

PCL2 antibody (mAb)

Catalog Nos: 61153, 61154

RRID: AB_2615020

Clone: 3A9

Isotype: IgG1

Application(s): WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 70 kDa

Background: PCL2 (Polycomb-like 2, MTF2, M96) is a member of the **Polycomb PRC2** protein complex and is important in cell fate decisions and may play a role in recruitment of PRC2 to target genes and histone methylation. **PCL2** regulates gene expression patterns controlling self-renewal and commitment in embryonic stem cells and plays a critical role in ESC differentiation and self-renewal. PCL2 contains two PHD fingers.

Immunogen: This PCL2 antibody was raised against a recombinant protein corresponding to amino acids 103 - 327 of mouse PCL2.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

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.