

Cathepsin S (human), (recombinant)

Highly active

Cathepsin S, a member of the papain family of lysosomal cysteine proteases, functions in the immune system.

Citations: 7

[View Online »](#)

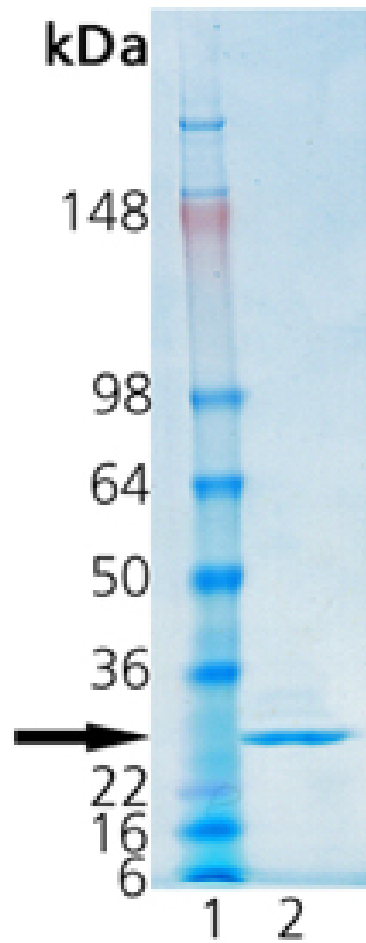
Ordering Information

[Order Online »](#)

BML-SE453-0010	10µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 0.5 μ g of Human Cathepsin S Protein

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application Notes

Useful to study enzyme kinetics, cleave target substrates, and screen for inhibitors.

Formulation

Liquid. In 100 mM MES, pH 6.5, containing 1 mM EDTA, 50 mM L-cysteine, 10 mM DTT, 0.5% Triton[®] X-100, 30% glycerol.

MW

26 kDa

Purity

≥95% (SDS-PAGE)

Purity Detail

Purified by multi-step chromatography.

Source

Produced in insect cells. Recombinant active cathepsin S cloned from human cDNA as full length proenzyme, auto-activated by incubation in pH 4.5 acetate buffer, and purified. Produced in a baculovirus expression system.

Specific Activity

≥20 U/mg protein. One unit will hydrolyze one μmole Z-Val-Val-Arg-AMC substrate (Prod. No. BML-P199, 40 μM) per minute at 37°C.

UniProt ID

P25774

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)