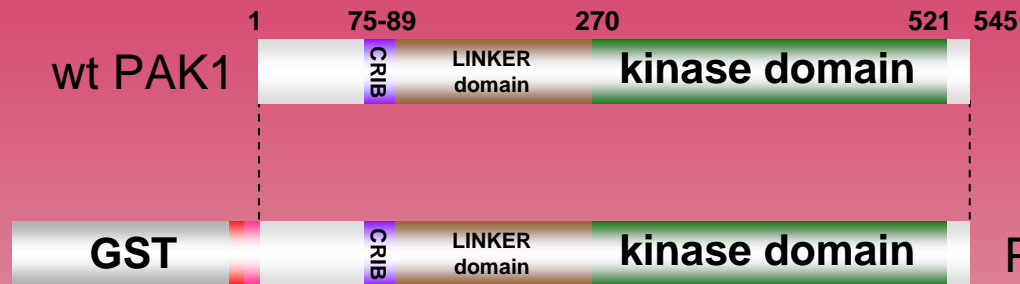


human **PAK1** kinase (GenBank Acc.No: NM_002576)



ProQinase PAK1 (Met₁-His₅₄₅)

805 aminoacids, 90.542 kDa

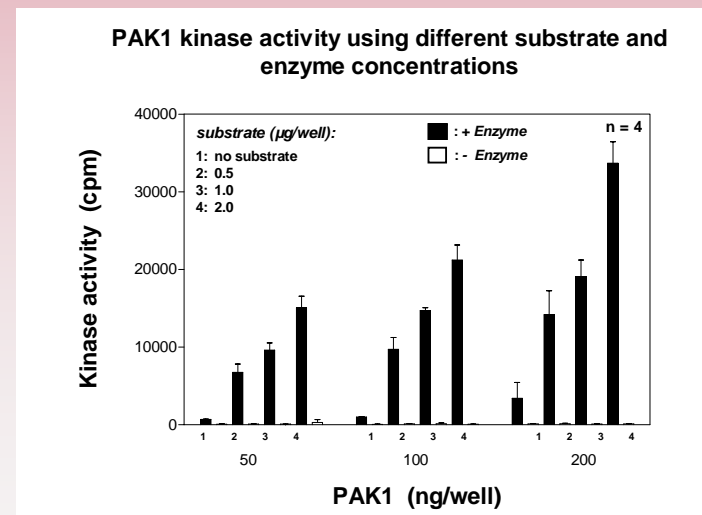
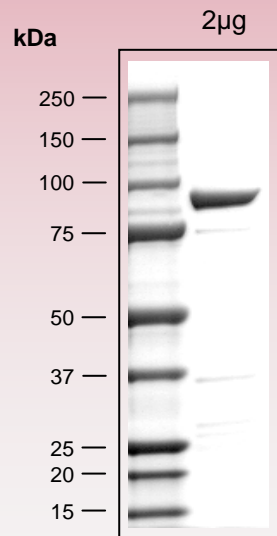
1-218 GST

224-229 HIS-tag

241-246 Thrombin cleavage site

261-805 PAK1 1-545

Representative lot: Coomassie stain & activity in PanQinase™ assay



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