

MOZ antibody (pAb)

Catalog No: 39867, 39868

Isotype: IgG

Application(s): ChIP, WB

Reactivity: Human

Purification: Affinity Purified

Host: Rabbit

Concentration: 1 µg/µl

Volume: 100 µl, 10 µl

Molecular Weight: 290 kDa

Background: MOZ (Monocytic leukemia zinc finger protein) / MYST3 (also known KAT6A) is a histone acetyltransferase specific for H3. It is part of the MOZ/MORF protein complex and is involved in the activation of transcription. Specifically, it is known to interact with and serve as a co-activator for RUNX1 and RUNX2. MOZ / MYST3 is found concentrated in sub-nuclear foci that are distinct from PML bodies. Chromosomal aberrations involving MYST3 may be a cause of acute myeloid leukemias. MYST3-NCoA2 has been shown to bind CBP and disrupt its transcriptional activation function.

Immunogen: This MOZ antibody was raised against a synthetic peptide within the N-terminal half of human MOZ (Monocytic leukemia zinc finger protein aka histone acetyltransferase MYST3).

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Validated Applications:

WB: 1:200 - 1:500 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.

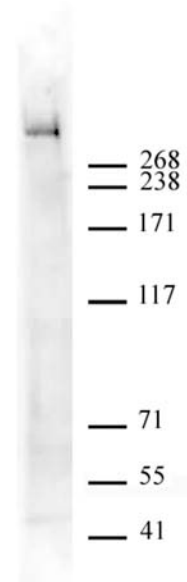
Published Applications:

ChIP

WB

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Upon receipt, unconjugated antibodies may be stored at -20°C for up to 2 years. Fluorophore- & enzyme-conjugated antibodies should be stored at 4°C. Fluorophore-conjugated antibodies should be protected from light. Keep reagents on ice when not in storage; to avoid repeated freeze/thaw cycles, we recommend aliquoting items that will be stored frozen into single-use fractions prior to freezing. This product is guaranteed for 6 months from date of receipt.

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



MOZ antibody (pAb) tested by Western blot.
Nuclear extract of K-562 cells (40 µg) probed with MOZ antibody at a dilution of 1:500.