

# human **GSK3-beta** kinase (GenBank Acc.No: L33801)



## ProQinase GSK3-beta (Met<sub>1</sub>-Thr<sub>420</sub>)

676 aminoacids, 76.312 kDa

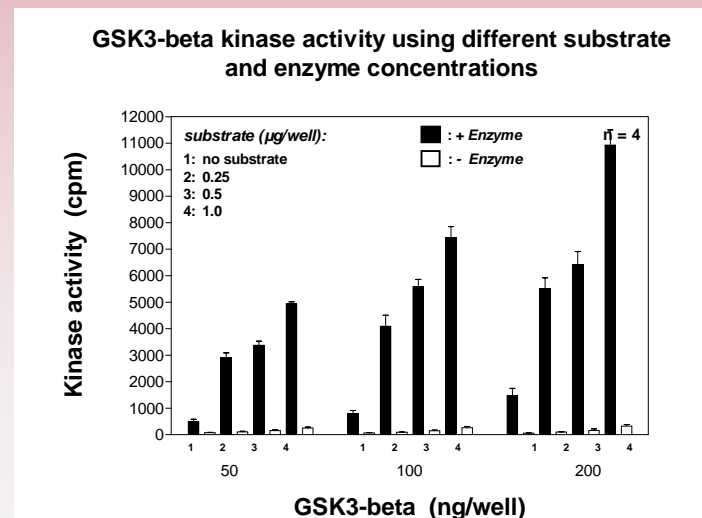
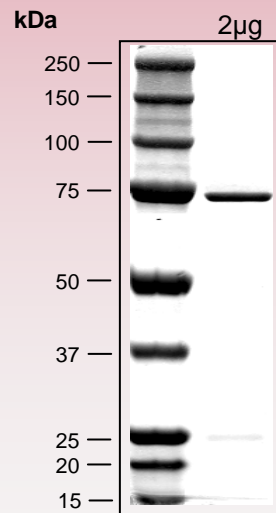
1-218 GST

224-229 HIS-tag

221-226 Thrombin cleavage site

257-676 GSK3-beta 1-420

## Representative lot: Coomassie stain & activity in PanQinase™ assay



Lot specific datasheet  
available on request.  
please contact:  
[info@proqinase.com](mailto:info@proqinase.com)