

Cathepsin B (human liver), (purified)

Highly pure lysosomal protease

Cathepsin B is a lysosomal protease (endopeptidase and dipeptidyl carboxypeptidase) that can also be secreted from the cell and is involved in protein degradation, apoptosis, arthritis, cancer, and Alzheimer's disease.

Citations: 16

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE198-0025	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.

Formulation Liquid. In 50mM sodium acetate, pH 5.0, containing 1mM EDTA.

MW ~27.5 kDa

Purity ≥95% (SDS-PAGE)

Source Purified from human liver.

UniProt ID P07858



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)