

Caspase-1 (human), (recombinant) (active)

Caspase-1 is the prototypical member of the caspase family of cysteine proteases. Caspase-1 exists in cells as an inactive 45kDa pro-enzyme. The pro-enzyme is matured by proteolysis to yield large (20kDa) and small (10kDa) subunits. The active enzyme is a heterotetramer consisting of two large and two small subunits. To date the mechanism of regulation of caspase-1 activation is complex and poorly understood. In THP-1 cells, a large proportion of the caspase-1 is present in the inactive pro-enzyme form.

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-056-U025	25U
ALX-201-056-U100	100U

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Interleukin 1 β converting enzyme, IL-1 β converting enzyme, ICE
Application Notes	Useful in studying enzyme regulation and inhibitors, determining target substrate, screening caspase inhibitors, or as a positive control in caspase assays. We recommend using 1 unit per assay for analyzing caspase activity.
Formulation	Lyophilized.
Quality Control	Routinely tested for its ability to cleave the caspase-1 colorimetric substrate Ac-YVAD- <i>p</i> NA (Prod. No. ALX-260-026).
Reconstitution	Reconstitute to 1U/ μ l in water
Source	Produced in <i>E. coli</i> . Contains an N-terminal His-tag.
Specific Activity	>5'000U/mg protein. One unit is defined as the amount of enzyme that cleaves 1nmol of the caspase substrate YVAD- <i>p</i> NA (<i>p</i> NA: <i>p</i> -nitroaniline) per hour at 37°C in a reaction solution containing 50mM HEPES, pH 7.2, 50mM sodium chloride, 0.1% CHAPS, 10mM EDTA, 5% glycerol and 10mM DTT.
UniProt ID	P29466

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)