

Cathepsin D recombinant monoclonal antibody (010)

This monoclonal antibody detects Cathepsin D (CTSD), a lysosomal protease involved in protein degradation and apoptosis.

Cathepsin D is a key aspartic protease in the lysosome, critical for protein catabolism and cellular homeostasis. It has been implicated in neurodegenerative diseases and cancer progression.

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

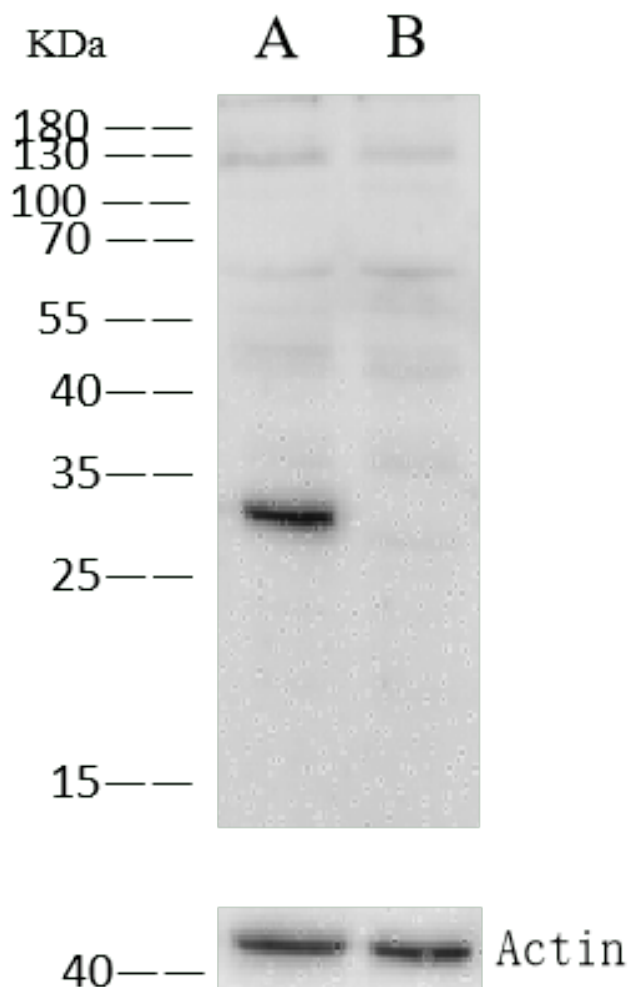
Ordering Information

[Order Online »](#)

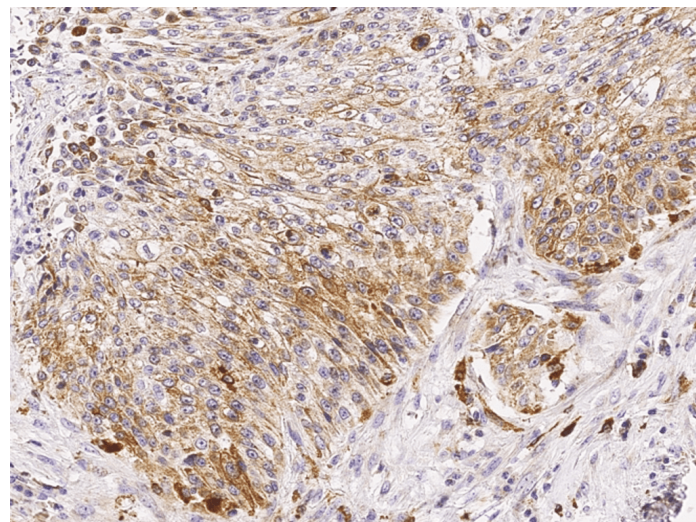
ENZ-ABS1070-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis using cathepsin D rabbit monoclonal antibody (ENZ-ABS1070) at 1:500 dilution. Secondary Antibody: Goat Anti-Rabbit IgG (H+L)/HRP at 1/10000 dilution. Lane A: 293T Whole Cell Lysate, Lane B: CTSD knockout 293T Whole Cell Lysate. Lysates/proteins at 30 µg per lane.



Immunochemical staining of human cathepsin D in human lung cancer with rabbit monoclonal antibody (ENZ-ABS1070, 1:300, formalin-fixed paraffin embedded sections)

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 CTSD

Application IHC (PS), WB

Clone 010

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

Host Rabbit

Immunogen Recombinant human Cathepsin D.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Western Blot (1:500-1:2'000)

Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:100-1:500)

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-
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-
fr@enzolifesciences.com](mailto:info-fr@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)