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 1β converting enzyme, IL-1β converting

enzyme, ICE

Application NotesUseful 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-pNA (Prod. No. ALX-260-

026).

Reconstitution Reconstitute to 1U/µI in water

Source Produced in *E. coli*. 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-pNA (pNA: p-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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com