

MLL / HRX antibody (mAb)

Catalog Nos: 39829, 39830

RRID: AB_2793358

Clone: N4.4

Isotype: IgG1

Application(s): ChIP, IP, WB

Reactivity: Human, Mouse

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 440 kDa

Background: MLL (myeloid/lymphoid or mixed-lineage leukemia (trithorax homolog, Drosophila)), also known as Histone-lysine N-methyltransferase **HRX**, is a SET-domain containing histone methyltransferase specific for lysine 4 of histone H3. It serves as a transcriptional activator for many target genes, including several HOX genes. MLL is a frequent target for translocations in acute myeloid leukemia (AML), acute lymphoblastic leukemia (ALL) and mixed lineage leukemia (MLL). A key feature of MLL fusion proteins is their ability to efficiently transform hematopoietic cells into leukemia stem cells. Leukemogenic MLL translocations encode MLL fusion proteins that have lost H3 Lys4 methyltransferase activity. MLL is found in a complex of proteins that includes ASH2L, WDR5, DPY30 and RBBP5.

Immunogen: This MLL / HRX antibody was raised against a recombinant protein corresponding to amino acids 161-356 of mouse MLL.

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:

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.