

PRODUCT INFORMATION



Cytokeratin 17 (C-Term) Rabbit Monoclonal Antibody (Clone RM351) Item No. 32275

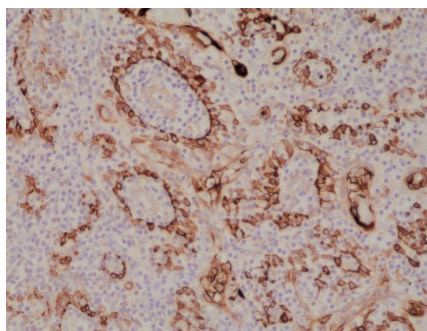
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	CK-17, Keratin-17, Keratin, Type II Cytoskeletal 17, KRT17
Immunogen:	Peptide from the C-terminal region of human cytokeratin 17
Cross Reactivity:	(+) Cytokeratin 17
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:	RM351
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:500-1:1,000 for IHC and 1:1,000-1:20,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



WB of HeLa cell lysates using Cytokeratin 17 (C-Term) Rabbit Monoclonal Antibody (Clone RM351) at a dilution of 1:10,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human tonsil tissue using Cytokeratin 17 (C-Term) Rabbit Monoclonal Antibody (Clone RM351) at a 1:1,000 dilution.

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, 01/31/2024

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 17 (CK-17) is an intermediate filament protein and a member of the cytokeratin protein family encoded by *KRT17* in humans.¹ It is a type I acidic cytokeratin expressed in the basal cells of epithelial appendages, such as hair follicles and sebaceous glands, and is localized to the cytoplasm. CK-17 expression can be induced in the epidermis under conditions of cellular stress, such as viral infection, injury, or psoriasis. CK-17 is also expressed in squamous, basal, and urothelial carcinomas, as well as adenocarcinomas, and is associated with poor prognosis in breast, cervical, oral squamous, and gastric carcinomas.^{1,2} Heterozygous dominant mutations in *KRT17* are associated with early onset alopecia.¹ Missense mutations in *KRT17* are associated with ectodermal dysplasias, including pachyonychia congenita and steatocystoma multiplex. Cayman's Cytokeratin 17 (C-Term) Rabbit Monoclonal Antibody (Clone RM351) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Yang, L., Zhang, S., and Wang, G. Keratin 17 in disease pathogenesis: From cancer to dermatoses. *J. Pathol.* **247(2)**, 158-165 (2019).
2. Kim, K., Lee, H.W., Chae, S.W., et al. Cytokeratin 17 expression is associated with poor prognosis in gallbladder adenocarcinoma. *Appl. Immunohistochem. Mol. Morphol.* **25(5)**, 346-350 (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