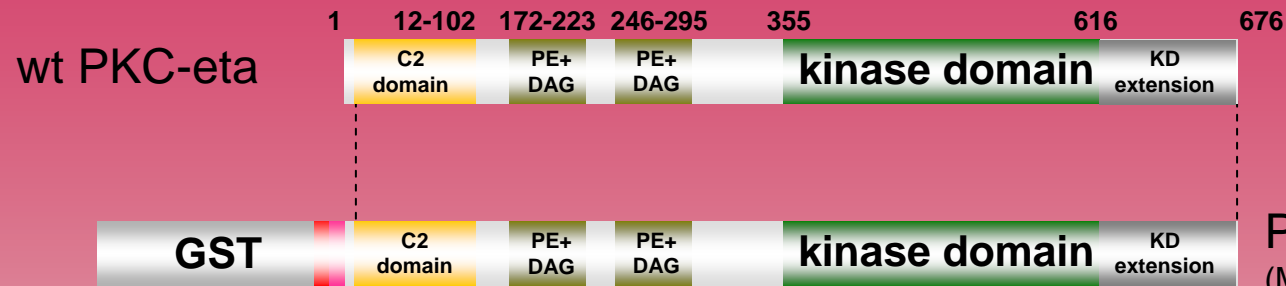
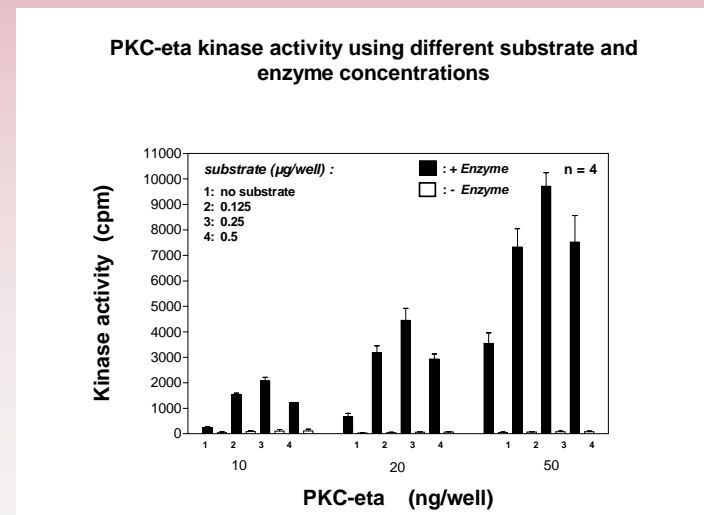
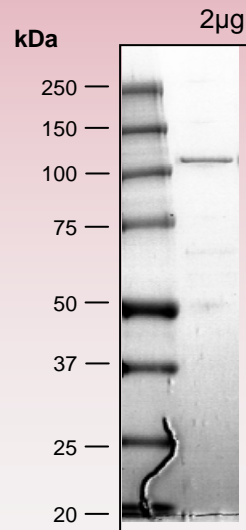


# murine PKC-eta kinase (GenBank Acc.No: D90242)



960 aminoacids, 109.613 kDa  
 1-218 GST  
 224-229 HIS-tag  
 241-246 Thrombin cleavage site  
 278-960 PKC-eta 1-683

## Representative lot: Coomassie stain & activity in PanQinase™ assay



Lot specific datasheet available on request.  
 please contact:  
[info@proqinase.com](mailto:info@proqinase.com)