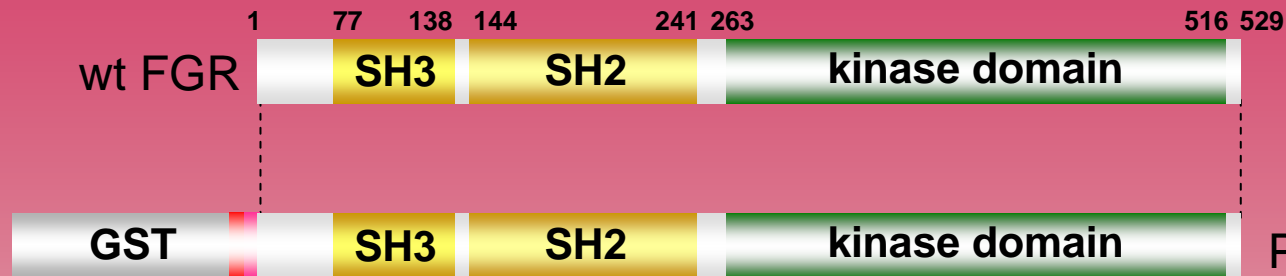


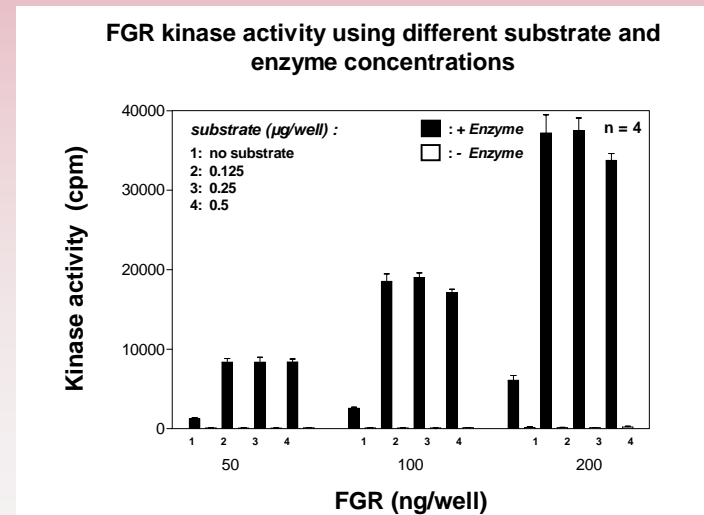
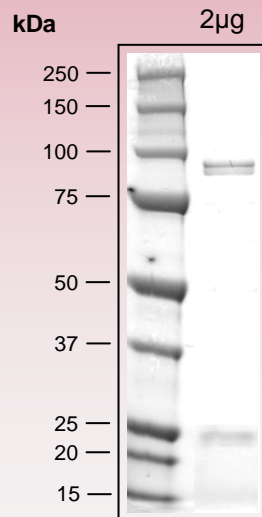
human FGR kinase (GenBank Acc.No: NM_005248)



ProQinase FGR (Met₁-Thr₅₂₉)

830 aminoacids, 93.934 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 262-830 FGR 1-529

Representative lot: Coomassie stain & activity in PanQinase™ assay



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