

Product Information



Caspase-9 Polyclonal Antibody

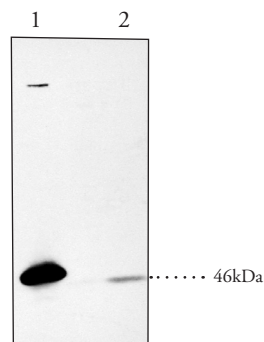
Catalog No. 160790

- Synonyms:** Apaf-3; ICE-LAP6; Mch6
Contents: This vial contains 100 µg of peptide affinity purified IgG in 200 µl PBS 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-9.¹⁻³
Application: This polyclonal antibody can be used for detection of caspase-9 by western blot at a dilution of 1:1000 to 1:2000. 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 Apaf3.¹⁻³ Caspase-9 and Apaf-1 bind to each other, which leads to caspase-9 activation.³ Caspase-9 is also activated by granzyme B and CPP32.^{1,2} Activated caspase-9 cleaves and activates caspase-3, one of the key proteases responsible for the proteolytic cleavage of many key proteins in apoptosis.³ 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.⁴ Caspase-9 is expressed in a variety of human tissues.^{1,2}

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).



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)

Related Products

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

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

WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.

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 under separate cover to the MSDS supervisor at 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, 12/11/2007