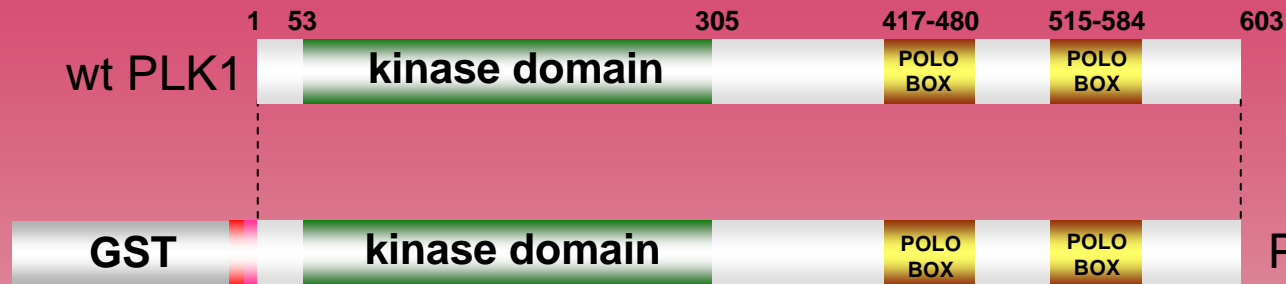


human **PLK1** kinase (GenBank Acc.No: NM_005030)



ProQinase PLK1 (Met₁-Ser₆₀₃)

873 aminoacids, 99.311 kDa

1-218

GST

224-229

HIS-tag

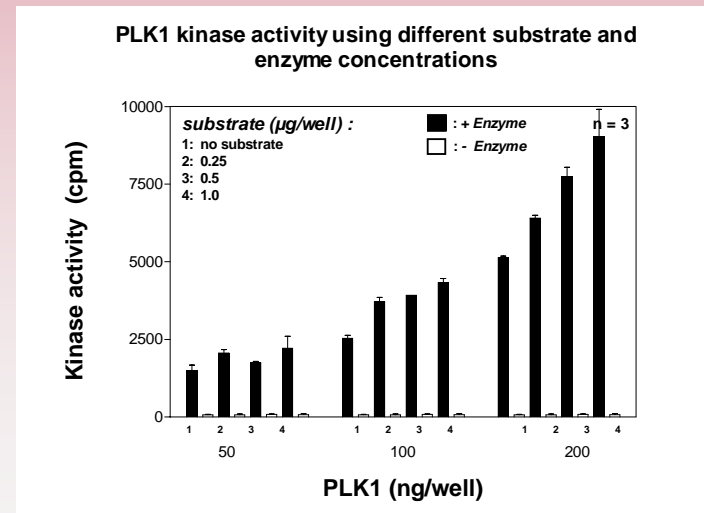
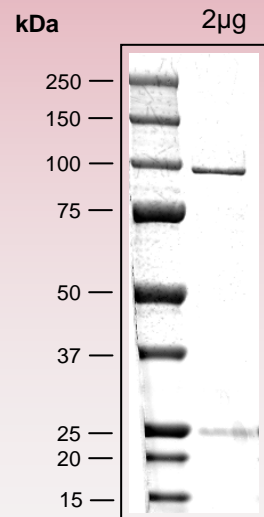
241-246

Thrombin cleavage site

271-873

PLK1 1-603

Representative lot: Coomassie stain & activity in PanQinase™ assay



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