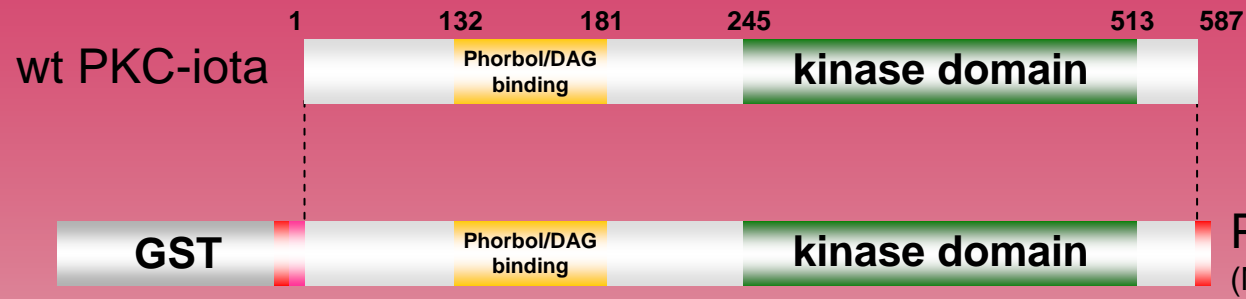
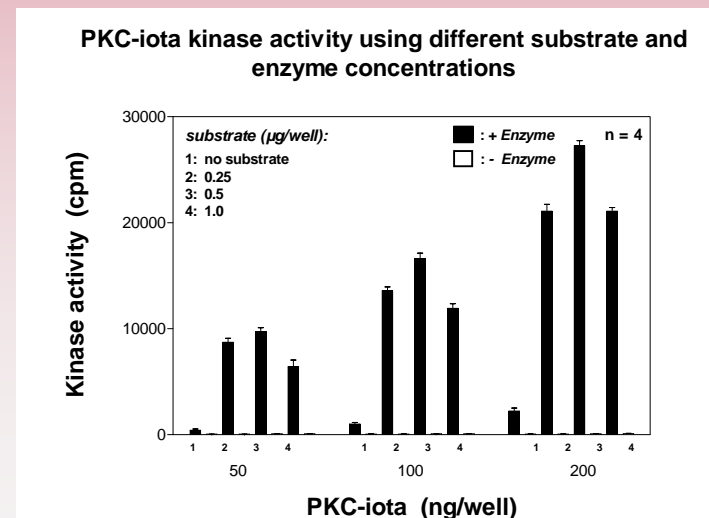
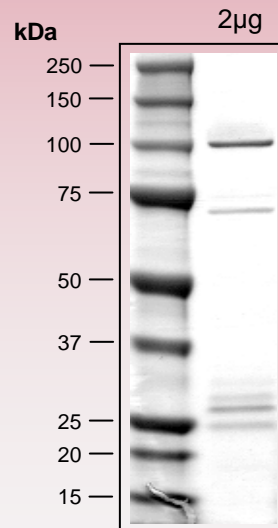


human PKC-iota kinase (GenBank Acc.No: NM_002740)



853 aminoacids, 98.027 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 261-847 PKC-iota 1-587
 848-853 HIS-tag

Representative lot: Coomassie stain & activity in PanQinase™ assay



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