

ERK1, Unactive

Full-length recombinant protein expressed in *E. coli* cells

Catalog # M29-14G

Lot # T1100-4

Product Description

Recombinant full-length human ERK1 was expressed in *E. coli* cells using an N-terminal GST tag. The gene accession number is [NM_002746](#).

Gene Aliases

PRKM3; P44ERK1; P44MAPK; MAPK3

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

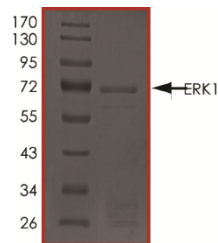
Scientific Background

ERK1 is a protein serine/threonine kinase that is a member of the extracellular signal-regulated kinases (ERKs) which are activated in response to numerous growth factors and cytokines (1). Activation of ERK1 requires both tyrosine and threonine phosphorylation that is mediated by MEK. ERK1 is ubiquitously distributed in tissues with the highest expression in heart, brain and spinal cord. Activated ERK1 translocates into the nucleus where it phosphorylates various transcription factors (e.g., Elk-1, c-Myc, c-Jun, c-Fos, and C/EBP beta).

References

1. Boulton, T G. et al: Purification and properties of extracellular signal-regulated kinase 1, an insulin-stimulated microtubule-associated protein 2 kinase. *Biochemistry*. 1991 Jan 8;30(1):278-86.

Purity



The purity was determined to be **>85%** by densitometry. Approx. MW **70kDa**.

ERK1, Unactive

Full-length recombinant protein expressed in *E. coli* cells

Catalog #	M29-14G
Lot #	T1100-4
Purity	>85%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.