## Mitochondrial Marker Monoclonal Mouse Antibody (MTC02)



## **Product Description**

This MAb recognizes a 60 kDa antigen associated with the mitochondria in human cells. It is a part of a new panel of reagents, which recognizes subcellular organelles or compartments of human cells. These markers may be useful in identification of these organelles in cells, tissues, and biochemical preparations. This MAb recognizes an antigen associated with the mitochondria in human cells only. It can be used to stain the mitochondria in cell or tissue preparations and can be used as a mitochondrial marker in subcellular fractions. It produces a spaghetti-like pattern in normal and malignant cells and may be used to stain mitochondria of cells in fixed or frozen tissue sections. It can also be used with paraformaldehyde fixed frozen tissue or cell preparations. This MAb is an excellent marker for human cells in xenographic model research. It reacts specifically with human cells, including neurons and embryonic stem cells.

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 0740, Anti-Mitochondrial Marker (MTC02)

## Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	0740		
Antibody reactivity (target)	Mitochondrial Marker		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	MTC02		
Isotype	lgG1, kappa		
Molecular weight	60 kDa		
Synonyms	Not Known		
Entrez gene ID	Not Known		
SwissProt	Not Known		
Unigene	ne Not Known		
Immunogen	Semi-purified mitochondrial preparation		
Antibody target cellular localization	Mitochondria		
Verified antibody applications	IHC (FFPE) (verified)		
Species reactivity	Human		
Antibody application notes  Higher concentration may be required for direct detection primary antibody conjugates than for indirect detection secondary antibody, Immunofluorescence: 0.5-1 ug/mL, not react with mouse or rat, others not known, Immunoh formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed is requires boiling tissue sections in 10 mM Tris with 1 mM pH 9.0, for 10-20 min followed by cooling at RT for 20 m Optimal dilution for a specific application should be determined by user			
Positive control	HeLa or HepG2 cells. Hepatic carcinoma.		
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		
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			
Antibody research areas	Organelle markers		
roduct origin  Product may contain either bovine serum albumin (BSA) fi bovine serum (Bos taurus), or recombinant BSA produced Chinese hamster ovary cells. Inquire for the specific lot.			

Email: btinfo@biotium.com

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 5, 2025 at 09:45:07 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/monoclonal-anti-mitochondrial-marker-mtc02/">https://biotium.com/product/monoclonal-anti-mitochondrial-marker-mtc02/</a>