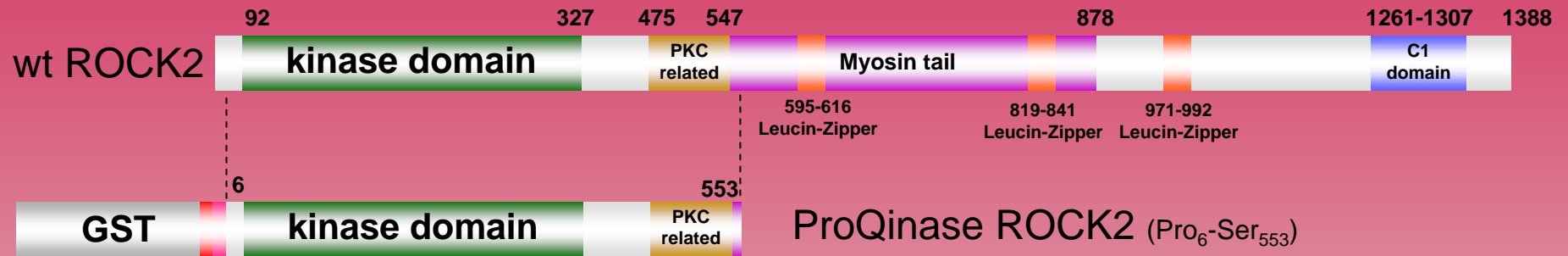


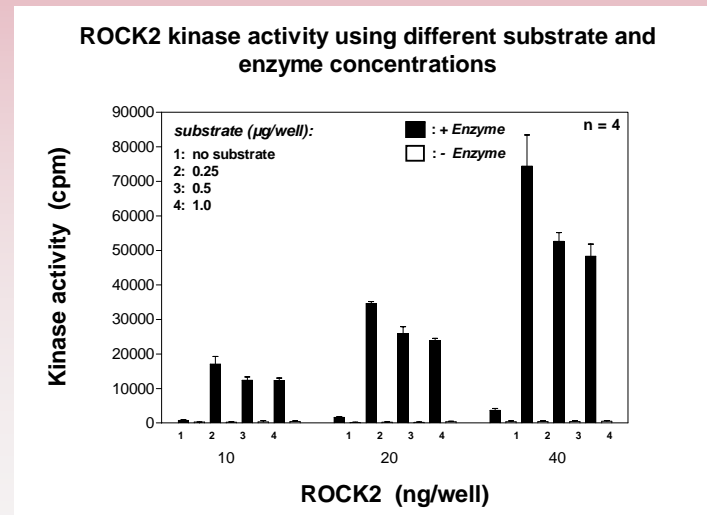
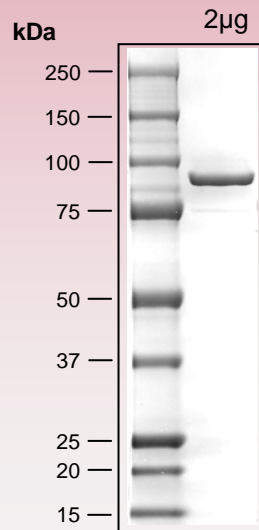
human **ROCK2** kinase (GenBank Acc.No: NM_004850)



ProQinase ROCK2 (Pro₆-Ser₅₅₃)

803 aminoacids, 92.279 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 256-803 ROCK2 6-553

Representative lot: Coomassie stain & activity in PanQinase™ assay



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