

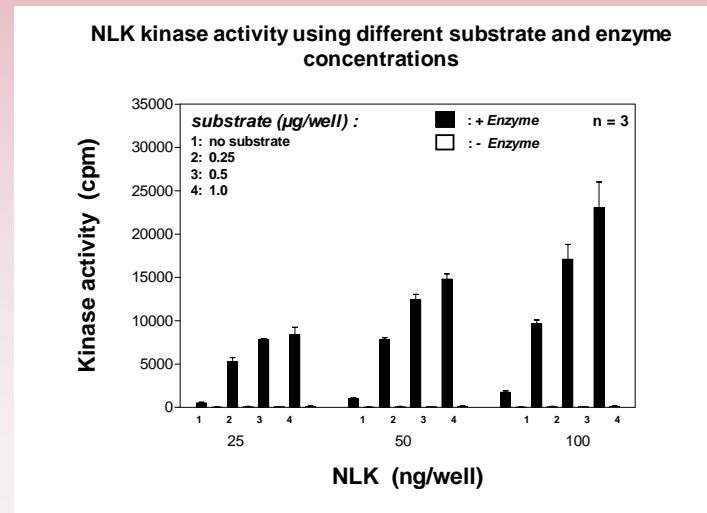
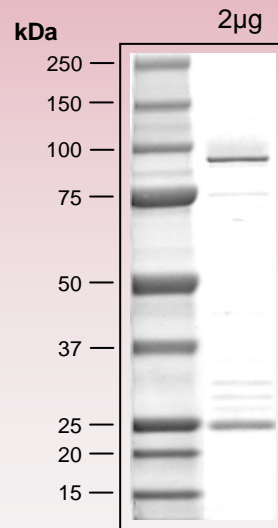
human **NLK** kinase (GenBank Acc.No: NM_016231)



ProQinase NLK (Met₁-Glu₅₁₅)

819 aminoacids, 91.818 kDa
 1-218 GST
 224-229 HIS-tag
 221-226 Thrombin cleavage site
 278-819 NLK 1-515

Representative lot: Coomassie stain & activity in PanQinase™ assay



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