

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-PDGF Receptor β (Tyr751) Assay Kit

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

Kit-Specificity

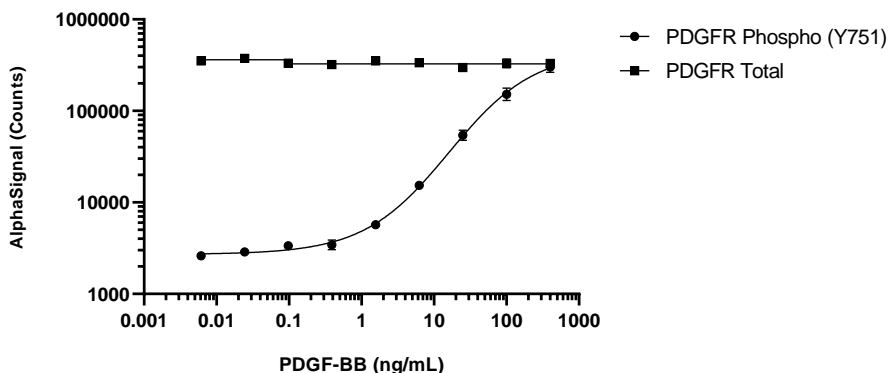
This assay kit contains antibodies which recognize the phospho-Tyr751 epitope, and a distal epitope, on PDGF Receptor β . The protein detected by this kit corresponds to UniProt ID P09619. These antibodies recognize PDGF Receptor β of human, rat and mouse origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from NIH/3T3 cells, cultured to 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. NIH/3T3 cells were seeded at 40K cells per well in a 96 well plate and incubated overnight. Cells were then treated for 5 mins with PDGF-BB at the indicated concentrations. Cells were lysed with 1X Lysis Buffer and assayed for Total or Phospho PDGFR β (Tyr751), using respective *SureFire Ultra* assay kits. Equivalent to approximately 4000 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>