

# EpCAM Recombinant monoclonal antibody (rVU-1D9)

Epithelial cellular adhesion molecule (EpCAM), is also identified as epithelial specific antigen (ESA) and the 40kDa transmembrane epithelial glycoprotein EGP40. EpCAM is expressed on the baso-lateral cell surface in most simple epithelia and a vast majority of carcinomas. EpCAM antibody is also useful in distinguishing serous carcinomas of the ovary from mesothelioma.

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

## Ordering Information

[Order Online »](#)

ENZ-ABS318-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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, IF, IHC (PS), WB

**Application Notes** The VU-1D9 antibody has been used to distinguish adenocarcinoma from pleural mesothelioma and hepatocellular carcinoma.

**Clone** rVU-1D9

**Formulation** Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.

**Gene/Protein Identifier** 4072 (Gene ID)

**Host** Mouse

**Immunogen** Small cell lung carcinoma cells.

**Isotype** IgG1k

**Purity Detail** Protein G affinity purified.

**Recommendation** Flow Cytometry (0.5-1 µg/106 cells) Immunofluorescence (1-2 µg/mL) Immunohistochemistry (formalin-fixed paraffin sections) (0.5-1 µg/mL for 30 min at RT) Western Blot (0.5-1 µg/mL) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Species Reactivity** Human

**UniProt ID** P16422



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

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

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

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

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

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

This antibody is covered by our **Worry-Free Guarantee**.