

Caspase-1 (human), (recombinant)

Highly active caspase essential for apoptosis.

First of the caspases described, caspase-1 was subsequently found to be homologous to Ced-3, the *C. elegans* caspase essential for apoptosis.

Citations: 15

[View Online »](#)

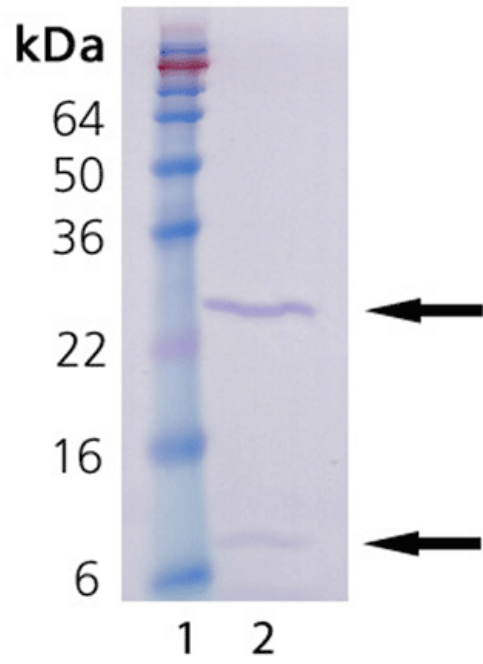
Ordering Information

[Order Online »](#)

BML-SE168-5000	5000U
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	After initial defrost, aliquot product into individual tubes and refreeze at -80°C. Avoid repeated freeze/thaw cycles. The enzyme is stable on ice for the time typically required to set up an experiment (30-60 min.), but may lose activity with prolonged storage on ice. It is recommended that thawing and dilution of the enzyme be done within as short a time as possible before start of the assay. The remaining, undiluted and unused enzyme should be refrozen quickly by, for example, snap-freezing in a dry ice ethanol bath or liquid nitrogen. The enzyme is stable to at least 4 freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Activity	100 U/μl
Alternative Name	Interleukin 1β converting enzyme, IL-1β converting enzyme, ICE
Application Notes	Useful tool to study enzyme regulation and kinetics, cleave target substrates, screen for inhibitors.
Formulation	Liquid. In 50mM HEPES, pH 7.4, containing 100mM sodium chloride, 0.5% CHAPS, 1mM EDTA, 10% glycerol and 10mM DTT.
MW	~20 + 10kDa
Purity	≥90% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> . cDNA encodes residues identical to Asn ¹²⁰ -His ⁴⁰⁴ (C-terminus) at Genbank Accession No. M87507, except for an Asp ³⁸¹ to Glu change, introduced to stabilize the enzyme against autoproteolysis.
Specific Activity	One U=1 pmol/min, using Ac-YVAD-pNA (200μM; Prod. No. ALX-260-026) as substrate, at 30°C.
UniProt ID	P29466



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)