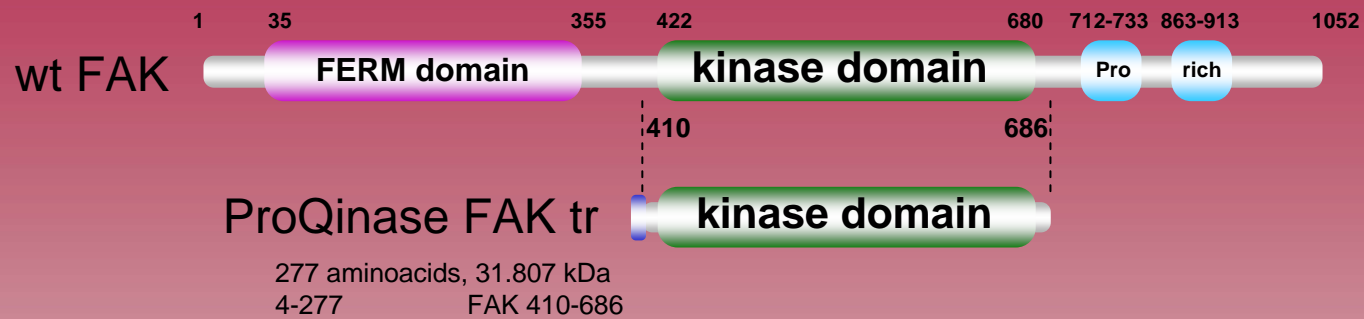
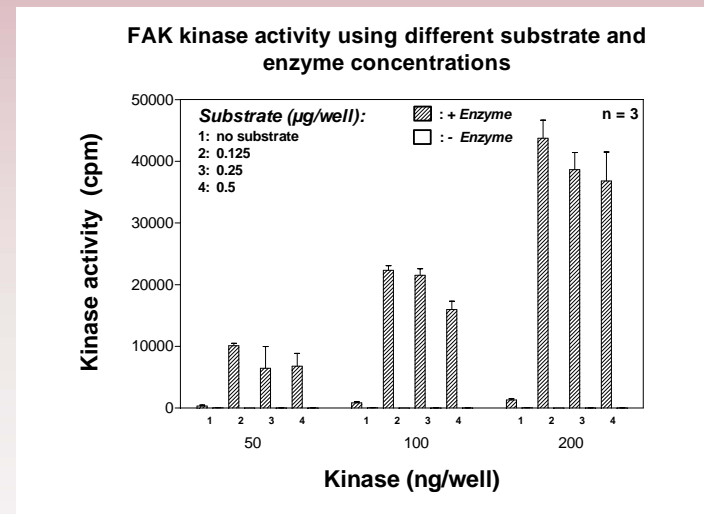
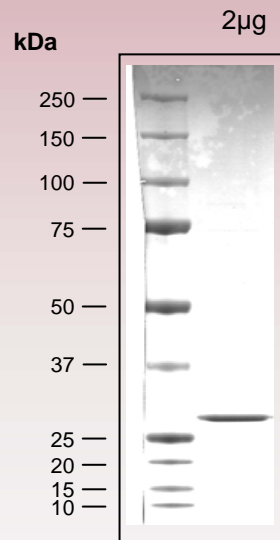


human **FAK tr** kinase (GenBank Acc.No: L13616)



Representative lot: Coomassie stain & activity in PanQinase™ assay



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