

[pSer55]Vimentin monoclonal antibody (4A4)

Vimentin is one of four type II intermediate filament (IF) proteins; Desmin, Glial fibrillary acidic protein (GFAP) and Peripherin being the others. Vimentin is the most widely distributed of all IF proteins. It is typically expressed in leukocytes, blood vessel endothelial cells, and mesenchymal cells such as fibroblasts. Vimentin filaments help support cellular membranes. Vimentin networks also may help keep the nucleus and other organelles in a defined place within the cell.

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

Citations: 5

[View Online »](#)

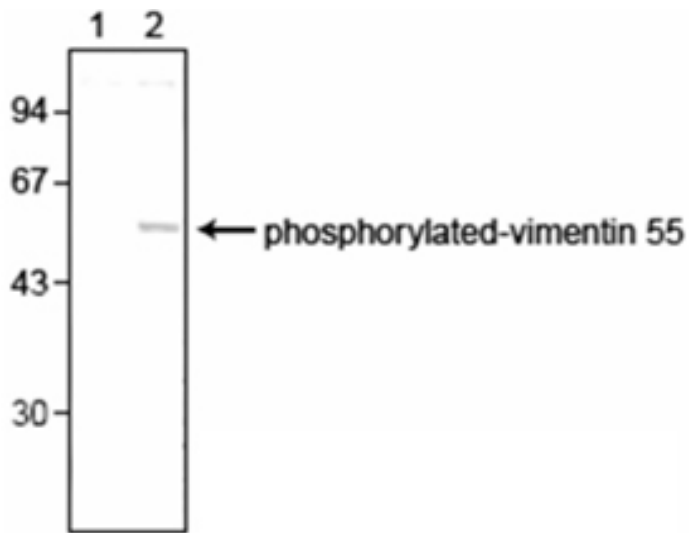
Ordering Information

[Order Online »](#)

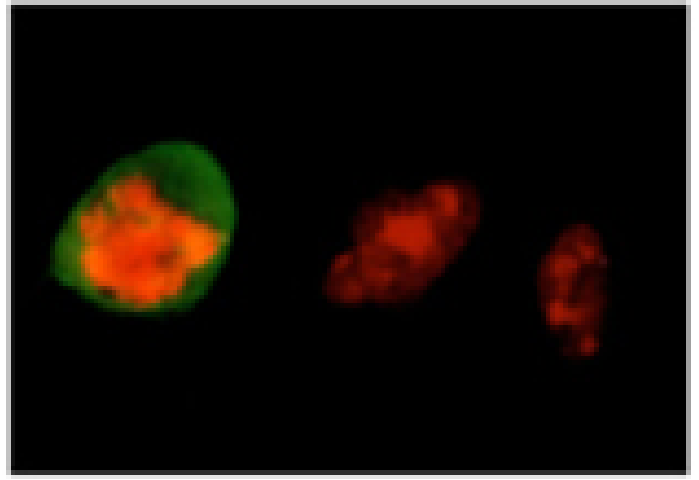
ADI-KAM-CC249-E	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of U251 interphase (1) and U251 M phase (2) cells, probed with Vimentin (pSer55) mAb (4A4).



Immunocytochemistry analysis of U251 cells using Vimentin (pSer55) mAb (4A4).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	VIM
Application	ELISA, ICC, WB
Application Notes	Detects a band of ~57kDa by Western blot.
Clone	4A4
Formulation	Liquid. In PBS containing 50% glycerol.
Host	Mouse
Immunogen	Synthetic peptide corresponding to a portion of mouse vimentin phosphorylated at Ser55.
Isotype	IgG2b
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	ELISA, Immunocytochemistry (1µg/ml)Western Blot (1-5µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	P20152

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)

Last modified: May 29, 2024