

## Suz12 antibody (mAb)

**Catalog Nos:** 39877, 39878

**RRID:** AB\_2793377

**Clone:** 2A09

**Isotype:** IgG2b

**Application(s):** ChIP, IHC, IP

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 95 kDa

**Background:** Suz12 (Suppressor-of-Zeste 12 protein, also designated as ChET9, JAZ1 protein) is a component of the Polycomb group complexes 2, 3 and 4 (PRC2, PRC3 and PRC4), which also contain the histone methyltransferase EZH2, EED (Embryonic Ectoderm Development protein) protein and the histone-binding proteins RbAP46 and RbAP48. PRC2 is responsible for di- and trimethylation of histone 3 at lysine 27 (Histone H3K27me2 and Histone H3K27me3), which leads to transcriptional repression of the target genes. Suz12 is a nuclear protein belonging to the VEFS family. Suz12 contains a zinc-finger binding motif and a VEFS box, which is required for the interaction between Suz12 and EZH2.

**Immunogen:** This Suz12 antibody was raised against a recombinant protein corresponding to full-length human Suz12.

**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. Sodium azide is highly toxic.

### Application Notes:

Applications Validated by Active Motif:

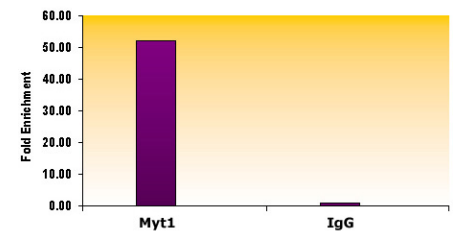
ChIP: 2.5 - 10 µg per ChIP

This antibody has been published, but has not been validated by Active Motif, for Western blot. For an antibody validated for use in Western blotting, see Catalog No. 39357.

This antibody is also available as an AbFlex® engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91163.

**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.



### Suz12 antibody (mAb) tested by ChIP.

Chromatin IP performed using the ChIP-IT® Express Kit (Catalog No. 53008) and HeLa chromatin ( $1.5 \times 10^6$  cell equivalents per ChIP) using 10 µg of Suz12 antibody or the equivalent amount of rabbit IgG as a negative control. Real time, quantitative PCR (RT-qPCR) was performed on DNA purified from each of the ChIP reactions using a primer pair specific for the Myt1 gene promoter. Data are presented as Fold Enrichment of the ChIP antibody signal versus the negative control IgG using the ddCT method.