

Mi-2 antibody (mAb)

Catalog Nos: 61463, 61464

RRID: AB_2793648

Clone: 4D8

Isotype: IgG2c

Application(s): IF, IP, WB

Reactivity: Drosophila

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 220 kDa

Background: **Mi-2** (ATP-dependent helicase Mi-2) is a member of the CHD family of ATP-dependent chromatin remodelers characterized by the presence of two tandemly arranged chromodomains. The Chd3/Chd4 subfamily contains a pair of PHD fingers preceding the chromodomains and these proteins exist in multisubunit complexes with nucleosome remodeling and deacetylation (NuRD) activity. *Drosophila* Mi-2 (dMI-2) is the orthologs of mammalian CHD4.

Immunogen: This antibody was raised against a synthetic peptide corresponding to amino acid residues 1641-1662 of *Drosophila* Mi-2.

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

Application Notes:

For Mi-2 beta (CHD4), we also offer AbFlex[®] Mi-2 beta (CHD4) Recombinant Antibody (rAb). For details, see Catalog No. 91275.

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.