

## Anti-JNK1

Rabbit Polyclonal Antibody

**Catalog # M33-63R**

Lot # D316-4

### Cited Applications

- Western blot (1:250)

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

### Specificity

Recognizes the JNK1 protein

### Cross Reactivity

- Western blot JNK1 from human, mouse, and rat cells

*JNK1 from other species may also be detectable*

### Host

Rabbit

### Immunogen

Protein Code: P45983

### Formulation

TBS pH 7.2, 50% glycerol

### Stability

Store at 4°C (add 0.1% NaN<sub>3</sub>) 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 optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

### Scientific Background

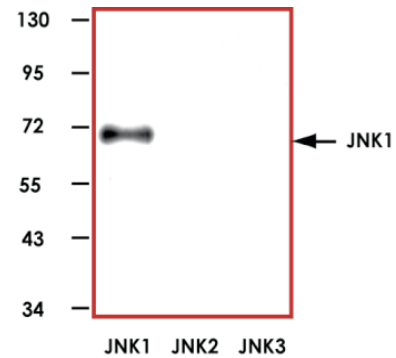
JNK1 is a member of the MAP kinase group that is activated by dual phosphorylation at thr and tyr residues during exposure to stress such as UV irradiation. JNK1 binds to the c-Jun transactivation domain and phosphorylates it on Ser-63 and Ser-73 (1). JNK1 has been shown to play an important role in 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. Derjard, 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.

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

### Sample Data



Representative western blot with Anti-JNK1 (1:250) using 20 ng of JNK1, JNK2, and JNK3 recombinant protein.

## Anti-JNK1

Rabbit Polyclonal Antibody

Catalog Number

M33-63R

Specific Lot Number

D316-4

Purification

Affinity purified by the peptide antigen

Concentration

0.25 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 optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.