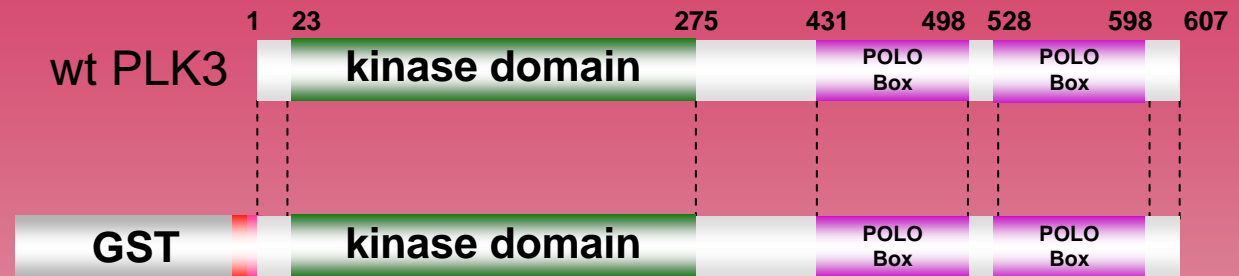


human **PLK3** kinase (GenBank Acc.No: U56998.1)



862 aminoacids, 97.131 kDa

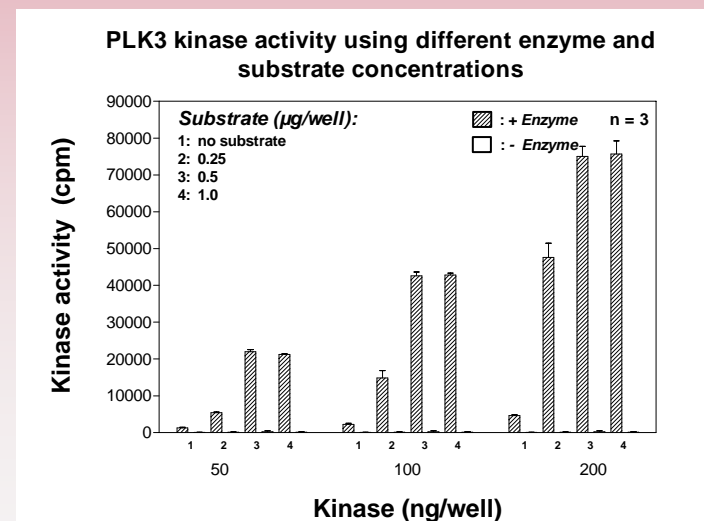
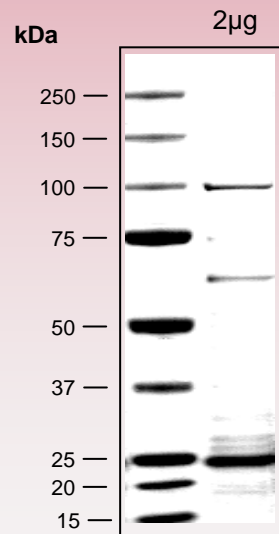
1-218 GST

224-229 HIS-tag

241-246 Thrombin cleavage site

256-862 PLK3 1-607

Representative lot: Coomassie stain & activity in PanQinase™ assay



Lot specific datasheet available on request.
 please contact:
info@proqinase.com