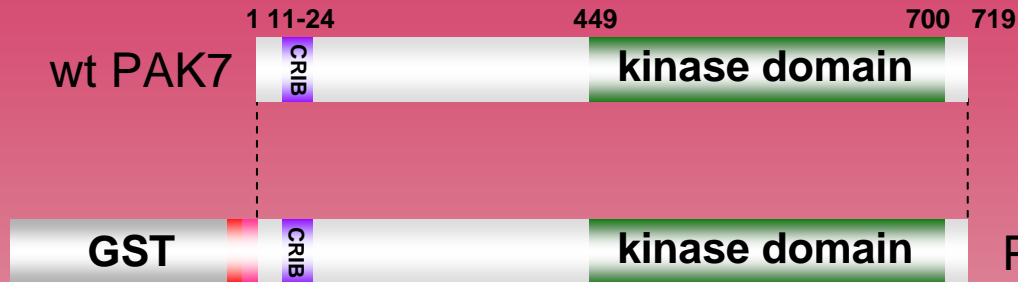


human **PAK7** kinase (GenBank Acc.No: NM_020341)

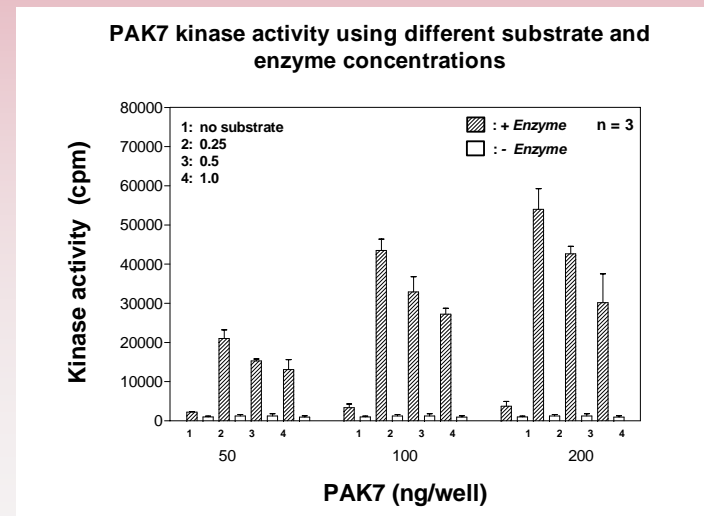
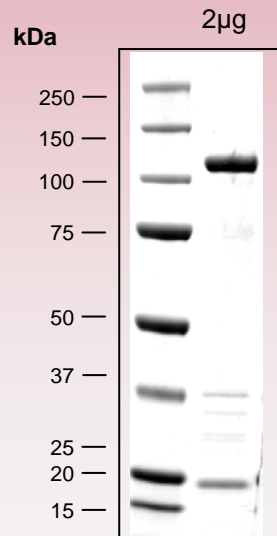


ProQinase PAK7 (Met₁-His₇₁₉)

979 aminoacids, 110.660 kDa

1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 261-979 PAK7 1-719

Representative lot: Coomassie stain & activity in PanQinase™ assay



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