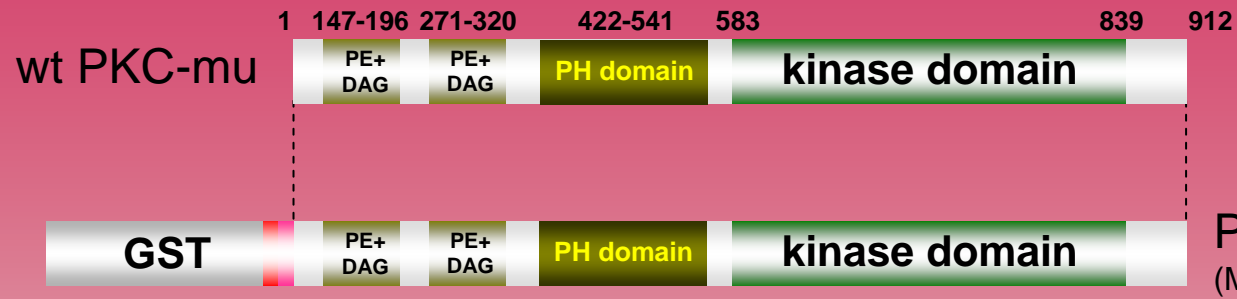
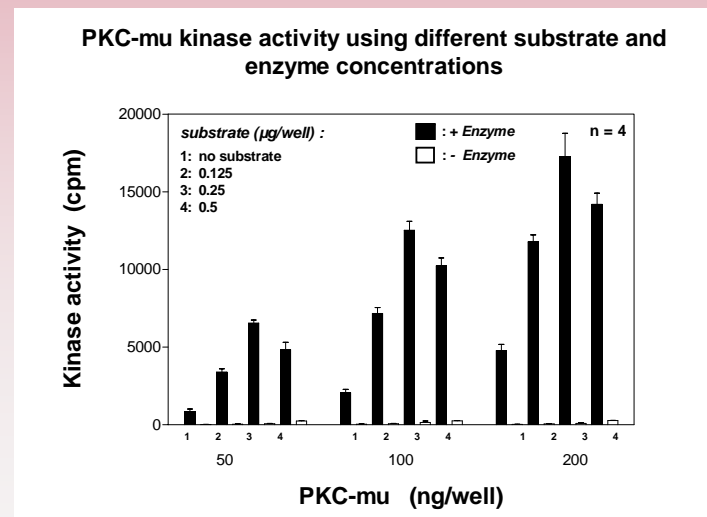
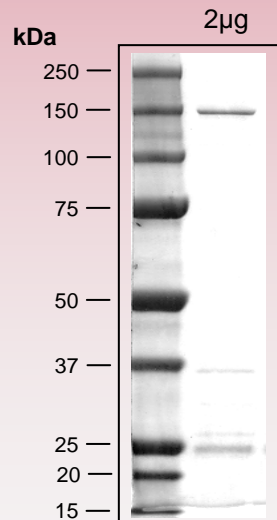


human PKC-mu kinase (GenBank Acc.No: NM_002742)



1246 aminoacids, 139.489 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 334-1246 PKC-mu 1-912

Representative lot: Coomassie stain & activity in PanQinase™ assay



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