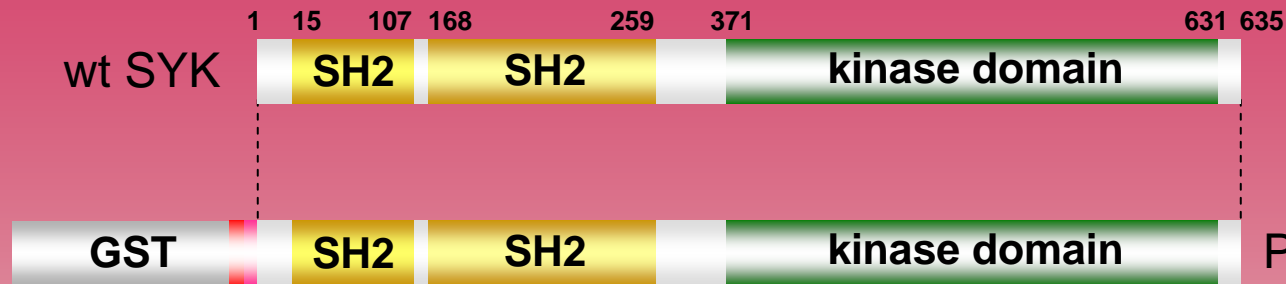


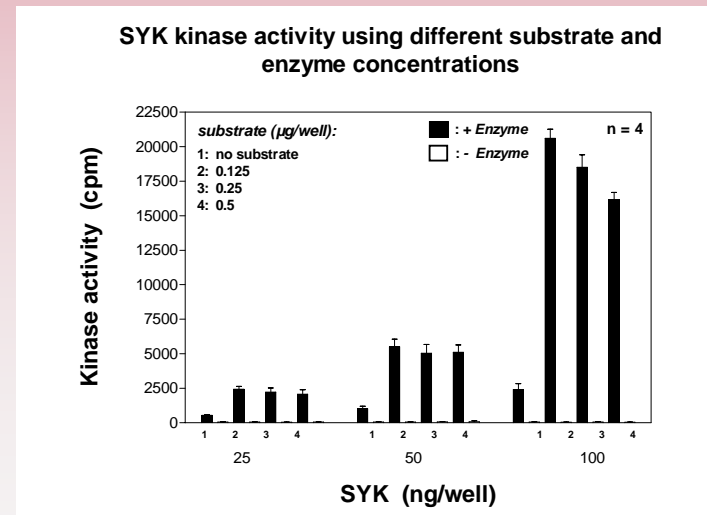
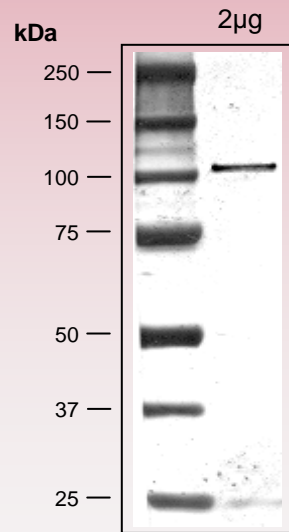
human SYK kinase (GenBank Acc.No: NM_003177)



ProQinase SYK (Met₁-Asn₆₃₅)

918 aminoacids, 104.408 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 284-918 SYK 1-635

Representative lot: Coomassie stain & activity in PanQinase™ assay



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