

# Cathepsin L (human liver), (purified)

Cathepsin L is a lysosomal protease that can also be secreted from the cell and is involved in protein degradation, arthritis, and cancer.

Citations: 6

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-SE201-0025	25µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -80°C

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application Notes** Useful for inhibitor screening, kinetic studies, and cleavage of substrates *in vitro* or in tissue culture.

**Purity** ≥95% by SDS-PAGE

**Source** Isolated from human liver. Negative for HbsAg, anti-HBc, anti-HCV, and anti-HIV-1 and -2.

**UniProt ID** P07711 (cathepsin L1)



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)