

# human p38-beta kinase (GenBank Acc.No: NM\_002751)

wt p38-beta



ProQinase p38-beta (Met<sub>1</sub>-Gln<sub>364</sub>)

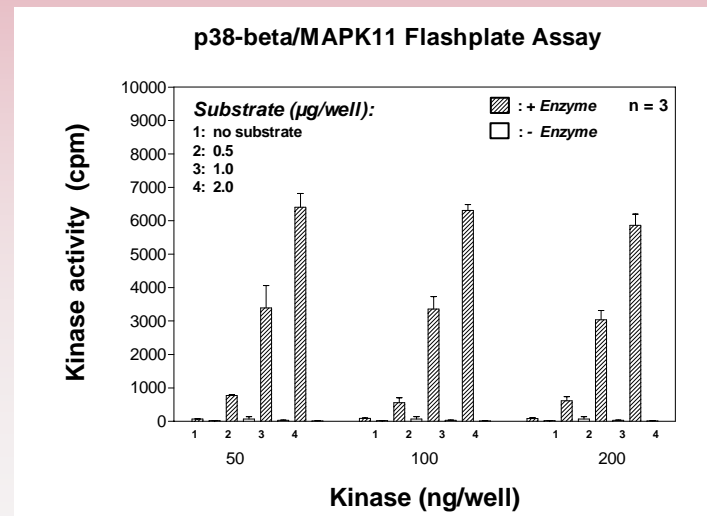
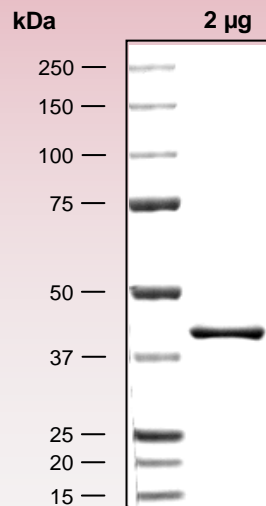


366 aminoacids, 41.485 kDa

3-366

p38-beta 1-364

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



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