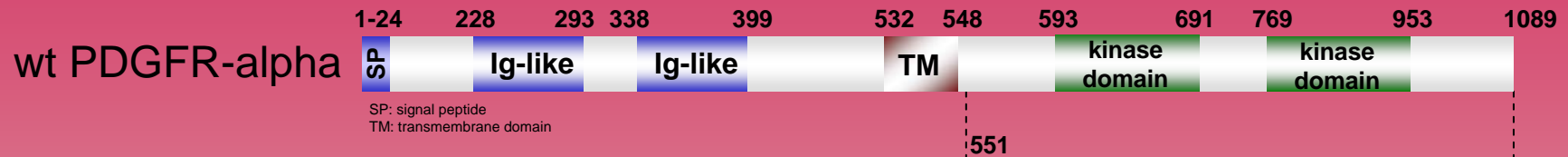


human PDGFR-alpha kinase (GenBank Acc.No: NM_006206)

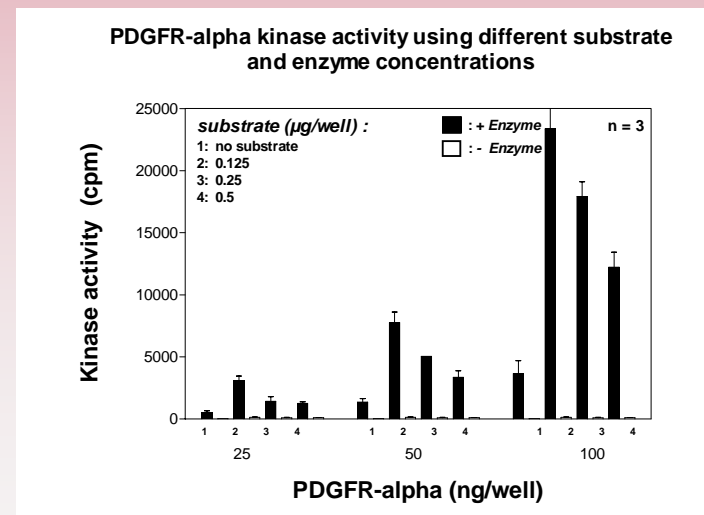
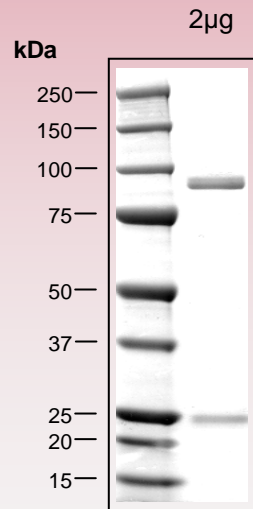


ProQinase PDGFR-alpha (Gln₅₅₁-Leu₁₀₈₉)

766 aminoacids, 87.851 kDa
 1-218 GST
 222-226 factor Xa cleavage site
 228-766 PDGFR-alpha 551-1089



Representative lot: Coomassie stain & activity in PanQinase™ assay



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