

## EZH2 antibody (pAb)

**Catalog Nos:** 61641, 61642

**RRID:** AB\_2793716

**Isotype:** IgG

**Application(s):** ChIP

**Reactivity:** Human, Mouse

**Quantities:** 100 µg, 10 µg

**Purification:** Protein A Chromatography

**Host:** Rabbit

**Concentration:** 1 µg/µl

**Molecular Weight:** 96 kDa

**Background:** EZH2 – (Enhancer of Zeste homolog 2) is a human homolog of the Drosophila Polycomb-group protein Enhancer of Zeste protein. It contains a SET domain that catalyzes the methylation of histone H3 at lysine 27 (Histone H3 Lys27 monomethylated, dimethylated and trimethylated). 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. EZH2 associates with the embryonic ectoderm development protein, the VAV1 oncoprotein, and the X-linked nuclear protein. EZH2 may play a role in the hematopoietic and central nervous systems. Deregulation of EZH2 is linked to a variety of cancers.

**Immunogen:** This antibody was raised against a recombinant fusion protein corresponding to amino acids 1-370 of mouse EZH2.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic. For your convenience, a sera version (Catalog No. 39901) of this antibody is also available.

### Application Notes:

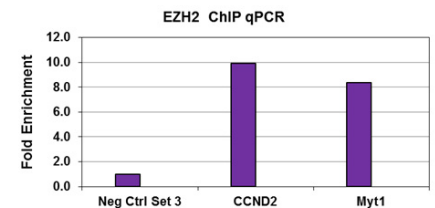
Applications Validated by Active Motif:

ChIP: 5 µg per ChIP

For EZH2, we also offer AbFlex<sup>®</sup> EZH2 Recombinant Antibody (rAb). For details, see Catalog No. 91159.

**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.



### EZH2 antibody (pAb) tested by ChIP.

Chromatin immunoprecipitation (ChIP) was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with 13 µg of chromatin from HeLa cells and 5 µg EZH2 antibody. ChIP DNA was used in qPCR with the control primer pairs or gene-specific primer pairs as indicated. Data are presented as Fold Enrichment of the ChIP antibody signal versus the negative control primer pair using the ddCT method.