

## CDT1 antibody (mAb)

---

**Catalog Nos:** 61039, 61040

**RRID:** AB\_2793483

**Clone:** ID8

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Hamster

**Concentration:** 1 µg/µl

**Molecular Weight:** 65 kDa

---

**Background:** CDT1 (chromatin licensing and DNA replication factor 1) is a key licensing factor which functions to allow DNA replication. It is recruited (along with CDC6) by the ORC (Origin Recognition Complex) to sites of DNA replication and helps form the pre-replicative complex (pre-RC). Its activity during the cell cycle is regulated by its association with the Geminin protein geminin, which both inhibits CDT1 activity during S phase in order to prevent unregulated DNA replication and also to inhibit CDT1 proteolysis.

**Immunogen:** This CDT1 antibody was raised against a recombinant protein corresponding to amino acids 177-557 of mouse CDT1.

**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:

**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.