

Vimentin (human) monoclonal antibody (65E)

Vimentin is a the type III intermediate filament (IF) protein expressed in a variety of mesenchymal and non-mesenchymal cells, including muscle cells, astrocytes, neurons, endothelial cells and fibroblasts. During development of the mammalian nervous system vimentin is transiently expressed in neuronal precursor cells, often with the type VI IF, nestin. Vimentin appears to play a role in many intracellular processes, recent examples from the literature include adhesion, glial scar formation, neuronal migration, and the normal development of Bergmann glia.

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

Ordering Information

[Order Online »](#)

BML-VG6100-0100	100µl
------------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC, WB

Clone 65E

Host Mouse

Immunogen Electrophoretically purified vimentin from human cultured fibroblasts.

Isotype IgG1

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-
be@enzolifesciences.com](mailto:info-be@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)