

AP-2 antibody (pAb)

Catalog No: 39001

RRID: AB_2793148

Isotype: IgG

Application(s): ChIP, EMSA, IHC, WB

Reactivity: Human

Volume: 100 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 50 and 40 kDa

Background: AP-2 (activator protein 2 or AP2) transcription factors activate transcription and regulate gene expression during embryogenesis. There are four different isoforms of AP-2 (AP-2 α , AP-2 β , AP-2 γ and AP-2 δ). The family members share a common structure in which the N-terminal region is involved with transcriptional activation and the C-terminal region is involved in DNA binding.

Immunogen: This AP-2 antibody was raised against a synthetic peptide corresponding to amino acid residues 120-134 of human AP-2.

Buffer: PBS containing 0.2% gelatin, 0.01% thimerosal and 0.1% sodium azide. Thimerosal and sodium azide are highly toxic.

Application Notes:

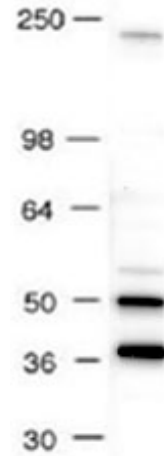
Applications Validated by Active Motif:

WB*: 1:500 - 1:1,000

*For optimal results, primary antibody incubations should be performed overnight at 4°C. 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.



AP-2 pAb tested by Western blot.

Detection of AP-2 by Western blot analysis. Activator protein 2 (AP-2) can be detected in JEG-3 nuclear extract using AP-2 rabbit polyclonal antibody at a 1:1,000 dilution.