

# Caspase-6 (human), (recombinant)

Highly, active caspase-6 for the study of  
enzyme kinetics or screening of inhibitors

Effector caspase responsible for the cleavage of nuclear lamins in apoptosis. Caspase-6 cleavages of various other protein targets have been implicated in the pathogenesis of neurodegenerative diseases, including Alzheimer's disease (b-amyloid precursor protein), Huntington's disease (huntingtin).

Citations: 11

[View Online »](#)

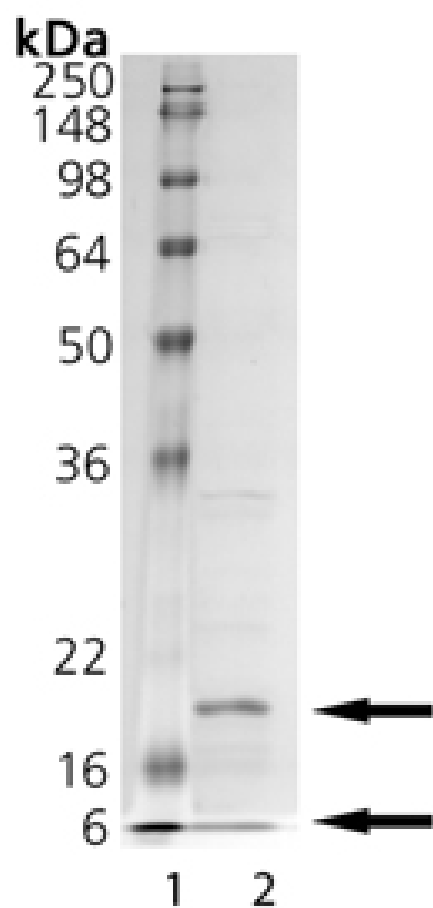
## Ordering Information

[Order Online »](#)

BML-SE170-5000	5000U
----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis – Lane 1: MW Marker, Lane 2: 1  $\mu$ g Caspase-6.

## Handling & Storage

Use/Stability	After initial defrost, aliquot product into individual tubes and refreeze the remaining, unused enzyme quickly by snap-freezing in a dry/ice ethanol bath or liquid nitrogen, if possible.
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	Mch2
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	18 + 11 kDa subunits
Purity	≥80% (SDS-PAGE)
Source	Produced in <i>E. coli</i> .
Specific Activity	One U=1 pmol/min at 30°C using Ac-VEID-pNA (200mM; Prod. No. BML-P437) as substrate.
UniProt ID	P55212

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)