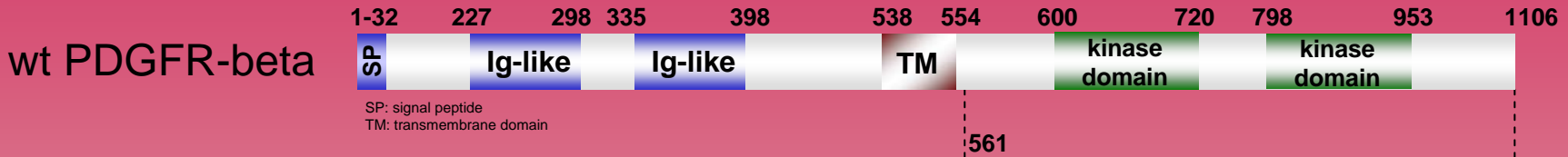


human PDGFR-beta kinase (GenBank Acc.No: NM_002609)

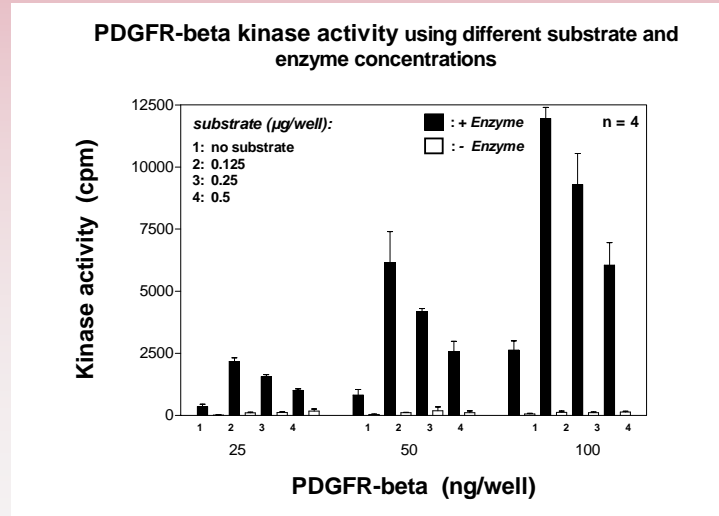
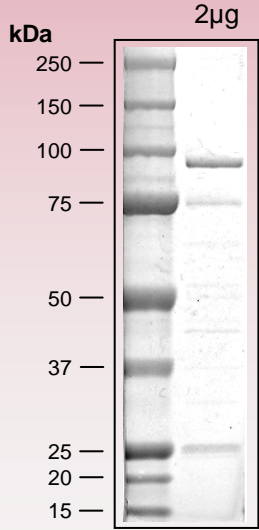


ProQinase PDGFR-beta (Arg₅₆₁-Leu₁₁₀₆)

777 aminoacids, 88.029 kDa
 1-218 GST
 222-226 factor Xa cleavage site
 232-777 PDGFR-beta 561-1106



Representative lot: Coomassie stain & activity in PanQinase™ assay



Lot specific datasheet available on request.
 please contact:
info@proqinase.com