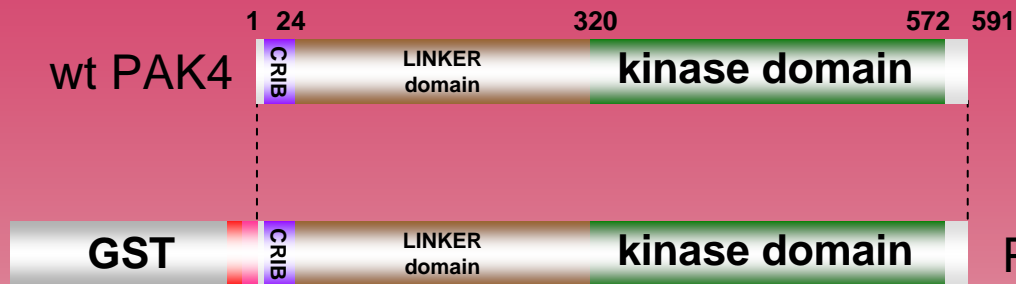


human **PAK4** kinase (GenBank Acc.No: NM_005884)



ProQinase PAK4 (Met₁-Arg₅₉₁)

851 aminoacids, 93.968 kDa

1-218

GST

224-229

HIS-tag

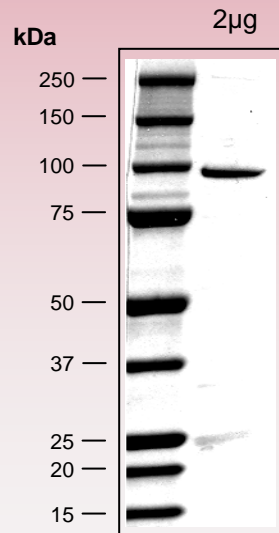
241-246

Thrombin cleavage site

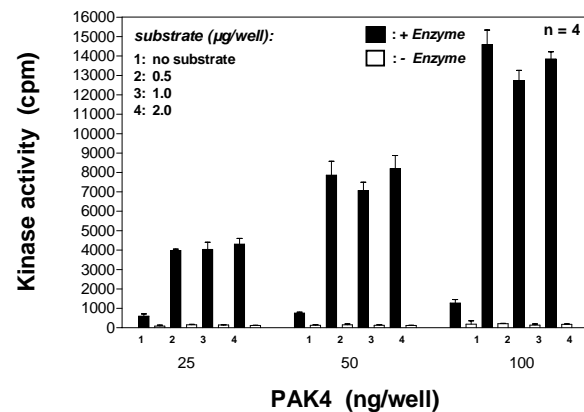
261-851

PAK4 1-591

Representative lot: Coomassie stain & activity in PanQinase™ assay



PAK4 kinase activity using different substrate and enzyme concentrations



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