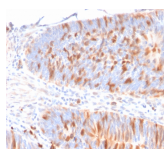


# MBD1 Monoclonal Mouse Antibody (CPTC-MBD1-1)



## Product Description

DNA methylation is the major modification of eukaryotic genomes and plays an essential role in mammalian development. Human proteins MECP2, MBD1, MBD2, MBD3, and MBD4 comprise a family of nuclear proteins related by the presence in each of a methyl-CpG binding domain (MBD). Each of these proteins, with the exception of MBD3, is capable of binding specifically to methylated DNA.

MECP2, MBD1 and MBD2 can also repress transcription from methylated gene promoters. MBD1 has also been predicted to be involved in the development and progression of tumors, and it has been shown that MBD1 is significantly upregulated in pancreatic cancer tissues when compared with surrounding normal tissues. Furthermore, MBD1 may serve as a potential therapeutic target for pancreatic cancer.

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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 2271, Anti-MBD1 (CPTC-MBD1-1)**

Call us : [800-304-5357](tel:800-304-5357)

## Product attributes

|                                       |                                                                                                                                                                                                                                     |
|---------------------------------------|-------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Antibody number                       | #2271                                                                                                                                                                                                                               |
| Antibody reactivity (target)          | MBD1                                                                                                                                                                                                                                |
| Antibody type                         | Primary                                                                                                                                                                                                                             |
| Host species                          | Mouse                                                                                                                                                                                                                               |
| Clonality                             | Monoclonal                                                                                                                                                                                                                          |
| Clone                                 | CPTC-MBD1-1                                                                                                                                                                                                                         |
| Isotype                               | IgG1, kappa                                                                                                                                                                                                                         |
| Molecular weight                      | ~80 kDa                                                                                                                                                                                                                             |
| Synonyms                              | CXXC-type zinc finger protein 3; CXXC3; MBD1; MECP1 COMPLEX; Methyl-CpG-binding domain protein 1 (MBD1); Protein containing methyl-CpG-binding domain 1 (PCM1); Regulator of fibroblast growth factor 2 (FGF 2) transcription (RFT) |
| Human gene symbol                     | MBD1                                                                                                                                                                                                                                |
| Entrez gene ID                        | 4152                                                                                                                                                                                                                                |
| SwissProt                             | Q9UIS9                                                                                                                                                                                                                              |
| Unigene                               | 405610                                                                                                                                                                                                                              |
| Immunogen                             | Recombinant human MBD1 protein                                                                                                                                                                                                      |
| Verified antibody applications        | IHC (FFPE) (verified)                                                                                                                                                                                                               |
| Antibody target cellular localization | Nucleus                                                                                                                                                                                                                             |
| Species reactivity                    | Human                                                                                                                                                                                                                               |
| Positive control                      | Human tonsil or colon (IHC).                                                                                                                                                                                                        |
| 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                                                                                                                     |
| 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                                   |
| Shelf life                            | Guaranteed for at least 24 months from date of receipt when stored as recommended                                                                                                                                                   |
| 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.                                                     |
| Tumor expression                      | Pancreatic cancer                                                                                                                                                                                                                   |

| Antibody # prefix | Conjugation        | Ex/Em (nm) | Laser line | Detection channel        | Dye Features                     |
|-------------------|--------------------|------------|------------|--------------------------|----------------------------------|
| BNC04             | CF®405S            | 404/431    | 405        | DAPI (microscopy), AF405 | <a href="#">CF®405S Features</a> |
| BNC88             | CF®488A            | 490/515    | 488        | GFP, FITC                | <a href="#">CF®488A Features</a> |
| BNC68             | CF®568             | 562/583    | 532, 561   | RFP, TRITC               | <a href="#">CF®568 Features</a>  |
| BNC94             | CF®594             | 593/614    | 561        | Texas Red®               | <a href="#">CF®594 Features</a>  |
| BNC40             | CF®640R            | 642/662    | 633-640    | Cy®5                     | <a href="#">CF®640R Features</a> |
| BNC47             | CF®647             | 650/665    | 633-640    | Cy®5                     | <a href="#">CF®647 Features</a>  |
| BNC74             | CF®740             | 742/767    | 633-685    | 775/50                   | <a href="#">CF®740 Features</a>  |
| 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 January 7, 2026 at 06:58:21 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/mbd1-monoclonal-mouse-antibody-cptc-mbd1-1/>