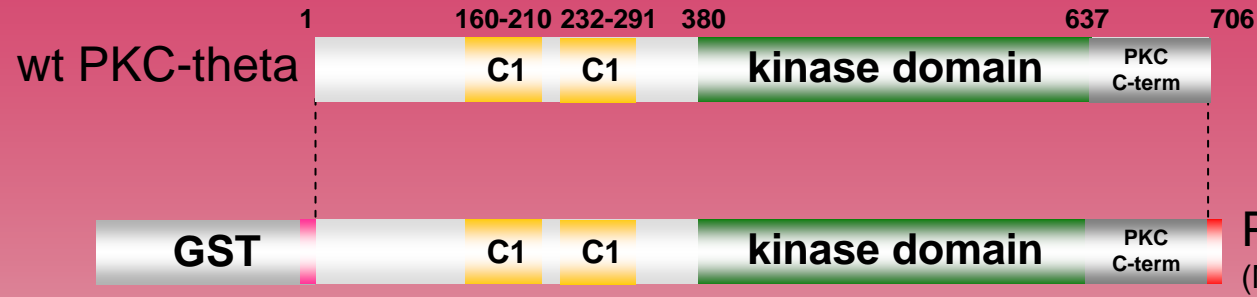


# human PKC-theta kinase (GenBank Acc.No: NM\_006257)



945 aminoacids, 109.677 kDa

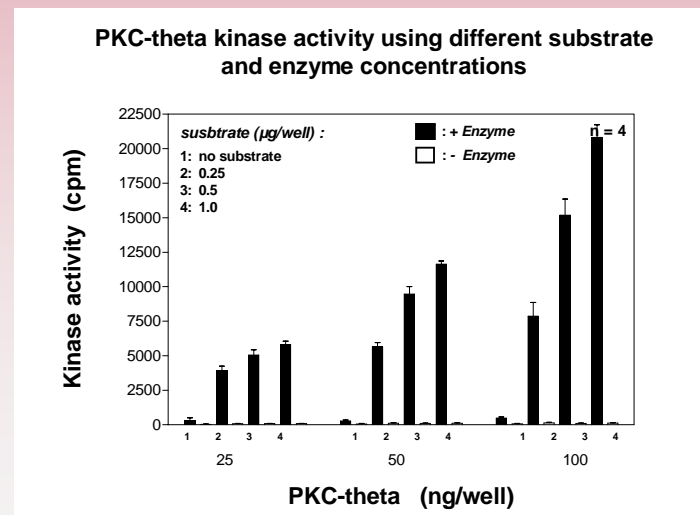
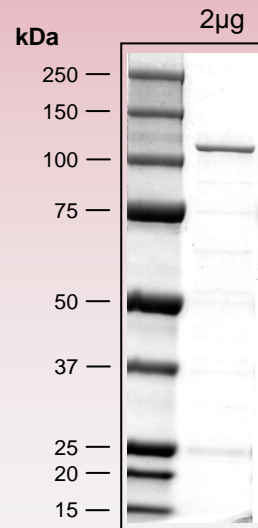
1-218 GST

222-225 factor Xa cleavage site

234-939 PKC-iota 1-587

940-845 HIS-tag

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



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