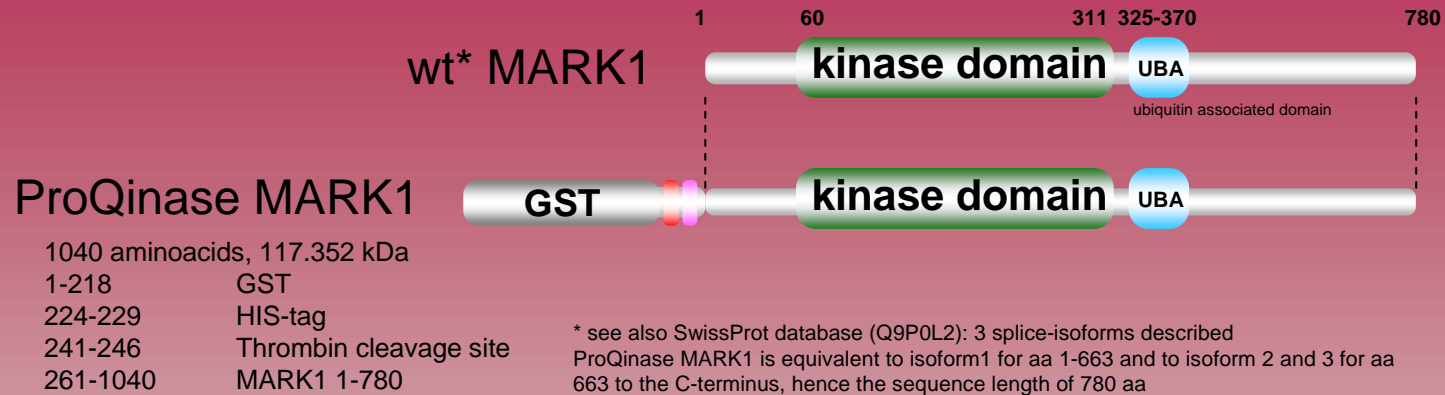
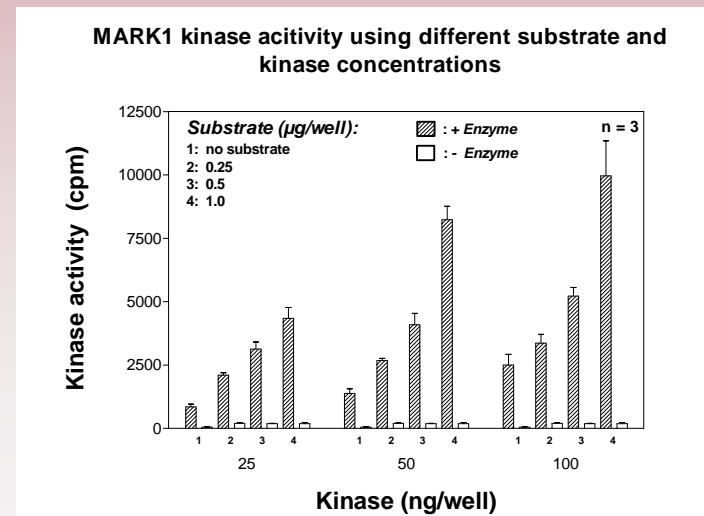
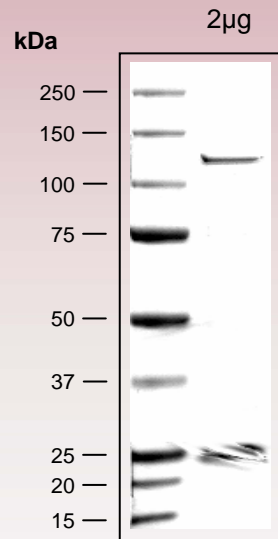


human **MARK1** kinase (GenBank Acc.No: NM_018650*)



Representative lot: Coomassie stain & activity in PanQinase™ assay



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