



Glowing products for science

## Neurofilament-H Recombinant Monoclonal Mouse Antibody (rRmdO-20) – Biotium Choice

A recombinant mouse monoclonal antibody that recognizes a nonphosphorylated epitope in the tail domain of Neurofilament-H. The antibody belongs to the Biotium Choice list of select antibodies that have been validated in-house for optimal performance.

### Product Description

Neurofilament-H Recombinant Monoclonal Mouse Antibody (rRmdO-20) is a recombinant mouse monoclonal antibody that recognizes a nonphosphorylated epitope in the tail domain of Neurofilament-H. Dephosphorylation of NF-H will increase immunoreactivity with this antibody. 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 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 and CF®790](#)
- Suitable for immunohistochemistry, immunofluorescence, and western

Neurofilaments, neuron-specific intermediate filament proteins, maintain neuronal caliber and consist mainly of NF-L, NF-M, and NF-H. Their expression is tissue- and development-specific. NF-H is crucial in mature axons and implicated in amyotrophic lateral sclerosis (ALS).

### 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](#)

## Neurofilament-H Recombinant Monoclonal Mouse Antibody (rRmdO-20) – Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
<a href="#">CF®488A</a>	490/516 nm	100 ug/mL	500 uL	<a href="#">P035-488A-500UL</a>	<a href="#">CF®488A Features</a>
<a href="#">CF®568</a>	562/584 nm	100 ug/mL	500 uL	<a href="#">P035-568-500UL</a>	<a href="#">CF®568 Features</a>
<a href="#">CF®594</a>	593/615 nm	100 ug/mL	500 uL	<a href="#">P035-594-500UL</a>	<a href="#">CF®594 Features</a>
<a href="#">CF®640R</a>	642/663 nm	100 ug/mL	500 uL	<a href="#">P035-640R-500UL</a>	<a href="#">CF®640R Features</a>
<a href="#">CF®647</a>	652/668 nm	100 ug/mL	500 uL	<a href="#">P035-647-500UL</a>	<a href="#">CF®647 Features</a>
<a href="#">CF®680R</a>	680/701 nm	100 ug/mL	500 uL	<a href="#">P035-680R-500UL</a>	<a href="#">CF®680R Features</a>
<a href="#">CF®740</a>	742/767 nm	100 ug/mL	500 uL	<a href="#">P035-740-500UL</a>	<a href="#">CF®740 Features</a>
<a href="#">CF®790</a>	783/808 nm	100 ug/mL	500 uL	<a href="#">P035-790-500UL</a>	<a href="#">CF®790 Features</a>



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

### Product attributes

Antibody Number	P035
Antibody reactivity (target)	Neurofilament H
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Primary, Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	rRmdO-20
Isotype	IgG1, kappa
Molecular weight	200 kDa
Synonyms	CMT2CC; I79_012512; KIAA0845; mKIAA0845; NEFH; NF-H; NF200; NFH
Human gene symbol	NEFH
Entrez gene ID	24587
SwissProt	P16884
Immunogen	Rat neurofilaments
Antibody target cellular localization	Cytoskeleton
Verified antibody applications	IF (verified), IHC (frozen) (verified)
Species reactivity	Bovine, Hamster, Human, Mollusc, Mouse, Rabbit, Rat
Expected antibody applications	IHC (published for clone), WB (published for clone)
Antibody application notes	Immunofluorescence (HeLa): 1-5 ug/mL Optimal concentration to be determined by end-user.
Positive control	Brain
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/conjugate formulation	Conjugates: PBS/2 mg/mL rBSA/0.05% azide
Antibody research areas	Neuroscience
Cell/tissue expression	Brain, Neurons
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line

This datasheet was generated on February 3, 2026 at 02:41:16 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/neurofilament-h-recombinant-monoclonal-mouse-antibody-rmdo-20-biotium-choice/>