

Phc1 antibody (mAb)

Catalog Nos: 39723, 39724

RRID: AB_2713961

Clone: 6-1-3

Isotype: IgG

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

Reactivity: Human, Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 150 kDa

Background: Phc1, polyhomeotic homolog 1 (also known as EDR1, early development regulatory protein 1) is a component of the polycomb group (PcG) multiprotein PRC1 complex. The PRC1 complex is required to maintain the stable repression of the Hox genes during development. This chromatin-associated protein complex acts via chromatin remodeling and histone modifications. Phc1 has been shown to interact with PHC2, BMI1, MAPKAPK2 and MCRS1.

Immunogen: This Phc1 antibody was raised against a recombinant mouse Phc1 protein fragment.

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:

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.