

Anti-JNK2

Mouse Monoclonal Antibody

Catalog # M34-60M

Lot # V336-2

Cited Applications

- Immunoprecipitation (2 µg)
- Western blot (1:1000 dilution)

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes only the JNK2 protein

Cross Reactivity

- Immunoprecipitation of human JNK2
- Western blot of human JNK2

JNK2 from other species may also be detectable

Host/Isotype/Clone#

Mouse, IgG2, 2G12

Immunogen

Human recombinant full length JNK2 protein

Formulation

10mM PBS pH 7.2

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. 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

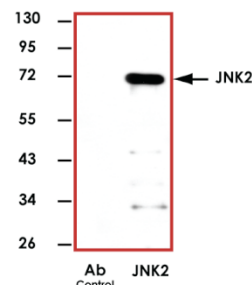
JNK2 is a member of the MAP kinase group that is activated by dual phosphorylation at thr and tyr residues during exposure to stress such as UV irradiation and treatment of cells with TNF α . JNK2 binds to the c-Jun transactivation domain and phosphorylates it on Ser-63 and Ser-73 (1). JNK2 has been 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

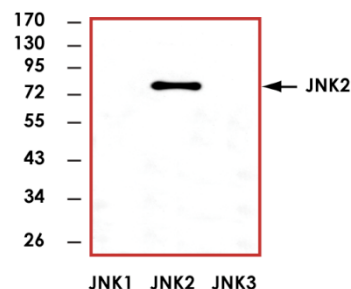
1. Sluss, H K. et al: Signal transduction by tumor necrosis factor mediated by JNK protein kinases. Mol Cell Biol. 1994 Dec;14(12):8376-84.
2. Vogel, V. et al: Cardiac hypertrophy in the Prague-hypertensive rat is associated with enhanced JNK2 but not ERK tissue activity. Kidney Blood Press Res. 2001;24(1):52-6.

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

Sample Data



Representative immunoprecipitation with Anti-JNK2 (2µg) using 100 ng of JNK2 recombinant protein.



Representative western blot with Anti-JNK2 (1:1000) using 20 ng of JNK1, JNK2, and JNK3 recombinant protein.

Anti-JNK2

Mouse Monoclonal Antibody

Catalog Number

M34-60M

Specific Lot Number

V336-2

Purification
Concentration

Affinity purified using Protein G
0.89 µg/µL

Stability

1 yr At -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody 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 ice packs.