

EZH2 Rabbit pAb

Catalog No: 39639

Host: Rabbit

Application: ChIP, WB

Reactivity: Human, Mouse

Purification: Affinity Purified

Volume: 100 μ l

Concentration: 1 μ g/ μ l

Molecular Weight: 96 kDa

Background: EZH2 is a human homolog of the Drosophila Polycomb-group protein Enhancer of Zeste protein, and contains a SET domain that catalyzes the methylation of histone H3 at lysine 27. Polycomb-group proteins repress gene expression by binding to chromatin and locally altering chromatin structure. EZH2, BMI-1 and SUZ12 are present in the PRC2 and PRC3 protein complexes that function as mediators of epigenetic transcriptional silencing. Deregulation of EZH2 is linked to a variety of cancers.

Immunogen: Recombinant fusion protein corresponding to amino acids 1-370 of mouse EZH2.

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.

Application Notes:

Recommended Dilution:

WB: 1:500 - 1:2,000

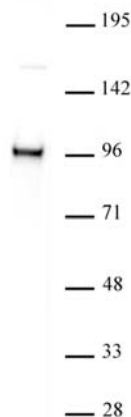
ChIP: 10 μ l per ChIP

References:

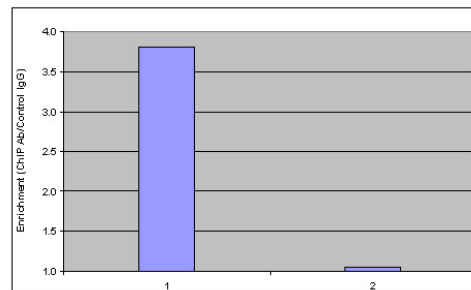
ChIP: van Dessel. et al (2010) Nucl Acids Res in press.

Storage and Guarantee: Antibodies in solution can be stored at -20°C for 2 years. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of arrival.

Application Key: ChIP = Chromatin Immunoprecipitation; FACS = Flow Cytometry; IF = Immunofluorescence; IHC = Immunohistochemistry; IP = Immunoprecipitation; WB = Western Blot



Western Blot: Nuclear extract of HeLa cells (20 μ g/lane) stained with 39639 (1:1,000 dilution).



ChIP performed on HeLa cell chromatin using 39639. qPCR was performed using primers of the human Myt1 gene. Lane 1: ChIP using 10 μ l of 39639. Lane 2: negative control rabbit IgG.