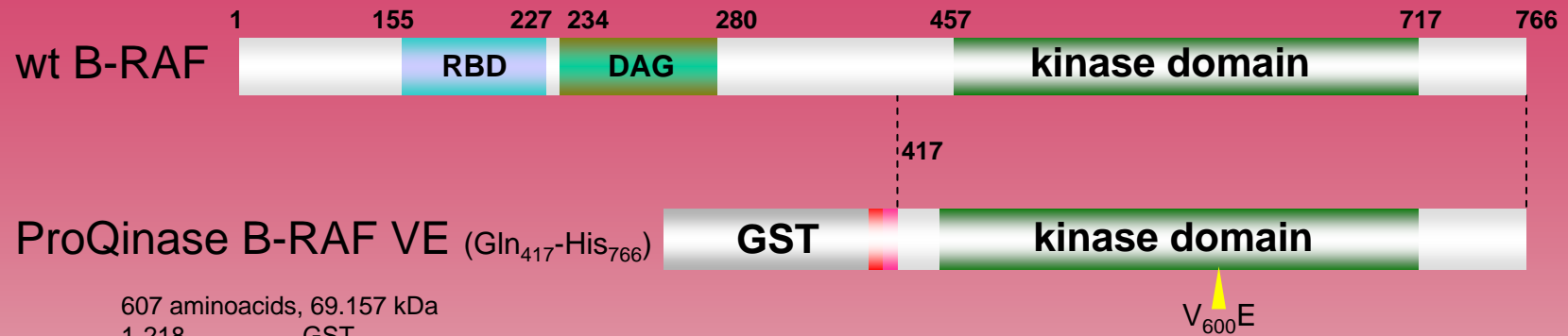
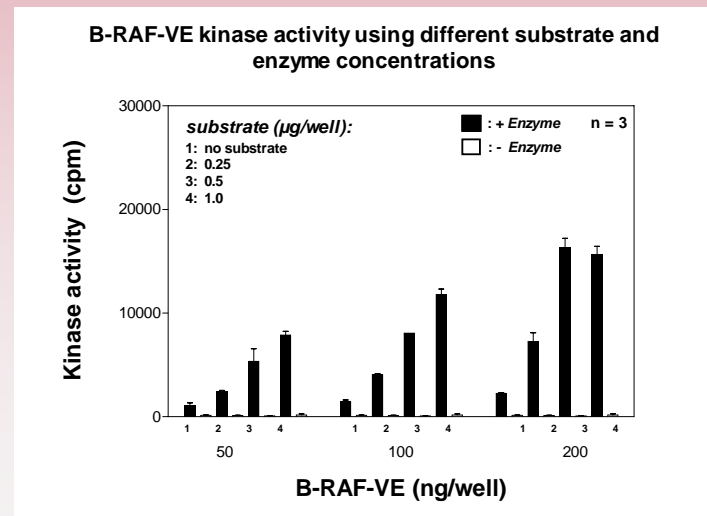
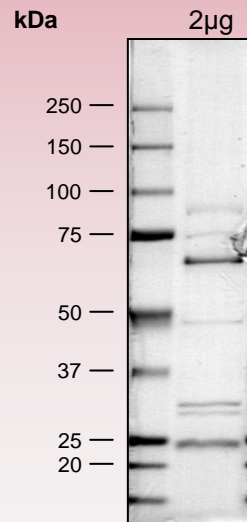


human **B-RAF VE** kinase (GenBank Acc.No: NM_004333.2)



607 aminoacids, 69.157 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 258-607 B-RAF 417-766

Representative lot: Coomassie stain & activity in PanQinase™ assay



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