

human WEE1 kinase (GenBank Acc.No: X62048)

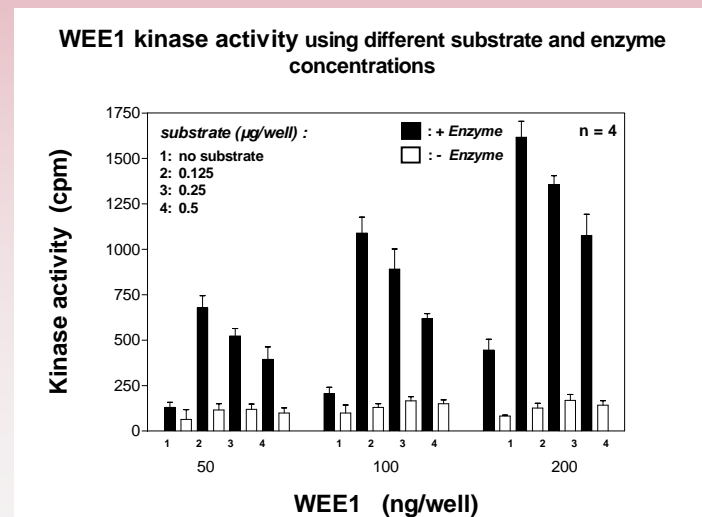
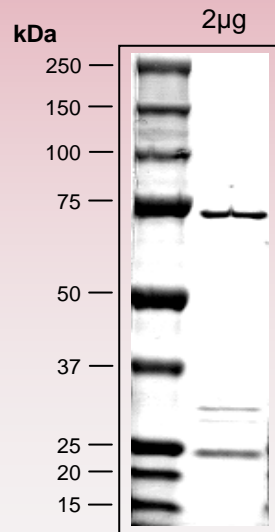


ProQinase WEE1 (Gln₂₅₀-Tyr₆₄₆)

655 aminoacids, 74.902 kDa
 1-218 GST
 224-228 HIS-tag
 241-246 Thrombin cleavage site
 258-655 WEE1 250-646



Representative lot: Coomassie stain & activity in PanQinase™ assay



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