

Ring1B antibody (mAb)

Catalog Nos: 39663, 39063, 39664

RRID: AB_2615006

Clone: 3-3

Isotype: IgG2b

Application(s): ChIP, ChIP-Seq, ICC, IF, IP, WB

Reactivity: Human, Monkey, Mouse

Quantities: 100 µg, 50 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 40 kDa

Background: Ring1B (also known as ring finger protein 2, RING2 or RNF2) is an essential component of the Polycomb group (PcG) multiprotein PRC1 complex. The PRC1 complex is important for the transcriptional repression of genes involved in development and cell proliferation, such as the *Hox* genes. Ring1B also acts as an E3 ubiquitin ligase on histone H2A of the PRC1 complex.

Immunogen: This Ring1B antibody was raised against the recombinant full-length mouse protein.

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:

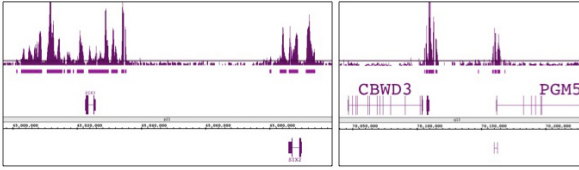
ChIP-Seq: 4 µl each

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.

Ring1B antibody (mAb) tested by ChIP-chip.

ChIP was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) with chromatin from 4.5 million human embryonic stem cells and 4 µl of antibody. ChIP DNA was amplified by WGA, labeled and hybridized to a human tiling array. The two images show regions of strong binding on chromosomes 2 and 9.



Ring1B antibody (mAb) tested by ChIP-Seq.

ChIP was performed using Ring1B antibody with 30 µg chromatin from the wild-type mouse ESC and 4 µg of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 19.5 million sequence tags were mapped to identify Ring1B binding sites. The image shows binding across a region of chromosome 2.

