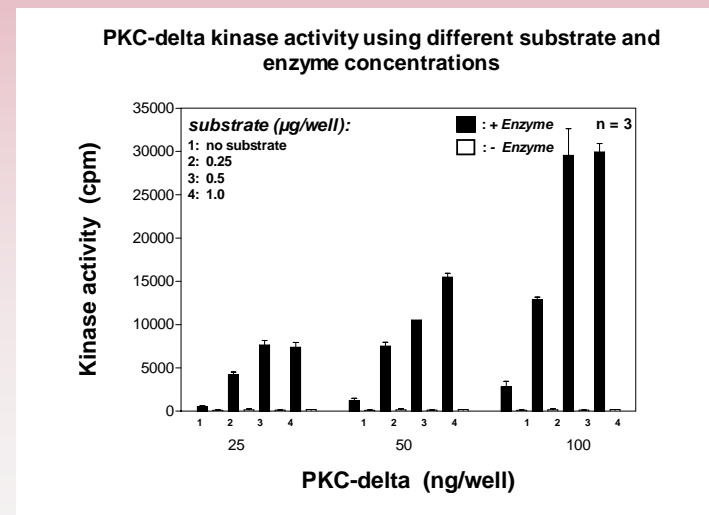
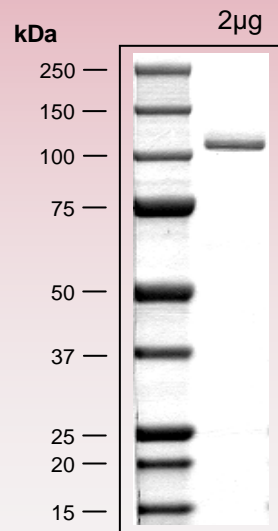


# human PKC-delta kinase (GenBank Acc.No: NM\_006254.3)



975 aminoacids, 111.765 kDa  
 1-218 GST  
 224-229 HIS-tag  
 241-246 Thrombin cleavage site  
 261-975 PKC-delta 1-675

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



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