

EpCAM recombinant monoclonal antibody (002)

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

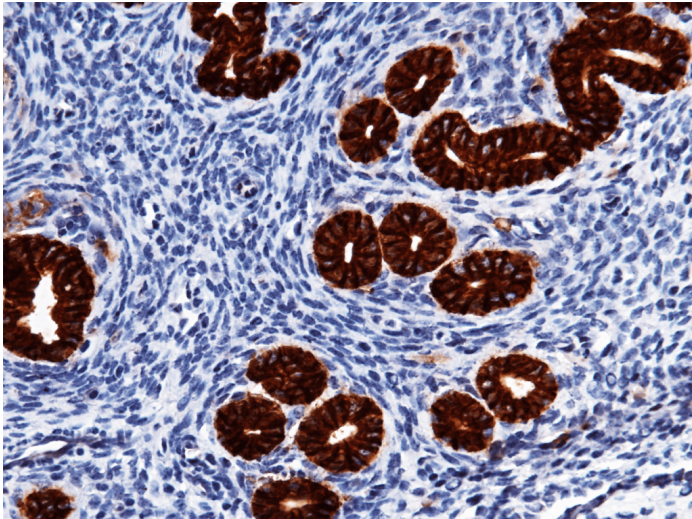
Ordering Information

[Order Online »](#)

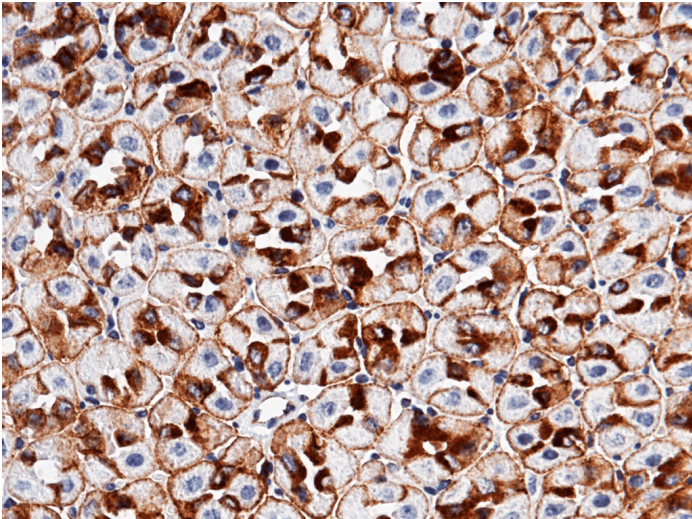
ENZ-ABS1021-0100	100µl
------------------	-------

Manuals, SDS & CofA

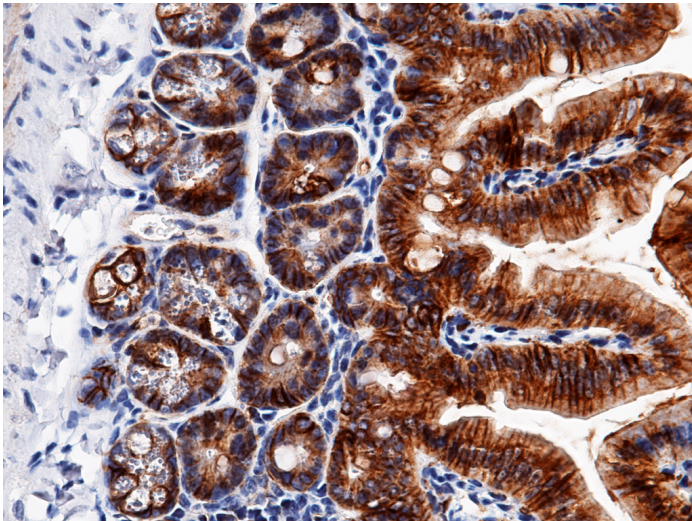
[View Online »](#)



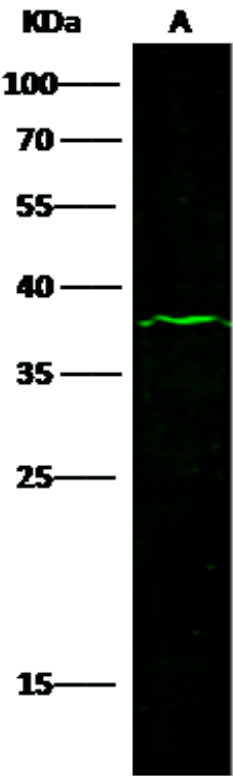
Immunochemical staining of mouse EpCAM in mouse uterus with rabbit monoclonal antibody at 1:1000 dilution, formalin-fixed paraffin embedded sections.



Immunochemical staining of mouse EpCAM in mouse stomach with rabbit monoclonal antibody at 1:1000 dilution, formalin-fixed paraffin embedded sections.



Immunochemical staining of mouse EpCAM in mouse intestine with rabbit monoclonal antibody at 1:1000 dilution, formalin-fixed paraffin embedded sections.



Anti-EPCAM rabbit monoclonal antibody at 1:500 dilution Lane A: 293T Whole Cell Lysate. Lysates/proteins at 30 µg per lane. Secondary Goat Anti-Rabbit IgG H&L (Dylight800) at 1/10000 dilution. Developed using the Odyssey technique. Performed under reducing conditions. Observed band size:38 kDa.

Handling & Storage

Handling Protected from light. Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Epithelial cell adhesion molecule, TROP-1, TACSTD1

Application ELISA, IHC (PS), WB

Clone 002

Formulation Liquid. 0.2 µm filtered solution in PBS. Preservative free.

Host Rabbit

Immunogen Recombinant mouse EpCAM / TROP-1 / TACSTD1 protein.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions Western Blot (1:500-1:1'000)Immunohistochemistry (paraffin sections) (1:500-1:2'000)ELISA (1:25'000-1:5'0000)ELISA (Detectino)(1:5'000-1:5'0000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Mouse

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



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

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

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

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

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

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

Last modified: September 17, 2024