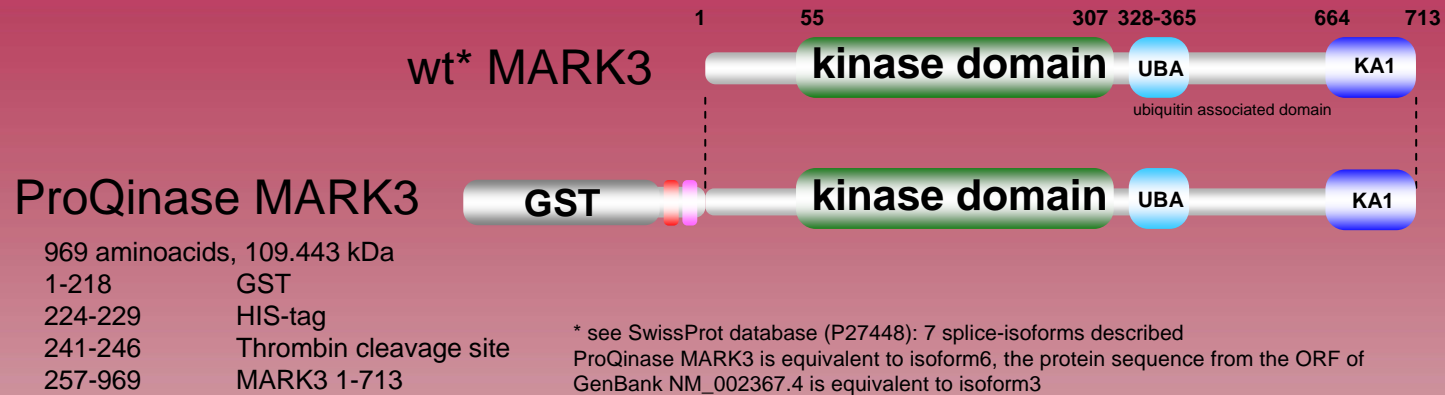
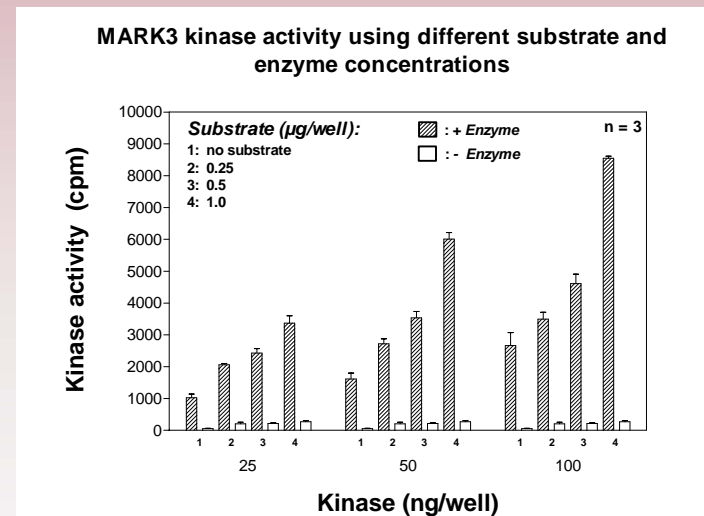
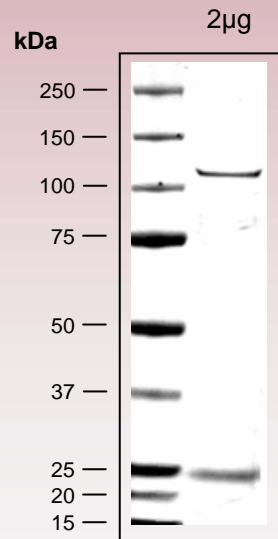


human **MARK3** kinase (GenBank Acc.No: NM_002367.4*, Swiss-Prot P27448)



Representative lot: Coomassie stain & activity in PanQinase™ assay



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