

## CDX1 antibody (mAb)

---

**Catalog Nos:** 61791, 61792

**RRID:** AB\_2793768

**Isotype:** IgG1

**Application(s):** ChIP, IF, IHC, WB

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein A Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

---

**Background:** The CDX family members (CDX1, CDX2, and CDX4) are Caudal-related transcription factors implicated in anterior-posterior patterning during embryonic development, regulation of cardiac differentiation in stem cells and down-regulation of genes associated with neuronal growth and vascular endothelial growth. The caudal-type homeobox 1 (CDX1) transcription factor controls enterocyte differentiation in the colon and is also central to the capacity of a colorectal cancer (CRC) cell line to differentiate. CDX1 is a negative marker of CRC stem cells. The expression of CDX1 is itself regulated by methylation at the CDX1 promoter region, and its expression is conversely correlated with expression of the polycomb complex protein BMI-1, which is involved in maintenance of quiescent stem cells and is expressed in cancer stem cells.

**Immunogen:** This antibody was raised against an N-terminal peptide (a.a. 30-47) of human CDX1 (NP\_001795).

**Buffer:** PBS containing 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.