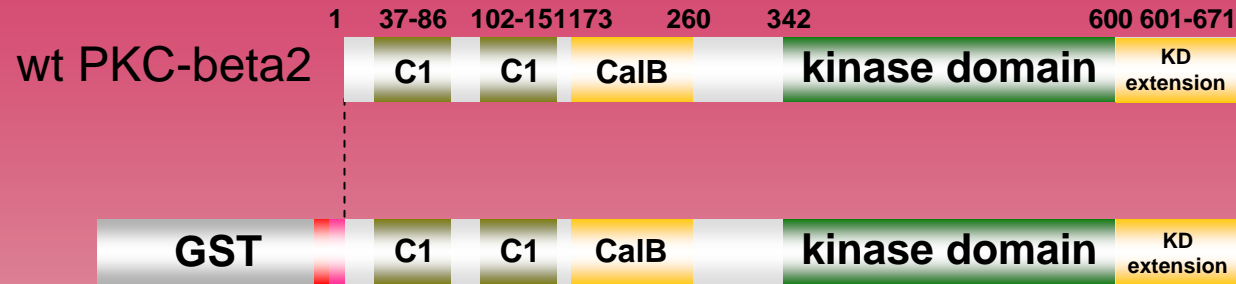


human PKC-beta2 kinase (GenBank Acc.No: X07109)



ProQinase PKC-beta2 (Met₁-Ser₆₇₃)

902 aminoacids, 103.660 kDa

1-218 GST

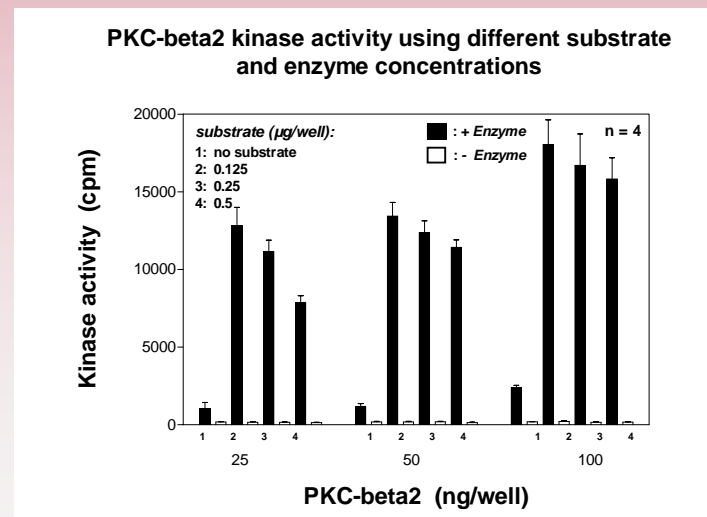
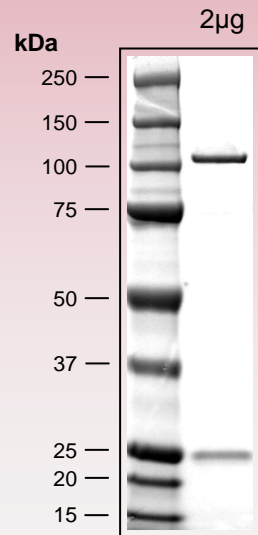
221-226

Thrombin cleavage site

230-902

PKC-beta2 1-673

Representative lot: Coomassie stain & activity in PanQinase™ assay



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