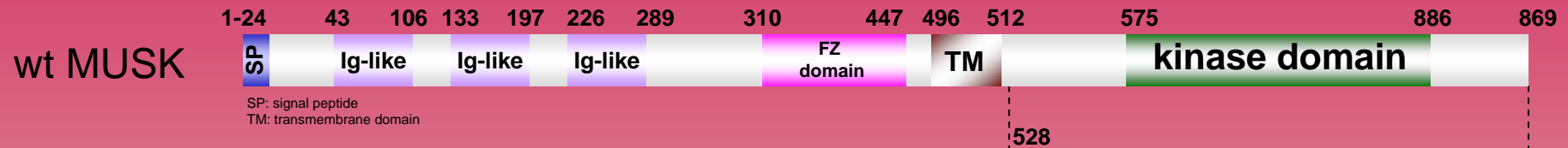


human MUSK kinase (GenBank Acc.No: NM_005592)



ProQinase MUSK (Arg₅₂₈-Val₈₆₉)

598 aminoacids, 68.345 kDa

1-218

GST

224-229

HIS-tag

241-246

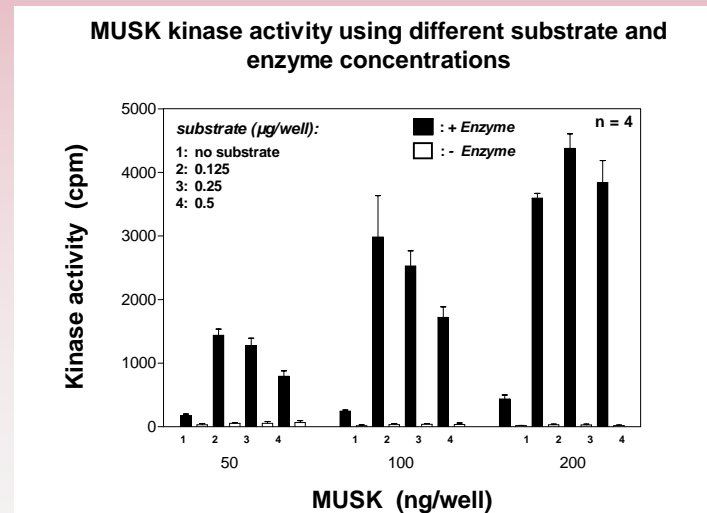
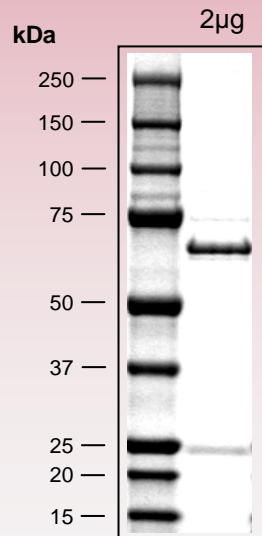
Thrombin cleavage site

258-598

MUSK 528-869



Representative lot: Coomassie stain & activity in PanQinase™ assay



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