

EpCAM Recombinant monoclonal antibody (EGP40/1555R)

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](#).

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	This antibody has been used to distinguish adenocarcinoma from pleural mesothelioma and hepatocellular carcinoma. It is also useful in distinguishing serous carcinomas of the ovary from mesothelioma.
Clone	EGP40/1555R
Formulation	Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide (around aa 20-60) from the N-terminus of human EpCAM.
Isotype	IgGk
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (0.5-1 µg/10 ⁶ cells in 0.1 mL)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

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)