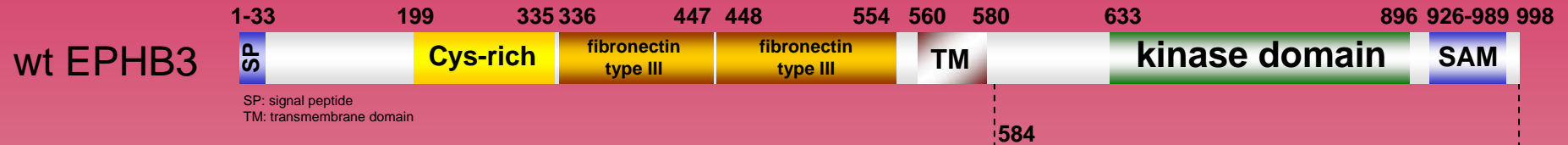


human **EPHB3** kinase (GenBank Acc.No: NM_004443)

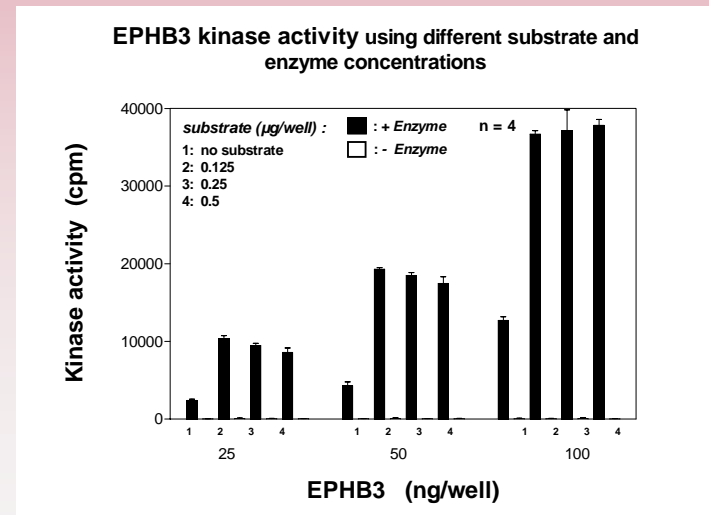
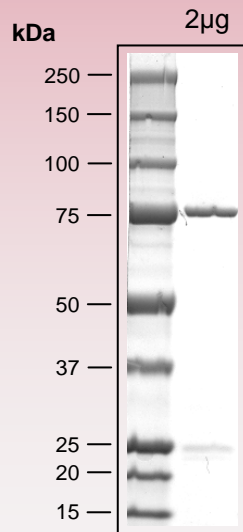


ProQinase EPHB3 (Gln₅₈₄-Val₉₉₈)

704 aminoacids, 80.148 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 280-704 EPHB2 584-998



Representative lot: Coomassie stain & activity in PanQinase™ assay



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