

JNK3, Unactive

Full-length recombinant protein expressed in E. coli cells

Catalog # M35-14G

Lot # K352-1

Product Description

Recombinant full-length human JNK3 was expressed in E.coli cells using an N-terminal GST tag. The gene accession number is [NM_002753](#).

Gene Aliases

MAPK10; JNK3A, MGC50974, PRKM10, p493F12, p54bSAPK; FLJ12099, FLJ33785

Formulation

Recombinant protein is stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 0.25mM DTT, 0.1mM EGTA, 0.1mM EDTA, 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

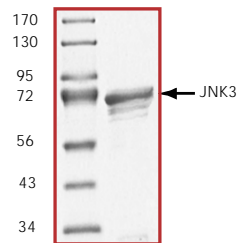
JNK3 is a member of the c-Jun N-terminal kinases (JNKs) which are part of the mitogen-activated protein (MAP) kinase family, and regulate signal transduction in response to environmental stress. JNK3 phosphorylates various transcription factors such as ATF2, Elk-1 and members of the Jun family (1). Activation and nuclear localization of JNK3, a neuronal-specific isoform of JNK, has been associated with hypoxic and ischemic damage of CA1 neurons in the hippocampus. Knockout mice lacking JNK3 showed reduced apoptosis of hippocampal neurons and reduced seizure induced by kainic acid, a glutamate-receptor agonist (2).

References

- Gupta, S. et al: Selective interaction of JNK protein kinase isoforms with transcription factors. EMBO J. 1996 Jun 3;15(11):2760-70.
- Yang, D D. et al: Absence of excitotoxicity-induced apoptosis in the hippocampus of mice lacking the Jnk3 gene. Nature. 1997 Oct 23;389(6653):865-70.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

Purity



The purity was determined to be **>75%** by densitometry.
Approx. MW **~71kDa**.

JNK3, Unactive

Full-length recombinant protein expressed in E. coli cells

Catalog Number M35-14G

Specific Lot Number K352-1

Purity	>75%
Concentration	0.2 $\mu\text{g}/\mu\text{l}$
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.