

human EPHB4 kinase (GenBank Acc.No: NM_004444.2)



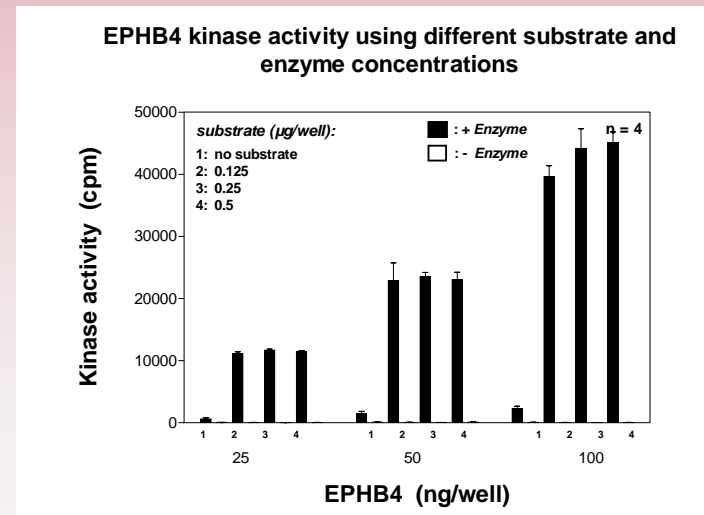
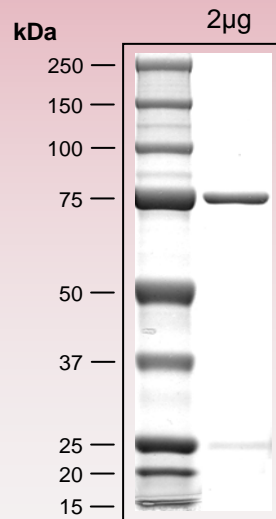
SP: signal peptide
TM: transmembrane domain

ProQinase EPHB4 (Leu₅₆₃-Tyr₉₈₇)

680 aminoacids, 76.794 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 256-680 EPHB4 563-987



Representative lot: Coomassie stain & activity in PanQinase™ assay



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