

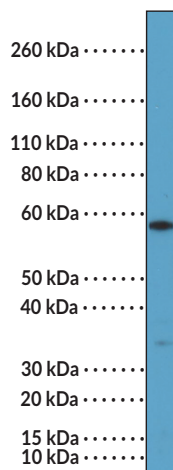
# PRODUCT INFORMATION

## Cytokeratin 10 (C-Term) Rabbit Monoclonal Antibody (Clone RM386) Item No. 32313

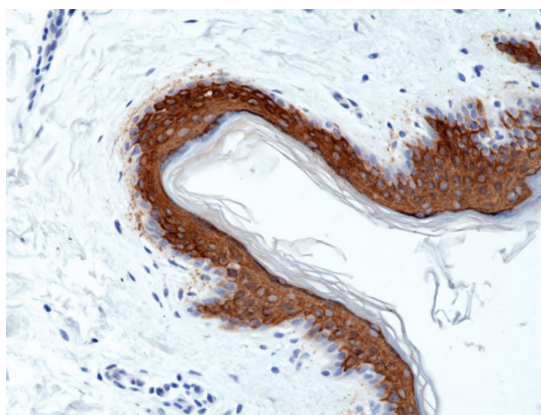
### Overview and Properties

<b>Contents:</b>	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
<b>Synonyms:</b>	CK-10, Keratin-10, Keratin, Type I Cytoskeletal 10, KRT10
<b>Immunogen:</b>	Peptide from the C-terminal region of human cytokeratin 10
<b>Cross Reactivity:</b>	(+) Cytokeratin 10
<b>Species Reactivity:</b>	(+) Human
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS, with 50% glycerol, 1% BSA, and 0.09% sodium azide
<b>Clone:</b>	RM386
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for IHC is 1:500-1:1,000 and 1:200-1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



WB of A431 cell lysate using Cytokeratin 10 (C-Term) Rabbit Monoclonal Antibody (Clone RM386) at a dilution of 1:200.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human skin tissue using Cytokeratin 10 (C-Term) Rabbit Monoclonal Antibody (Clone RM386) at a dilution of 1:1,000.

**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, 02/22/2024

**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

---

Cytokeratin 10 (CK-10) is a type I acidic epithelial intermediate filament protein.<sup>1</sup> It is composed of a central rod containing four  $\alpha$ -helical domains, which are important for self-assembly, and non-helical head and tail domains at the N- and C-termini, respectively.<sup>2</sup> CK-10 expression is induced upon keratinocyte differentiation, increasing during keratinocyte migration from the basal layer to the suprabasal layer of the epidermis.<sup>1</sup> It dimerizes with the type II epithelial intermediate filament protein CK-1 *via* heptad repeats in the central rod domain to form a network of filament bundles throughout the cytoplasm.<sup>3</sup> Increased tumor CK10 expression has been found in patients with hepatocellular carcinoma (HCC) and is associated with poor prognosis.<sup>4</sup> CK10 mutations have been found in patients with epidermolysis bullosa simplex (EBS), a disorder characterized by skin blistering.<sup>3</sup> Cayman's Cytokeratin 10 (C-Term) Rabbit Monoclonal Antibody (Clone RM386) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

## References

---

1. Moll, R., Divo, M., and Langbein, L. The human keratins: biology and pathology. *Histochem. Cell Biol.* **129(6)**, 705-733 (2008).
2. Gu, L.-H. and Coulombe, P.A. Keratin function in skin epithelia: A broadening palette with surprising shades. *Curr. Opin. Cell Biol.* **19(1)**, 13-23 (2007).
3. Chan, Y.M., Yu, Q.C., LeBlanc-Straceski, J., *et al.* Mutations in the non-helical linker segment L1-2 of keratin 5 in patients with Weber-Cockayne epidermolysis bullosa simplex. *J. Cell Sci.* **107(Pt 4)**, 765-774 (1994).
4. Yang, X.-R., Xu, Y., Shi, G.-M., *et al.* Cytokeratin 10 and cytokeratin 19: Predictive markers for poor prognosis in hepatocellular carcinoma patients after curative resection. *Clin. Cancer Res.* **14(12)**, 3850-3859 (2008).

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