

# human PBK kinase (GenBank Acc.No: NM\_018492)



## ProQinase PBK (Met<sub>1</sub>-Val<sub>322</sub>)

596 aminoacids, 67.521 kDa

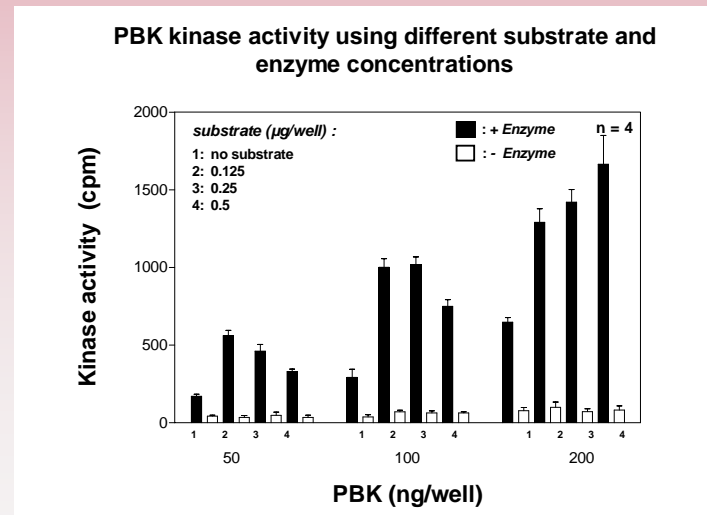
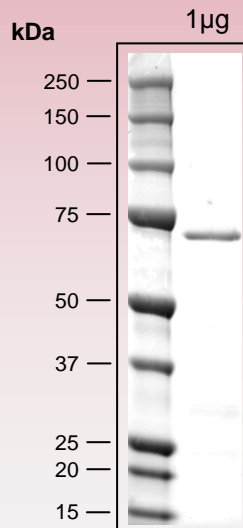
1-218 GST

224-229 HIS-tag

241-246 Thrombin cleavage site

270-596 PBK 1-322

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



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