

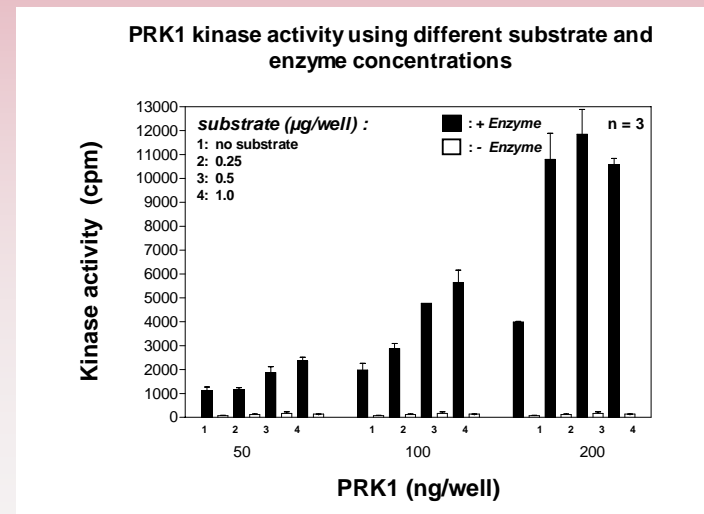
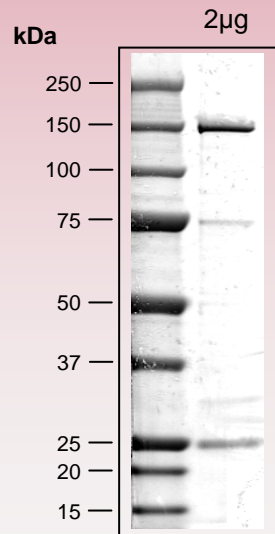
human PRK1 kinase (GenBank Acc.No: NM_002759)



ProQinase PRK1 (Met₁-Cys₉₄₂)

1230 aminoacids, 136.910 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 278-1219 PRK1 1-942

Representative lot: Coomassie stain & activity in PanQinase™ assay



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