

Alpha SureFire® Ultra™ Multiplex

Human and Mouse p-PDGF Receptor β (Tyr751) + Total PDGF Receptor β Detection Kit

Product number: MPSU-PTPDGF-K500, MPSU-PTPDGF-K10K,
MPSU-PTPDGF-K50K, MPSU-PTPDGF-K-HV



Kit specificity:

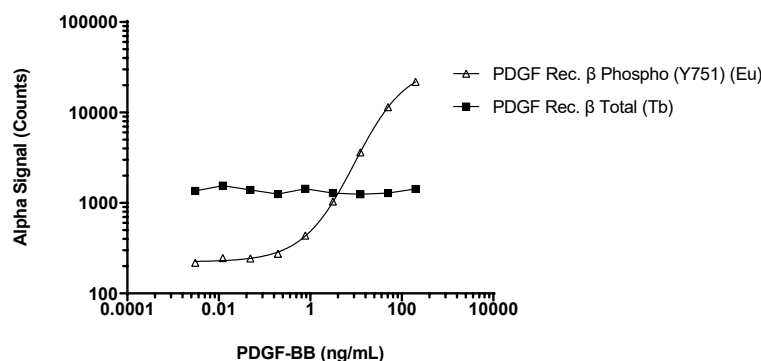
This assay kit contains antibodies which recognize the phospho-Tyr751 epitope and a distal epitope on PDGFR β , plus antibodies to detect Total Platelet-derived growth factor (PDGF) Receptor β . The protein detected by this kit corresponds to UniProt ID P09619. PDGF Receptor β is also known as CD140 antigen-like family member B. These antibodies recognize PDGF Receptor β 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 medium, treated with 50 ng/mL PDGF-BB for 5 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 20K cells/well in a 96 well plate and incubated overnight. Cells were treated with PDGF-BB at the indicated concentrations for 5 minutes. Cells were lysed with Lysis Buffer and assayed for Phospho (Tyr751) and Total PDGF Receptor β in the same wells of an assay plate using the *SureFire Ultra* Multiplex kit. Equivalent to approximately 1,000 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"