

# EpCAM Recombinant monoclonal antibody (EGP40/1556R)

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
Clone	EGP40/1556R
Formulation	Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
Host	Rabbit
Immunogen	Recombinant human EpCAM.
Isotype	IgGκ
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (0.5-1 µg/10 <sup>6</sup> 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 <a href="#">Worry-Free Guarantee</a>

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)