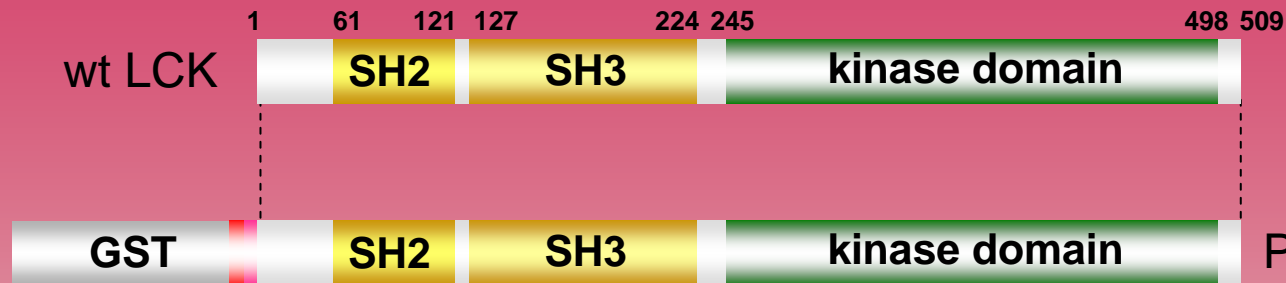


# human LCK kinase (GenBank Acc.No: NM\_005356)



## ProQinase LCK (Met<sub>1</sub>-Pro<sub>509</sub>)

813 aminoacids, 92.792 kDa

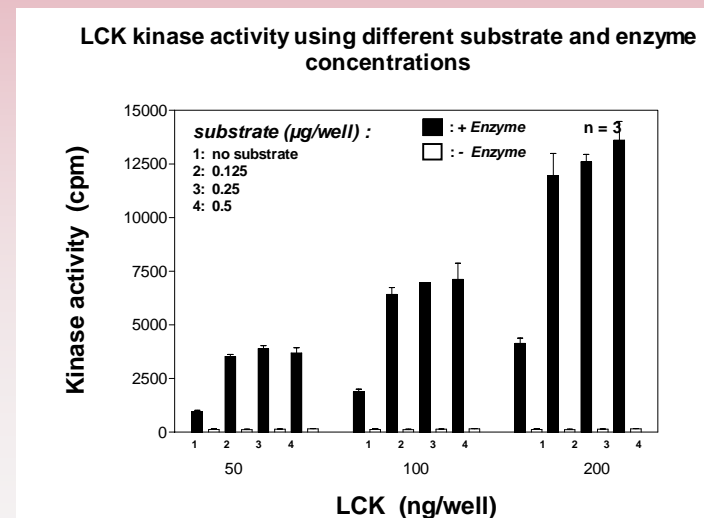
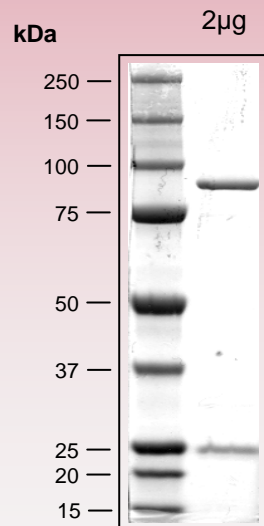
1-218 GST

224-229 HIS-tag

241-246 Thrombin cleavage site

278-786 LCK 1-509

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



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