

# Clathrin, Heavy Chain Monoclonal Mouse Antibody (CHC/1432)

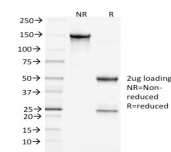
## Product Description

Recognizes protein of 192 kDa, which is identified as Clathrin Heavy Chain. Clathrin is composed of three heavy chains and three light chains, which associate non-covalently to form a triskelion structure. Clathrin heavy chain (HC) is composed of a terminal globular domain, a distal segment and a proximal segment containing a light chain-binding site. The proximal segment of the Clathrin HC protein is essential for interactions between Clathrin heavy chains and light chains, which result in the formation of the triskelion structure.

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 1432, Anti-Clathrin, Heavy Chain (CHC/1432)**



Call us : [800-304-5357](tel:800-304-5357) Email: [techsupport@biotium.com](mailto:techsupport@biotium.com)

### Product attributes

|                                       |   |
|---------------------------------------|---|
| Antibody number                       | #1432   |
| Antibody reactivity (target)          | Clathrin, Heavy Chain   |
| Antibody type                         | Primary   |
| Host species                          | Mouse   |
| Clonality                             | Monoclonal  |
| Clone                                 | CHC/1432  |
| Isotype                               | IgG1, kappa   |
| Molecular weight                      | 192 kDa   |
| Synonyms                              | Clathrin heavy chain; Clathrin heavy chain 1; Clathrin heavy chain 2; Clathrin heavy chain on chromosome 17 (CHC17 or CHC-17); CLH22; CLTC; CLTCL; CLTCL1; CLTCL2; CLTD   |
| Human gene symbol                     | CLTC  |
| Entrez gene ID                        | 1213  |
| SwissProt                             | P53675; Q00610  |
| Unigene                               | 491351  |
| Immunogen                             | Recombinant full-length human Clathrin Heavy Chain protein  |
| Verified antibody applications        | IF (verified)   |
| Antibody target cellular localization | Cytoplasmic, Lysosomes, Vesicular   |
| Species reactivity                    | Bird, Cow, Dog, Green monkey, Human, Mouse, Pig, Rat, Xenopus laevis  |
| Antibody application notes            | Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL; Optimal dilution for a specific application should be determined. |
| Positive control                      | HeLa cells, Placenta or Prostate 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   |
| 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.   |

| 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 September 17, 2025 at 09:57:08 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/clathrin-heavy-chain-monoclonal-mouse-antibody-chc-1432/>