

# 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

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

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Application Notes</b>	Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors.
<b>Formulation</b>	Liquid. In 20mM NaOAc, pH 5.0, containing 2.5mM EDTA, 250mM sodium chloride and 20mM L-cysteine.
<b>MW</b>	~39kDa
<b>Purity</b>	≥80% (SDS-PAGE)
<b>Purity Detail</b>	Purified by multi-step chromatography.
<b>Source</b>	Produced in insect cells (secreted) as untagged proenzyme; activated during purification. Produced in a baculovirus expression system.
<b>Specific Activity</b>	≥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.
<b>UniProt ID</b>	Q9UBX1



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)