



Glowing products for science

Nucleolin Recombinant Monoclonal Mouse Antibody (2507.NCL) - Biotium Choice

A recombinant mouse monoclonal antibody that recognizes nucleolin (NCL). The antibody belongs to the Biotium Choice list of select antibodies that have been validated in-house for optimal performance.

Product Description

Nucleolin Recombinant Monoclonal Mouse Antibody (2507.NCL) is a recombinant mouse monoclonal antibody that recognizes nucleolin (NCL). This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for fluorescence microscopy. 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 immunofluorescence and western

Recognizes Nucleolin (NCL), the major nucleolar phosphoprotein of growing eukaryotic cells. NCL is located mainly in dense fibrillar regions of the nucleolus. It is found associated with intranucleolar chromatin and pre-ribosomal particles. Human NCL gene consists of 14 exons with 13 introns and spans approximately 11 kb. It has a predicted molecular weight of ~76 kDa with an apparent MW of 100-110 kDa, attributed to phosphorylation of its N-terminal domain. It induces chromatin decondensation by binding to histone H1. It is thought to play a role in pre-rRNA transcription and ribosome assembly.

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](#)

Nucleolin Recombinant Monoclonal Mouse Antibody (2507.NCL) - Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	100 ug/mL	500 uL	P039-488A-500UL	CF®488A Features
CF®568	562/584 nm	100 ug/mL	500 uL	P039-568-500UL	CF®568 Features
CF®594	593/615 nm	100 ug/mL	500 uL	P039-594-500UL	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	500 uL	P039-640R-500UL	CF®640R Features
CF®647	652/668 nm	100 ug/mL	500 uL	P039-647-500UL	CF®647 Features
CF®680R	680/701 nm	100 ug/mL	500 uL	P039-680R-500UL	CF®680R Features
CF®740	742/767 nm	100 ug/mL	500 uL	P039-740-500UL	CF®740 Features

This datasheet was generated on January 8, 2026 at 04:08:02 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/nucleolin-recombinant-monoclonal-mouse-antibody-2507-ncl-biotium-choice/>



Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody Number	P039
Antibody reactivity (target)	Nucleolin
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Primary, Primary
Host species	Rat
Clonality	Recombinant Monoclonal
Clone	2507.NCL
Isotype	IgG1, lambda
Molecular weight	76 kDa (predicted), 100-110 kDa (observed)
Synonyms	NCL; Nucl; Nucleolin; Protein C23
Human gene symbol	NCL
Entrez gene ID	4691
SwissProt	P19338
Immunogen	Lysate of SU-DHL-1 Nuclei
Antibody target cellular localization	Cytoplasm, Nucleus, Nucleoli
Verified antibody applications	IF (verified), WB (verified)
Species reactivity	Human
Expected antibody applications	IF (verified), WB (verified)
Antibody application notes	Immunofluorescence: 1-5 ug/mL; Optimal concentration to be determined by end-user.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Store at 2 to 8 °C or below, Protect fluorescent conjugates 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)
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line
Antibody/conjugate formulation	Organelle markers
Cell/tissue expression	All human cells
Filter by antibody application	IF (verified), WB (verified)