

CA-125 monoclonal antibody

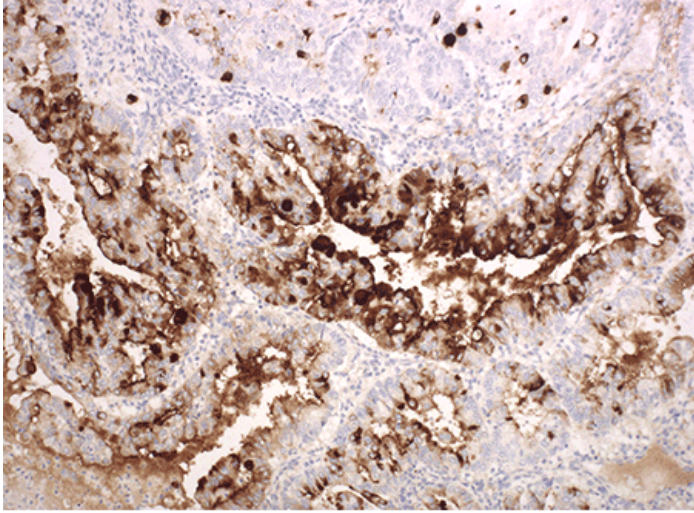
Monoclonal antibody to CA-125 for use with IHC applications

CA-125 is a key marker for ovarian carcinomas. It reacts with a majority of epithelial ovarian neoplasms of various types including serous, endometrioid, clear cell and undifferentiated but is also found in other neoplasms.

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

Manuals, SDS & CofA

[View Online »](#)



Formalin-fixed paraffin-embedded endometrial tissue stained using Enzo's CA-125 antibody (Prod. No. ENZ-ABS775). IHC staining was performed on an automated Leica BOND III.

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Cancer antigen 125, Mucin-16
Application	IHC
Application Notes	<p>The CA-125 antibody can be used to qualitatively identify the presence of CA-125 protein in FFPE tissue sections using IHC test methods.</p> <p>For IHC applications: Dilute the concentrate as appropriate in antibody diluent, and then apply 150 µL to the slide, as indicated in the user's IHC protocol.</p>
Formulation	Liquid. In Tris buffer, pH 7.5, containing 1% BSA and <0.1% sodium azide.
Host	Mouse
Immunogen	Epitope group B1
Isotype	IgG
Purity Detail	Affinity-purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (1:50-1:200)Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q8WXI7

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)