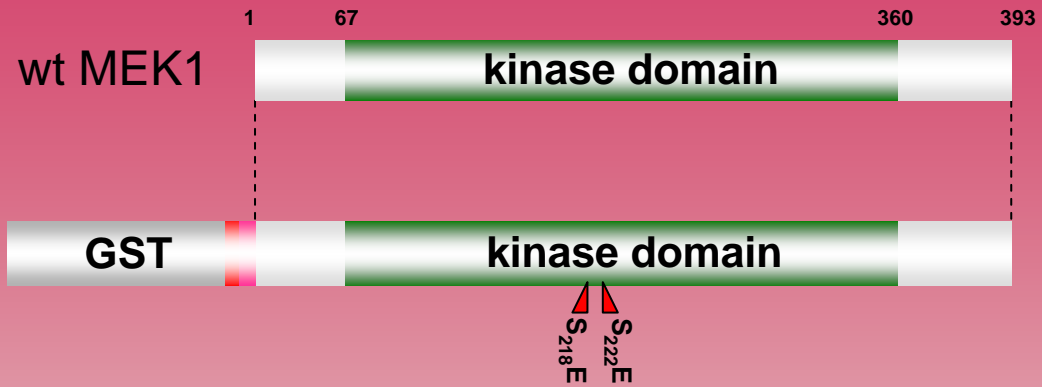
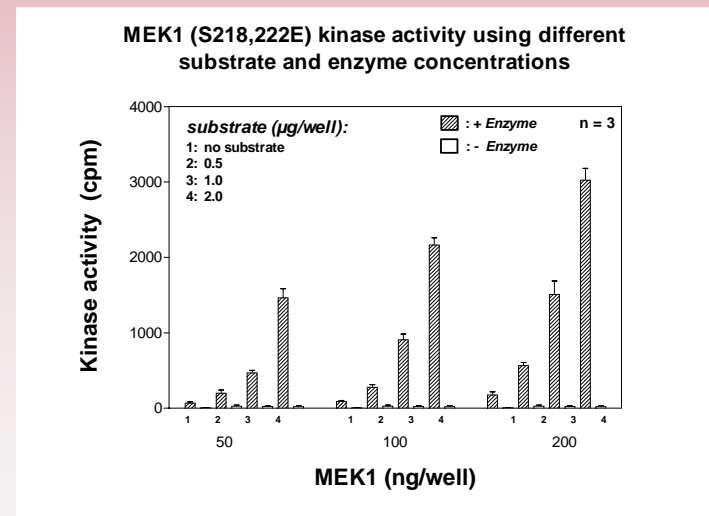
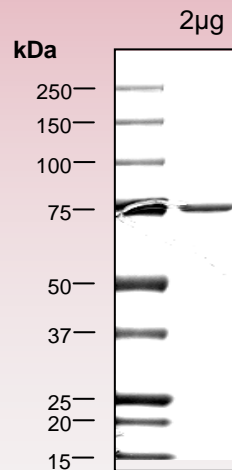


human MEK1 SESE kinase (S_{218,222}E; GenBank Acc.No: NM_002755)



Representative lot: Coomassie stain & activity in PanQinase™ assay



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