

Product Information

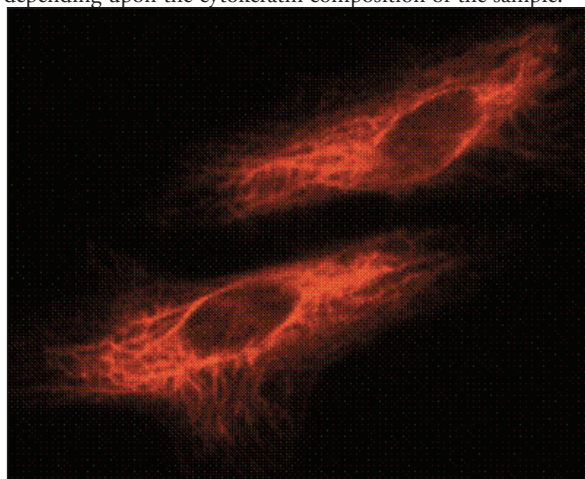


Cytokeratin Monoclonal Antibody (Clone C-11)

Item No. 10004600

Contents:	This vial contains 100 µg affinity-purified IgG ₁ in 500 µl PBS, pH 7.4.
Synonym:	pan-Cytokeratin
Host:	Mouse
Cross Reactivity:	(+) human cytokeratins 4, 5, 6, 8, 10, 13, and 18; other species not tested
Stability:	≥1 year at 4°C
Applications:	Western blot, flow cytometry, immunocytochemistry, and immunohistochemistry (paraffin sections). The recommended starting dilution for WB, flow cytometry, and ICC is 1:200 (1 µg/ml) and 1:200 - 1:50 (1-4 µg/ml) for IHC (paraffin sections). Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

This pan-cytokeratin antibody detects cytokeratin isoforms 4, 5, 6, 8, 10, 13, and 18.¹ Cytokeratins make up a large family of proteins classified as intermediate filament polypeptides.^{1,2} They are subdivided into type I (acidic) and type II (basic/neutral) proteins.³ Under normal conditions, these proteins are found in epithelial cells as type I/type II pairs.^{1,4} These pairs provide cellular support as well as contributing to more dynamic processes. They undergo extensive post translational modifications, such as phosphorylation, and are important for cancer diagnostics.^{4,5} This antibody recognizes human cytokeratins 4, 5, 6, 8, 10, 13, and 18.¹ Accordingly the molecular migration is isoform dependent and expected to range from 40 to 70 kDa depending upon the cytokeratin composition of the sample.



Human cervical carcinoma cells stained with Cytokeratin Monoclonal Antibody (1 µg/ml) followed by goat anti-mouse IgG-Cy3.

References

1. Bártek, J., Vojtěšek, B., Stašková, Z., *et al.* *Journal of Pathology* **164**, 215-224 (1991).
2. Magin, T.M., Vijayaraj, P., and Leube, R.E. *Exp. Cell Res.* **313**, 2021-2032 (2007).
3. Gu, L.-H. and Coulombe, P.A. *Curr. Opin. Cell Biol.* **19**, 13-23 (2007).
4. Ditzel, H.J., Strik, M.C.M., Larsen, M.K., *et al.* *J. Biol. Chem.* **277**(24), 21712-21722 (2002).
5. Plowman, J.E. *Journal of Chromatography B* **849**, 181-189 (2007).

Related Products

Cytokeratin Monoclonal APC Antibody (Clone C-11) - Item No. 10568 • Goat Anti-Mouse IgG HRP - Item No. 10004302 • Goat Anti-Mouse (IgG+IgM) FITC - Item No. 10006617 • 20S Proteasome Assay Kit - Item No. 10008041 • Apoptotic Blebs Assay Kit - Item No. 10010750

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, 03/04/2011

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