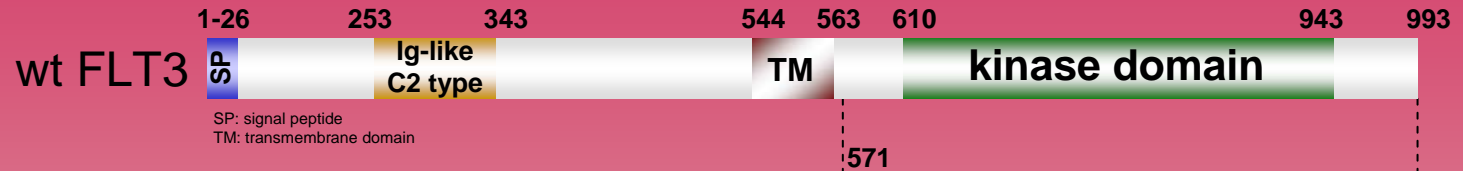


human **FLT3** kinase (GenBank Acc.No: NM_004119)

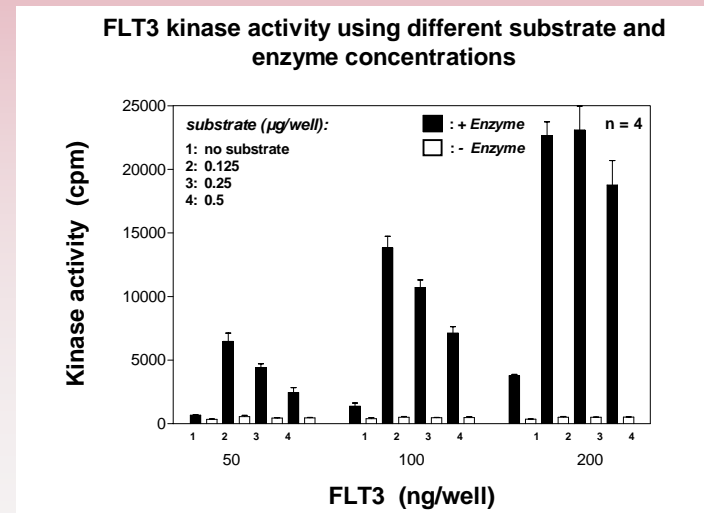
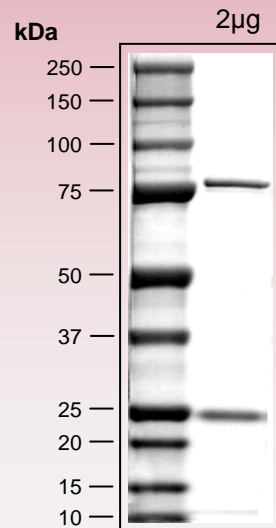


ProQinase FLT3 (Arg₅₇₁-Ser₉₉₃)

714 aminoacids, 82.076 kDa
 1-218 GST
 224-229 HIS-tag
 241-246 Thrombin cleavage site
 280-714 FLT3 571-993



Representative lot: Coomassie stain & activity in PanQinase™ assay



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