

Product Information

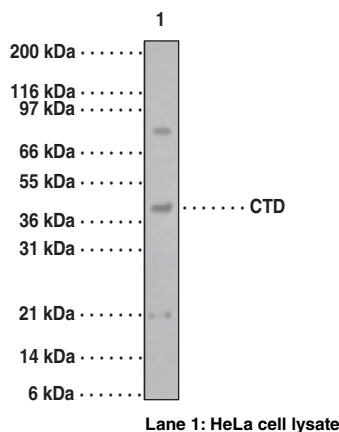


Caspase-9 (carboxyl-terminal divergent) Polyclonal Antibody

Item No. 13915

Contents:	This vial contains IgG 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
Synonym:	CTD
Antigen:	Mixture of synthetic peptides corresponding to rat caspase-9 amino acids 345-359 and 366-383
Host:	Rabbit
Cross Reactivity:	(+) Human, mouse, and rat caspase-9
Storage:	≤6 months at 4°C; ≥6 months at -20°C
Applications:	Western blot (WB); the recommended starting dilution for WB is 1:250-1:1,000

Caspase-9 and Apaf-1 bind to each other, which leads to caspase activation. Activated caspase-9 cleaves and activates caspase-3, one of the key proteases responsible for the proteolytic cleavage of many proteins in apoptosis. Caspase-9 plays a central role in cell death induced by a wide variety of apoptosis inducers including TNFα, TRAIL, anti-CD95, FADD, and TRADD. Caspase-9 is expressed in a variety of human tissues. Recently, a novel isoform of rat caspase-9 has been identified in which the C-terminus of full-length caspase-9 is replaced with an alternative peptide sequence. This protein, called Caspase-9 (CTD) (where CTD is carboxyl-terminal divergent) is expressed in multiple tissues, with the relative highest levels observed in ovary and heart. The variant C-terminus of CTD in rat is derived from an alternative exon of the rat caspase-9 gene. CTD was found primarily in the cytoplasm and was not detected in the nucleus. CTD proenzyme is not processed in the cells and lacks apoptotic activity. The CTD-transfected cells are resistant to caspase induction by cytochrome c, suggesting the CTD acts as a dominant-negative variant.



Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13915

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, 12/19/2011

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