

## BMI-1 antibody (mAb)

**Catalog Nos:** 61089, 61090

**RRID:** AB\_2793501

**Clone:** DC9

**Isotype:** IgG1, k

**Application(s):** ICC, IF, WB

**Reactivity:** Human, Mouse

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 40 kDa

**Background:** BMI-1 is a member of the Polycomb PRC1 complex that is recruited to transcriptionally repressed genes subsequent to histone H3 lysine 27 methylation in order to maintain repression. In the PRC1 complex, BMI-1 is required to stimulate the E3 ubiquitin-protein ligase activity of RNF2/RING2, resulting in the ubiquitylation of histone H2A at lysine 119. BMI-1 acts as an oncogene and cooperates with c-myc in the initiation of lymphoma. High levels of expression of BMI-1 are observed in metastatic melanoma, and BMI-1 has been implicated in several other cancers. BMI-1 is required for hematopoietic stem cell self renewal.

**Immunogen:** This BMI-1 antibody was raised against a recombinant protein corresponding to full-length mouse BMI-1.

**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:

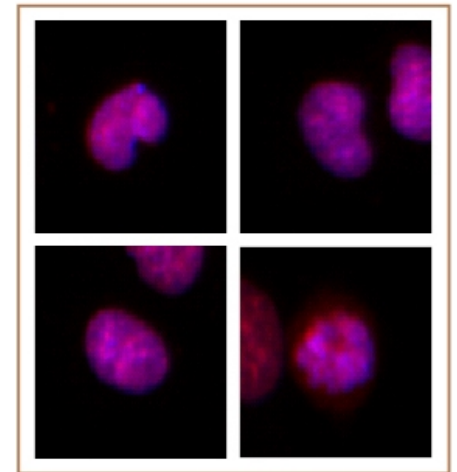
Applications Validated by Active Motif:

ICC/IF: 0.5 µg/ml dilution

For ChIP-Seq analysis of BMI-1, we offer AbFlex® BMI-1 Recombinant Antibody (rAb). For details, see Catalog No. 91195.

**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.



### BMI-1 antibody (mAb) tested by Immunofluorescence.

Formaldehyde fixed HeLa cells stained with BMI-1 antibody at a 0.5 µg/ml dilution.