

Vimentin Recombinant Monoclonal Mouse Antibody (rV9) – Biotium Choice



A recombinant mouse monoclonal antibody that recognizes Vimentin. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.

Product Description

Vimentin Recombinant Monoclonal Mouse Antibody (rV9) is a recombinant mouse monoclonal antibody that recognizes Vimentin. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance. The antibody is available conjugated to a selection of CF® Dyes and HRP.

- Available in 7 bright and photostable CF® Dyes, including [near-IR CF®740](#), and HRP.
- Suitable for immunohistochemistry, immunofluorescence, and western

Vimentin is a developmentally regulated intermediate filament protein expressed in mesenchymal cells and early cell development, where it supports intracellular protein transport and is later replaced by tissue-specific IFPs. It also plays a key role in steroid synthesis by serving as a storage network for cholesterol-rich lipid droplets. Phosphorylation of vimentin leads to filament breakdown and cytoskeletal rearrangement, enabling lipid droplet-mitochondria interactions that promote cholesterol transport and increase steroid production.

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](#)

Vimentin Recombinant Monoclonal Mouse Antibody (rV9) – Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	100 ug/mL	500 uL	P044-488A-500UL	CF®488A Features
CF®568	562/584 nm	100 ug/mL	500 uL	P044-568-500UL	CF®568 Features
CF®594	593/615 nm	100 ug/mL	500 uL	P044-594-500UL	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	500 uL	P044-640R-500UL	CF®640R Features
CF®647	652/668 nm	100 ug/mL	500 uL	P044-647-500UL	CF®647 Features
CF®680R	680/701 nm	100 ug/mL	500 uL	P044-680R-500UL	CF®680R Features
CF®740	742/767 nm	100 ug/mL	500 uL	P044-740-500UL	CF®740 Features
CF®790	783/808 nm	100 ug/mL	500 uL	P044-790-500UL	CF®790 Features
HRP	N/A	100 ug/mL	500 uL	P044-HRP-500UL	

Product attributes

Antibody number	P044
Antibody reactivity (target)	Vimentin
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice, Primary
Host species	Mouse
Clonality	Monoclonal, Recombinant Monoclonal
Clone	rV9
Isotype	IgG1, kappa
Molecular weight of antigen	~54 kDa
Synonyms	VIM; CTRCT30; HEL113
Human gene symbol	VIM
SwissProt	P08670
Antibody target cellular localization	Cytoskeleton
Verified antibody applications	IF (verified), IHC (frozen) (verified)
Species reactivity	C. Elegans, Cat, Chicken, Cow, Dog, Gerbil, Hamster, Horse, Human, Mouse, Non-human primates, Pig, Rabbit, Rat
Antibody application notes	Immunofluorescence: 1-5 ug/mL; Optimal concentration to be determined by end-user
Positive control	HeLa cells.
Shipping condition	Room temperature
Storage Conditions	Store CF® Dye conjugates at 2 °C to 8 °C, Store HRP conjugates at -10 °C to -35 °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, HRP conjugates: PBS/50% glycerol/2 mg/mL rBSA
Antibody research areas	Cancer
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line
Expected antibody applications	IHC (FFPE) (published for clone), WB (published for clone)

This datasheet was generated on June 7, 2026 at 08:25:01 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/vimentin-recombinant-monoclonal-mouse-antibody-rv9-biotium-choice/>