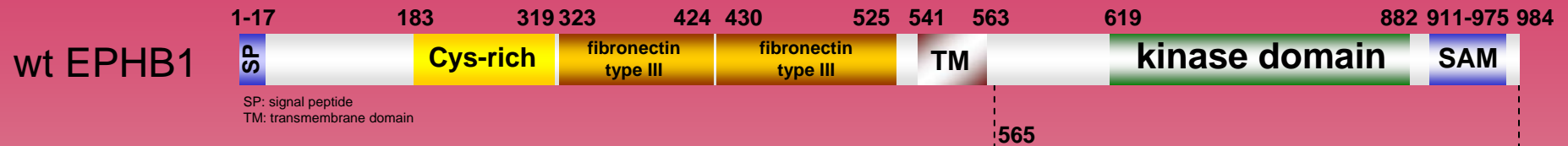


human EPHB1 kinase (GenBank Acc.No: NM_004441.2)

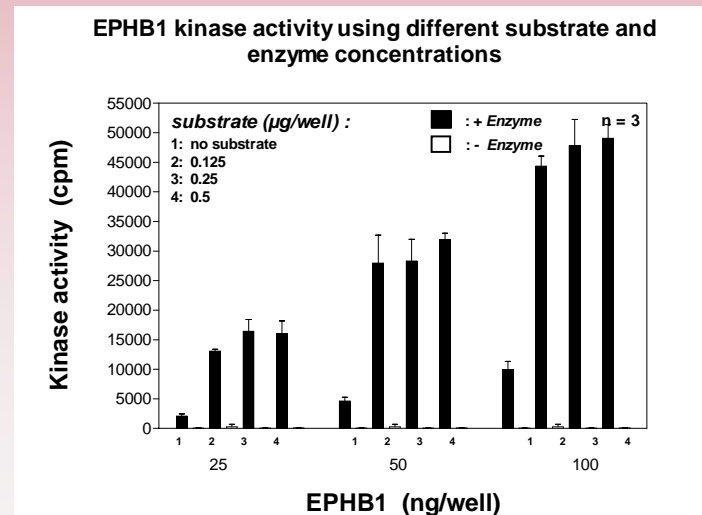
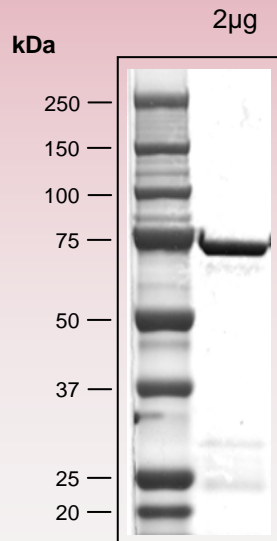


ProQinase EPHB1 (Arg₅₆₅-Ala₉₈₄)

675 aminoacids, 76.906 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 256-675 EPHB1 565-984



Representative lot: Coomassie stain & activity in PanQinase™ assay



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