

Product Information

Erk5(MAPK7)[inactive mutant]

Product Number : **04-145-11**

Product description

Human Erk5, catalytic domain [1-398 amino acids and K84R of accession number NP_002740] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system. GST-Erk5 was purified by using glutathione sepharose chromatography.

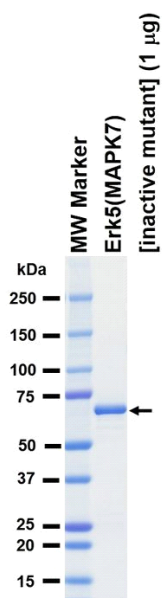
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

Store at -80C.
Avoid repeating freeze-thaws.

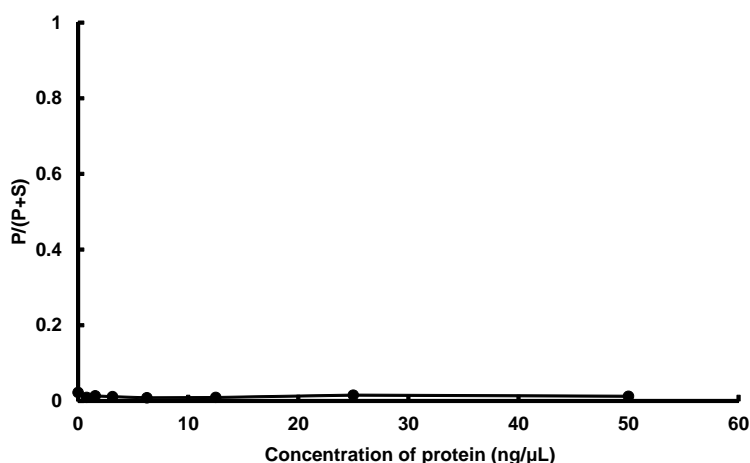
SDS-PAGE



Purity: 93 %

The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was measured by off-chip mobility shift assay. The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by LabChip™3000.

Substrate : EGFR-derived peptide

ATP : 1000 µM