

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® Dves.

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 LeapTM 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 <u>Astral Leap™ tandem dyes</u> for expanding multiplexing
- New antibody clones and dye conjugates continuously in development

View our full catalog of Biotium Choice antibodies

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				
<u>CF®568</u>	562/584 nm	100 ug/mL	200 uL	P023-568-200UL	CF®568 Features
1 mL	P023-568-1ML				
<u>CF®594</u>	593/615 nm	100 ug/mL	200 uL	P023-594-200UL	CF®594 Features
1 mL	P023-594-1ML				
<u>CF®640R</u>	642/663 nm	100 ug/mL	200 uL	P023-640R-200UL	CF®640R Features
1 mL	P023-640R-1ML				
<u>CF®647</u>	652/668 nm	100 ug/mL	200 uL	P023-647-200UL	CF®647 Features
1 mL	P023-647-1ML				
<u>CF®740</u>	742/767 nm	100 ug/mL	200 uL	P023-740-200UL	CF®740 Features
1 mL	P023-740-1ML				
<u>CF®750</u>	755/779 nm	100 ug/mL	200 uL	P023-750-200UL	CF®750 Features
1 mL	P023-750-1ML				

Call us: 800-304-5357

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.		