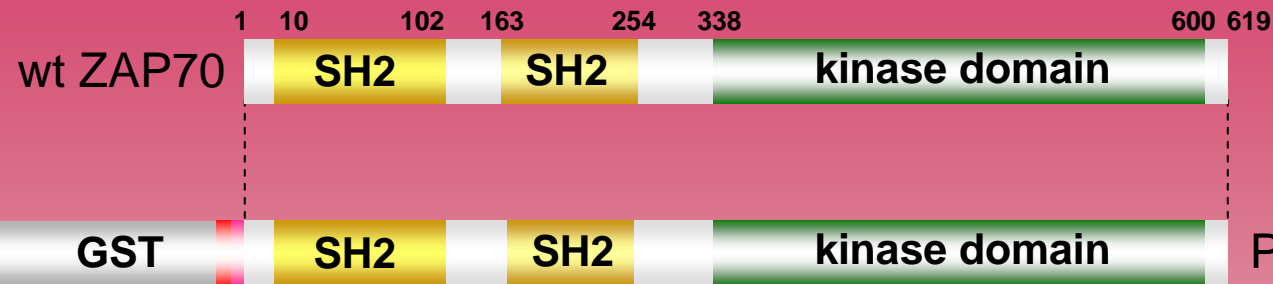


human ZAP70 kinase (GenBank Acc.No: NM_001079)



ProQinase ZAP70 (Met₁-Ala₆₁₉)

877 aminoacids, 99.496 kDa

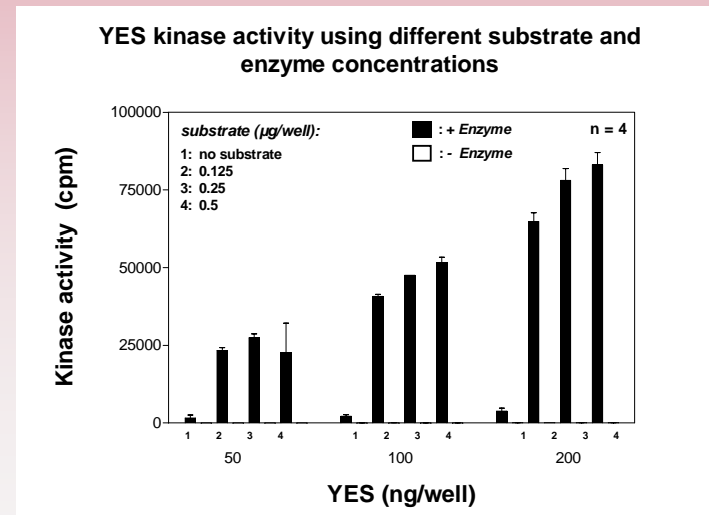
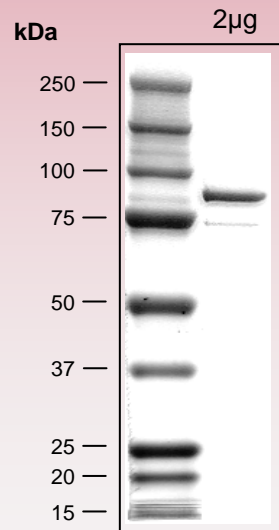
1-218 GST

224-229 HIS-tag

241-246 Thrombin cleavage site

259-877 ZAP70 1-619

Representative lot: Coomassie stain & activity in PanQinase™ assay



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