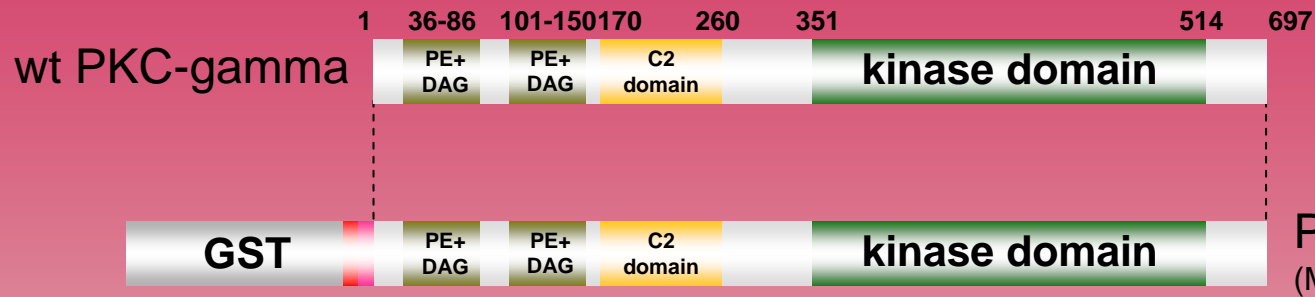
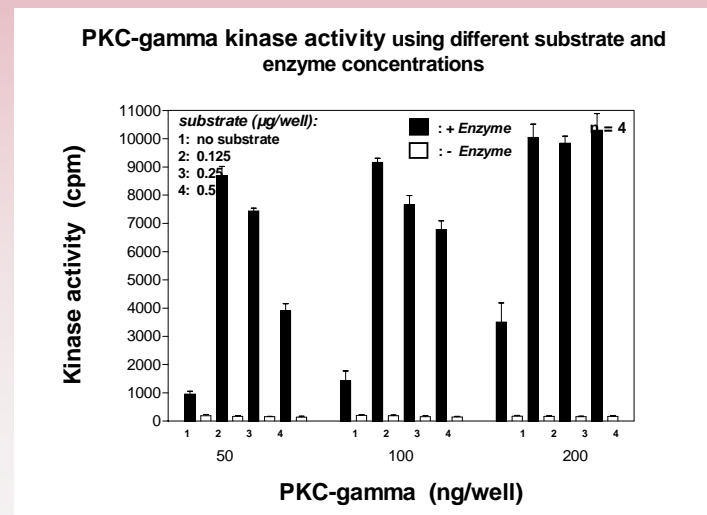
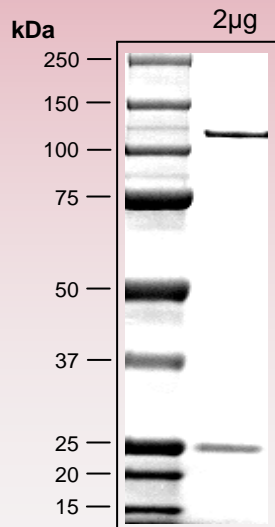


human PKC-gamma kinase (GenBank Acc.No: NM_002739.3)



993 aminoacids, 112.289 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 270-966 PKC-gamma 1-697

Representative lot: Coomassie stain & activity in PanQinase™ assay



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