

CHD3 antibody (mAb)

Catalog Nos: 61469, 61470

RRID: AB_2793650

Clone: 7A11

Isotype: IgG1

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

Reactivity: Drosophila

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 103 kDa

Background: CHD3 (chromodomain helicase DNA binding protein 3) is a component of the CHD family of ATP-dependent chromatin remodelers and forms histone deacetylase NuRD complexes which participate in the remodeling of chromatin by deacetylating histones.

Immunogen: This antibody was raised against a synthetic peptide corresponding to amino acid residues 865-885 of Drosophila CHD3.

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

Application Notes:

This antibody is also available as an AbFlex[®] engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91273.

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.