

## MLL / HRX antibody (pAb)

**Catalog Nos:** 61295, 61296

**RRID:** AB\_2793585

**Isotype:** IgG

**Application(s):** ChIP, ChIP-Seq, ICC, IF, IP, WB

**Reactivity:** Human, Mouse

**Quantities:** 100 µg, 10 µg

**Purification:** Protein A Chromatography

**Host:** Rabbit

**Concentration:** 1 µg/µl

**Molecular Weight:** 180 kDa

**Background:** MLL (myeloid/lymphoid or mixed-lineage leukemia (trithorax homolog, Drosophila)), also known as Histone-lysine N-methyltransferase **HRX**, is a histone methyltransferase that plays an essential role in early development and hematopoiesis. Catalytic subunit of the MLL1 /MLL complex, a multiprotein complex that mediates both methylation of 'Lys-4' of histone H3 (H3K4me) complex and acetylation of 'Lys-16' of histone H4 (H4K16ac). In the MLL1/MLL complex, it specifically mediates H3K4me, a specific tag for epigenetic transcriptional activation. Has weak methyltransferase activity by itself, and requires other component of the MLL1/MLL complex to obtain full methyltransferase activity. Has no activity toward histone H3 phosphorylated on 'Thr-3', less activity toward H3 dimethylated on 'Arg-8' or 'Lys-9', while it has higher activity toward H3 acetylated on 'Lys-9'. Required for transcriptional activation of HOXA9. Promotes PPP1R15A-induced apoptosis.

**Immunogen:** This MLL / HRX antibody was raised against a recombinant protein corresponding to amino acids 2829-2883 of human MLL.

**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:

ChIP: 5 - 10 µg per ChIP

ChIP-Seq: 5 - 10 µg each

ICC/IF: 0.5 µg/ml dilution

WB: 0.5 - 2 µg per ChIP

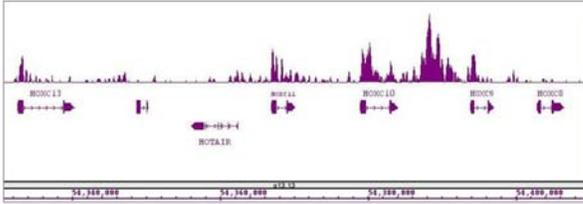
ChIP-Seq validation was performed by Active Motif's Epigenetics Services; the complete data set is available in the UCSC Genome Browser by clicking [here](#).

**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.

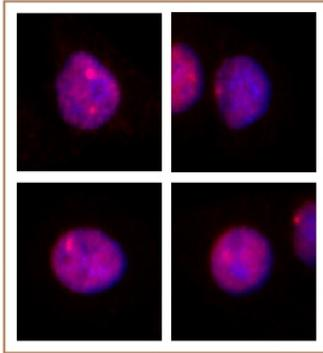
**MLL / HRX antibody (pAb) tested by ChIP-Seq.**

ChIP was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with 30 ug of chromatin from HeLa cells and 4 µg of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 20 million sequence tags were mapped to identify MLL / HRX binding sites. The image shows binding across a region of chromosome 12. You can view the complete data set in the UCSC Genome Browser, starting at this specific location, here.



**MLL / HRX antibody (pAb) tested by Immunofluorescence.**

Formaldehyde fixed HeLa cells stained with MLL / HRX antibody at a 0.5 µg/ml dilution.



**MLL/HRX antibody (pAb) tested by Western blot.**

HeLa whole cell extract extract (20 µg) probed with MLL/HRX antibody (pAb) at 2 µg/ml.

