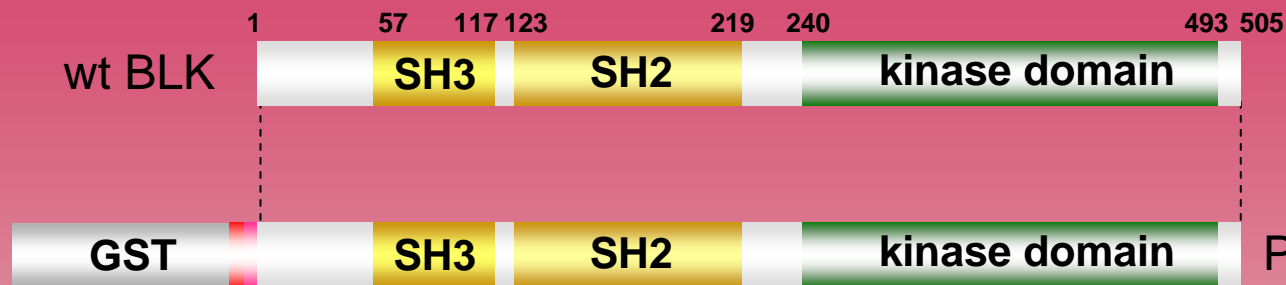


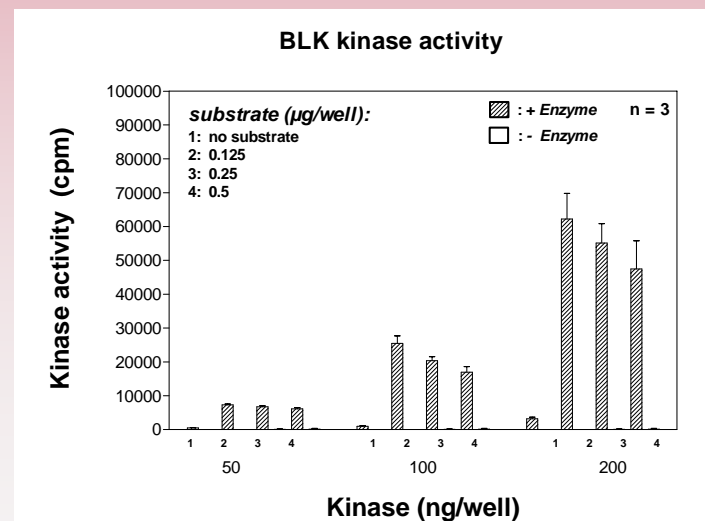
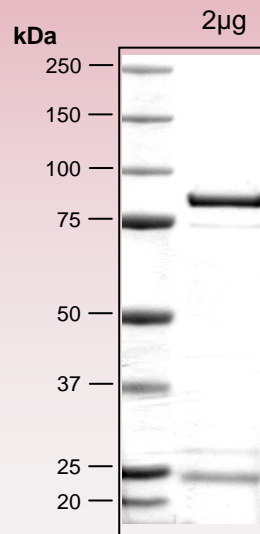
human **BLK** soluble tyrosine kinase (GenBank Acc.No: NM_001715.2)



ProQinase BLK (Met₁-Pro₅₀₅)

760 aminoacids, 87.102 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 256-760 BLK 1-505

Representative lot: Coomassie stain & activity in PanQinase™ assay



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