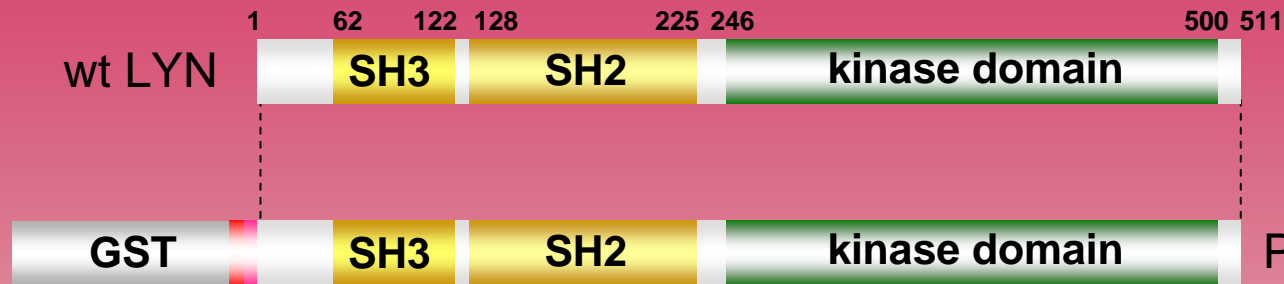


human LYN kinase (GenBank Acc.No: NM_002350)



ProQinase LYN (Met₁-Pro₅₁₁)

772 aminoacids, 88.470 kDa

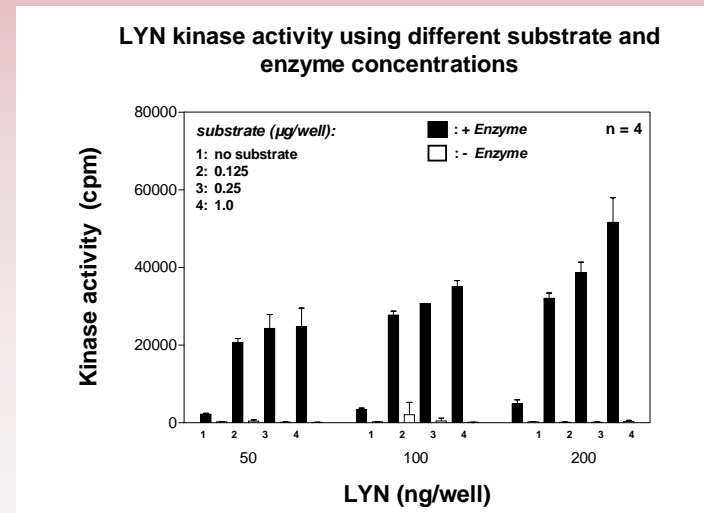
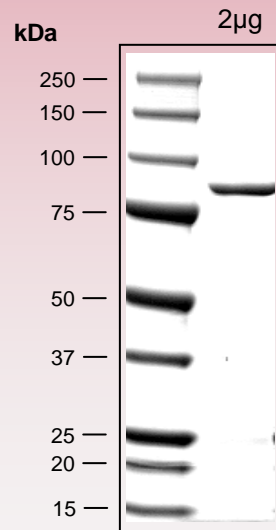
1-218 GST

224-229 HIS-tag

241-246 Thrombin cleavage site

261-772 LYN 1-511

Representative lot: Coomassie stain & activity in PanQinase™ assay



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