

JNK2, Unactive

Full-length recombinant protein expressed in E. coli cells

Catalog # M34-14G

Lot # K064-1

Product Description

Recombinant full-length rat JNK2 was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM_017322](#).

Gene Aliases

SAPK; MAPK9

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.

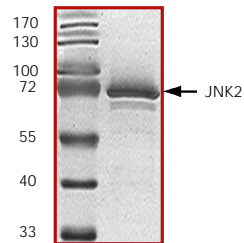
Scientific Background

JNK2 is a member of the MAP kinase group that is activated by dual phosphorylation at thr and tyr residues during exposure to stress such as UV irradiation and treatment of cells with $\text{TNF}\alpha$. JNK2 binds to the c-Jun transactivation domain and phosphorylates it on Ser-63 and Ser-73 (1). JNK2 has been shown to play an important role in disease processes. JNK2 plays a role in ventricular hypertrophy and is thought to be involved in hypertensive cardiac disease (2).

References

1. Sluss, H K. et al: Signal transduction by tumor necrosis factor mediated by JNK protein kinases. *Mol Cell Biol.* 1994 Dec;14(12):8376-84.
2. Vogel, V. et al: Cardiac hypertrophy in the Prague-hypertensive rat is associated with enhanced JNK2 but not ERK tissue activity. *Kidney Blood Press Res.* 2001;24(1):52-6.

Purity



The purity was determined to be **>90%** by densitometry.
Approx. MW **~74kDa**.

JNK2, Unactive

Full-length recombinant protein expressed in E. coli cells

Catalog Number M34-14G

Specific Lot Number K064-1

Purity >90%

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.

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

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.