

JNK1, Unactive

Full-length recombinant protein expressed in E. coli cells

Catalog # M33-14G

Lot # K089-2

Product Description

Recombinant full-length mouse JNK1 was expressed in E.coli cells using an N-terminal GST tag. The gene accession number is [NM_016700](#).

Gene Aliases

JNK; JNK1; PRKM8; SAPK1; JNK1A2; JNK21B1/2

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.

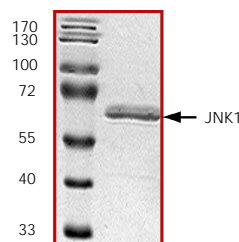
Scientific Background

JNK1 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. JNK1 binds to the c-Jun transactivation domain and phosphorylates it on Ser-63 and Ser-73 (1). JNK1 has been shown to play an important role in disease processes. Activation of JNK1 results in defects in myotube viability and integrity leading to dystrophic myofiber destruction (2). JNK1 activity is also abnormally elevated in obesity and removal of JNK1 results in decreased adiposity and significantly improved insulin sensitivity.

References

1. Derijard, B. et al: JNK1: a protein kinase stimulated by UV light and Ha-Ras that binds and phosphorylates the c-Jun activation domain. *Cell*. 1994 Mar 25;76(6):1025-37.
2. Kolodziejczyk, S M. et al: Activation of JNK1 contributes to dystrophic muscle pathogenesis. *Curr Biol*. 2001 Aug 21;11(16):1278-82.

Purity



The purity was determined to be **>90%** by densitometry.
Approx. MW **~70kDa**.

JNK1, Unactive

Full-length recombinant protein expressed in E. coli cells

Catalog Number M33-14G

Specific Lot Number K089-2

Purity	>90%
Concentration	0.2 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.

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.