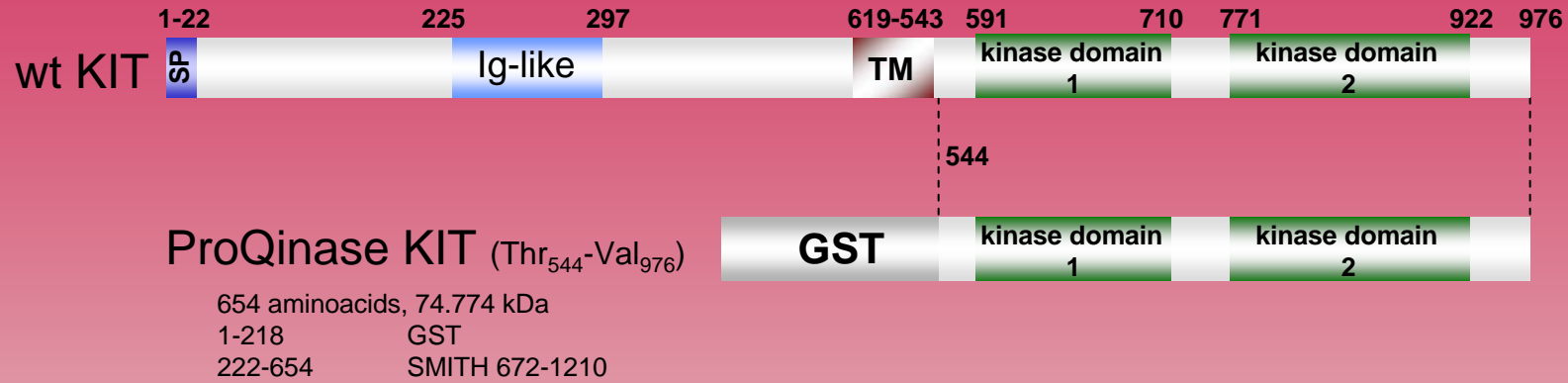
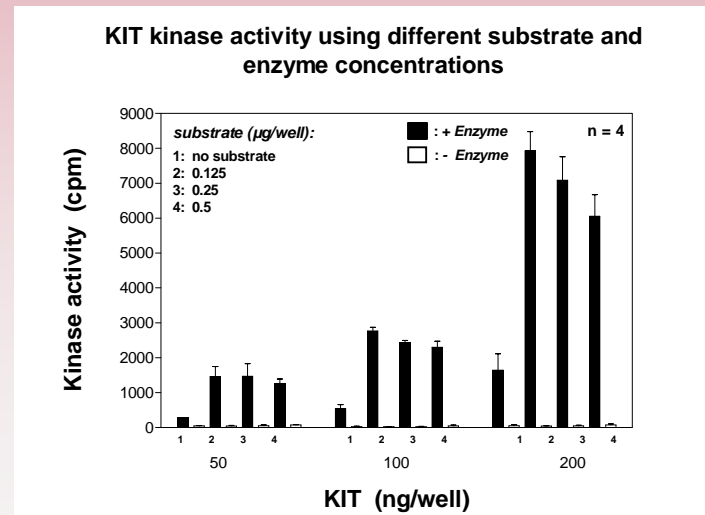
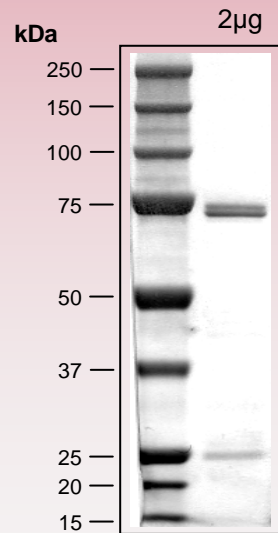


human **KIT** kinase (GenBank Acc.No: NM_000222)



Representative lot: Coomassie stain & activity in PanQinase™ assay



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