



Glowing products for science

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

Biotin Recombinant Monoclonal Mouse Antibody (rBN-34) – Biotium Choice

A recombinant mouse monoclonal antibody that recognizes free biotin and biocytin. The antibody belongs to the Biotium Choice list of select antibodies that have been validated in-house for optimal performance.



Product Description

Biotin Recombinant Monoclonal Mouse Antibody (rBN-34) is a recombinant mouse monoclonal antibody that recognizes free biotin and biocytin. 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 purified, or conjugated to a selection of CF® Dyes.

- Available in 10 bright and photostable CF® Dyes
- Suitable for flow cytometry, immunohistochemistry, immunofluorescence, and western

See our full selection of [anti-tag and anti-hapten antibody conjugates](#).

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 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](#)

Product attributes

Antibody Number	P033
Antibody reactivity (target)	Biotin
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	rBN-34
Synonyms	Bios II, Coenzyme R, D-Biotin, Vitamin B7, Vitamin H
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, 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)
Antibody research areas	Protein tag
Filter by antibody application	Flow, IHC, WB, IF (verified)
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line

Biotin Recombinant Monoclonal Mouse Antibody (rBN-34) -Biotium-Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF®500R	500/500 nm	100 ug/mL	1 mL	P033-500R-WML	CF®500R Features

This datasheet was generated on September 9, 2025 at 08:38:37 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-recombinant-monoclonal-mouse-antibody-rbn-34-biotium-choice/>