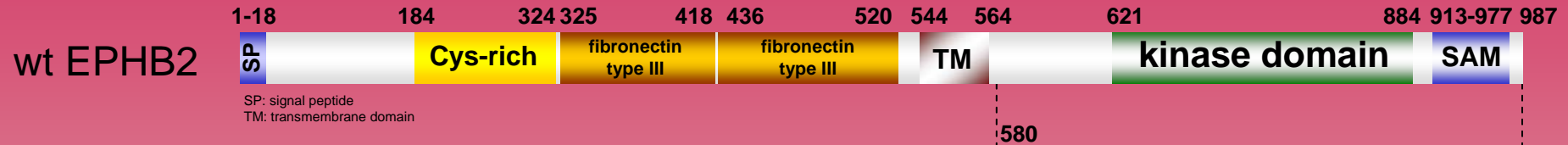


# human EPHB2 kinase (GenBank Acc.No: NM\_004442)

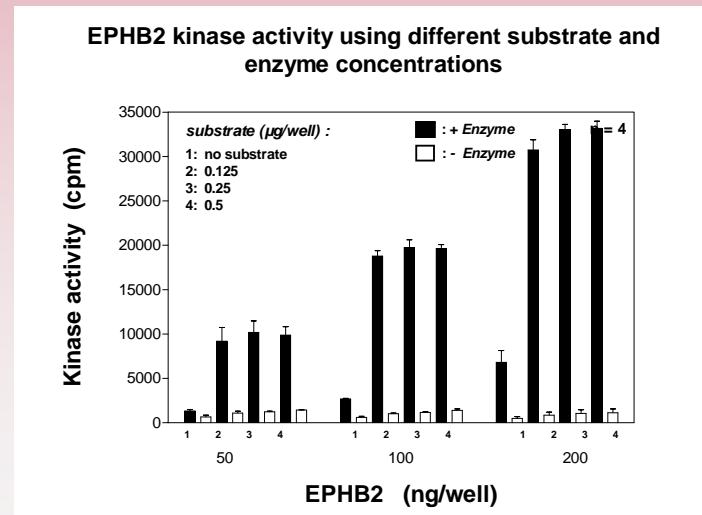
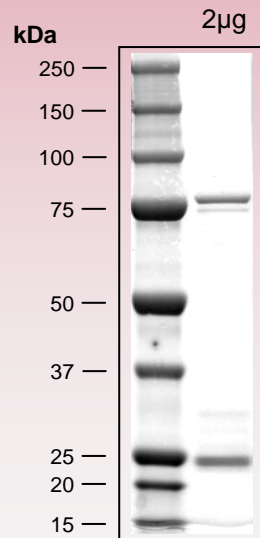


## ProQinase EPHB2(Gly<sub>580</sub>-Val<sub>987</sub>)

697 aminoacids, 79.283 kDa  
 1-218 GST  
 224-229 HIS-tag  
 241-246 Thrombin cleavage site  
 280-697 EPHB2 580-987



### Representative lot: Coomassie stain & activity in PanQinase™ assay



Lot specific datasheet available on request.  
 please contact:  
[info@proqinase.com](mailto:info@proqinase.com)