

human **MAPKAPK5** kinase (GenBank Acc.No: NM_139078.1)

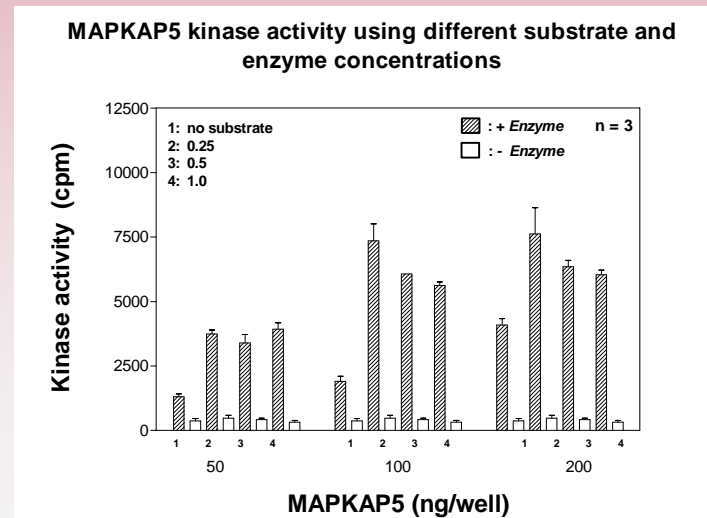
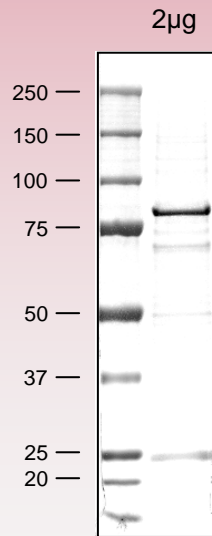


ProQinase MAPKAPK5 (Met₁-Gln₄₇₃)

733 aminoacids, 84.116 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 260-733 MAPKAPK5 1-473



Representative lot: Coomassie stain & activity in PanQinase™ assay



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