

human **TYRO3** kinase (GenBank Acc.No: NM_006293.2)



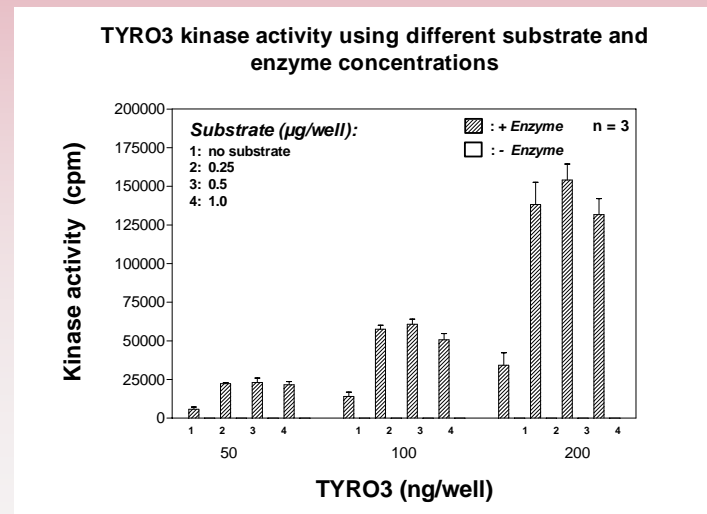
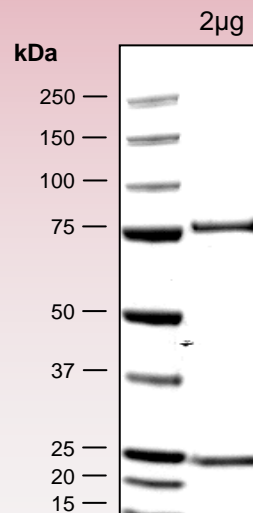
ProQinase TYRO3 (Arg₄₅₂-Cys₈₉₀)

694 aminoacids, 78.384 kDa

1-218 GST
224-229 HIS-tag
241-246 Thrombin cleavage site
256-694 TYRO3 452-890



Representative lot: Coomassie stain & activity in PanQinase™ assay



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