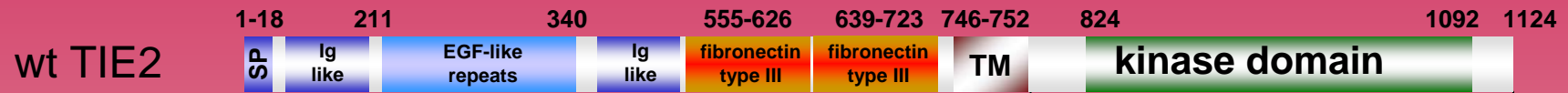


human TIE2 kinase (GenBank Acc.No: NM_000459)



ProQinase TIE2 (Gln₇₇₁-Ala₁₁₂₄)

581 aminoacids, 66.882 kDa

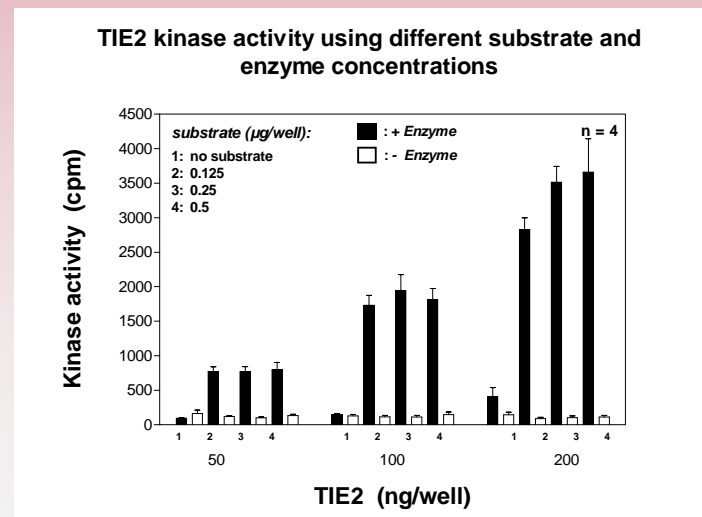
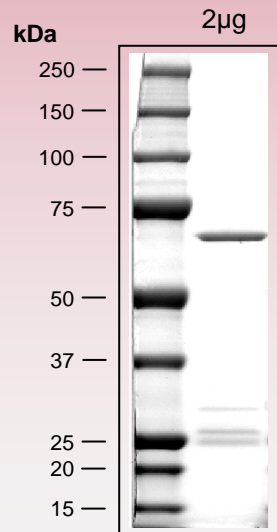
1-218 GST

222-225 factor Xa cleavage site

228-581 TIE2 771-1124



Representative lot: Coomassie stain & activity in PanQinase™ assay



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