

Product Information

Erk2(MAPK1)[inactive mutant]

 Product Number : **04-143-11**

Product description

Full-length human Erk2 [1-360(end) amino acids and K54R of accession number NP_002736.3] was expressed as N-terminal GST-fusion protein (69 kDa) using E.coli expression system. GST-Erk2 was purified by using glutathione sepharose chromatography.

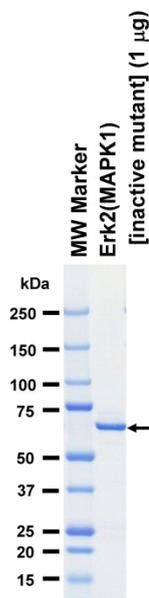
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,
 1 mM DTT, 10% glycerol, pH 7.5

Storage and Handling:

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

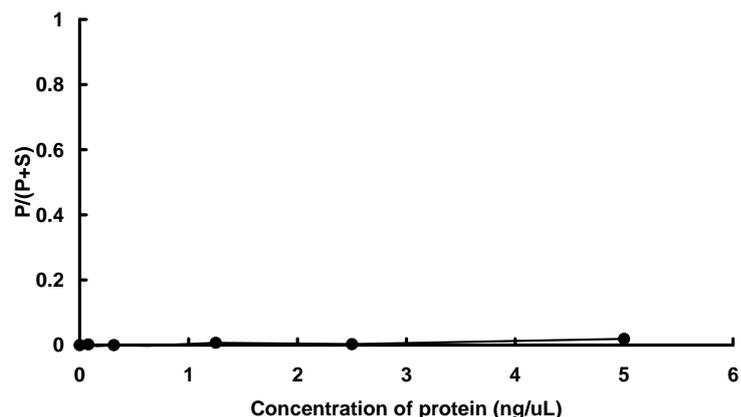
SDS-PAGE



Purity: 96 %

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 : Modified Erktide

ATP : 100 µM