

PRODUCT INFORMATION



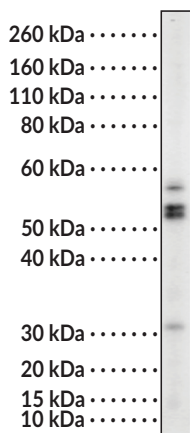
WT1 Rabbit Monoclonal Antibody (Clone RM516)

Item No. 42782

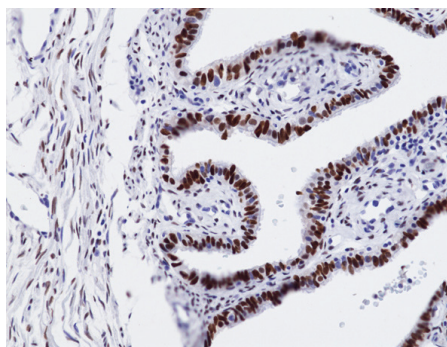
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonym:	Wilms Tumor Protein 1
Immunogen:	A peptide corresponding to residues near the N-terminus of human WT1
Cross Reactivity:	(+) WT1
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:	RM504
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:200 for IHC and 1:500-1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



WB of K562 cell lysates using WT1 Rabbit Monoclonal Antibody (Clone RM516) at a 1:500 dilution.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human fallopian tube tissue using GDFP-15 Rabbit Monoclonal Antibody (Clone RM504) at a dilution of 1:100 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.

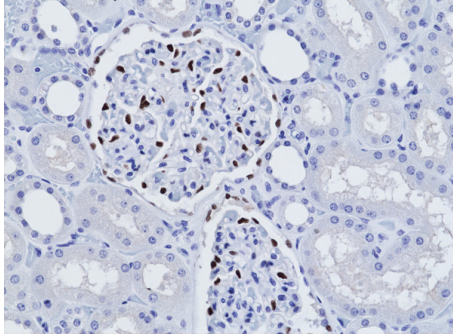
WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/02/2025

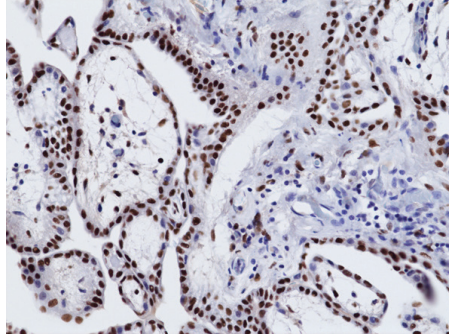
CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

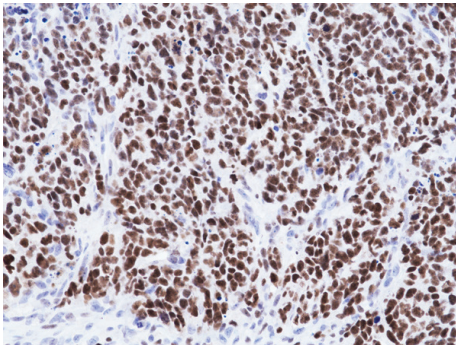
Images



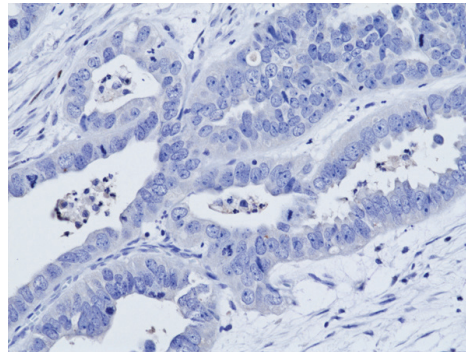
Immunohistochemical staining of formalin-fixed and paraffin-embedded human normal kidney tissue using WT1 Rabbit Monoclonal Antibody (Clone RM516) at a dilution of 1:100 for one hour at room temperature.



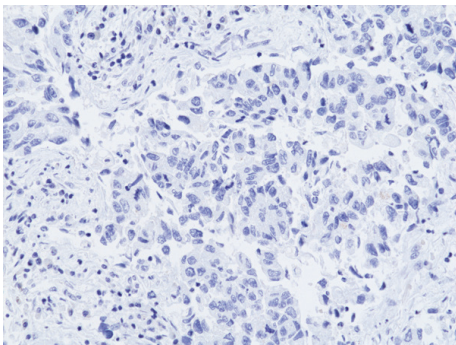
Immunohistochemical staining of formalin-fixed and paraffin-embedded human mesothelioma tissue using WT1 Rabbit Monoclonal Antibody (Clone RM516) at a dilution of 1:100 for one hour at room temperature.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human ovarian serous adenocarcinoma (positive control verified) tissue using WT1 Rabbit Monoclonal Antibody (Clone RM516) at a dilution of 1:100 for one hour at room temperature.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human ovarian serous adenocarcinoma (negative control verified) tissue using WT1 Rabbit Monoclonal Antibody (Clone RM516) at a dilution of 1:100 for one hour at room temperature.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human lung cancer (as negative control) tissue using WT1 Rabbit Monoclonal Antibody (Clone RM516) at a dilution of 1:100 for one hour at room temperature.

PRODUCT INFORMATION



Description

Wilms tumor protein 1 (WT1) is a transcription factor and member of the SP1 family of transcription factors.¹ It is composed of an N-terminal dimerization domain, a transcriptional repressor domain, a transactivation domain, and a C-terminal DNA/RNA-binding domain containing four zinc-finger motifs.^{1,2} There are many isoforms of WT1 formed by various processes.¹ The most common isoforms contain or lack a lysine, threonine, and serine sequence (+KTS or -KTS) in the zinc-finger region, contain or lack 17 amino acids encoded by exon 5 (+17aa or -17aa), lack the KTS sequence but contain the 17-amino acid sequence (-KTS/+17aa), or contain both sequences (+KTS/+17aa).^{1,3} WT1 is expressed in, and crucial for the development of, the genitourinary system, gonads, and adrenal glands.^{2,3} WT1 binds to DNA and acts as a transcriptional repressor or activator in a context-dependent manner to modulate the expression of its targets, including *NR5A1/SF1*, which is involved in adrenal gland and gonadal development.^{1,3} It also binds to and modulates mRNA splicing and stability.¹ Certain mutations in *WT1* disrupt differentiation of metanephric mesenchyme cells and are associated with Wilms tumor, a renal cancer primarily affecting children.⁴ Other *WT1* mutations are associated with non-cancer pathologies, such as Frasier syndrome and Denys-Drash syndrome, which include genitourinary and gonadal development disruptions, as well as heart and central and peripheral nervous system disorders.^{3,4} Cayman's WT1 Rabbit Monoclonal Antibody (Clone RM516) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Torban, E. and Goodyer, P. Wilms' tumor gene 1: Lessons from the interface between kidney development and cancer. *Am. J. Physiol. Renal. Physiol.* **326(1)**, F3-F19 (2024).
2. Brown, K.W. and Malik, K.T. The molecular biology of Wilms' tumour. *Expert. Rev. Mol. Med.* 1-16 (2001).
3. Ferrari, M.T.M., Elias, F.M., Gomes, N.L.R.G., et al. WT1: A single gene associated with multiple and severe phenotypes. *Endocrinol. Metab.* **13**, 100143 (2023).
4. Morrison, A.A., Viney, R.L., Saleem, M.A., et al. New insights into the function of the Wilms tumor suppressor gene *WT1* in podocytes. *Am. J. Physiol. Renal. Physiol.* **295(1)**, F12-17 (2008).

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