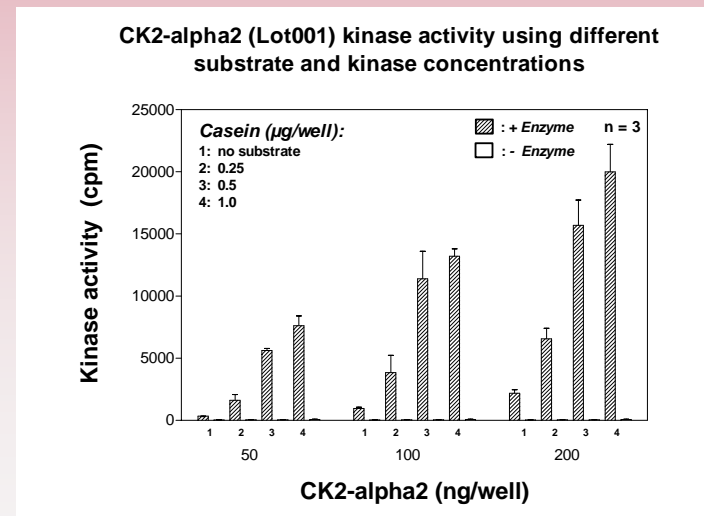
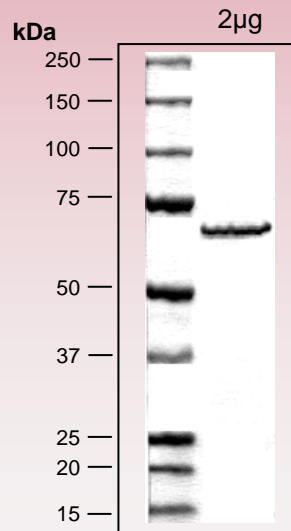


human **CK2-alpha2** kinase (GenBank Acc.No: NM_001896.2)



610 aminoacids, 71.108 kDa
1-218 GST
260-610 CK2-alpha1 1-350

Representative lot: Coomassie stain & activity in PanQinase™ assay



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