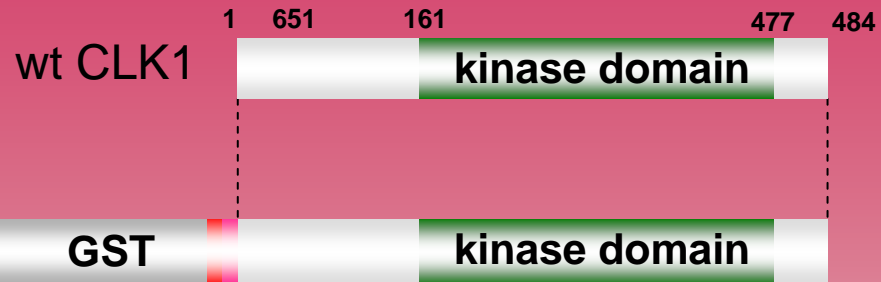


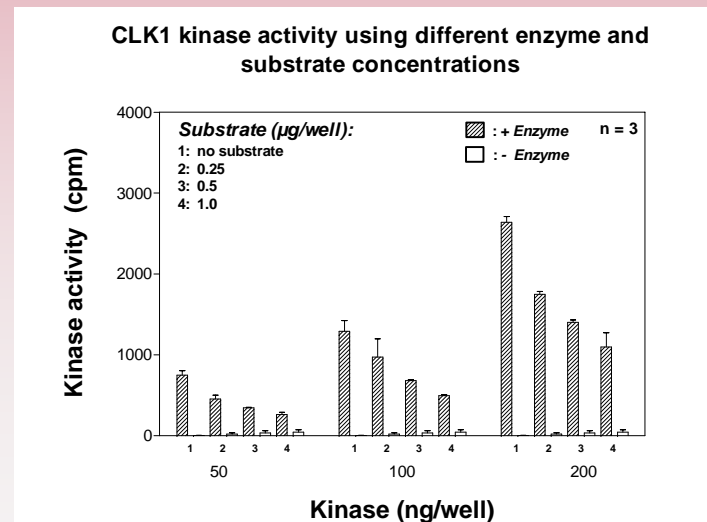
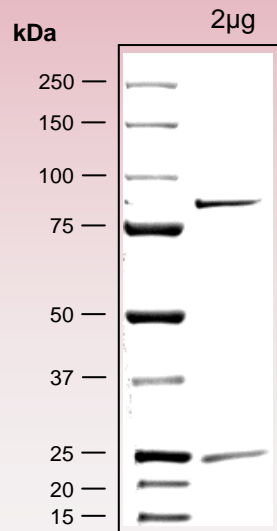
human **CLK1** kinase (GenBank Acc.No: NM_004071.2)



ProQinase CLK1 (Met₁-Ile₄₈₄)

739 aminoacids, 86.695 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 256-739 CLK1 1-484

Representative lot: Coomassie stain & activity in PanQinase™ assay



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