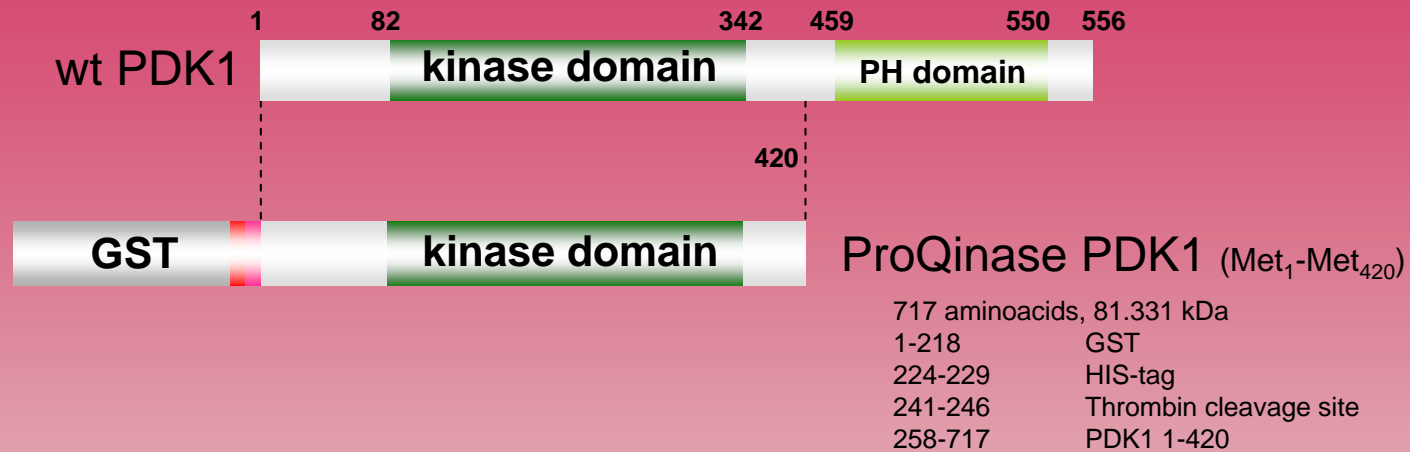
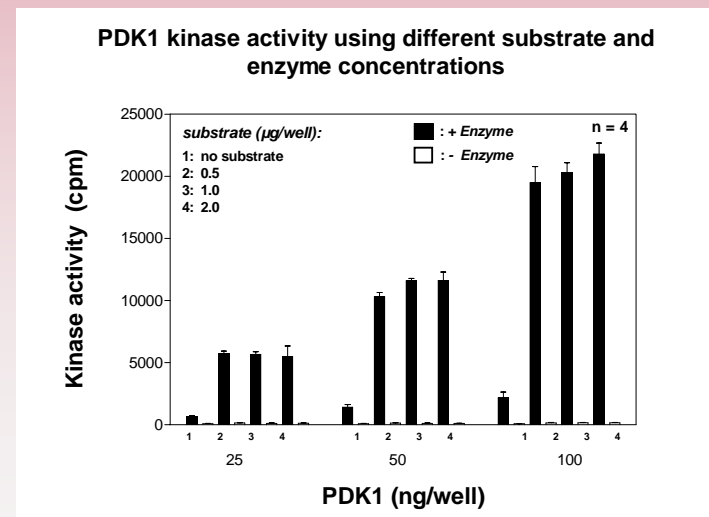
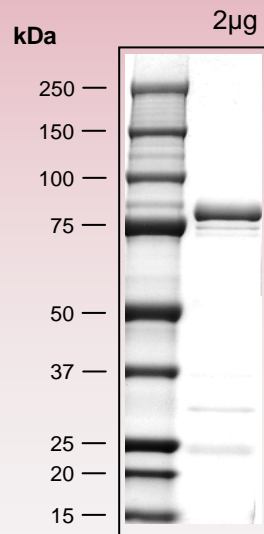


# human PDK1 kinase (GenBank Acc.No: NM\_002613)



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



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