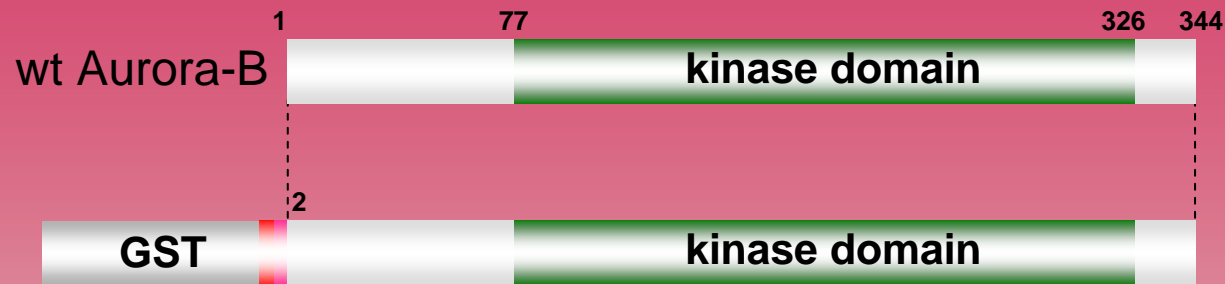


human **Aurora-B** kinase (GenBank Acc.No: NM_004217)



ProQinase Aurora-B (Ala₂-Ala₃₄₄)

613 aminoacids, 70.073 kDa

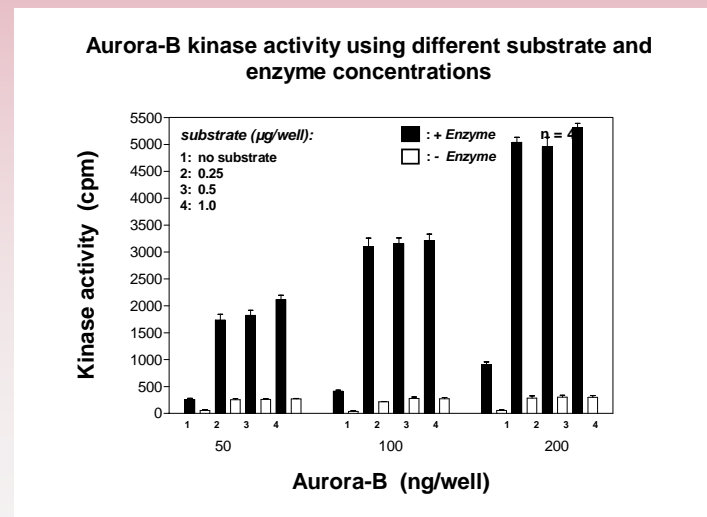
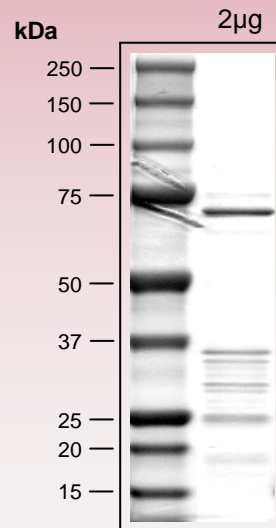
1-218 GST

224-229 HIS-tag

241-246 Thrombin cleavage site

271-613 Aurora-B 2-344

Representative lot: Coomassie stain & activity in PanQinase™ assay



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