

# Caspase-7 (human), (recombinant)

Caspase-7 acts as an effector caspase in apoptosis and has substrate preferences similar to caspase-3. It displays an affinity for ADP-ribose polymers and automodification of PARP-1 (ADP-ribosylation) makes it a better caspase-7 substrate.

Citations: 6

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                |       |
|----------------|-------|
| BML-SE177-5000 | 5000U |
|----------------|-------|

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

|                   |                                                                                                                        |
|-------------------|------------------------------------------------------------------------------------------------------------------------|
| Use/Stability     | After initial defrost, aliquot product into individual tubes and refreeze at -80°C. Avoid repeated freeze/thaw cycles. |
| Long Term Storage | -80°C                                                                                                                  |
| Shipping          | Dry Ice                                                                                                                |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                   |                                                                                                                   |
|-------------------|-------------------------------------------------------------------------------------------------------------------|
| Activity          | 100 U/μl                                                                                                          |
| Alternative Name  | Mch3, CMH-1                                                                                                       |
| Application Notes | Useful tool to study enzyme regulation and kinetics, cleave target substrates, screen for inhibitors.             |
| Formulation       | Liquid. In 50mM HEPES, pH 7.4, containing 100mM sodium chloride, 0.5% CHAPS, 1mM EDTA, 10% glycerol and 10mM DTT. |
| MW                | ~20 + 12kDa                                                                                                       |
| Purity            | ≥95% (SDS-PAGE)                                                                                                   |
| Source            | Produced in <i>E. coli</i> . Human caspase-7 (aa 24-303).<br>Genbank Accession No. U40281.                        |
| Specific Activity | One U=1 pmol/min at 30°C using Ac-DEVD-pNA (200μM; Prod. No. ALX-260-033) as substrate.                           |
| UniProt ID        | P55210                                                                                                            |

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-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)