

PRODUCT INFORMATION



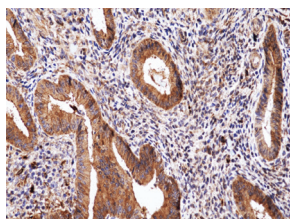
PTEN Rabbit Monoclonal Antibody (Clone RM511)

Item No. 42778

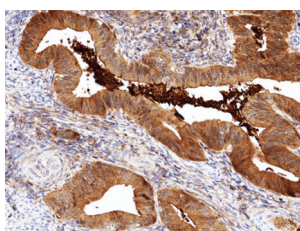
Overview and Properties

Contents:	This vial contains 100 µg of protein A-affinity purified monoclonal antibody.
Synonyms:	MMAC1, Mutated in Multiple Advanced Cancers 1, Phosphatase and Tensin Homolog, TEP1
Immunogen:	A peptide corresponding to residues at the C-terminus of human PTEN
Cross Reactivity:	(+) PTEN
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:	RM511
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry; the recommended starting concentration is 1:100-1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

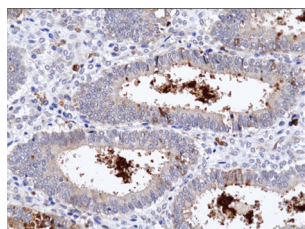
Images



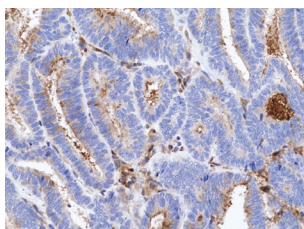
Immunohistochemical staining of formalin-fixed and paraffin-embedded human endometrium (proliferative) tissue section using PTEN Rabbit Monoclonal Antibody (Clone RM511) at a 1:100 dilution for one hour at room temperature.



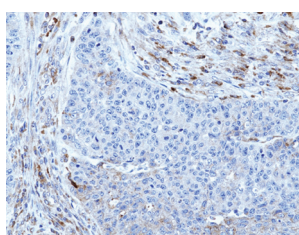
Immunohistochemical staining of formalin-fixed and paraffin-embedded human endometrium (secretory) tissue section using PTEN Rabbit Monoclonal Antibody (Clone RM511) at a 1:100 dilution for one hour at room temperature.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human endometrium neoplasia (Stage IA) tissue section using PTEN Rabbit Monoclonal Antibody (Clone RM511) at a 1:100 dilution for one hour at room temperature.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human endometrium adenocarcinoma (Stage IB) tissue section using PTEN Rabbit Monoclonal Antibody (Clone RM511) at a 1:100 dilution for one hour at room temperature.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human endometrium adenocarcinoma (Stage III) tissue section using PTEN Rabbit Monoclonal Antibody (Clone RM511) 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, 02/03/2025

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

Phosphatase and tensin homolog (PTEN) is a tumor suppressor with roles in cell division, cell migration, and apoptosis.¹⁻⁴ It is ubiquitously expressed and localized to the cytosol and nucleus where it dephosphorylates phosphatidylinositol-(3,4,5)-triphosphate (PIP₃) to phosphatidylinositol-(3,4)-diphosphate (PIP₂) to antagonize PI3K signaling and regulates spliceosome assembly and pre-mRNA splicing, respectively.² PTEN inhibits cell cycle progression and induces apoptosis, as well as inhibits cell adhesion and migration *in vitro*.⁵ PTEN deletions have been found in various cancers, including glioblastomas and prostate, breast, and kidney cells. Germline mutations in PTEN induce a variety of syndromes, including Cowden, Bannayan-Riley-Ruvalcaba, Proteus, and Proteus-like syndromes, which are characterized by increased risk of cancers, developmental delays, autism spectrum disorders, and macrocephaly.⁴ Cayman's PTEN Rabbit Monoclonal Antibody (Clone RM511) can be used for immunohistochemistry (IHC).

References

1. Di Cristofano, A. and Pandolfi, P.P. The multiple roles of PTEN in tumor suppression. *Cell* **100**(4), 387-390 (2000).
2. Liang, H. and Yin, Y. Multiple roles of PTEN isoforms PTEN α and PTEN β in cellular activities and tumor development. *Sci. China Life Sci.* **62**(12), 1722-1724 (2019).
3. Boossani, C.S., Gunasekar, P., and Agrawal, D.K. An update on PTEN modulators - a patent review. *Expert Opin. Ther. Pat.* **29**(11), 881-889 (2019).
4. Nunes-Santos, C.J., Uzel, G., and Rosenzweig, S.D. PI3K pathway defects leading to immunodeficiency and immune dysregulation. *J. Allergy Clin. Immunol.* **143**(5), 1676-1687 (2019).
5. Dahia, P.L.M. PTEN, a unique tumor suppressor gene. *Endocr. Relat. Cancer* **7**(2), 115-129 (2000).

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