

## Anti-phospho-ATF2 (Ser112)

Rabbit Polyclonal Antibody

**Catalog # A10-565R**

Lot # J1178-17

### Cited Applications

IHC

*Suggested Dilutions:*

IHC: 1:50-1:100

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

### Specificity

Recognizes the ATF2 protein phosphorylated at serine 112

### Cross Reactivity

Human and Mouse

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser112

### Formulation

PBS (pH 7.4) 150mM NaCl, 0.02% sodium azide and 50% glycerol.

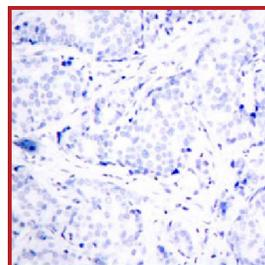
### Scientific Background

The activating transcription factor ATF2 (also known CREBP1) is a member of the ATF/CREB family of leucine zipper transcription factors. These bind both AP-1 and CRE DNA response elements (1). ATF2 has two unique characteristics: it mediates adenovirus E1A-induced trans-activation and forms a heterodimer with c-Jun (2). ATF2 is a target of the JNK signal transduction pathway (3). ATF2 is phosphorylated by JNK at two closely spaced threonine residues located in the NH<sub>2</sub>-terminal activation domain. The replacement of these phosphorylation sites with alanine inhibited the transcriptional activity of ATF2. Furthermore, ATF2 is a target for p38 MAPK phosphorylation.

### References

1. Maekawa, T. et al: Leucine zipper structure of the protein CRE-BP1 binding to the cyclic AMP response element in brain. EMBO J 1989 8:2023-2028.
2. 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.
3. Gupta, S. et al: Transcription factor ATF2 regulation by the JNK signal transduction pathway. Science, 1995 Jan 20;267(5196):389-93.

### Sample Data



Immunohistochemistry image of paraffin-embedded human breast carcinoma tissue using the anti-phospho-ATF (Ser112) antibody.

## Anti-phospho-ATF2 (Ser112)

Rabbit Polyclonal Antibody

Catalog Number

A10-565R

Specific Lot Number

J1178-17

Purification

Affinity chromatography

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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**