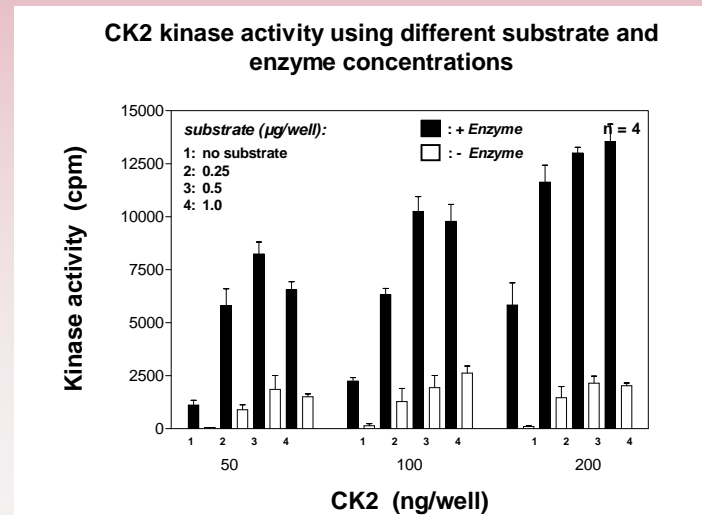
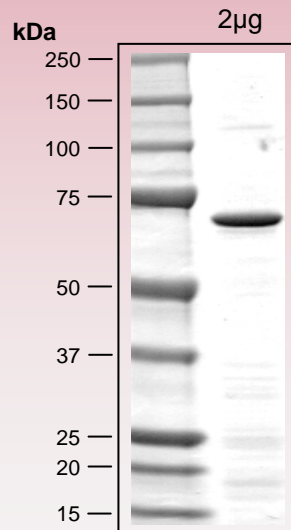


human **CK2-alpha1** kinase (GenBank Acc.No: NM_001895)



612 aminoacids, 70.924 kDa
 1-218 GST
 222-612 CK2-alpha1 1-391

Representative lot: Coomassie stain & activity in PanQinase™ assay



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