

Technical Data Sheet

Caution: Research use only.

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

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

Kit-Specificity Information

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 Entrez-Gene ID 5159. These antibodies recognize PDGF Receptor β of human, rat and mouse origin. Other species should be tested on a case-by-case basis.

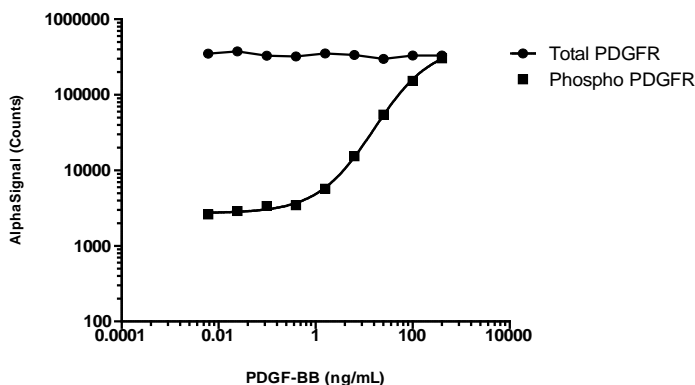
Control Lysate Information

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 50ng/mL PDGF-BB for 5min and lysed in 5mL of 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. NIH/3T3 cells were seeded at 2×10^4 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.

PDGF-BB Dose Response - NIH/3T3 cells



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>

TGRSU284.0