



Glowing products for science

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

Nuclear Membrane Recombinant Monoclonal Mouse Antibody (2406.NM) - Biotium Choice



This nuclear membrane antibody (clone 2406.NM) is a validated antibody that recognizes an antigen associated with the nuclear membrane expressed in human cells. Validated for IF.

Product Description

Nuclear Membrane Recombinant Monoclonal Mouse Antibody (clone 2406.NM) recognizes an antigen associated with the nuclear membrane expressed in human cells. This antibody belongs to the Biotium Choice list of select antibodies that have been formulated and validated in-house for optimal performance. This antibody has been validated for immunofluorescence microscopy and is available conjugated to CF® Dyes.

The nuclear envelope or nuclear membrane is the double membrane of the cell nucleus. This antibody stains the rims of cell nuclei in fixed and permeabilized cells or frozen or paraffin tissue sections.

Biotium Choice Antibody Features

- Robust and validated clones against common targets
- Developed and optimized for flow cytometry 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](#)

Learn more about our innovative reagents for [immunofluorescence microscopy](#), such as [NucSpot® Nuclear Stains](#) for bright and nuclear-specific staining in a wide color selection, and [CytoLiner™ Fixed Cell Membrane Stains](#) for robust membrane staining in formaldehyde-fixed cells.

Discover Sharper Signals and Unmatched Panel Flexibility with Biotium Choice Antibodies Powered by CF® Dyes and Astral Leap™ Tandem Dyes

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise. Labeled with our advanced CF® Dyes and Astral Leap™ tandem dyes, they are our top-recommended antibodies for flow cytometry and other applications.

Biotium Choice Antibody Features

- Robust and validated clones against common targets
- Developed and optimized for flow cytometry and other applications
- Conjugated to bright, photostable [CF® Dyes](#) for superior signal and clarity
- Also available with [Astral Leap™ tandem dyes](#) for expanding multiplexing
- New antibody clones and dye conjugates continuously in development

[View our full catalog of Biotium Choice antibodies](#)

Product attributes

Antibody number	#P023
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Primary, Primary
Clonality	Recombinant Monoclonal
Host species	Mouse
Clone	2406.NM
Isotype	IgG1, kappa
Antibody reactivity (target)	Nuclear Membrane
Synonyms	Nuclear Envelope, Perinuclear Envelope, Nuclear Membrane, Nucleolemma, Karyotheca
Species reactivity	Human
Antibody target cellular localization	Nuclear membrane
Cell/tissue expression	All cells
Verified antibody applications	IF (verified)
Positive control	HeLa or other human cell line
Antibody application notes	Does not react with rat, Immunofluorescence (PFA- or MeOH-fixed): 1-5 ug/mL
Antibody research areas	Organelle markers
Antibody/conjugate formulation	Conjugates: PBS/50% glycerol/2 mg/mL rBSA/ 0.05% azide
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Storage Conditions	Store conjugates at 2 °C to 8 °C
Regulatory status	For research use only (RUO)
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.

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	100 ug/mL	200 uL	P023-488A-200UL	CF®488A Features
			1 mL	P023-488A-1ML	
CF®568	562/584 nm	100 ug/mL	200 uL	P023-568-200UL	CF®568 Features
			1 mL	P023-568-1ML	
CF®594	593/615 nm	100 ug/mL	200 uL	P023-594-200UL	CF®594 Features
			1 mL	P023-594-1ML	
CF®640R	642/663 nm	100 ug/mL	200 uL	P023-640R-200UL	CF®640R Features
			1 mL	P023-640R-1ML	
CF®647	652/668 nm	100 ug/mL	200 uL	P023-647-200UL	CF®647 Features
			1 mL	P023-647-1ML	
CF®740	742/767 nm	100 ug/mL	200 uL	P023-740-200UL	CF®740 Features
			1 mL	P023-740-1ML	
CF®750	755/779 nm	100 ug/mL	200 uL	P023-750-200UL	CF®750 Features
			1 mL	P023-750-1ML	

This datasheet was generated on August 2, 2025 at 05:17:13 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/nuclear-membrane-recombinant-monoclonal-mouse-antibody-2406-nm-biotium-choice/>