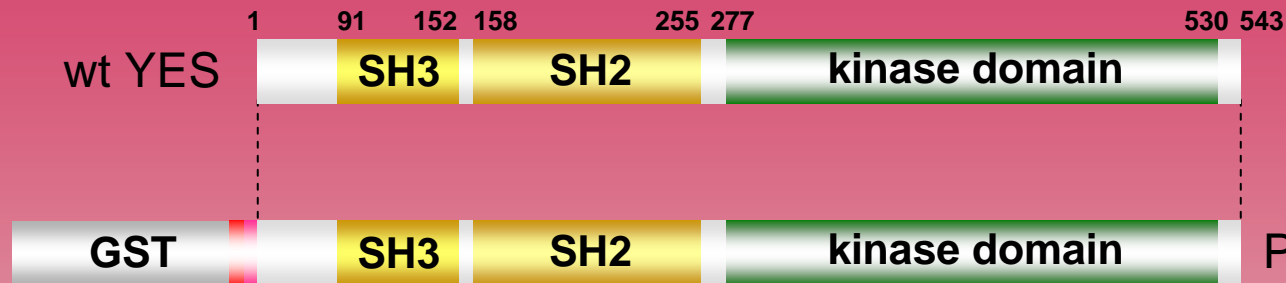


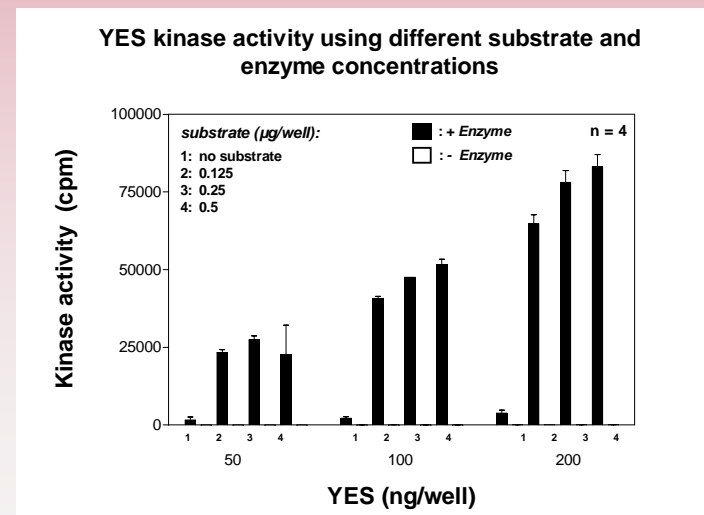
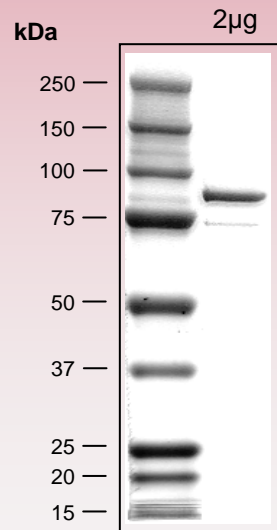
human YES kinase (GenBank Acc.No: NM_005433)



ProQinase YES (Met₁-Leu₅₄₃)

803 aminoacids, 90.997 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 261-803 YES 1-543

Representative lot: Coomassie stain & activity in PanQinase™ assay



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