

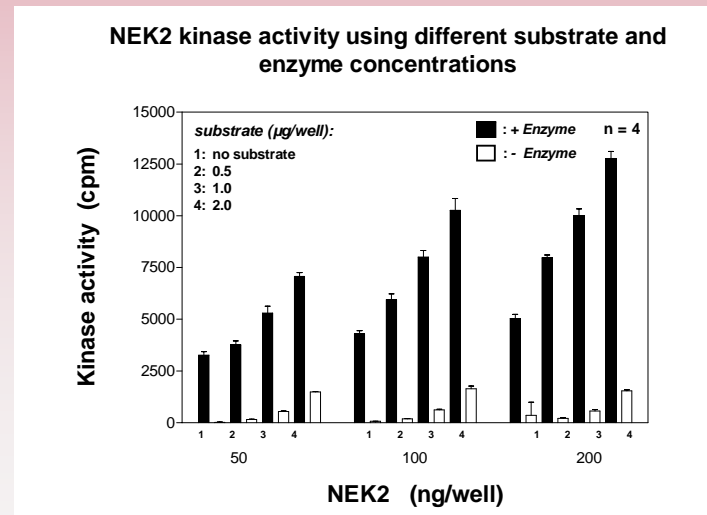
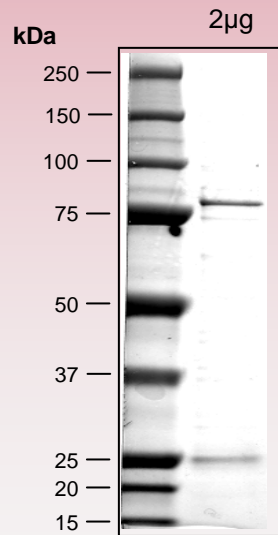
human **NEK2** kinase (GenBank Acc.No: NM_002492)



ProQinase NEK2 (Met₁-Arg₄₄₅)

715 aminoacids, 82.687 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 271-7156 NEK2 1-445

Representative lot: Coomassie stain & activity in PanQinase™ assay



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