

Anti-c-JUN

Rabbit Polyclonal Antibody

Catalog # C05-63R

Lot # M191-6

Cited Applications

- Western blot (1:1000)

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

Specificity

Recognizes the c-JUN protein

Cross Reactivity

- Western blot human and mouse c-JUN

c-JUN from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 71-75 of human c-JUN

Formulation

PBS, pH7.4, 0.02%NaN₃, 50% glycerol

Stability

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

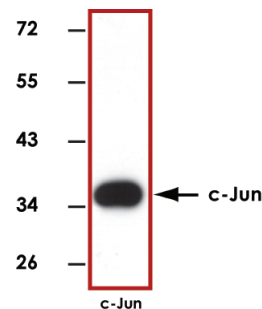
Scientific Background

c-Jun is a proto-oncogene that forms a complex with c-Fos which regulates transcription from promoters containing AP-1 activation elements (1). c-Jun has specific DNA binding activity and following in vitro translation, c-Jun binds as a homodimer to the AP-1 DNA site. The transactivating function of c-Jun is acutely regulated by a wide variety of cellular signals via modulation of phosphorylation of two serines (63 and 73). Jun N-terminal domain kinases (JNKs), are responsible for mediating S63/73 phosphorylation on c-Jun in response to a variety of cellular stimuli including TNF- α , heat stress and u.v. light (2).

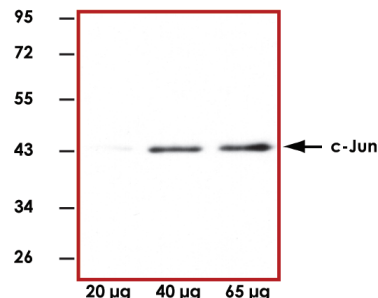
References

1. Halazonetis, T D. et al: c-Jun dimerizes with itself and with c-Fos, forming complexes of different DNA binding affinity. Cell. 1988 Dec 2; 55(5):917-24.
2. Dai, T. et al: Stress-activated protein kinases bind directly to the delta domain of c-Jun in resting cells: implications for repression of c-Jun function. Oncogene. 1995 Mar 2; 10(5):849-55.

Sample Data



Representative Western blot with Anti-c-Jun (1:1000) using 15ng of GST-tagged human recombinant c-JUN.



Representative Western blot with Anti-c-JUN (1:1000) using 20µg, 40µg and 65µg of 3T6 cell lysate.

Anti-c-JUN

Rabbit Polyclonal Antibody

Catalog Number

C05-63R

Specific Lot Number

M191-6

Purification

Affinity chromatography

Concentration

1.0 µg/µL

Stability

1yr 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.

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.