Caspase-3 (human), (recombinant) (active)

Highly active caspase essential for apoptosis

Caspase-3 is a member of the interleukin-1β converting enzyme (ICE) family of cysteine proteases. It exists in cells as an inactive 32kDa proenzyme. During apoptosis procaspase-3 is processed at aspartate residues by self-proteolysis and/or cleavage by upstream caspases, such as caspase-6, caspase-8 and granzyme B. The processed form of caspase-3 consists of large (17kDa) and small (11kDa) subunits which associate to form the active enzyme. Active caspase-3 has been shown involved in the proteolysis of several important molecules, such as poly(ADP-ribose) polymerase (PARP), the sterol regulatory element binding proteins (SREBPs), focal adhesion kinase (FAK) and others.

Citations: 17

View Online »

Ordering Information

Order Online »

ALX-201-059-U025	25U
ALX-201-059-U100	100U

Manuals, SDS & CofA

View Online »

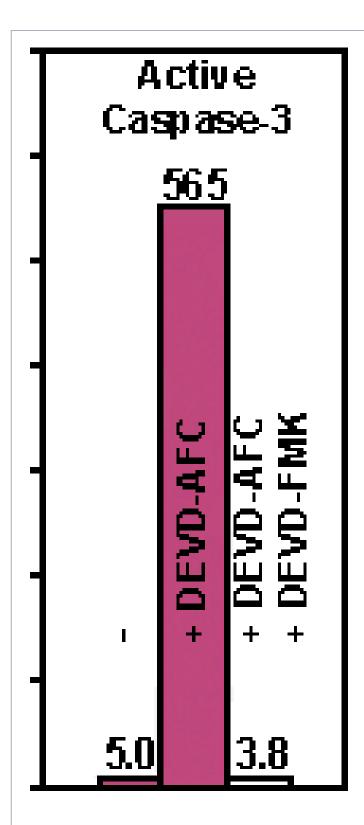


Figure: Active human caspase was expressed in *E. coli* and purified. The activity of recombinant caspase-3 was determined by cleaving AFC conjugates of DEVD. The cleavage activity was effectively inhibited by the corresponding peptide inhibitor as indicated.

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 CPP32, Yama, Apopain

Application NotesUseful in studying enzyme regulation, 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.

For a complete caspase-3 assay protocol, please refer to Caspase-3 Fluorometric or Colorimetric Assay Kits (Prod.

No. ALX-850-216 or ALX-850-215).

Formulation Lyophilized.

Quality Control Routinely tested for its ability to cleave the caspase-3

substrates Ac-DEVD-pNA (Prod. No. ALX-260-033) and

Ac-DEVD-AFC (Prod. No. ALX-260-032).

Reconstitution Reconstitute to 1U/μl in water.

Source Produced in E. coli. Contains an N-terminal His-tag.

Specific Activity ≥15,000 U/mg protein. One unit is defined as the amount

of enzyme that cleaves 1nmol of the caspase substrate DEVD-pNA (Prod. No. ALX-260-033) per hour at 37°C in a reaction solution containing 50mM HEPES, pH 7.2, 50mM NaCl, 0.1% CHAPS, 10mM EDTA, 5% glycerol and 10mM

DTT.

UniProt ID P42574



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