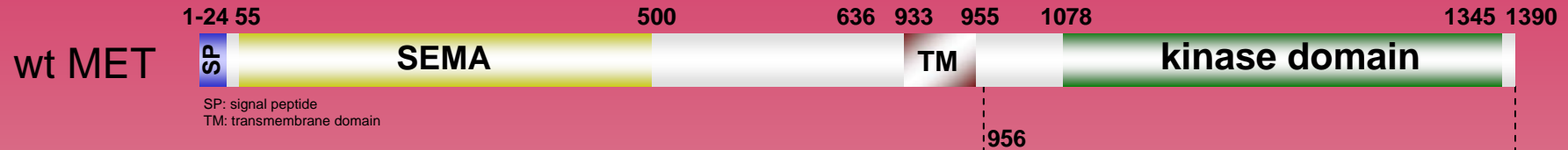


human MET kinase (GenBank Acc.No: NM_000245)



ProQinase MET (Lys₉₅₆-Ser₁₃₉₀)

693 aminoacids, 78.788 kDa

1-218 GST

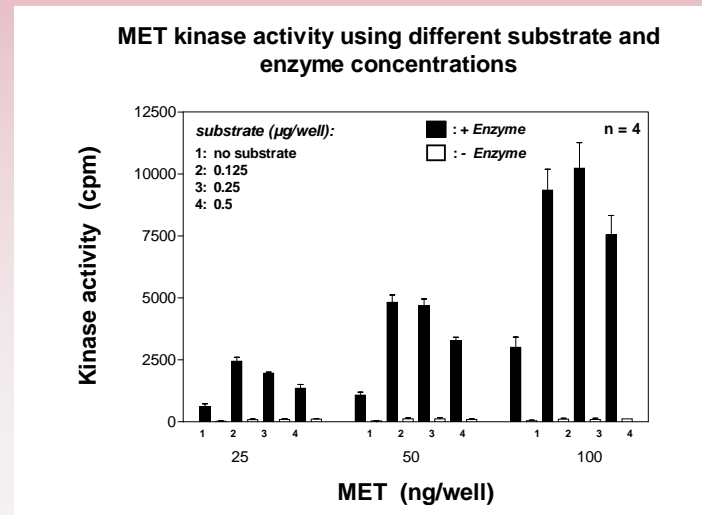
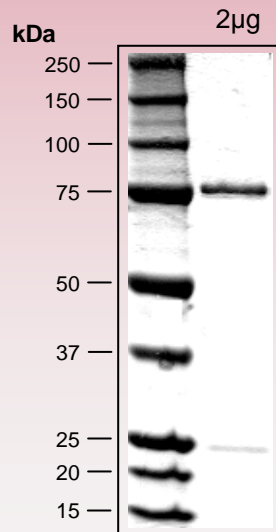
224-229 HIS-tag

241-246 Thrombin cleavage site

259-693 MET 956-1390



Representative lot: Coomassie stain & activity in PanQinase™ assay



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