

## ATF-2 antibody (pAb)

**Catalog No:** 40977**RRID:** AB\_2614939**Application(s):** WB**Reactivity:** Human**Quantity:** 100 µg**Purification:** Affinity Purified**Host:** Rabbit**Isotype:** IgG**Concentration:** 0.5 µg/µl**Molecular Weight:** 60 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 antibody was raised against a synthetic peptide corresponding to amino acid residues 13-27 of human 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. Store at 4°C for short term. 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 pAb tested by Western blot.**

ATF-2 detection by Western blot. The analysis was performed using 293 nuclear extract and ATF-2 pAb at a 2 µg/ml dilution.