

MLL1/KMT2A CUTANA™ CUT&RUN Antibody



EpiCypher®

Catalog No. 13-2004

Lot No. 21013001-38

Pack Size 100 µL

Type Monoclonal [BL-175-7E8] **Target Size** 431 kDa

Host Rabbit **Format** Pur. IgG

Reactivity Human

Applications CUT&RUN, WB, IP

Product Description:

This antibody meets EpiCypher's "CUTANA Compatible" criteria for performance in Cleavage Under Targets and Release Using Nuclease (CUT&RUN) and/or Cleavage Under Targets and Tagmentation (CUT&Tag) approaches to genomic mapping. Every lot of a CUTANA Compatible antibody is tested in the indicated CUTANA approach using EpiCypher optimized protocols (EpiCypher.com/resources/protocols) and determined to yield peaks that show a genomic distribution pattern consistent with reported function(s) of the target protein. MLL1 antibody produces CUT&RUN peaks above background (Figure 1) that overlap with H3K4me3 (Figures 1-2), consistent with its known role as the catalytic subunit of the MLL1/MLL histone methyltransferase complex.

Immunogen:

A synthetic peptide corresponding to human MLL1 amino acids 720 to 780.

Formulation:

Purified recombinant monoclonal antibody (1.0 mg/mL) in borate buffered saline (BBS) pH 8.2 with 0.09% sodium azide.

Storage and Stability:

Stable for 1 year at 4°C from date of receipt.

Application Notes:

Recommended Dilutions:

CUT&RUN: 0.5 µg

WB: 1:1,000

IP: 10 µL / 0.5 mg lysate

References:

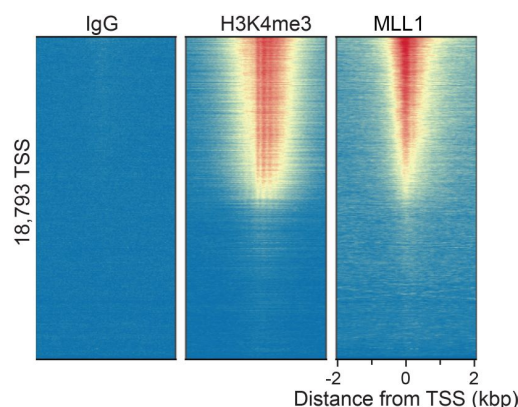


Figure 1: MLL1 enrichment at annotated transcription start sites (TSSs) in CUT&RUN. CUT&RUN was performed using 500,000 K562 cells with MLL1 and control antibodies (0.5 µg each; IgG negative control, EpiCypher 13-0042; H3K4me3 positive control, EpiCypher 13-0041). Sequencing reads were aligned to annotated TSSs (+/- 2 kbp) of 18,793 genes. High, medium, and low signal is ranked by intensity (top to bottom) and reflected by red, yellow, and blue colors, respectively. All rows aligned relative to MLL1.

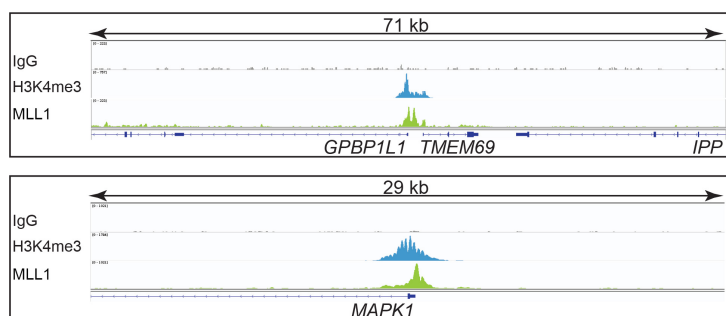


Figure 2: MLL1 CUT&RUN peaks and functional overlap. Two representative gene loci from the CUT&RUN data in Figure 1 are shown. MLL1 peaks overlap with H3K4me3 CUT&RUN peaks (EpiCypher 13-0041), consistent with the reported function of MLL1 as an H3K4 histone methyltransferase. Images were generated using the Integrative Genomics Viewer (Broad Institute).

This product is provided for commercial sale under license from Bethyl Laboratories, Inc.

