

EpCAM (human) monoclonal antibody (VU-1D9)

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Epithelial cell adhesion molecule, Epithelial specific antigen, ESA, CD326
Application	Flow Cytometry, FUNC, IHC (FS), IHC (PS), IP, WB
Application Notes	Functional Application: Blocking.
Clone	VU-1D9
Formulation	Liquid. In PBS.
Host	Mouse
Immunogen	Small cell lung carcinoma cells.
Isotype	IgG1
Purity Detail	Protein A-affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections, microwave treatment not necessary)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Specificity	Recognizes human EpCAM (epithelial specific antigen), a 40 kDa epithelial protein expressed on baso-lateral cell surface in many epithelial tissues (but absent from mesothelial tissue). Differentiates between mesothelioma and carcinoma.

UniProt ID

P16422

Worry-free Guarantee

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

Last modified: May 29, 2024



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)