

Cathepsin D (human liver), (purified)

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

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE199-0025	25µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquotes and store at -80°C.
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 <i>in vitro</i> or in tissue culture.
Formulation	Lyophilized with phosphate buffer.
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute just before first use with original volume of water.
Source	Isolated from human liver.
UniProt ID	P07339

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)