

CID antibody (pAb)

Catalog Nos: 39713, 39714

RRID: AB_2793317

Isotype: Serum

Application(s): WB

Reactivity: *Drosophila*

Volumes: 100 µl, 10 µl

Purification: None

Host: Rabbit

Molecular Weight: 30 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 CID antibody was raised against recombinant CID protein from *Drosophila melanogaster*.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

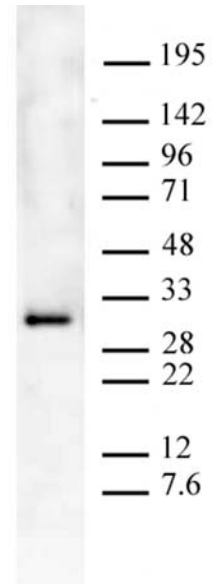
Validated Applications:

WB: 1:500 dilution

For an antibody against CID that works in immunofluorescence, use Catalog No. 39719.

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.



CID pAb tested by Western blot.

Drosophila Schneider L2 cells nuclear extract (10 µg) probed with CID pAb (1:500 dilution).