

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex p-PDGF Receptor $\beta$ (Tyr751) + Total PDGF Receptor $\beta$ Assay Kit

**Products: MPSU-PTPDGF-K500, MPSU-PTPDGF-K10K, MPSU-PTPDGF-K50K,  
MPSU-PTPDGF-K-HV**

### Kit-Specificity Information

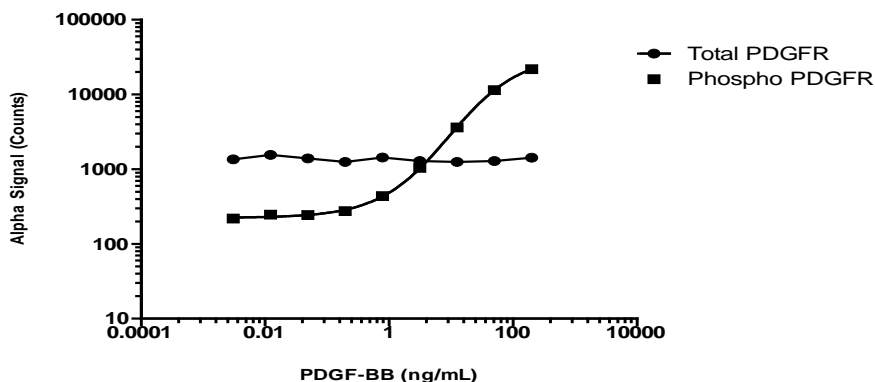
This assay kit contains antibodies which recognize the phospho-Tyr751 epitope, and a distal epitope, on PDGF Receptor  $\beta$  for phospho PDGF Receptor  $\beta$  analysis, plus antibodies to measure the total levels of PDGF Receptor  $\beta$ . The protein detected by this kit corresponds to Entrez Gene ID 5159. These antibodies recognize PDGF Receptor  $\beta$  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  $2 \times 10^4$  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 100 $\mu$ L of 1X Lysis buffer and samples were assayed for both total and phospho (Tyr751) PDGFR  $\beta$ , as detailed in the MPSU manual.



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>