
EHMT2 Polyclonal Antibody

(Catalog # A-3019)

Background

EHMT2 (Euchromatic Histone Lysine Methyltransferase 2) is a histone methyltransferase that specifically mono- and dimethylates Lys-9 of histone H3 (H3K9me1 and H3K9me2, respectively) in euchromatin. H3K9me is a specific tag for epigenetic transcriptional repression by recruiting HP1 proteins to methylated histones. This gene also mediates monomethylation of Lys-56 of histone H3 (H3K56me1) in G1 phase, promoting interaction between histone H3 and PCNA and regulating DNA replication. Also weakly methylates Lys-27 of histone H3 (H3K27me). Also required for DNA methylation, the histone methyltransferase activity is not required for DNA methylation, suggesting that these 2 activities function independently. Most likely targeted to histone H3 by different DNA-binding proteins like E2F6, MGA, MAX and/or DP1. May also methylate histone H1. In addition to the histone methyltransferase activity, also methylates non-histone proteins: mediates dimethylation of Lys-373 of p53/TP53. Also methylates CDYL, WIZ, ACIN1, DNMT1, HDAC1, ERCC6, KLF12 and itself.

Description

EHMT2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human

Isotype

IgG

Uniprot ID

Q96KQ7

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 877-1176 of human EHMT2 (NP_079532.5).

Storage

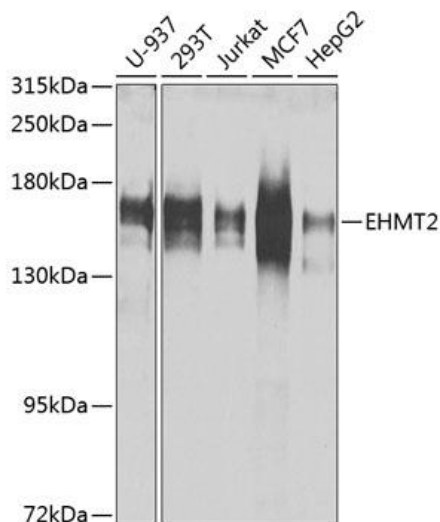
Shipped at 4°C. Upon receipt, store at -20°C. Avoid repeated freeze.

Alternative Names

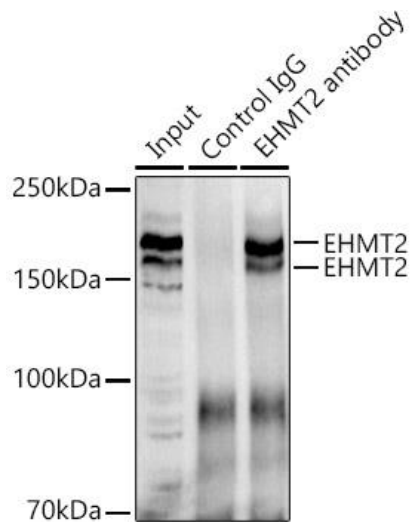
EHMT2; BAT8; C6orf30; G9A; GAT8; KMT1C; NG36; histone-lysine N-methyltransferase EHMT2

Application

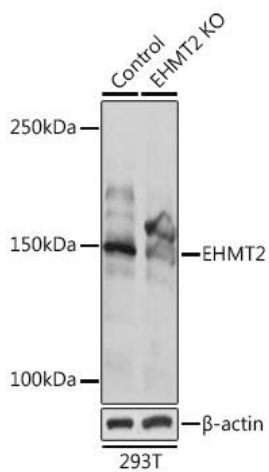
WB, IP, ELISA; Recommended dilution: WB 1:500 - 1:1000, IP 0.5 µg-4 µg antibody for 200 µg-400 µg extracts of whole cells



Western blot analysis of various lysates using EHMT2 Polyclonal Antibody at 1:1000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25µg per lane. Blocking buffer: 3% nonfat dry milk in TBST.



Immunoprecipitation analysis of 300 µg extracts of 293T cells using 3µg EHMT2 Polyclonal Antibody. Western blot was performed from the immunoprecipitate using EHMT2 Polyclonal Antibody at a dilution of 1:1000



Western blot analysis of lysates from wild type (WT) and EHMT2 Rabbit pAb knockout (KO) 293T cells, using EHMT2 Polyclonal Antibody at 1:1000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25µg per lane. Blocking buffer: 3% nonfat dry milk in TBST.