

# E-Cadherin recombinant monoclonal antibody (032)

The anti-E-Cadherin (CDH1/CD324) antibody is specifically designed to detect E-Cadherin, a calcium-dependent cell adhesion protein critical for maintaining epithelial tissue structure and integrity.

E-Cadherin (epithelial cadherin), also known as CDH1 or CD324, is a calcium-dependent cell adhesion molecule crucial for maintaining epithelial tissue integrity and polarity. It belongs to the cadherin superfamily and mediates cell-cell adhesion through homophilic interactions, playing a central role in tissue architecture, cellular communication, and signaling. E-Cadherin is a key regulator of epithelial-to-mesenchymal transition (EMT), a process critical during embryonic development, wound healing, and cancer metastasis. Loss or dysfunction of E-Cadherin disrupts cell adhesion, contributing to increased cell motility and invasion, often correlating with poor prognosis in cancers such as breast, gastric, and ovarian carcinomas.

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

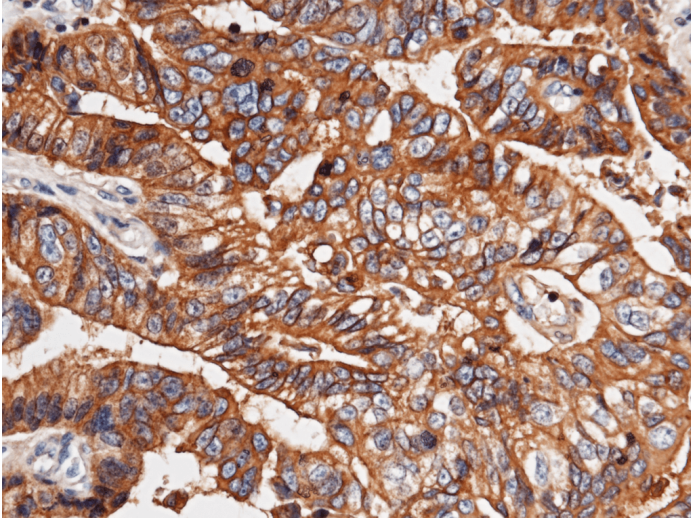
## Ordering Information

[Order Online »](#)

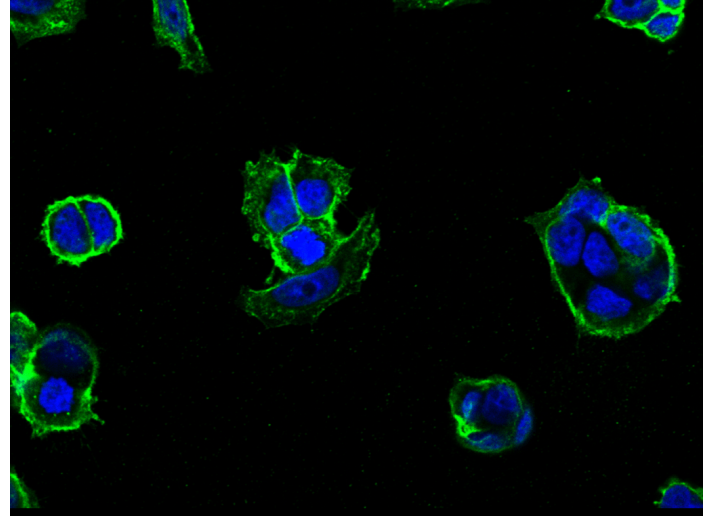
ENZ-ABS1046-0100	100µl
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Immunohistochemical staining of human E-cadherin in human gastric cancer with rabbit monoclonal Antibody (1:200, formalin-fixed paraffin embedded sections). Positive staining was localized to membrane of epithelial cells.



Immunofluorescence staining of CD324 in MCF7 cells. Cells were fixed with 4% PFA, permeabilized with 0.3% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit anti-human CD324 monoclonal antibody (1:60) at 4°C overnight. Then cells were stained with the Alexa Fluor® 488-conjugated Goat Anti-rabbit IgG secondary antibody (green) and counterstained with DAPI (blue).

# Handling & Storage

Handling	Avoid freeze-thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	CDH1, Cadherin-1, E-Cad, Uvomorulin, CAM 120/80, Epithelial cadherin, CD324
Application	ICC, IF, IHC (PS)
Clone	032
Formulation	Liquid. 0.2 µm filtered solution in PBS. Carrier-free. Preservative-free.
Host	Rabbit
Immunogen	Recombinant human E-cadherin.
Isotype	IgG
Purity Detail	Protein A purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections) (1:100-1:500) Immunocytochemistry/Immunofluorescence (1:20-1:100)  Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .

Last modified: January 9, 2025



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)