

[pSer71]Vimentin monoclonal antibody (TM71)

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](#).

Ordering Information

[Order Online »](#)

ADI-KAM-CC250-E	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

kDa

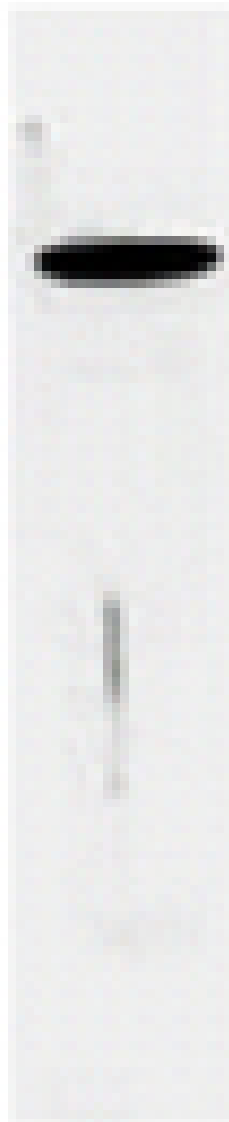
97 -

66 -

43 -

30 -

22 -



Western blot analysis of phosphorylated vimentin (Ser71) expression in U251 using [pSer71]Vimentin, mAb (TM71).



Immunocytochemistry analysis of phosphorylated vimentin (Ser71) on 4% PFA fixed U251 cells with [pSer71]Vimentin, mAb (TM71).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	VIM
Application	ICC, WB
Application Notes	Detects a band of ~58kDa by Western blot.
Clone	TM71
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol.
Host	Rat
Immunogen	Synthetic peptide corresponding to aa 66-77 of human Vimentin phosphorylated at Ser71.
Isotype	IgG2a
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (1µg/ml)Western Blot (1µg/ml for chemiluminescence detection system)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	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)