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.

statum 8

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. They are supplied in PBS, 2 mg/mL rBSA, and 0.05% sodium azide.

- Available in 7 bright and photostable CF® Dyes, including near-IR CF®740
- 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

Product attributes

| i ioddci attiibates | | | |
|---------------------------------------|--|--|--|
| Antibody number | P044 | | |
| Antibody reactivity (target) | Vimentin | | |
| Biotium Choice Antibody | Primary | | |
| Antibody type | Primary | | |
| Host species | Mouse | | |
| Clonality | Recombinant Monoclonal | | |
| Clone | rV9 | | |
| Isotype | IgG1, kappa | | |
| Molecular weight | ~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 conjugates at 2 °C 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 | 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) | | |
| | | | |

Call us: 800-304-5357 Email: btinfo@biotium.com

Vimentin Recombinant Monoclonal Mouse Antibody (rV9) - Biotium Choice

| Conjugation | Ex/Em | Conc. | Size | Catalog No. | Dye Features | | |
|--|---|---|--------|-----------------|---------------------|--|--|
| CF@501R This datasheet was genera | 562/565 nm ted on November 6, 2025 at 0 | 100 ug/ml 2:18:33 PM. Visit product pa | 500 uL | P044-58675500UL | CF®669 FF Eatangses | | |
| Product link: https://biotium.com/product/vimentin-recombinant-monoclonal-mouse-antibody-ry9-biotium-choice/ | | | | | | | |