

Kat6a antibody (mAb)

Catalog No: 61327

Clone: 2MOZ-2D4

Isotype: IgG1k

Application(s): ChIP, IP, WB

Reactivity: Human, Mouse

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Quantity: 100 µg

Molecular Weight: 290 kDa

Background: **Kat6a** (K(lysine) acetyltransferase 6a, MOZ, MYST3) is a histone acetyltransferase that acetylates lysine residues in histone H3 and histone H4 (in vitro). Component of the MOZ/MORF complex which has a histone H3 acetyltransferase activity. May act as a transcriptional coactivator for RUNX1 and RUNX2.

Immunogen: This antibody was raised against a peptide corresponding to amino acid residue(s) 856-869 of mouse Kat6a.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 32 mM glycine, 0.07 mM EDTA, 30% glycerol, and 0.035% sodium azide.

Application Notes:

Published Applications:

ChIP

IP

WB

See references for more information. Individual optimization may be required.

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.