

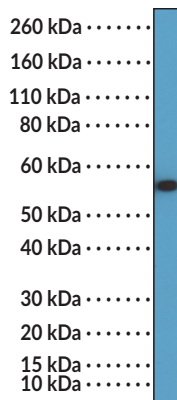
PRODUCT INFORMATION

Cytokeratin 7 (C-Term) Rabbit Monoclonal Antibody (Clone RM416) Item No. 32337

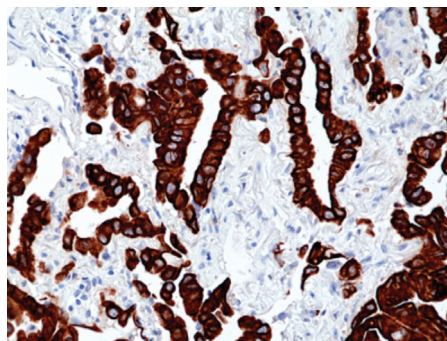
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	CK-7, Keratin-7, Keratin, Type II Cytoskeletal 7, KRT7
Immunogen:	Peptide from the C-terminal region of human cytokeratin 7
Cross Reactivity:	(+) Cytokeratin 7
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:	RM416
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:200 for IHC and 1:1,000-1:10,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



WB of HeLa cell lysate using Cytokeratin 7 (C-Term) Rabbit Monoclonal Antibody (Clone RM416) at a dilution of 1:10,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human lung adenocarcinoma tissue using Cytokeratin 7 (C-Term) Rabbit Monoclonal Antibody (Clone RM416) at a dilution of 1:100.

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, 01/13/2021

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 7 (CK-7) is an intermediate filament protein and a member of the cytokeratin protein family.^{1,2} It is a type II neutral-basic cytokeratin expressed in ductal, glandular, and transitional epithelial tissues, as well as endothelial cells, but not stratified squamous epithelia, and is localized to the cytoplasm.^{2,3} CK-7 is found in various tumors, including ductal, lobular, and mucinous carcinomas, as well as ovarian, but not colorectal, adenocarcinomas, malignant mesotheliomas, and transitional cell carcinomas.³ It is also found in chromophobe renal cell carcinomas but not oncocytoma and conventional carcinomas and has been used as a biomarker for differentiation of tumor origin.⁴ Hepatocyte expression of CK-7 is positively correlated with fibrosis stage in patients with hepatitis B and hepatitis C.⁵ Cayman's Cytokeratin 7 (C-Term) Rabbit Monoclonal Antibody (Clone RM416) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Weng, Y.-R., Cui, Y., and Fang, J.-Y. Biological functions of cytokeratin 18 in cancer. *Mol. Cancer Res.* **10(4)**, 485-493 (2012).
2. Rabban, J.T., Soslow, R.A., and Zaloudek, C.Z. Immunohistology of the female genital tract. *Diagnostic immunohistochemistry*. Dabbs, D.J., editor, 3rd edition, W.B. Saunders (2010).
3. Campbell, F. and Herrington, C.S. Application of cytokeratin 7 and 20 immunohistochemistry to diagnostic pathology. *Curr. Diagn. Pathol.* **7(2)**, 113-122 (2001).
4. Mathers, M.E., Pollock, A.M., Marsh, C., et al. Cytokeratin 7: A useful adjunct in the diagnosis of chromophobe renal cell carcinoma. *Histopathology* **40(6)**, 563-567 (2002).
5. Matsukuma, S., Takeo, H., Kono, T., et al. Aberrant cytokeratin 7 expression of centrilobular hepatocytes: A clinicopathological study. *Histopathology* **61(5)**, 857-862 (2012).

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