

PRODUCT INFORMATION



Caspase-7 (human, recombinant)

Item No. 43692

Overview and Properties

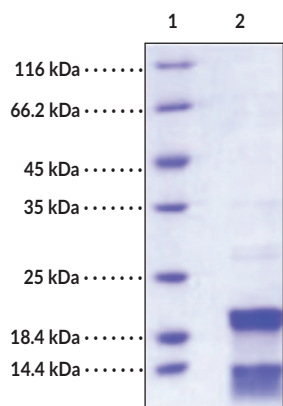
Synonyms: CASP-7, Apoptotic Protease Mch-3, CMH-1, ICE-like Apoptotic Protease 3, ICE-LAP3
Source: Recombinant human C-terminal His-tagged caspase-7 expressed in *E. coli*
Amino Acids: Met1; 24-303
Uniprot No.: P55210
Molecular Weight: 35 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: ≥90% estimated by SDS-PAGE
Supplied in: Lyophilized from sterile 20 mM HEPES, pH 7.5, with 100 mM sodium chloride, 1 mM EDTA, 0.10% sucrose, and 0.1% CHAPS

Protein

Concentration: *batch specific* mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers
Lane 2: Caspase-7

SDS-PAGE Analysis of Caspase-7. This protein has a calculated molecular weight of 35 kDa.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/23/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Caspase-7 is a cysteinyl aspartic protease and a member of the caspase family of proteases.¹ It is expressed as a proprotein that upon pro-apoptotic signaling is cleaved into two subunits, p20 and p11, which form an active homodimer that recognizes and cleaves proteins containing the amino acid sequence DEVD.^{1,2} Caspase-7 is ubiquitously expressed but is found at low levels in the brain and localizes to the cytoplasm.³ It is an effector caspase and targets the same protein substrates as caspase-3 but is less effective at inducing apoptosis.^{4,5} However, it can induce apoptosis in response to specific signaling, such as granzyme B-induced poly(ADP-ribose) polymerase (PARP) cleavage.^{3,4} Caspase-7 cleaves and activates acid sphingomyelinase to repair gasdermin D (GSDMD) and perforin pores in the plasma membrane and prevents non-apoptotic forms of cell death.⁶ Increased intratumoral levels of cleaved caspase-7 are positively associated with increased overall survival in patients with breast cancer.⁷ Cayman's Caspase-7 (human, recombinant) protein was synthesized from a DNA sequence encoding the mature form of human caspase-7 (Ala24-Gln303) with an N-terminal translation-initiating methionine (Met1). The expressed protein consists of 313 amino acids, has a calculated molecular weight of 35 kDa, and a predicted N-terminus of Met1.

References

1. Lamkanfi, M., and Kanneganti, T.D. Caspase-7: A protease involved in apoptosis and inflammation. *Int. J. Biochem. Cell Biol.* **42(1)**, 21-24 (2010).
2. Talanian, R.V., Quinlan, C., Trautz, S., *et al.* Substrate specificities of caspase family proteases. *J. Biol. Chem.* **272(15)**, 9677-9682 (1997).
3. Cohen, G.M. Caspases: The executioners of apoptosis. *Biochem. J.* **326(Pt 1)**, 1-16 (1997).
4. Slee, E.A., Adrain, C., and Martin, S.J. Executioner caspase-3, -6, and -7 perform distinct, non-redundant roles during the demolition phase of apoptosis. *J. Biol. Chem.* **276(10)**, 7320-7326 (2001).
5. Walsh, J.G., Cullen, S.P., Sheridan, C., *et al.* Executioner caspase-3 and caspase-7 are functionally distinct proteases. *Proc. Natl. Acad. Sci. USA* **105(35)**, 12815-12819 (2008).
6. Nozaki, K., Maltez, V.I., Rayamajhi, M., *et al.* Caspase-7 activates ASM to repair gasdermin and perforin pores. *Nature* **606(7916)**, 960-967 (2022).
7. Lindner, A.U., Lucantoni, F., Varešlija, D., *et al.* Low cleaved caspase-7 levels indicate unfavourable outcome across all breast cancers. *J. Mol. Med. (Berl)* **96(10)**, 1025-1037 (2018).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM