

AEBP2 antibody (pAb)

Catalog Nos: 61539, 61540

RRID: AB_2614969

Isotype: IgG

Application(s): WB

Reactivity: Mouse

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 52 kDa

Background: **AEBP2** (AE binding protein 2) is a DNA-binding transcriptional repressor and component of the PRC2 complex, which methylates 'Lys-27' residues of Histone H3.

Immunogen: This antibody was raised against a peptide within the C-terminal region of mouse AEBP2.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

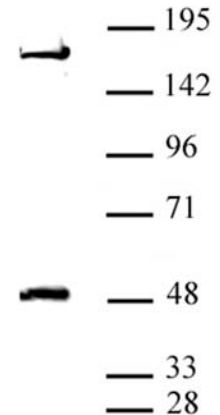
WB*: 1:1,000 - 1:5,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

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.



AEBP2 antibody (pAb) tested by Western blot.

AEBP2 detection by Western blot. The analysis was performed using 20 μ g P19 nuclear cell extract and AEBP2 antibody at a 1:2,500 dilution. An unspecified band is detected around 160 kDa.