

human VEGF-R1 kinase (GenBank Acc.No: NM_002019)



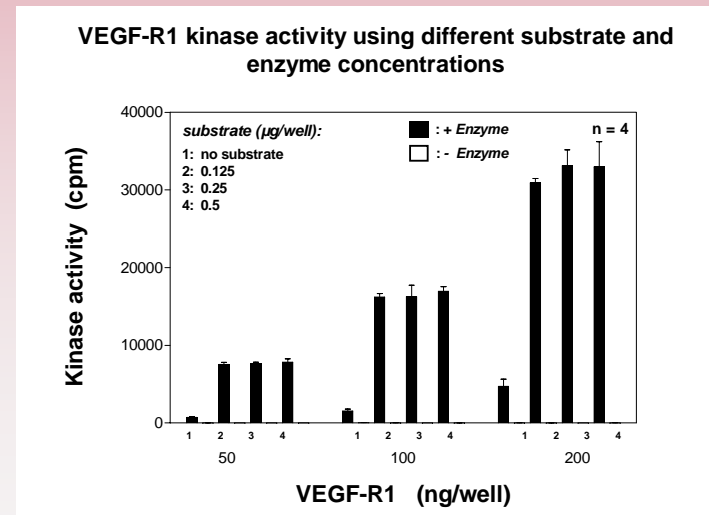
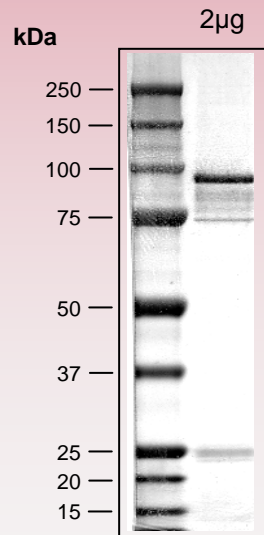
SP: signal peptide
TM: transmembrane domain

ProQinase VEGF-R1 (Lys₇₈₄-Ile₁₃₃₈)

783 aminoacids, 89.357 kDa
1-218 GST
222-225 Factor Xa cleavage site
229-783 VEGF-R1 784-1338



Representative lot: Coomassie stain & activity in PanQinase™ assay



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