Applications Key: ChIP: Chromatin immunoprecipitation; CUT&RUN: Cleavage Under Targets and Release Using Nuclease; CUT&Tag: Cleavage Under Targets and Tagmentation; E: ELISA; FACS: Flow cytometry; ICC: Immunocytochemistry; IF: Immunofluorescence; IHC: Immunohistochemistry; IP: Immunoprecipitation; L: Luminex; WB: Western Blot. **Reactivity Key:** B: Bovine; Ce: C. elegans; Ch: Chicken; Dm: Drosophila; Eu: Eukaryote; H: Human; M: Mouse; Ma: Mammal; R: Rat; Sc: S.cerevesiae; Sp: S. pombe; WR: Wide Range (predicted); X: Xenopus; Z: Zebrafish

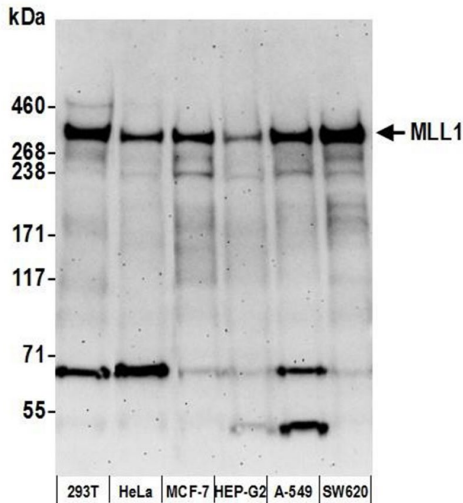


Figure 3: Western blot detection of human MLL1. Whole cell lysates were isolated from HEK293T, HeLa, MCF-7, HEP-G2, A-549, and SW620 cells using NETN lysis buffer. Fifty micrograms (50 µg) of lysate were loaded onto SDS-PAGE gel and analyzed under standard western blot conditions using MLL1 antibody (1:1,000 dilution).

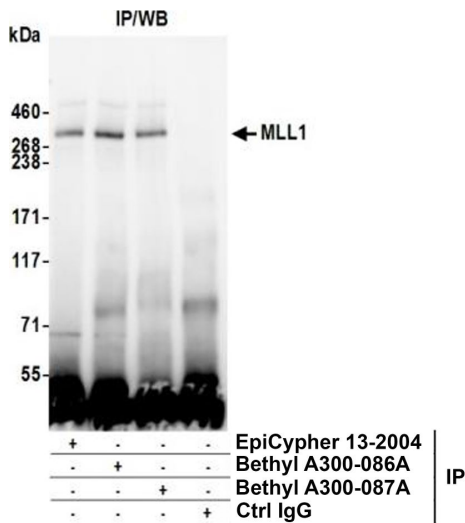


Figure 4: Immunoprecipitation of human MLL1. EpiCypher MLL1 antibody (10 µL) was used to immunoprecipitate whole cell lysates isolated from HEK293T cells using NETN lysis buffer (0.5 mg per IP). A negative control IgG antibody and positive control antibodies to various MLL1 epitopes (Bethyl Laboratories) were also used to demonstrate specificity of the IP. Immunoprecipitates were loaded onto SDS-PAGE gel (20% of IP loaded) and probed via western blot with EpiCypher MLL1 antibody (1:1,000 dilution).

This product is provided for commercial sale under license from Bethyl Laboratories, Inc.

Applications Key: ChIP: Chromatin immunoprecipitation; CUT&RUN: Cleavage Under Targets and Release Using Nuclease; CUT&Tag: Cleavage Under Targets and Tagmentation; E: ELISA; FACS: Flow cytometry; ICC: Immunocytochemistry; IF: Immunofluorescence; IHC: Immunohistochemistry; IP: Immunoprecipitation; L: Luminex; WB: Western Blot. **Reactivity Key:** B: Bovine; Ce: *C. elegans*; Ch: Chicken; Dm: *Drosophila*; Eu: Eukaryote; H: Human; M: Mouse; Ma: Mammal; R: Rat; Sc: *S. cerevisiae*; Sp: *S. pombe*; WR: Wide Range (predicted); X: Xenopus; Z: Zebrafish