Cdc34 homolog Monoclonal Mouse Antibody (CPTC-CDC34-2)



Product Description

Cell cycle events are regulated by the sequential activation and deactivation of cyclin dependent kinases (Cdks) and by the proteolysis of cyclins. The cell division cycle (Cdc) genes are required at various points in the cell cycle. Cdc25A, Cdc25B and Cdc25C protein tyrosine phosphatases function as mitotic activators by dephosphorylating Cdc2 p34 on regulatory tyrosine residues. Cdc6 is the human homolog of Saccharomyces cerevisiae Cdc6, which is involved in the initiation of DNA replication. Cdc37 appears to facilitate Cdk4/cyclin D1 complex formation and has been shown to form a stable complex with HSP 90. Cdc34, Cdc27 and Cdc16 function as ubiquitinconjugating enzymes. Cdc34 is thought to be the structural and functional homolog of Saccharomyces cerevisiae Cdc34, which is essential for the G1 to S phase transition. Cdc16 and Cdc27 are components of the APC (anaphase promoting complex) which ubiquitinates cyclin B, resulting in cyclin B/Cdk complex degradation.

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 2267, Anti-Cell Division Cycle 34 homolog (CPTC-CDC34-2)

Product attributes

Call us: 800-304-5357

| Fiduuci attiibutes | | | |
|--|--|--|--|
| Antibody number | #2267 | | |
| Antibody reactivity (target) | ell Division Cycle 34 homolog | | |
| Antibody type | Primary | | |
| Host species | Mouse | | |
| Clonality | Monoclonal | | |
| Clone | CPTC-CDC34-2 | | |
| Isotype | IgG1, kappa | | |
| Molecular weight | 34 kDa | | |
| Synonyms | Cdc34; cell division cycle 34; UBC3; UBCH3; UBE2R1; Ubiquitin carrier protein; Ubiquitin conjugating enzyme E2 32 kDa complementing; Ubiquitin protein ligase R1; Ubiquitin conjugating enzyme E2-CDC34 | | |
| Human gene symbol | CDC34 | | |
| Entrez gene ID | 997 | | |
| SwissProt | P49427 | | |
| Unigene | 514997 | | |
| Immunogen | Recombinant human full-length CDC34 protein | | |
| Verified antibody applications | IHC (FFPE) (verified) | | |
| Antibody target cellular localization | Nucleus & cytoplasm | | |
| Species reactivity | Human | | |
| Positive control | 293, K562, Jurkat cell lysates. Human small intestine. | | |
| Shipping condition | Room temperature | | |
| Storage Conditions | Store at 2 to 8 $^{\circ}$ C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 $^{\circ}$ 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/o.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide | | |
| Validated in protein array | Monospecific | | |
| 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. | | |

| 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 November 22, 2025 at 01:33:30 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cdc34-homolog-monoclonal-mouse-antibody-cptc-cdc34-2/