

PRODUCT INFORMATION

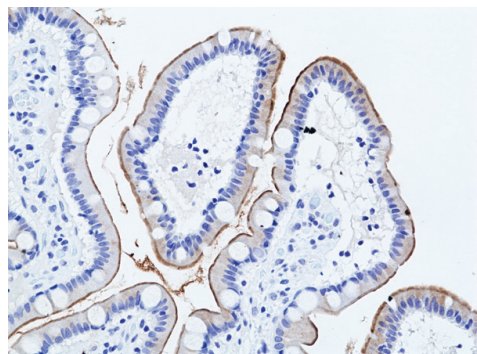
MUC17 Rabbit Monoclonal Antibody (Clone RM517)

Item No. 42783

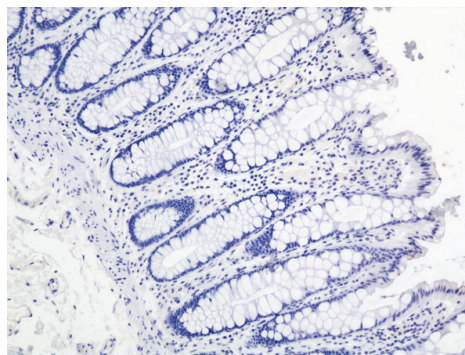
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified antibody.
Synonym:	Mucin 17
Immunogen:	Peptide corresponding to residues near the C-terminus of human MUC17
Cross Reactivity:	(+) MUC17
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Clone:	RM517
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC); the recommended starting dilution is 1:100-1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Immunohistochemical staining of formalin-fixed and paraffin-embedded human small intestine tissue sections using MUC17 Rabbit Monoclonal Antibody (Clone RM517) at a 1:100 dilution for one hour at room temperature.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human normal colon tissue sections using MUC17 Rabbit Monoclonal Antibody (Clone RM517) at a 1:100 dilution for one hour at room temperature.

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, 03/19/2025

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

Mucin 17 (MUC17) is a member of the mucin family.¹ It is composed of an N-terminal tandem repeat domain, which contains O- and N-linked glycosylation sites, two EGF-like domains, a SEA module, a transmembrane region, and a short C-terminal cytoplasmic domain.^{1,2} MUC17 is alternatively expressed as a membrane-bound or secreted protein.² Membrane-bound MUC17 is predominantly expressed in the small intestine and colon by absorptive epithelial cells.^{1,3} The levels of MUC17 are increased in tumor tissue from patients with colon adenocarcinoma.³ Cayman's MUC17 Rabbit Monoclonal Antibody (Clone RM517) can be used for immunohistochemistry (IHC).

References

1. Gum, J.R., Jr., Crawley, S.C., Hicks, J.W., *et al.* MUC17, a novel membrane-tethered mucin. *Biochem. Biophys. Res. Commun.* **291**(3), 466-475 (2002).
2. Moniaux, N., Junker, W.M., Singh, A.P., *et al.* Characterization of human mucin MUC17. Complete coding sequence and organization. *J. Biol. Chem.* **281**(33), 23676-23685 (2006).
3. Krishn, S.R., Kaur, S., Smith, L.M., *et al.* Mucins and associated glycan signatures in colon adenoma-carcinoma sequence: Prospective pathological implication(s) for early diagnosis of colon cancer. *Cancer Lett.* **374**(2), 304-314 (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