

## CREB antibody (mAb)

**Catalog No:** 40961

**RRID:** AB\_2614962

**Clone:** 82B514

**Isotype:** IgG1

**Application(s):** WB

**Reactivity:** Human

**Quantity:** 100 µg

**Purification:** Affinity Purified

**Host:** Mouse

**Concentration:** 0.5 µg/µl

**Molecular Weight:** 43 kDa

**Background:** CREB (cAMP response element binding) transcription factor binds as a homodimer to specific DNA sequences called cAMP response elements (CRE) and regulates transcription. CREB belongs to the leucine zipper family of DNA binding proteins. CREB induces transcription in response to hormonal stimulation.

**Immunogen:** This CREB antibody was raised against a synthetic peptide corresponding to human CREB.

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

**Application Notes:**

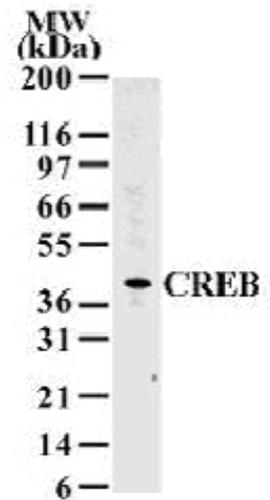
Applications Validated by Active Motif:

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.



**CREB mAb tested by Western blot.**

Detection of CREB by Western blot analysis. A protein band of approximate molecular weight of 43 kDa was detected using HeLa whole-cell extract and CREB mAb at a dilution of 2 µg/ml.