

Caspase-3 (human), (recombinant)

The major effector caspase of apoptosis, caspase-3 is responsible for cleavage of multiple cellular targets.

Citations: 15

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE169-5000	5000U
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid 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	CPP32, Yama, Apopain
Application Notes	Can be used to study enzyme regulation and kinetics, cleave target substrates and screen for inhibitors.
Formulation	Liquid. In 50mM HEPES, pH 7.4, 100mM sodium chloride, 0.5% CHAPS, 1mM EDTA, 10% glycerol and 10mM dithiothreitol.
MW	17 + 12 kDa
Purity	≥95% (SDS-PAGE)
Source	Produced in <i>E. coli</i> . Caspase-3 (CPP32; Yama; apopain) from human cDNA. The enzyme was cleaved and activated from the proenzyme
Specific Activity	One U=1 pmol/min using Ac-DEVD-pNA (200 μM; Prod. No. ALX-260-033) as substrate at 30°C
UniProt ID	P42574

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