

BMI-1 (IN) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat BMI-1 (Polycomb Group RING Finger Protein 4, RNF51)

Instruction Manual

Catalog Number	PK-AB718-3755
Synonyms	BMI-1 Antibody: Polycomb group RING finger protein 4, RNF51
Description	The transcriptional repressor BMI-1 was first identified as a proto-oncogene frequently activated by Moloney murine leukemia proviral insertions in mice and cooperating with c-myc in the generation of mouse lymphomas. BMI-1 is involved in segment specification, cell growth and maintenance, transcriptional regulation, and chromatin modification. A major target of BMI-1 is the ink4a locus which encodes tumor suppressor proteins p16 and p19Arf, which are important in tumor progression and thought to be critical in cell proliferation and senescence. Recent studies have also shown that BMI-1 is required for the maintenance of adult normal and leukemic stem cells, suggesting that BMI-1 could be an attractive therapeutic target for stem cell proliferation and renewal as well as for anti-cancer strategies.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	BMI-1 antibody was raised against a peptide corresponding to 15 amino acids near the center of human BMI-1. Antigen: Human BMI-1 (Intermediate Domain) Peptide (Cat. No. PK-AB718-3755P).
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, ICC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. BMI-1 antibody can be used for detection of BMI-1 by Western blot at 0.5 to 2 µg/mL. Antibody can also be used for immunocytochemistry starting at 10 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Alkema MJ, Wiegant J, Raap AK, et al. Characterization and chromosomal localization of the human proto-oncogene BMI-1. <i>Hum. Mol. Genet.</i> 1993; 2:1597-603. Jacobs JJ, Kieboom K, Marino S, et al. The oncogene and polycomb-group gene bmi-1 regulates cell proliferation and senescence through the ink4a locus. <i>Nature</i> 1999; 397:164-8. Lessard J and Sauvageau G. BMI-1 determines the proliferative capacity of normal and leukaemic stem cells. <i>Nature</i> 2003; 255-60.
Images	Available upon request.
Related Products	BMI-1 Peptide; Cat. No.: PK-AB718-3755P K562 Cell Lysate (CT); Cat. No.: PK-AB718-1204

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.