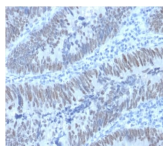


CDX2 / Caudal Type Homeobox 2 Monoclonal Mouse Antibody (PCRP-CDX2-1A3)



Product Description

The specificity of this monoclonal antibody to its intended target was validated by HuProt™Array, containing more than 19,000, full-length human proteins. The intestine-specific transcription factors CDX1 and CDX2 are important for directing intestinal development, differentiation, proliferation and maintenance of the intestinal phenotype. CDX2 protein expression has been seen in GI carcinomas. Anti-CDX2 has been useful to establish GI origin of metastatic adenocarcinomas and carcinoids and is especially useful to distinguish metastatic colorectal adenocarcinoma from lung adenocarcinoma. However, mucinous carcinomas of the ovary also express CDX2 protein. It limits the usefulness of this marker in the distinction of metastatic colorectal adenocarcinoma from mucinous carcinoma of the ovary.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 1920, Anti-CDX2 / Caudal Type Homeobox 2 (PCRP-CDX2-1A3)

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #1920 |
| Antibody reactivity (target) | CDX2 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | PCRP-CDX2-1A3 |
| Isotype | IgG1 |
| Molecular weight | 40 kDa |
| Synonyms | Caudal type homeobox 2; Caudal type homeobox transcription factor 2; Caudal-type homeobox protein 2; CDX2 |
| Human gene symbol | CDX2 |
| Entrez gene ID | 1045 |
| SwissProt | Q99626 |
| Unigene | 174249 |
| Immunogen | Recombinant full-length human CDX2 protein |
| Verified antibody applications | IHC (FFPE) (verified) |
| Antibody target cellular localization | Nucleus |
| Species reactivity | Human |
| Antibody application notes | Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunohistology (formalin): 0.5-1 ug/mL for 30 minutes at RT. Immunoprecipitation 2-4 ug/mg protein. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes. Western blotting 0.5-1.0 ug/mL. Optimal dilution for a specific application should be determined by user |
| Positive control | HT29 cells or Colon Carcinoma |
| Shipping condition | Room temperature |
| Storage Conditions | Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C |
| Shelf life | Guaranteed for at least 24 months from date of receipt when stored as recommended |
| Regulatory status | For research use only (RUO) |
| Antibody/conjugate formulation | Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide |
| Validated in protein array | Monospecific |
| Antibody research areas | Cancer, Developmental biology, Transcription factors |
| Product origin | Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. |
| Cell/tissue expression | Intestine |
| Tumor expression | Colorectal cancer, Ovarian cancer |

| Antibody # prefix | Conjugation | Ex/Em (nm) | Laser line | Detection channel | Dye Features |
|-------------------|--------------------|------------|------------|--------------------------|----------------------------------|
| BNC04 | CF®405S | 404/431 | 405 | DAPI (microscopy), AF405 | CF®405S Features |
| BNC88 | CF®488A | 490/515 | 488 | GFP, FITC | CF®488A Features |
| BNC68 | CF®568 | 562/583 | 532, 561 | RFP, TRITC | CF®568 Features |
| BNC94 | CF®594 | 593/614 | 561 | Texas Red® | CF®594 Features |
| BNC40 | CF®640R | 642/662 | 633-640 | Cy®5 | CF®640R Features |
| BNC47 | CF®647 | 650/665 | 633-640 | Cy®5 | CF®647 Features |
| BNC74 | CF®740 | 742/767 | 633-685 | 775/50 | CF®740 Features |
| BNCB | Biotin | N/A | N/A | N/A | |
| BNUB | Purified | N/A | N/A | N/A | |
| BNUM | Purified, BSA-free | N/A | N/A | N/A | |

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on August 28, 2025 at 03:37:07 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cdx2-caudal-type-homeobox-2-monoclonal-mouse-antibody-pcrp-cdx2-1a3/>