

Clathrin, Heavy Chain Monoclonal Mouse Antibody (CLTC/1431)

Product Description

This antibody 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 to inquire about stock status and lead times before placing your order. **Catalog number key for antibody number 1431, Anti-Clathrin, Heavy Chain (CLTC/1431)**

Product attributes

Antibody number	#1431
Antibody reactivity (target)	Clathrin, Heavy Chain
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CLTC/1431
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	Purified recombinant N-terminal fragment of human Clathrin Heavy Chain (exact sequence is proprietary)
Antibody target cellular localization	Cytoplasmic, Lysosomes, Vesicular
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
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
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
Antibody research areas	Membrane trafficking

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF®488A	490/515	488	GFP, FITC
BNC68	CF®568	562/583	532, 561	RFP, TRITC
BNC94	CF®594	593/614	561	Texas Red®
BNC40	CF®640R	642/662	633-640	Cy®5
BNC47	CF®647	650/665	633-640	Cy®5
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

Dye Features

[CF®405S Features](#)

[CF®488A Features](#)

[CF®568 Features](#)

[CF®594 Features](#)

[CF®640R Features](#)

[CF®647 Features](#)

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, LI-COR Bioscience.