

ATF-2 phospho Thr71 antibody (mAb)

Catalog No: 40975

RRID: AB_2793467

Clone: 103C411

Isotype: IgG1

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Concentration: 0.5 µg/µl

Molecular Weight: 61 kDa

Background: ATF-2 (activating transcription factor 2) transcription factor is a member of the leucine-zipper family of proteins and was previously known as cAMP responsive element binding protein 2 (or CREB2). ATF 2 is a transcriptional activator that binds the cAMP-responsive element (CRE) and interacts with c-Jun to stimulate CRE transcription. This protein is also capable of acetylating Histone H2B and Histone H4 *in vitro* as a histone acetyltransferase (HAT).

Immunogen: This ATF-2 phospho Thr71 antibody was raised against a synthetic peptide containing phosphorylated threonine at position 71 corresponding to human Phosphorylated-ATF-2.

Buffer: PBS containing 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

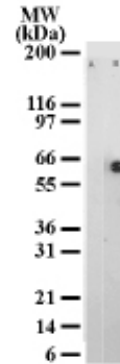
Applications Validated by Active Motif:

WB: 1 - 2 µg/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



ATF-2 phospho Thr71 mAb tested by Western blot.

ATF-2 (phospho) detection by Western blot. Detection of ATF-2 phosphorylation in HeLa cell extract with ATF-2 phospho Thr71 mAb.

Lane A: Untreated HeLa nuclear extract;
Lane B: HeLa nuclear extract treated with TNF-α for 1 hour.