

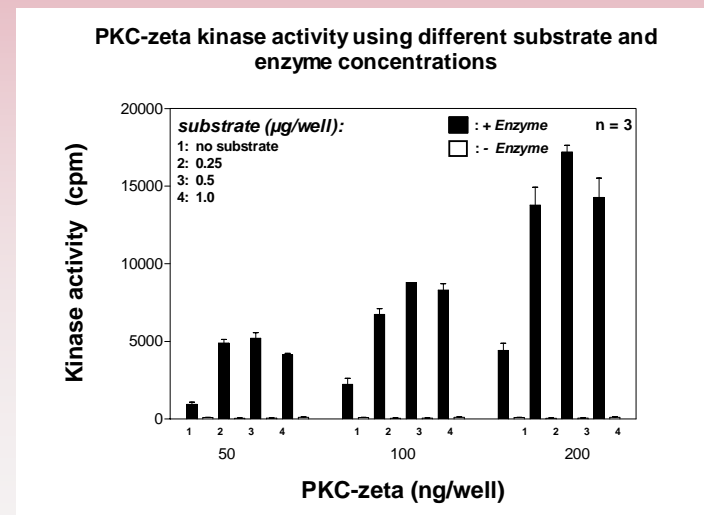
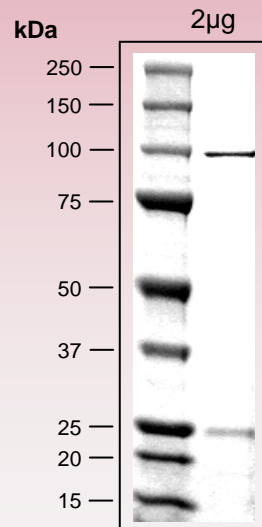
human PKC-zeta kinase (GenBank Acc.No: Z15108)



ProQinase PKC-zeta
(Gly₆-Val₅₈₄)

843 aminoacids, 96.654 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 265-939 PKC-iota 6-584

Representative lot: Coomassie stain & activity in PanQinase™ assay



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