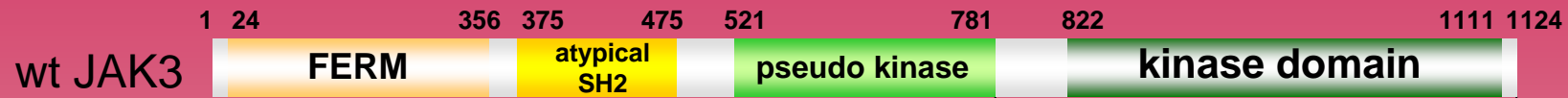


human **JAK3** kinase (GenBank Acc.No: NM_000215)



ProQinase JAK3 (Ile₇₈₁-Ser₁₁₂₄)

601 aminoacids, 68.379 kDa

1-218 GST

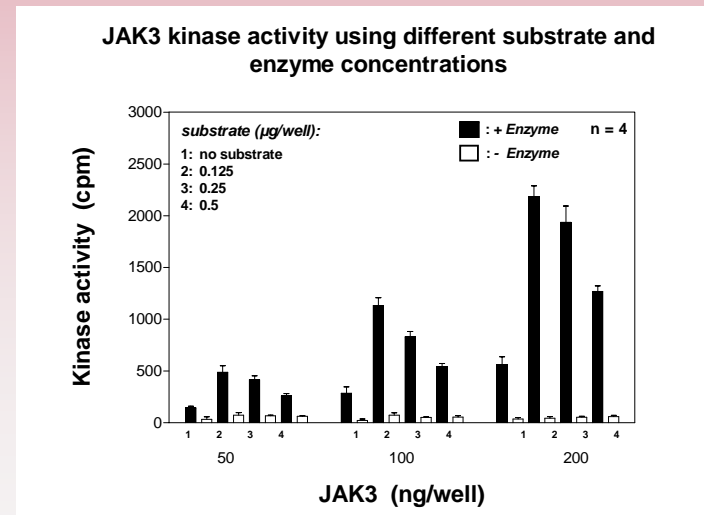
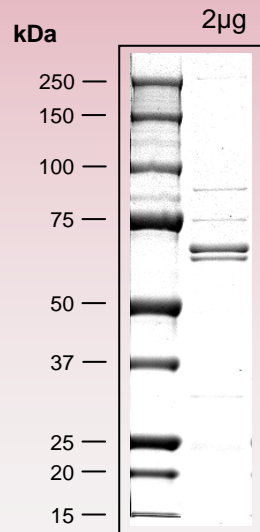
224-229 HIS-tag

241-246 Thrombin cleavage site

258-601 JAK3 781-1124



Representative lot: Coomassie stain & activity in PanQinase™ assay



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