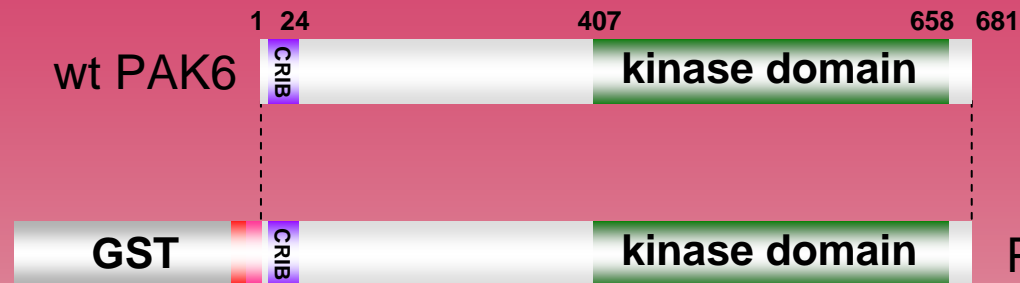


human **PAK6** kinase (GenBank Acc.No: NM_020168)



ProQinase PAK6 (Met₁-Cys₆₈₁)

941 aminoacids, 104.780 kDa

1-218

GST

224-229

HIS-tag

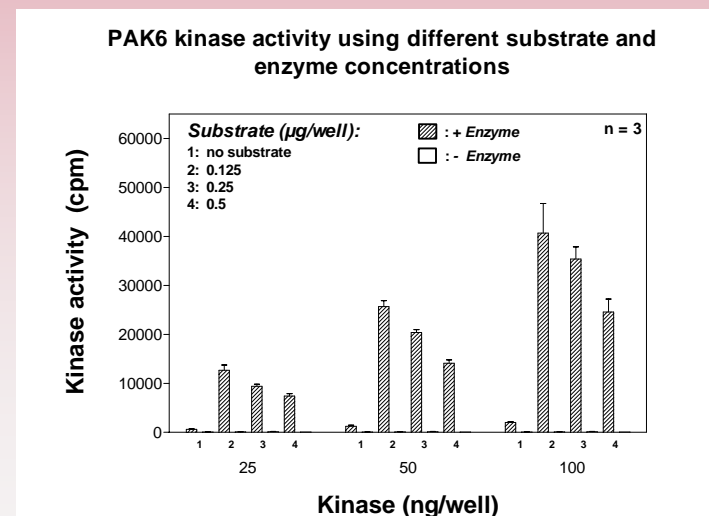
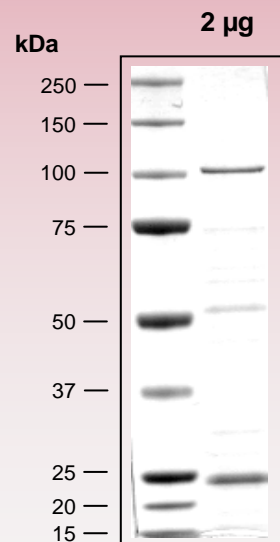
241-246

Thrombin cleavage site

261-941

PAK6 1-681

Representative lot: Coomassie stain & activity in PanQinase™ assay



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