

Anti-ATF2

Rabbit Polyclonal Antibody

Catalog # A10-63R

Lot # M191-5

Cited Applications

- Western blot (1:1000)

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

Specificity

Recognizes the ATF2 protein

Cross Reactivity

- Western blot human and mouse ATF2

ATF2 from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 60-64 of human ATF2

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.

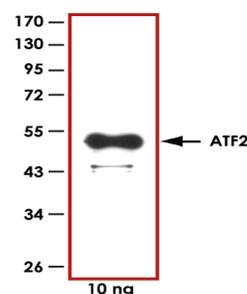
Scientific Background

Activating transcription factor (ATF)-2, also designated CRE-BP1, is a member of the ATF/cyclic AMP-responsive element binding protein family of transcription factors. ATF2 has two unique characteristics: it mediates the adenovirus E1A-induced transactivation and forms a heterodimer with c-Jun (1). ATF2 is a target of the JNK signal transduction pathway (2). ATF2 is phosphorylated by JNK on two closely spaced threonine residues within the NH₂-terminal activation domain. The replacement of these phosphorylation sites with alanine inhibited the transcriptional activity of ATF2. Furthermore, p38 MAPK can also phosphorylate ATF2.

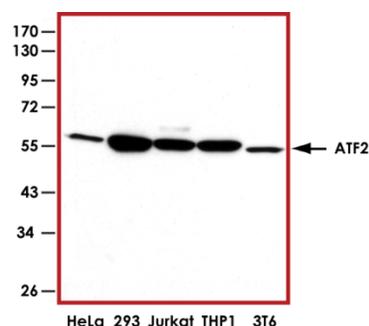
References

1. Nomura, N. et al: Isolation and characterization of a novel member of the gene family encoding the cAMP response element-binding protein CRE-BP1. *J Biol Chem.* 1993 Feb 25; 268(6):4259-66.
2. Gupta, S. et al: Transcription factor ATF2 regulation by the JNK signal transduction pathway. *Science.* 1995 Jan 20; 267(5196):389-93.

Sample Data



Representative Western blot with Anti-ATF2 (1:1000) using 10ng of GST-tagged human recombinant ATF2.



Representative Western blot with Anti-ATF2 (1:1000) using 30 µg of HeLa, 293, Jurkat, THP1 and 3T6 cell lysates.

Anti-ATF2

Rabbit Polyclonal Antibody

Catalog Number

A10-63R

Specific Lot Number

M191-5

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.