

human **ABL1** kinase (GenBank Acc.No: NM_005157)



ProQinase ABL1 (Pro₁₁₈-Ser₅₃₅)

645 aminoacids, 74.070 kDa

1-218

GST

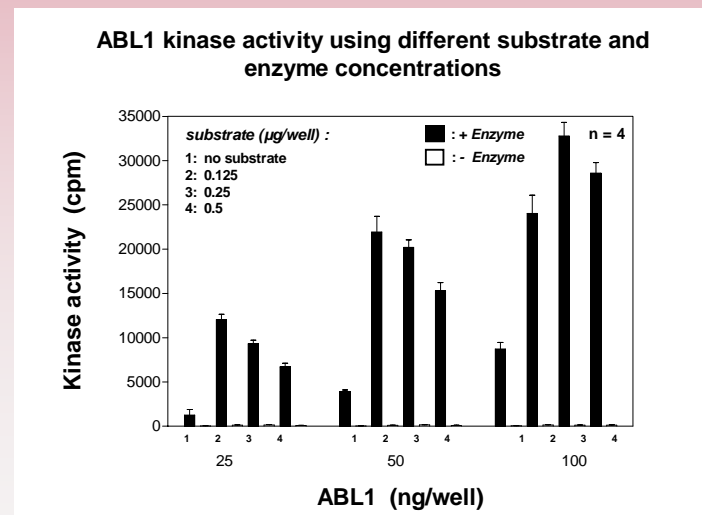
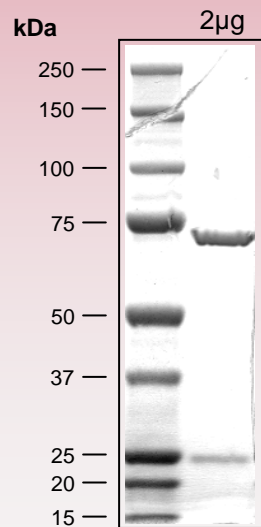
222-226

Factor Xa cleavage site

228-645

ABL1 118-535

Representative lot: Coomassie stain & activity in PanQinase™ assay



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