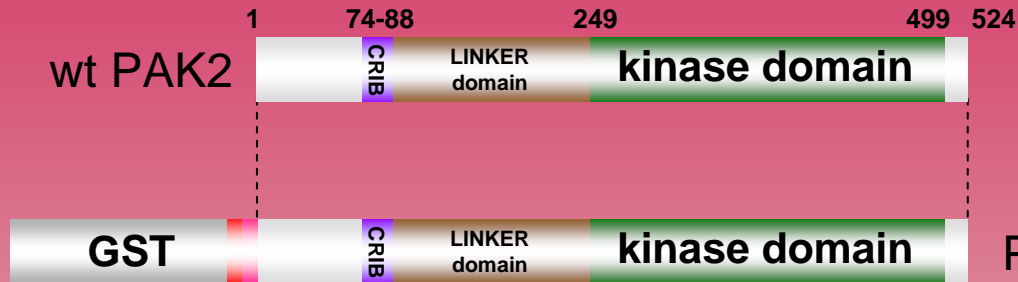


human **PAK2** kinase (GenBank Acc.No: NM_002577)



ProQinase PAK2 (Asp₃-Arg₅₂₄)

784 aminoacids, 87.954 kDa

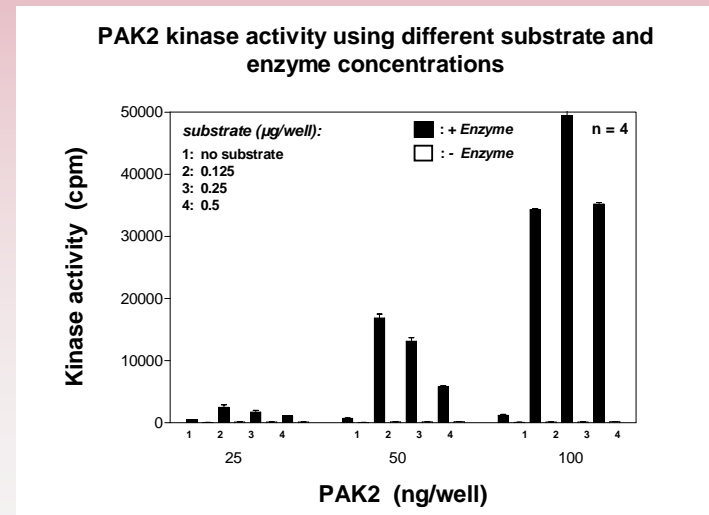
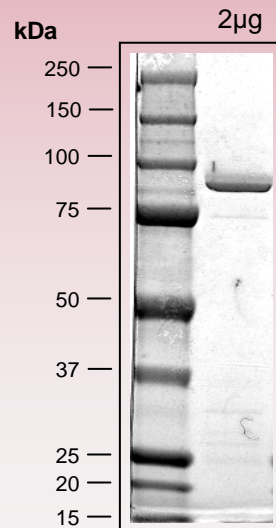
1-218 GST

224-229 HIS-tag

241-246 Thrombin cleavage site

263-784 PAK2 3-524

Representative lot: Coomassie stain & activity in PanQinase™ assay



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