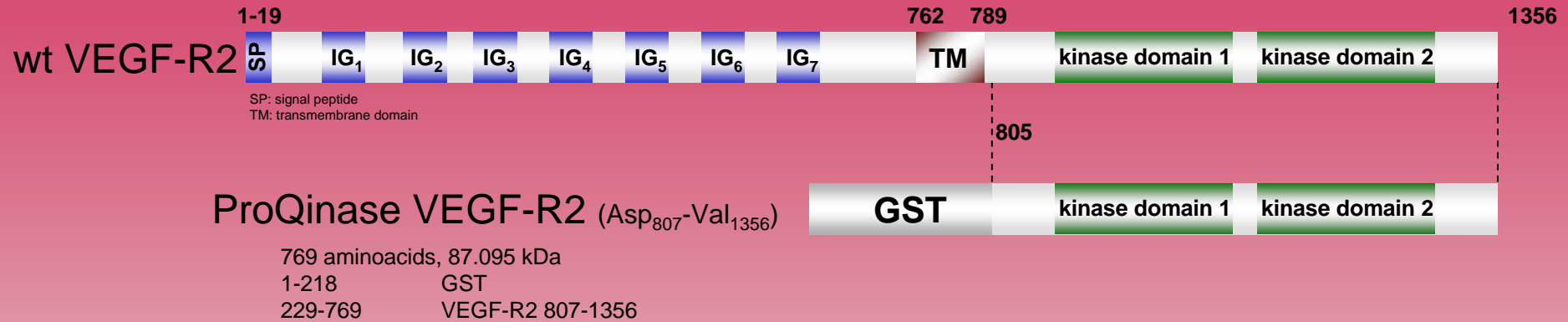
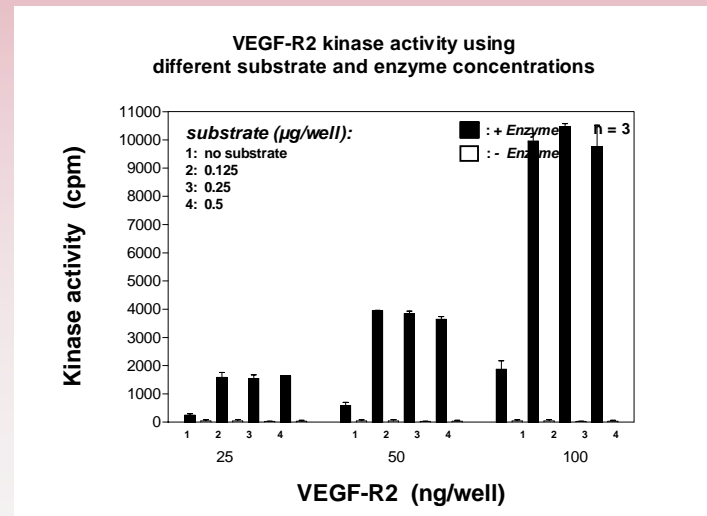
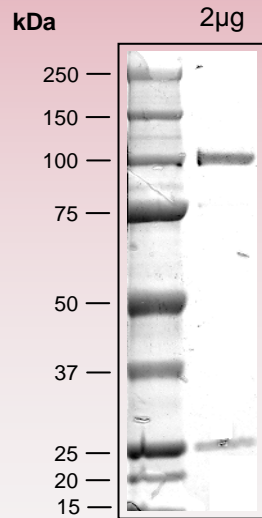


# human VEGF-R2 kinase (GenBank Acc.No: NM\_002253)



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



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