

# Caspase-3 (human), (recombinant)

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

Citations: 14

[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 *E. coli*. 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)