

PRODUCT INFORMATION



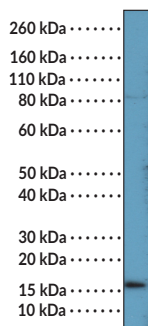
UP2 Rabbit Monoclonal Antibody (Clone RM507)

Item No. 42858

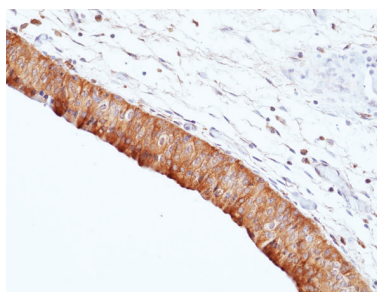
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified antibody.
Synonyms:	UPII, Uroplakin II, Uroplakin-2
Immunogen:	Peptide corresponding to residues near the C-terminus of human UP2
Cross Reactivity:	(+) UP2
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:	RM507
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:200 for IHC and 1:100-1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

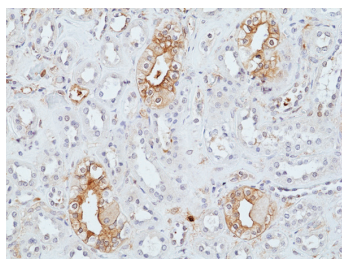
Images



WB of rhesus macaque bladder tissue lysate using UP2 Rabbit Monoclonal Antibody (Clone RM507) at a 1:100 dilution.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human ureter tissue using UP2 Rabbit Monoclonal Antibody (Clone RM507) at a dilution of 1:200.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human kidney tissue using UP2 Rabbit Monoclonal Antibody (Clone RM507) 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, 06/10/2025

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

Uroplakin-2 (UP2) is a member of the uroplakin family.¹ It is expressed as a proprotein composed of an N-terminal signal peptide and prodomain, which are cleaved in the endoplasmic reticulum (ER), and a C-terminal domain, which contains the mature peptide sequence, which is composed of a serine- and threonine-rich extracellular domain, a transmembrane domain, and a short cytoplasmic tail.^{1,2} It is expressed in urothelial umbrella cells and forms protective plaques with UP1a, UP1b, and UP3 in the urinary tract.¹ Antibodies targeting UP2 have been used in diagnostic panels for identifying cancers that originate in urothelial tissue.³ Cayman's UP2 Rabbit Monoclonal Antibody (Clone RM507) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Matuszewski, M.K., Tupikowski, K., Dołowy, Ł., *et al.* Uroplakins and their potential applications in urology. *Cent. European J. Urol.* **69(3)**, 252-257 (2016).
2. Lin, J.H., Wu, X.R., Kreibich, G., *et al.* Precursor sequence, processing, and urothelium-specific expression of a major 15-kDa protein subunit of asymmetric unit membrane. *J. Biol. Chem.* **269(3)**, 1775-1784 (1994).
3. Hoang, L.L., Tacha, D., Bremer, R.E., *et al.* Uroplakin II (UPII), GATA3, and p40 are highly sensitive markers for the differential diagnosis of invasive urothelial carcinoma. *Appl. Immunohistochem. Mol. Morphol.* **23(10)**, 711-716 (2015).

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