

# Product Information

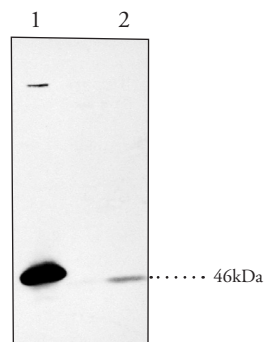


## Caspase-9 Polyclonal Antibody

Item No. 160790

- Synonyms:** Apaf-3; ICE-LAP6; Mch6
- Contents:** This vial contains 50 µg of peptide affinity purified IgG in 500 µl PBS, pH 7.2, containing 0.02% azide.
- Host:** Rabbit
- Antigen:** The anti-caspase-9 polyclonal antibody was raised against a peptide corresponding to amino acids 299 to 318 of human caspase-9.<sup>1-3</sup>
- Application:** This polyclonal antibody can be used for detection of caspase-9 by western blot at a dilution of 1:200. A whole cell lysate from HeLa cells can be used as a positive control and a 46 kDa band should be detected.
- Stability:** ≥1 year at 4°C

Apoptosis is related to many diseases and is induced by a family of cell death receptors and their ligands. Cell death signals are transduced by death domain containing adapter molecules and members of the caspase family of proteases. A novel member in the caspase family was recently identified and designated ICE-LAP6, Mch6, and Apaf3.<sup>1-3</sup> Caspase-9 and Apaf-1 bind to each other, which leads to caspase-9 activation.<sup>3</sup> Caspase-9 is also activated by granzyme B and CPP32.<sup>1,2</sup> Activated caspase-9 cleaves and activates caspase-3, one of the key proteases responsible for the proteolytic cleavage of many key proteins in apoptosis.<sup>3</sup> Caspase-9 plays a central role in cell death induced by a wide variety of apoptosis inducers including TNFα, TRAIL, anti-CD-95, FADD, and TRADD.<sup>4</sup> Caspase-9 is expressed in a variety of human tissues.<sup>1,2</sup>



Lane 1: Caspase 9 transfected 293 T-cell lysate (20 µg)  
Lane 2: 293 T-cell lysate (50 µg)  
Lane 3: HeLa cell lysate (50 µg)

### References

1. Srinivasula, S.M., Fernandes-Alnemri, T., Zangrilli, J., *et al.* The Ced-3/interleukin 1β converting enzyme-like homolog Mch6 and the lamin-cleaving enzyme Mch2α are substrates for the apoptotic mediator CPP32. *J. Biol. Chem.* **271**, 27099-27106 (1996).
2. Duan, H., Orth, K., Chinnaiyan, A.M., *et al.* ICE-LAP6, a novel member of the ICE/Ced-3 gene family, is activated by the cytotoxic T cell protease granzyme B. *J. Biol. Chem.* **271**, 16720-16724 (1996).
3. Li, P., Nijhawan, D., Budihardjo, I., *et al.* Cytochrome c and dATP-dependent formation of Apaf-1/caspase-9 complex initiates an apoptotic protease cascade. *Cell* **91**, 479-489 (1997).
4. Pan, G., O'Rourke, K., and Dixit, V.M. Caspase-9, Bcl-XL, and Apaf-1 form a ternary complex. *J. Biol. Chem.* **273**, 5841-5845 (1998).

### Related Products

Apaf-1 Polyclonal Antibody - Item No. 160780 • Caspase-10 Polyclonal Antibody - Item No. 160795

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 06/22/2011

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

custserv@caymanchem.com

#### Web

www.caymanchem.com