

Product Information

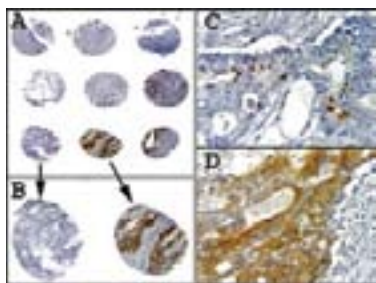


Caspase-1 Monoclonal Antibody (Clone 14F468)

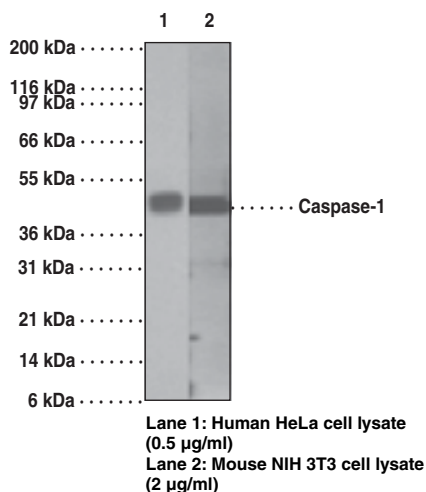
Item No. 13907

Contents:	This vial contains 100 µg of protein G-purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
Synonym:	IL-1β Converting Enzyme
Antigen:	Synthetic peptide corresponding to human caspase-1 amino acids 371-390 (RKVRFSEQPDGRAQMPTTE)
Host:	Mouse, Clone 14F468
Isotype:	IgG _{1κ}
Cross Reactivity:	(+) Human and mouse caspase-1; predicted to react with rat caspase-1
Storage:	≤6 months at 4°C; ≥6 months at -20°C
Applications:	Immunohistochemistry (paraffin-embedded sections) and western blot (WB); the recommended starting concentration for western blot is 0.5-2 µg/ml. The antibody will recognize full-length caspase-1 and cleaved caspase-1 forms that retain amino acids 371-390 of the caspase-1 protein.

Caspases are a family of cysteine proteases that are key mediators of programmed cell death or apoptosis.¹ The precursor form of all caspases is composed of a prodomain and large and small catalytic subunits. The active forms of caspases are generated by several stimuli including ligand-receptor interactions, growth factor deprivation, and inhibitors of cellular functions. All known caspases require cleavage adjacent to aspartates to liberate one large and one small subunit, which associate into a tetramer to form the active enzyme. Caspase-1 is similar to the cell death gene CED-3 of *C. elegans* and regulates multiple proinflammatory cytokines, including interleukin-1β and interferon-γ-inducing factor.



Immunohistochemical analysis of Caspase-1 using Caspase-1 monoclonal antibody. Formalin-fixed, paraffin-embedded sections of human colon cancer stained for Caspase-1 expression using a DAB chromogen and Hematoxylin counterstain. Nine sections, from different human patients were stained. Differential staining was observed (top right panel). Two of these sections are shown at higher magnifications (arrows and A, B). Only rare staining in A is observed. In contrast, abundant staining is observed in B.



Reference

1. Cerretti, D.P., Kozlosky, C.J., Mosley, B., *et al.* Molecular cloning of the interleukin-1β converting enzyme. *Science* 256, 97-100 (1992).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13907

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