

human VEGF-R3 kinase (GenBank Acc.No: NM_002020)



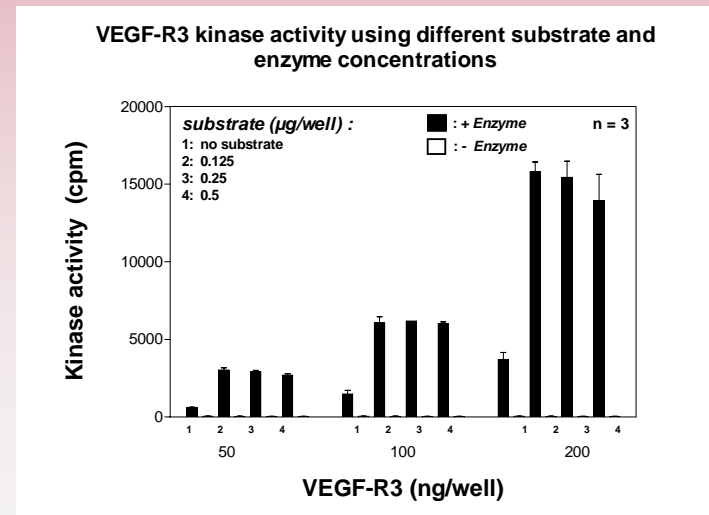
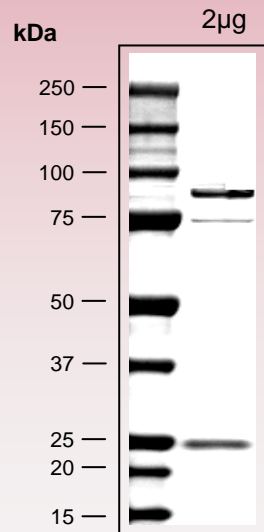
SP: signal peptide
TM: transmembrane domain

ProQinase VEGF-R3 (Asn₇₉₉-Arg₁₂₉₈)

756 aminoacids, 85.565 kDa
 1-218 GST
 224-228 HIS-tag
 241-246 Thrombin cleavage site
 257-783 VEGF-R3 799-1298



Representative lot: Coomassie stain & activity in PanQinase™ assay



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