# 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**

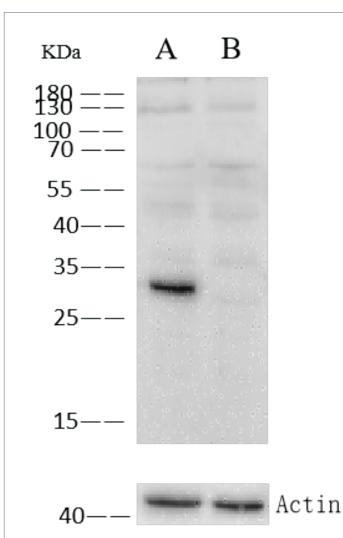
**Order Online** »

**ENZ-ABS1070-0100** 

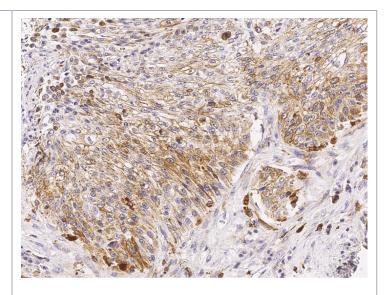
100µl

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 konckout 293T Whole Cell Lysate. Lysates/proteins at 30  $\mu$ g 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.

lgG lgG

Purity Detail Protein A purified.

**Recommendation Dilutions/Conditions**Western Blot (1:500-1:2'000)

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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com