EpCAM Recombinant monoclonal antibody (rMOC-31)

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-ABS317-0100

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 NotesThis antibody has been used to distinguish

adenocarcinoma from pleural mesothelioma and hepatocellular carcinoma. This antibody is also useful in

distinguishing serous carcinomas of the ovary from

mesothelioma.

Clone rMOC-31

Formulation Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced)

and 0.05% sodium azide.

Host Mouse

Immunogen Neuraminidase treated GLS-1 human small cell lung

carcinoma cells.

lgG1κ

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (0.5-1 μg/106 cells in 0.1

mL)Immunofluorescence (1-2 μ g/mL)Western blot (0.5-1 μ g/mL)Immunohistochemistry (formalin-fixed paraffin sections) (0.5-1 μ g/mL for 30 min at RT)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined

individually for each application.

Species Reactivity Human

Specificity

Binding epitope of this antibody is located in the first EGFlike repeat domain (EGF1) between amino acids 27-59 of Ep-CAM.

UniProt ID

P16422

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.