

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

Blottum

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

Call us: 800-304-5357

Antibody Number P033 Antibody reactivity (target) Biotin Biotium Choice Antibody Primary Host species Mouse Clonality Recombinant Monoclonal Clone rBN-34
Biotium Choice Antibody Primary Host species Mouse Clonality Recombinant Monoclonal Clone rBN-34
Host species Mouse Clonality Recombinant Monoclonal Clone rBN-34
Clonality Recombinant Monoclonal Clone rBN-34
Clone rBN-34
151101
0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0
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 Flow, IHC, WB, IF (verified) application
Secondary/tag antibody Flow cytometry, IHC, IF (cells or tissue sections), Western blot applications
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®488A	490/516 nm	100 ug/mL	1 mL	P033-488A-1ML	CF®488A Features
<u>CF®568</u>	562/584 nm	100 ug/mL	1 mL	P033-568-1ML	CF®568 Features
CF®594	593/615 nm	100 ug/mL	1 mL	P033-594-1ML	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	1 mL	P033-640R-1ML	CF®640R Features
CF®647	652/668 nm	100 ug/mL	1 mL	P033-647-1ML	CF®647 Features
<u>CF®660R</u>	662/682 nm	100 ug/mL	1 mL	P033-660R-1ML	CF®660R Features
CF®680R	680/701 nm	100 ug/mL	1 mL	P033-680R-1ML	CF®680R Features
<u>CF®740</u>	742/767 nm	100 ug/mL	1 mL	P033-740-1ML	CF®740 Features
CF®750	755/779 nm	100 ug/mL	1 mL	P033-750-1ML	CF®750 Features
<u>CF®770</u>	770/797 nm	100 ug/mL	1 mL	P033-770-1ML	CF®770 Features

This datasheet was generated on November 23, 2025 at 03:16:45 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/