

Product Information

Erk5(MAPK7)[inactive]

 Product Number : **04-146-10**

Product description

Human Erk5, catalytic domain [1-398 amino acids of accession number NP_002740.2] was expressed as N-terminal GST-fusion protein (72 kDa) using E.coli expression system. GST-Erk5 was purified by using glutathione sepharose chromatography.

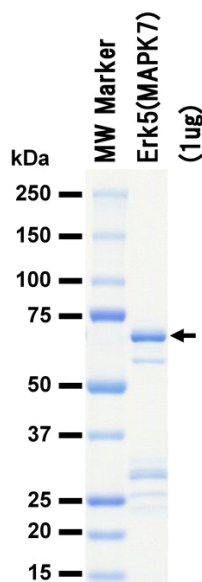
Storage buffer:

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.

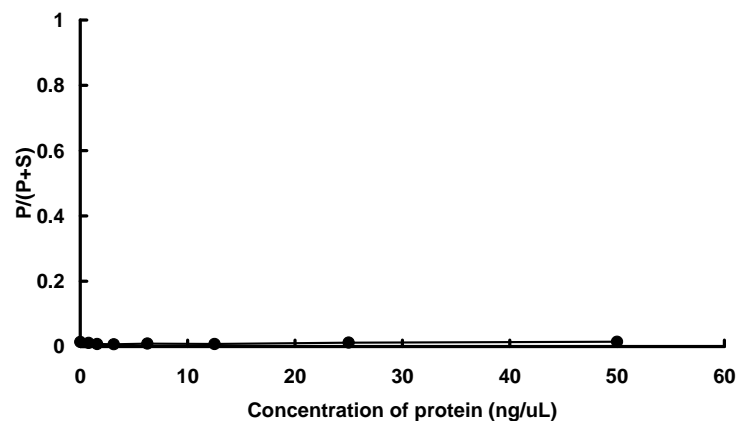
SDS-PAGE



Purity: 55 %

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

Activity data



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

Substrate : EGFR-derived peptide

 ATP : 1000 μ M