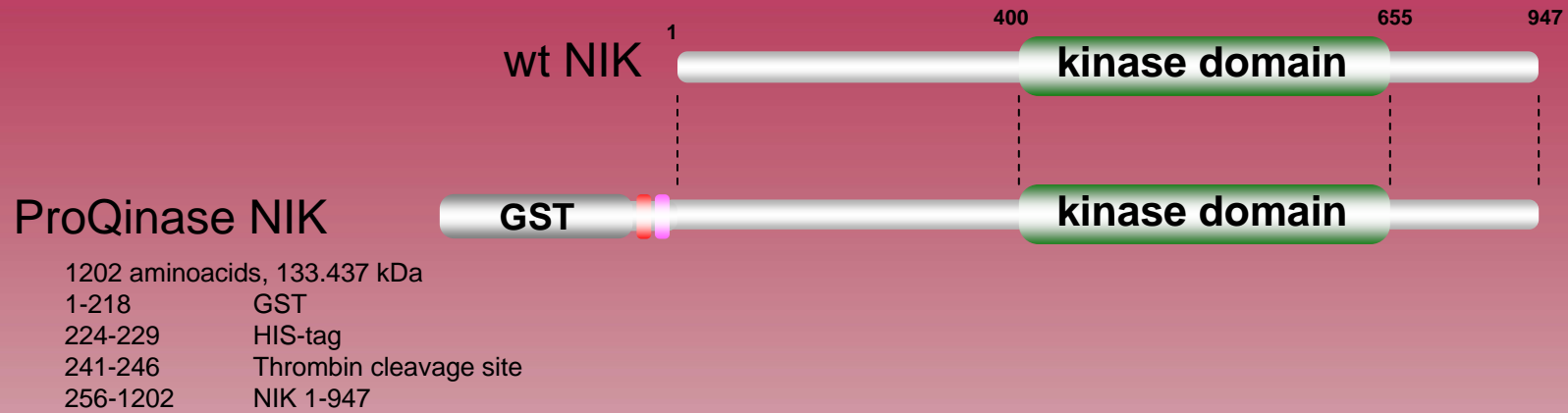
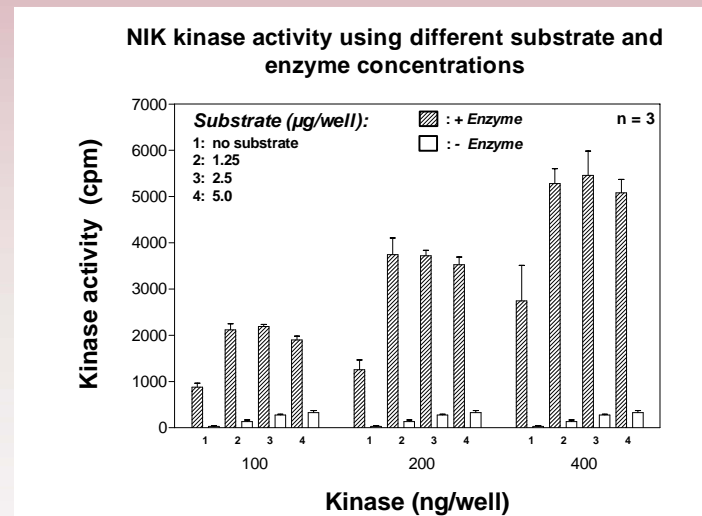
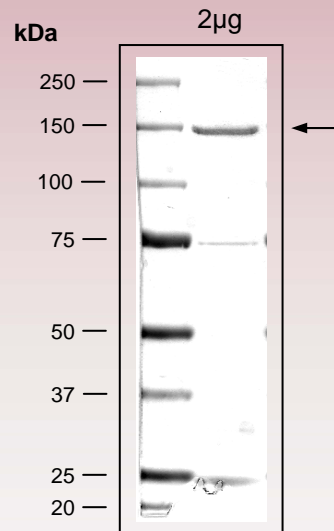


human **NIK** kinase (GenBank Acc.No: NM_003954)



Representative lot: Coomassie stain & activity in PanQinase™ assay



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