

PRODUCT INFORMATION



Cytokeratin 19 (C-Term) Rabbit Monoclonal Antibody (Clone RM364) Item No. 32289

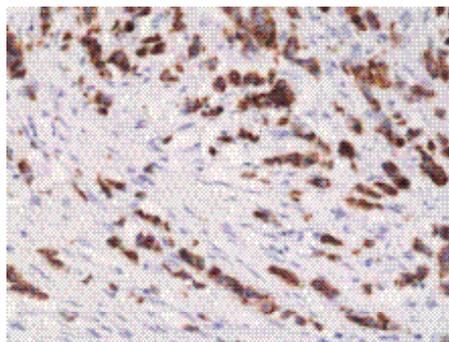
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	CK-19, Keratin, Type I Cytoskeletal 19, Keratin-19, KRT19
Immunogen:	Peptide from the C-terminal region of human cytokeratin 19
Cross Reactivity:	(+) Cytokeratin 19
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM364
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:500-1:1,000 for IHC and 1:1,000-1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



WB of MCF-7 cell lysate using Cytokeratin 19 (C-Term) Rabbit Monoclonal Antibody (Clone RM364) at a dilution of 1:1,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human breast cancer tissue using Cytokeratin 19 (C-Term) Rabbit Monoclonal Antibody (Clone RM364) 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, 02/15/2024

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 19 is a type I epithelial intermediate filament protein.¹ It is composed of a central rod containing four α -helical domains, which are important for self-assembly, and non-helical head and tail domains at the N- and C-termini, respectively. Cytokeratin 19 is expressed in bile duct epithelial cells and non-keratinized stratified squamous epithelial cells, as well as other epithelial cells.² Cytokeratin 19 is associated with a variety of cancers, including salivary gland, lung, and thyroid papillary carcinomas and various gastrointestinal cancers.³ The presence of cytokeratin 19 in tumor tissue from patients with hepatocellular carcinoma is associated with reduced recurrence-free and overall survival.⁴ Cayman's Cytokeratin 19 (C-Term) Rabbit Monoclonal Antibody (Clone RM364) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. 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).
2. Chu, P.G. and Weiss, L.M. Keratin expression in human tissues and neoplasms. *Histopathology* **40(5)**, 403-439 (2002).
3. Caviglia, G.P., Ciruolo, M., Olivero, A., *et al.* Prognostic role of serum cytokeratin-19 fragment (CYFRA 21-1) in patients with hepatocellular carcinoma. *Cancers* **12(10)**, 2776 (2020).
4. Kawai, T., Yasuchika, K., Ishii, T., *et al.* Keratin 19, a cancer stem cell marker in human hepatocellular carcinoma. *Clin. Cancer Res.* **21(13)**, 3081-3091 (2015).

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