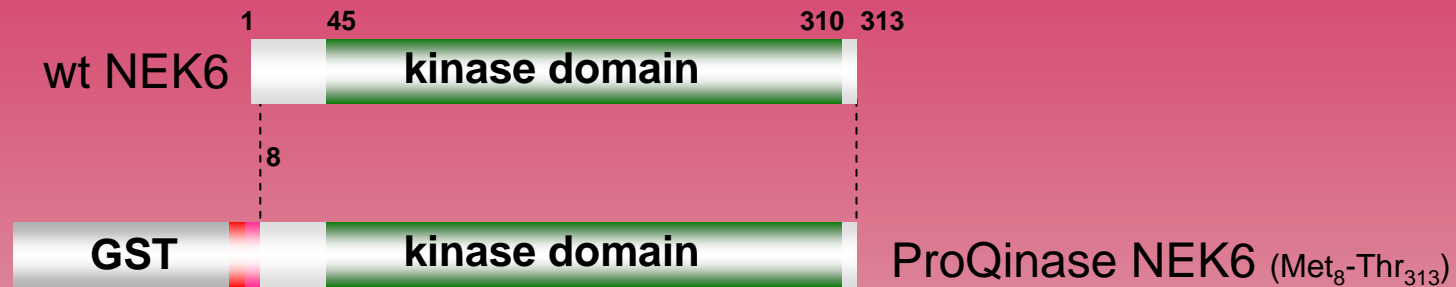
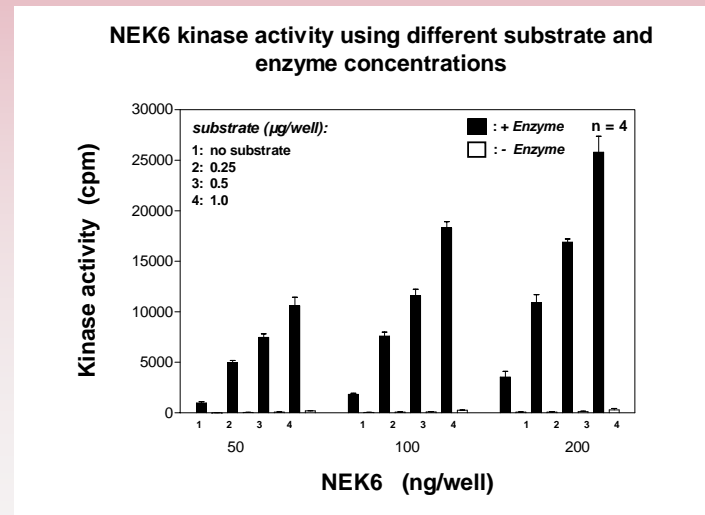
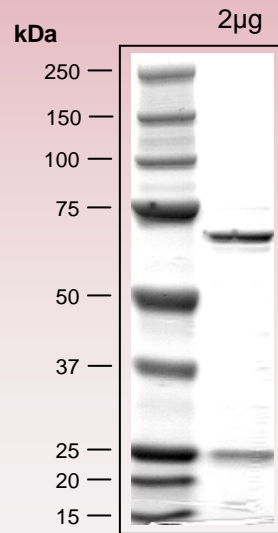


human **NEK6** kinase (GenBank Acc.No: NM_014397)



594 aminoacids, 68.041 kDa
 1-218 GST
 224-229 HIS-tag
 221-226 Thrombin cleavage site
 278-594 NEK6 8-313

Representative lot: Coomassie stain & activity in PanQinase™ assay



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