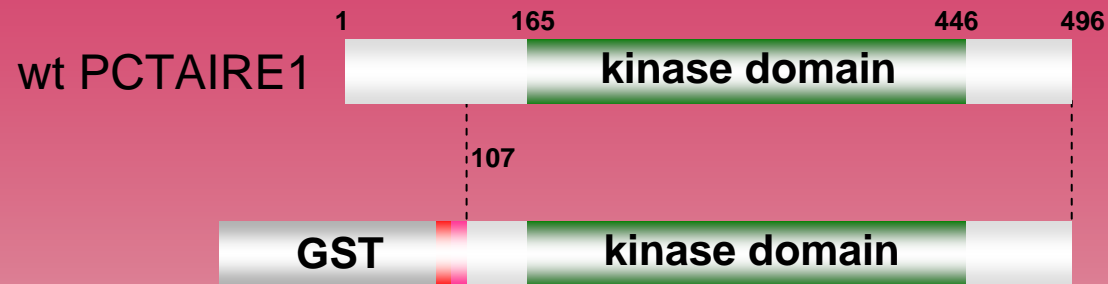


# human PCTAIRE1 kinase (GenBank Acc.No: NM\_006201)



ProQinase PCTAIRE1 (Arg<sub>107</sub>-Phe<sub>496</sub>)

654 aminoacids, 74.926 kDa

1-218

GST

224-229

HIS-tag

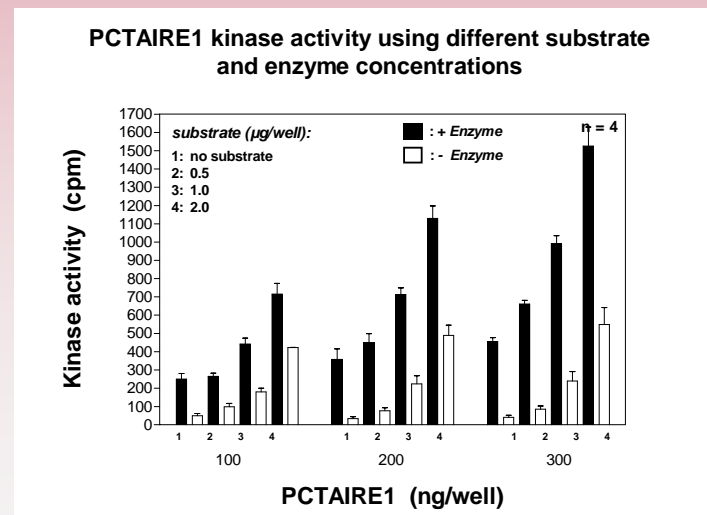
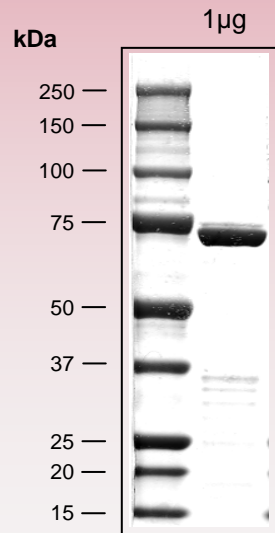
241-246

Thrombin cleavage site

263-654

PCTAIRE1 107-496

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



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