

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](#).

Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| ENZ-ABS365-0100 | 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 |
| 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 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 |



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)