

# Mitochondrial Marker Recombinant Monoclonal Mouse Antibody (2509.MTC)

## - Biotium Choice

A recombinant mouse monoclonal antibody that recognizes a 60 kDa antigen associated with the mitochondria in cells. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.

# Biothum

## **Product Description**

Mitochondrial Marker (2509.MTC) is a recombinant mouse monoclonal antibody that recognizes a 60 kDa antigen associated with the mitochondria in cells. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance. The antibody is available conjugated to a selection of CF® Dyes. They are supplied in PBS, 2 mg/mL rBSA, and 0.05% sodium azide.

- Available in 7 bright and photostable CF® Dyes, including near-IR CF®740
- Suitable for immunohistochemistry, immunofluorescence, and western

This mitochondrial marker antibody selectively stains mitochondria in cell or tissue preparations and subcellular fractionation studies.

# Discover Sharper Signals and Unmatched Panel Flexibility with Biotium Choice Antibodies – Powered by CF® Dyes

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise. Labeled with our advanced CF® Dyes, they are our top-recommended antibodies for immunofluorescence and other applications.

#### **Biotium Choice Antibody Features**

- Robust and validated clones against common targets
- Developed and optimized for immunofluorescence and other applications
- Conjugated to bright, photostable CF® Dyes for superior signal and clarity
- New antibody clones and dye conjugates continuously in development

View our full catalog of Biotium Choice antibodies

#### Product attributes

Call us: 800-304-5357

Antibody number	P045			
Antibody reactivity (target)	Mitochondrial Marker			
<b>Biotium Choice Antibody</b>	Primary			
Antibody type	Primary			
Host species	Mouse			
Clonality	Recombinant Monoclonal			
Clone	2509.MTC			
Isotype	IgG1, kappa			
Molecular weight	60 kDa			
Antibody target cellular localization	Mitochondria			
Verified antibody applications	IF (verified)			
Species reactivity	Human			
Antibody application notes	Immunofluorescence: 1-5 ug/mL; Optimal concentration to be determined by end-user			
Positive control	HeLa cells.			
Shipping condition	Room temperature			
Storage Conditions	Store conjugates at 2 °C to 8 °C, Protect fluorescent conjugate from light			
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: PBS/2 mg/mL rBSA/0.05% azide			
Antibody research areas	Cancer, Neuroscience			
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line			
Expected antibody applications	Flow (intracellular), IHC (frozen), IHC (FFPE), WB			

### Mitochondrial Marker Recombinant Monoclonal Mouse Antibody (2509.MTC) - Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	100 ug/mL	500 uL	P045-488A-500UL	CF®488A Features
CF®568	562/584 nm	100 ug/mL	500 uL	P045-568-500UL	CF®568 Features
CF®594	593/615 nm	100 ug/mL	500 uL	P045-594-500UL	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	500 uL	P045-640R-500UL	CF®640R Features
CF®647	652/668 nm	100 ug/mL	500 uL	P045-647-500UL	CF®647 Features
CF®680R	680/701 nm	100 ug/mL	500 uL	P045-680R-500UL	CF®680R Features
CF®740	742/767 nm	100 ug/mL	500 uL	P045-740-500UL	CF®740 Features

This datasheet was generated on November 24, 2025 at 11:34:08 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/mitochondrial-marker-recombinant-monoclonal-mouse-antibody-2509-mtc-biotium-choice/">https://biotium.com/product/mitochondrial-marker-recombinant-monoclonal-mouse-antibody-2509-mtc-biotium-choice/</a>