

# Caspase-8 (human), (recombinant) (active)

Caspase-8 is a member of the interleukin-1 $\beta$  converting enzyme (ICE) family of cysteine proteases. Same as other caspases, caspase-7 also exists in cells as an inactive proenzyme. During apoptosis procaspase-8 is processed at aspartate residues by self-proteolysis and/or cleavage by another caspase. The processed form of caspase-8 consists of large and small subunits which associate to form the active enzyme.

Citations: 6

[View Online »](#)

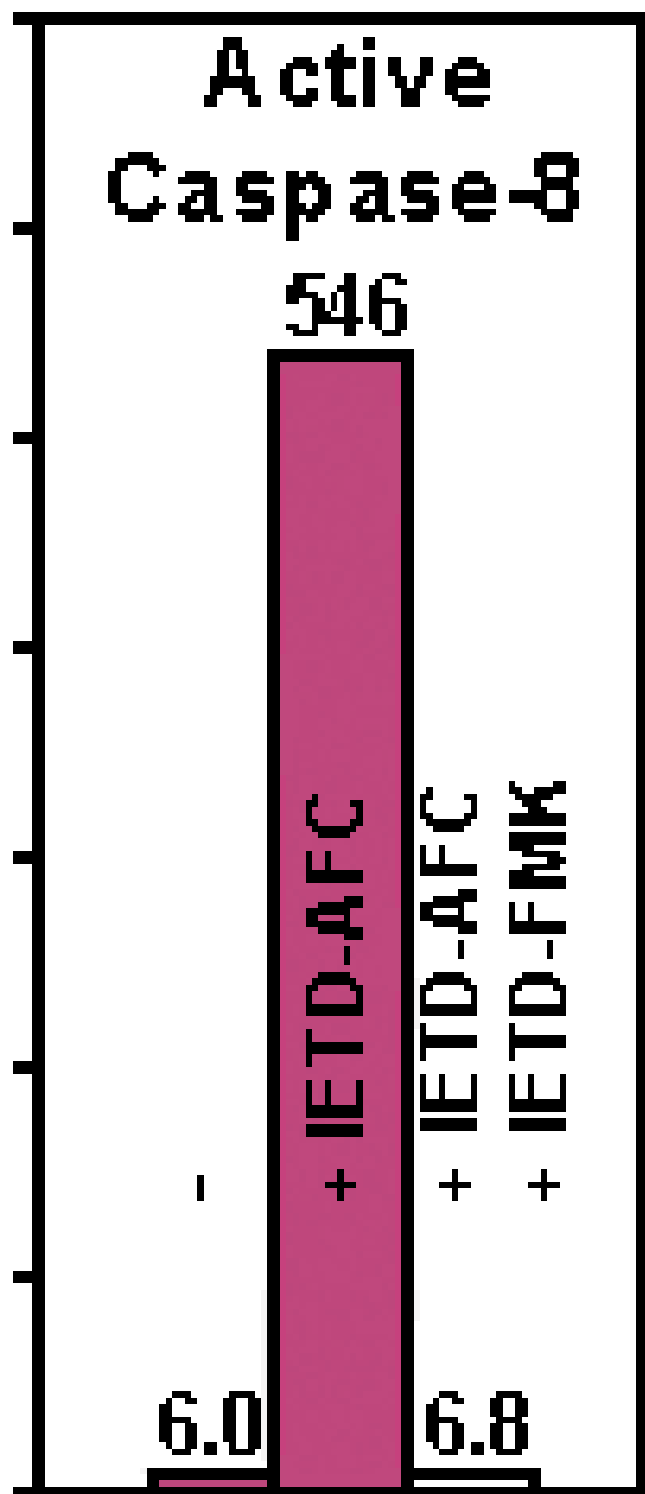
## Ordering Information

[Order Online »](#)

ALX-201-062-U100	100U
------------------	------

## Manuals, SDS & CofA

[View Online »](#)



**Figure:** Active human caspase was expressed in *E. coli* and purified. The activity of recombinant caspase-8 was determined by cleaving AFC conjugates of IETD. 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	FLICE, Mch5, MACH
Application Notes	Useful in screening caspase inhibitors, studying enzyme regulation, determining specificity of caspase substrates or as positive control in caspase activity assays. We recommend using 1U per assay for analyzing caspase activity.
Formulation	Lyophilized.
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute to 1U/μl with PBS containing 15% glycerol.
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 IETD-pNA 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	Q14790

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)