

Caspase-8 (human), (recombinant)

Highly active caspase-8 for the study of
enzyme regulation and kinetics

Caspase-8 is an initiator caspase activated in the death-inducing signaling complexes (DISCs) of the TNF receptor superfamily (e.g. TRAIL, FasL, TNF α receptors).

Citations: 8

[View Online »](#)

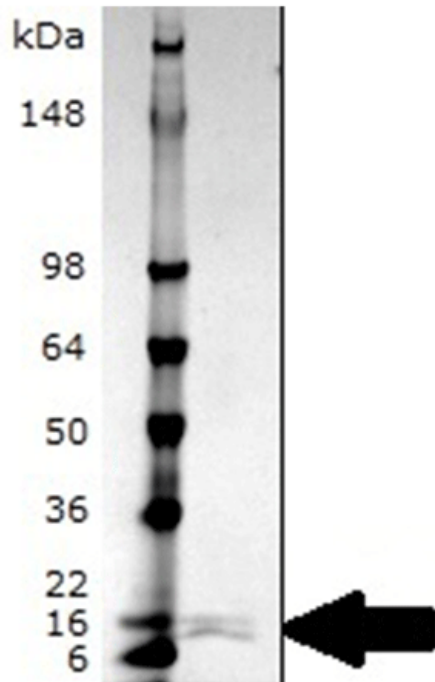
Ordering Information

[Order Online »](#)

BML-SE172-5000	5000U
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS PAGE Analysis: Lane 1: MW Marker, Lane 2: 1 μ g (Prod. No. BML-SE172).

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. Avoid repeated freeze/defrost cycles. NOTE: When stored under the above conditions, this enzyme is stable at the concentration supplied, in its current storage buffer. Procedures such as dilution of the enzyme followed by refreezing, could lead to loss of activity.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Activity	100 U/μl
Alternative Name	FLICE, Mch5, MACH
Application Notes	Useful tool to study enzyme regulation and kinetics, cleave target substrates, screen for inhibitors.
Formulation	Liquid. In 50mM HEPES, pH 7.4, 100mM sodium chloride, 0.5% CHAPS, 1mM EDTA, 10% glycerol and 10mM DTT.
MW	18 + 11 kDa
Purity	≥90% (SDS-PAGE)
Source	Produced in <i>E. coli</i>
Specific Activity	One U=1 pmol/min using Ac-IETD-pNA (200mM; Prod. No. ALX-260-045) as substrate, at 30°C.
UniProt ID	Q14790



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)