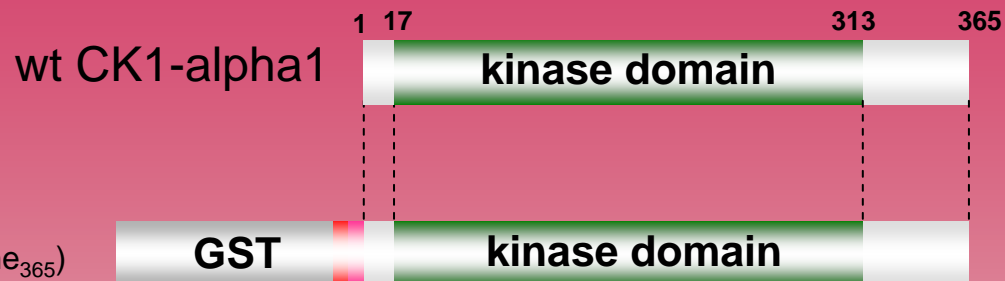


human **CK1-alpha1** kinase (GenBank Acc.No: DQ082865.1)



621 aminoacids, 71.434 kDa

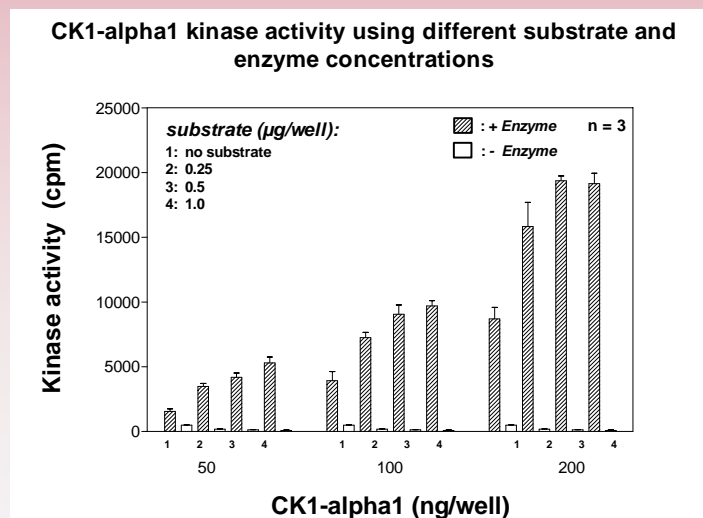
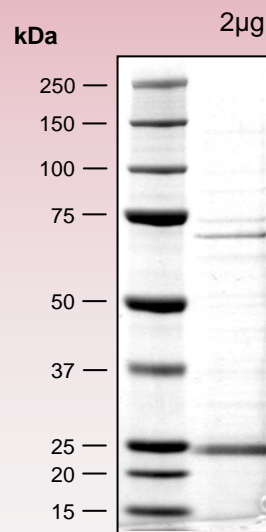
1-218 GST

224-229 HIS-tag

241-246 Thrombin cleavage site

257-621 CK1-alpha1 1-365

Representative lot: Coomassie stain & activity in PanQinase™ assay



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