

Vimentin monoclonal antibody

Monoclonal antibody to Vimentin for use with IHC applications

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| ENZ-ABS694-1000 | 1ml |
| ENZ-ABS694-0100 | 100µl |

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC

Application Notes For IHC applications: Dilute the concentrate as appropriate in antibody diluent, and then apply 150 µL to the slide, as indicated in the user's IHC protocol.

Formulation Liquid. In Tris buffer, pH 7.5, containing 1% BSA and <0.1% sodium azide.

Host Mouse

Immunogen Recombinant vimentin.

Isotype IgG

Purity Detail Affinity-purified.

Recommendation Dilutions/Conditions Immunohistochemistry (1:50-1:100) Optimal conditions must be determined individually for each application.

Species Reactivity Human

UniProt ID P08670

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)