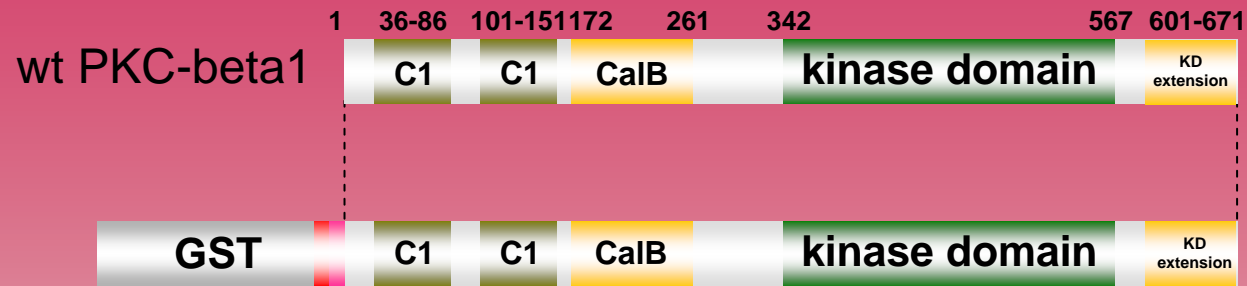


human PKC-beta1 kinase (GenBank Acc.No: NM_002738)



ProQinase PKC-beta1 (Met₁-Val₆₇₁)

900 aminoacids, 103.517 kDa

1-218 GST

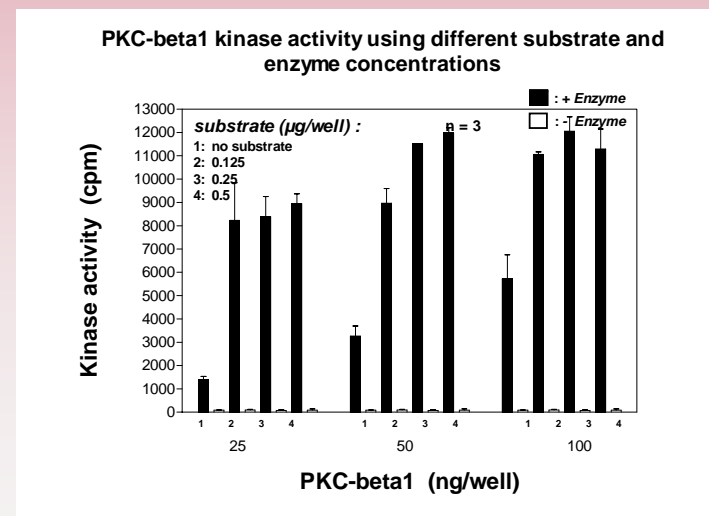
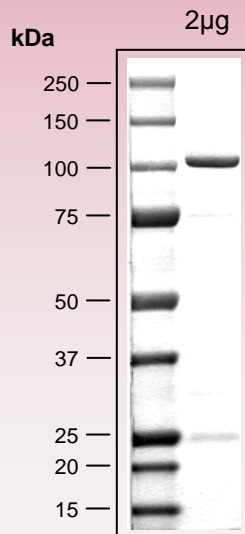
221-226

Thrombin cleavage site

230-900

PKC-beta1 1-671

Representative lot: Coomassie stain & activity in PanQinase™ assay



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