

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 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](#)

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

Product attributes

Antibody type	Biotium Choice, Primary, Tag Antibody
Antibody Number	P033
Antibody reactivity (target)	Biotin
Biotium Choice Antibody	Primary, Tag Antibody
Host species	Mouse
Clonality	Monoclonal, 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
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

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
CF@568	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
CF@660R	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
CF@740	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
CF@770	770/797 nm	100 ug/mL	1 mL	P033-770-1ML	CF@770 Features
CF@790	783/808 nm	100 ug/mL	1 mL	P033-790-1ML	CF@790 Features

This datasheet was generated on June 7, 2026 at 08:36:40 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/>