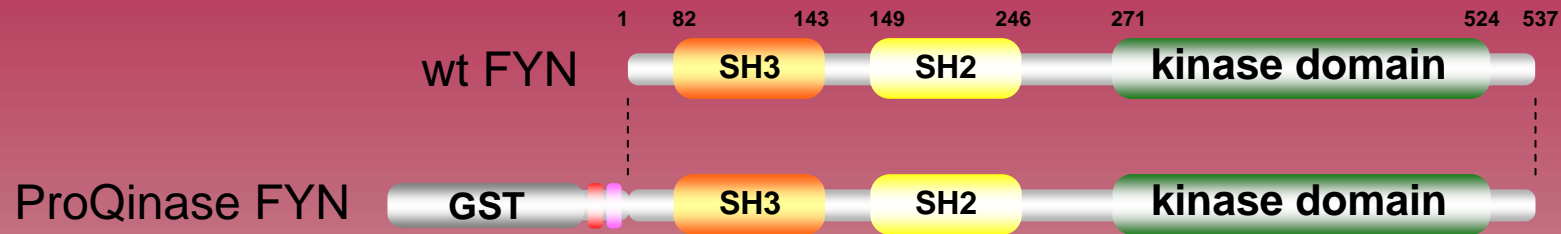
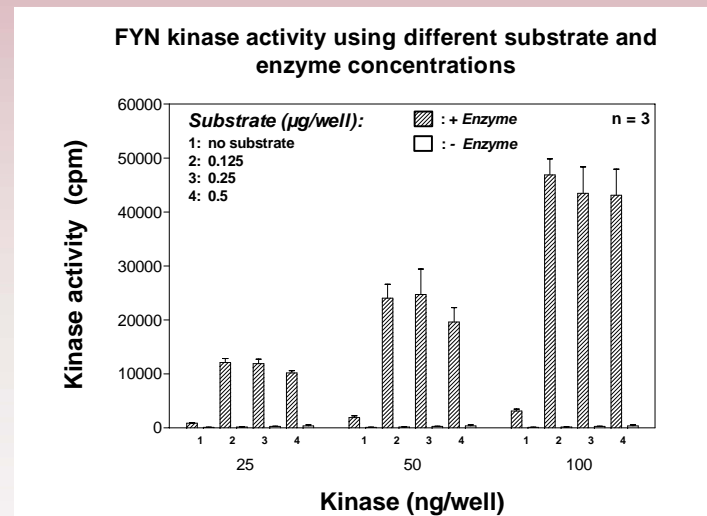
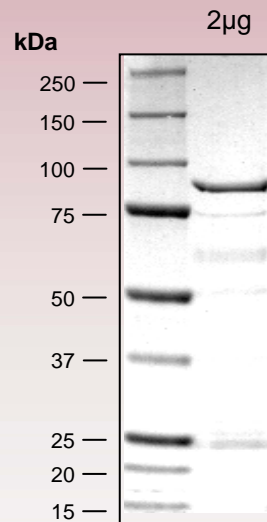


human FYN kinase (GenBank Acc.No: NM_002037)



794 aminoacids, 90.358 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 258-794 FYN 1-537

Representative lot: Coomassie stain & activity in PanQinase™ assay



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