

# Caspase-4 (human), (recombinant) (active)

Caspase-4 is a member of the caspase family of cysteine proteases. It exists in cells as an inactive pro-enzyme. The pro-enzyme is matured by proteolysis to yield large and small subunits. The active enzyme is a heterotetramer consisting of two large and two small subunits. The expressed caspase-4 spontaneously undergoes auto-processing to yield subunits characteristic of the native enzyme. The active caspase-4 is routinely tested for its ability to enzymatically cleave the substrate WEHD-AFC or WEHD-pNA.

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-201-093A-U025	25U
ALX-201-093A-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	ICH-2, Protease TX, ICERel-II
Application Notes	Useful in screening caspase inhibitors, studying enzyme regulation and kinetics, determining target substrate or as a positive control in caspase activity assays. We recommend using 1 unit per assay for analyzing caspase activity.
Formulation	Lyophilized.
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with PBS containing 15% glycerol to 1U/μl.
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 WEHD-pNA per hour at 37°C in reaction solution containing 50mM HEPES, pH 7.2, 50mM NaCl, 0.1% CHAPS, 10mM EDTA, 5% glycerol and 10mM DTT.
UniProt ID	P49662



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-  
be@enzolifesciences.com](mailto:info-be@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)