

## Recombinant ERK1 protein

---

**Catalog No:** 31152

**Lot No:** 17207002

**Expressed In:** *E. coli*

**Quantity:** 10 µg

**Concentration:** 0.5 µg/µl

**Source:** Human

**Buffer Contents:** 10 µg of Recombinant ERK1 protein in 50 mM Tris-HCl, 150 mM NaCl, 1 mM DTT, 50% glycerol pH 8.5.

**Protein Details:** Recombinant human ERK1 protein was expressed in *E. coli* with an N-terminal His-Tag. The accession number is NM\_002746. This highly active form was produced by phosphorylation of the purified protein *in vitro* with MEK1 and is suitable for labeling ERK1 substrates. This kinase is phosphorylated at Thr202. Purified by chromatography to > 95% by SDS-PAGE.

**Application Notes:** Recombinant ERK1 is suitable for kinase assays. The molecular weight of the protein is ~44.508 kDa. The activity of the protein is ~ 257,100 units/mg with 1 unit defined as the amount of enzyme that will catalyze the transfer of 1 pmol phosphate to myelin basic protein per minute at 30°C.

Recommended Kinase Reaction Buffer: 20 mM Tris-HCL, 25 mM β-glycerophosphate, 5 mM EGTA, 1 mM sodium orthovanadate and 1 mM DTT, pH 7.5.

Kinase activity may vary depending on the substrate and reaction conditions used.

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.