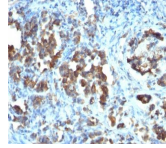


Cdc20 Monoclonal Mouse Antibody (CDC20/1102)



Product Description

Cyclins, regulatory subunits that associate with kinases, control many of the important steps in cell cycle progression. The Cdc2 protein kinase (p34Cdc2) exhibits protein kinase activity in vitro and exists in a complex with both cyclin B and a protein homologous to p13SUC1. Cdc2 kinase is the active subunit of the M phase promoting factor (MPF) and the M phase-specific Histone H1 kinase. The p34Cdc2/cyclin B complex is required for the G2 to M transition. An additional cell cycle-dependent protein kinase, termed p55cdc, exhibits a high degree of homology with the *S. cerevisiae* proteins Cdc20 and Cdc4. The p55cdc transcript is readily detectable in a variety of cultured cell lines in growth phase, but disappears when cell growth is chemically arrested. 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 1102, Anti-Cdc20 (CDC20/1102)**

Product attributes

| | |
|--|--|
| Antibody number | #1102 |
| Antibody reactivity (target) | Cdc20 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | CDC20/1102 |
| Isotype | IgG1, kappa |
| Molecular weight | 55 kDa |
| Synonyms | CDC20, CDC20A, p55CDC, P55CDC-LSB |
| Human gene symbol | CDC20 |
| Entrez gene ID | 991 |
| SwissProt | Q12834 |
| Unigene | 524947 |
| Immunogen | Recombinant human Cdc20 protein |
| Antibody target cellular localization | Cytoplasmic |
| Verified antibody applications | IHC (FFPE) (verified) |
| Species reactivity | Human |
| Antibody application notes | Immunohistochemistry formalin-fixed 2-4 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Optimal dilution for a specific application should be determined by user |
| Positive control | Ramos or HeLa cells. Tonsil or gastric 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 buffer 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 |
| Antibody research areas | Cell cycle |

Catalog Number Key for Primary Antibody Conjugates

| Antibody # prefix | Conjugation | Ex/Em (nm) | Laser line | Detection channel | Special applications/notes |
|-------------------|----------------------|-------------------|------------|----------------------------------|---|
| BNC04 | CF@405S | 404/431 | 405 | DAPI (microscopy), AF405 | SIM |
| BNC05 | CF@405M | 408/452 | 405 | DAPI (microscopy), Pacific Blue™ | STORM, STED, SIM, 2-Photon, TIRF |
| BNC06 | CF@405L | 395/545 | 405 | Pacific Orange™ | Spectral flow/imaging |
| BNC88 | CF@488A | 490/515 | 488 | GFP, FITC | STORM, STED, 2-photon, TIRF |
| BNC14 | CF@514 | 516/548 | 488 | GFP, FITC | Spectral flow/imaging |
| BNC43 | CF@543 | 541/560 | 532 | RFP, TRITC | |
| BNC55 | CF@555 | 555/565 | 532, 561 | RFP, TRITC | STORM |
| BNC68 | CF@568 | 562/583 | 532, 561 | RFP, TRITC | STORM, STED, SIM, TIRF |
| BNC94 | CF@594 | 593/614 | 561 | Texas Red® | STORM, STED, 2-photon |
| BNC40 | CF@640R | 642/662 | 633-640 | Cy@5 | SIM, TIRF, and FLImP |
| BNC47 | CF@647 | 650/665 | 633-640 | Cy@5 | STORM |
| BNC60 | CF@660C | 667/685 | 633-640 | Cy@5 | Spectral flow/imaging, STORM |
| BNC61 | CF@660R | 663/682 | 633-640 | Cy@5 | Highly photostable |
| BNC80 | CF@680 | 681/698 | 633-640 | Cy@5.5, IR@700 | STORM, 3D SMLM, LI-COR® Odyssey® |
| BNC81 | CF@680R | 680/701 | 633-640 | Cy@5.5, IR@700 | 2-photon, STED, SMT, LI-COR® Odyssey®; Highly photostable |
| BNC00 | CF@700 | 695/720 | 633-640 | AF700 | Spectral flow/imaging |
| BNC70 | CF@770 | 770/797 | 785, 808 | IR@800 | LI-COR® Odyssey® |
| BNCR | R-PE (PE) | 496, 546, 565/578 | 488-561 | PE | |
| BNCA | APC | 650/660 | 594-633 | APC | |
| BNCP | PerCP | 482/677 | 488-561 | PerCP, PE-Cy@5 | |
| BNCB | Biotin | N/A | N/A | N/A | |
| BNCAP | Alkaline Phosphatase | N/A | N/A | N/A | |
| BNCH | HRP | 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 GE Healthcare; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.