

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 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 |
| CF®790 | 783/808 nm | 100 ug/mL | 500 uL | P039-790-500UL | CF®790 Features |

Product attributes

| | |
|---------------------------------------|--|
| Antibody Number | P039 |
| Antibody reactivity (target) | Nucleolin |
| Biotium Choice Antibody | Primary |
| Antibody type | Biotium Choice, Primary |
| Host species | Rat |
| Clonality | Monoclonal, Recombinant Monoclonal |
| Clone | 2507.NCL |
| Isotype | IgG1, lambda |
| Molecular weight of antigen | 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 |

This datasheet was generated on June 7, 2026 at 08:41:53 PM. 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/>