

PRODUCT INFORMATION



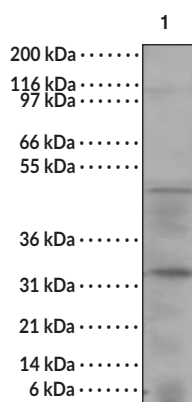
Caspase-14 Monoclonal Antibody (Clone 70A1426)

Item No. 13916

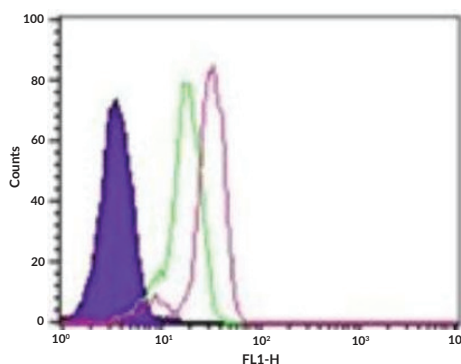
Overview and Properties

Contents: This vial contains 200 µl antibody solution with 0.09% sodium azide.
Synonyms: CASP-14, MICE, Mini-interleukin-1 Converting Enzyme
Immunogen: A synthetic peptide corresponding to amino acids 2-18 of mouse caspase-14
Cross Reactivity: (+) Caspase-14
Species Reactivity: (+) Human, mouse
Storage: -20°C (as supplied)
Stability: ≥6 months
Storage Buffer: Ascites with PBS
Clone: 70A1426
Host: Mouse
Isotype: IgG1κ
Applications: Flow cytometry (FC) (intracellular), Immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution is 1:500/10⁶ cells for FC (intracellular) and 0.1-4 µg/ml for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: Mouse spleen lysate (1:1,000)



Intracellular flow cytometric analysis of caspase-14 in 10⁶ 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.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. 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
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/28/2025

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Caspase-14 is a cysteinyl aspartic protease.¹⁻³ It is expressed during embryonic development and in differentiating keratinocytes and localizes to the cytoplasm and nucleus.^{2,3} Caspase-14 is produced as a zymogen containing a large and small subunit and is activated *via* proteolytic cleavage by kallikrein 7 (KLK7).⁴ It is involved in keratinocyte differentiation, cornification, UVB protection, and profilaggrin degradation.¹ A SNP in *CASP14*, the gene encoding caspase-14, is associated with ichthyosis.⁵ Cayman's Caspase-14 Monoclonal Antibody (Clone 70A1426) can be used for intracellular flow cytometry (FC), immunohistochemistry (IHC), and Western blot (WB) applications. The antibody recognizes caspase-14 from human and mouse samples.

References

1. Sahoo, G., Samal, D., Khandayataray, P., *et al.* A review on caspases: Key regulators of biological activities and apoptosis. *Mol. Neurobiol.* **60**(10), 5805-5837 (2023).
2. Hu, S., Snipas, S.J., Vincenz, C., *et al.* Caspase-14 is a novel developmentally regulated protease. *J. Biol. Chem.* **273**(45), 29648-29653 (1998).
3. Lippens, S., Kockx, M., Knaapen, M., *et al.* Epidermal differentiation does not involve the pro-apoptotic executioner caspases, but is associated with caspase-14 induction and processing. *Cell Death Differ.* **7**(12), 1218-1224 (2000).
4. Yamamoto, M., Miyai, M., Matsumoto, Y., *et al.* Kallikrein-related peptidase-7 regulates caspase-14 maturation during keratinocyte terminal differentiation by generating an intermediate form. *J. Biol. Chem.* **287**(39), 32825-32834 (2012).
5. Kirchmeier, P., Zimmer, A., Bouadjar, B., *et al.* Whole-exome-sequencing reveals small deletions in *CASP14* in patients with autosomal recessive inherited ichthyosis. *Acta Derm. Venereol.* **97**(1), 102-104 (2017).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM