

# Product Information

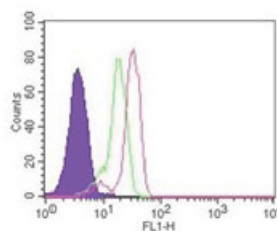
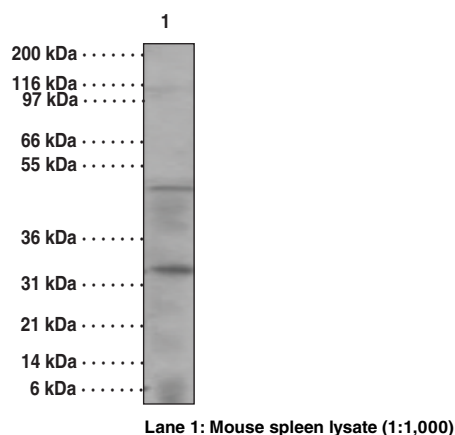


## Caspase-14 Monoclonal Antibody (Clone 70A1426)

Item No. 13916

<b>Contents:</b>	This vial contains 200 µl antibody solution containing 0.09% sodium azide.
<b>Antigen:</b>	Synthetic peptide corresponding to mouse caspase-14 amino acids 2-18 (ESEMSPQLQEERYDM)
<b>GenBank</b>	
<b>Accession No.:</b>	NP_033939
<b>Host:</b>	Mouse, clone 70A1426
<b>Cross Reactivity:</b>	(+) Human and mouse caspase-14
<b>Storage:</b>	≥6 months at -20°C
<b>Applications:</b>	Flow cytometry (FC) (intracellular) and western blot (WB); the recommended starting dilution for FC (intracellular) is 1:500/10 <sup>6</sup> cells and 1:500-1:2,000 for WB.

Caspases are a family of cysteine proteases that are key mediators of programmed cell death or apoptosis.<sup>1</sup> 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 an a2b2 tetramer to form the active enzyme. Caspase-14 has a conserved active site pentapeptide, QACRG. However, no proteolytic or biological activities have been identified so far. Its high expression in embryonic tissues and limited expression in adult tissues suggests that it may have some role during ontogenesis.



Intracellular flow cytometric analysis of caspase-14 in 10<sup>6</sup> Jurkat cells probed with 1:500 dilution of caspase-14 antibody. Shaded histogram represents cells without antibody; green represents isotype control antibody; red represents caspase-14 antibody.

### Reference

1. Cohen, G.M. Caspases: The executioners of apoptosis. *Biochem. J.* **326**, 1-16 (1997).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/13916](http://www.caymanchem.com/catalog/13916)

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

#### 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 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, 08/07/2013

### 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](http://www.caymanchem.com)