

Phc2 antibody (mAb)

Catalog Nos: 39661, 39662

Isotype: IgG

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

Reactivity: Human, Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 90 kDa

Background: Phc2, polyhomeotic homolog 2 (also known as EDR2, early development regulatory protein 2) 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. Phc2 has been shown to interact with PHC1, BMI1, MAPKAPK2 and MCRS1.

Immunogen: This Phc2 antibody was raised against a recombinant fragment of human Phc2.

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:

Published Applications:

ChIP

IP

ICC/IF

WB

See references for more information. Individual optimization may be required.

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.