate information:

Positive Control Lysate: Prepared from NIH/3T3 cells, cultured to confluence in T175 flasks in 10% FBS containing media, then treated with 50 ng/mL PDGF-BB for 10 minutes and lysed with 5 mL of Lysis Buffer.

### Representative data

Data obtained with a 2-plate, 2-incubation protocol. NIH/3T3 cells were seeded at 40K/well in a 96 well plate and incubated overnight. Cells were treated with PDGF-BB at the indicated concentrations for 10 minutes. Cells were lysed in 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.



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"