

# human EGF-R kinase (GenBank Acc.No: NM\_005228)



SP: signal peptide  
TM: transmembrane domain

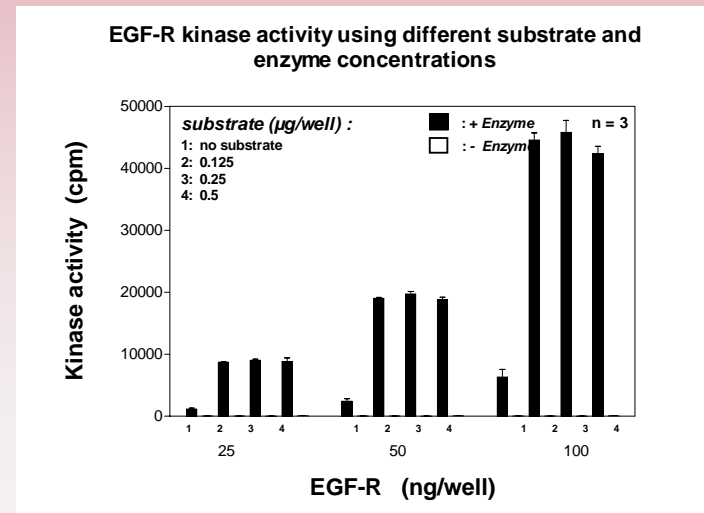
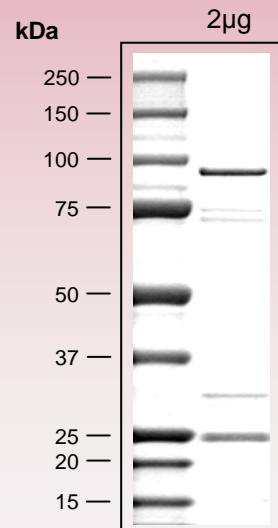
## ProQinase EGF-R (His<sub>672</sub>-Ala<sub>1210</sub>)

801 aminoacids, 91.002 kDa

1-218	GST
224-229	HIS-tag
241-246	Thrombin cleavage site
263-801	EGF-R 672-1210



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



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