

CID antibody (mAb)

Catalog Nos: 61735, 61736

RRID: AB_2793749

Clone: 4F8

Isotype: IgG2a

Application(s): ICC, IF

Reactivity: Drosophila

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 35 kDa

Background: CID (Centromer Identifier) is the *Drosophila* histone H3 variant which replaces canonical H3 in centromeric chromatin at the inner plate of the kinetochore (similar to mammalian CENP-A). CID is required for recruitment and assembly of kinetochore proteins, mitotic progression and chromosome segregation. It may serve as an epigenetic mark that propagates centromere identity through DNA replication and cell division.

Immunogen: This antibody was raised against a recombinant protein corresponding to the N-terminus of drosophila CID.

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.