

Cathepsin F (human), (recombinant)

Highly active

Recombinant glycosylated procathepsin F cloned from human cDNA (NM_003793), expressed in insect cells, and purified as the active form of the enzyme. Cathepsin F, a member of the papain family of lysosomal cysteine proteases, acts upon proteins such as MHC Class II-associated invariant chain and ApoB-100. It is inhibited by the cysteine proteinase inhibitor E-64 (Prod. No. BML-EI105) and cystatin C (Prod. No. BML-SE479). It is a lysosomal protease expressed in numerous tissues, notably brain and macrophages. It functions in the immune system and is implicated in disease states such as atherosclerosis, cancer, and angiogenesis.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE568-0005	5µg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored as supplied at -80°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application Notes	Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors.
Formulation	Liquid. In 20mM NaOAc, pH 5.0, containing 2.5mM EDTA, 250mM sodium chloride and 20mM L-cysteine.
MW	~39kDa
Purity	≥80% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in insect cells (secreted) as untagged proenzyme; activated during purification. Produced in a baculovirus expression system.
Specific Activity	≥1 U/mg. One unit hydrolyzes 1 μmol Z-Phe-Arg-AMC substrate (OMNICATHEPSIN [®] Fluorogenic Substrate, Prod. No. BML-P139) per min. at 37°C.
UniProt ID	Q9UBX1

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-
fr@enzolifesciences.com](mailto:info-fr@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)