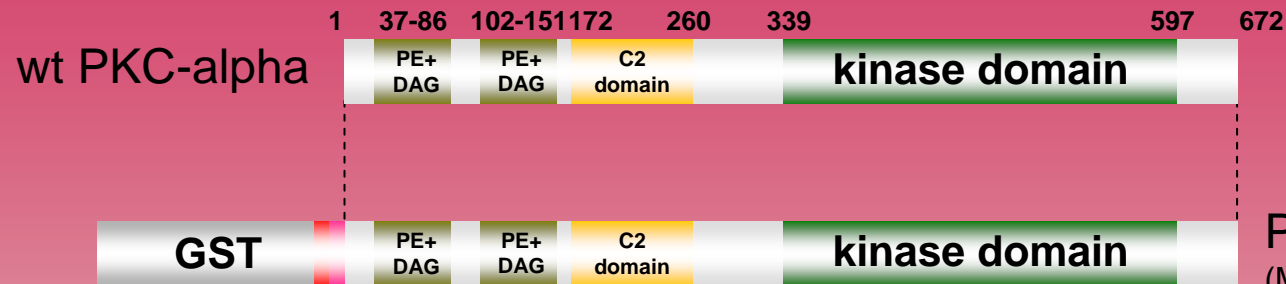
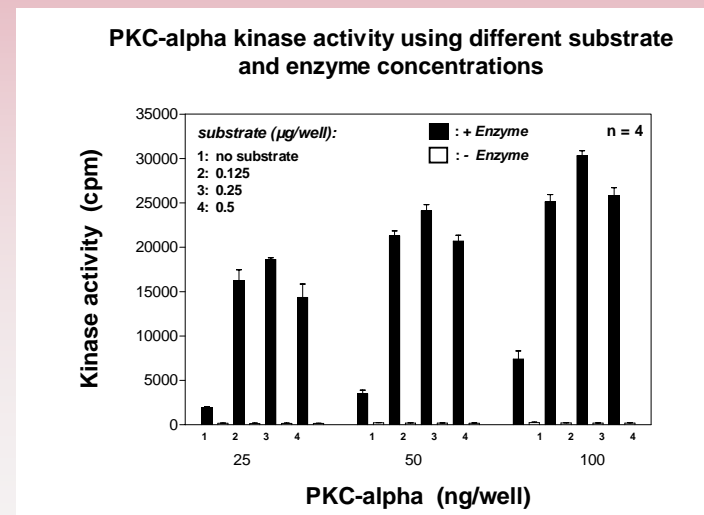
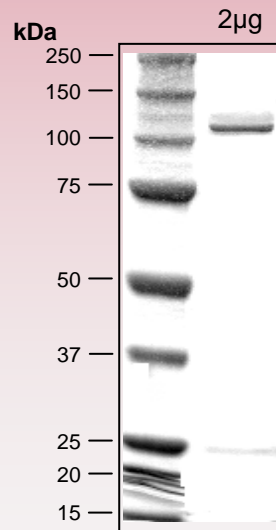


human PKC-alpha kinase (GenBank Acc.No: NM_002737)



962 aminoacids, 109.972 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 279-950 PKC-alpha 1-672

Representative lot: Coomassie stain & activity in PanQinase™ assay



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