

PRODUCT INFORMATION

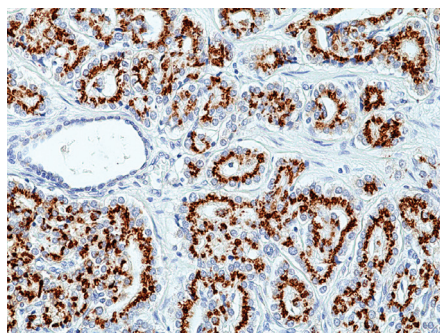
Prostein (C-Term) Rabbit Monoclonal Antibody (Clone RM426)

Item No. 35318

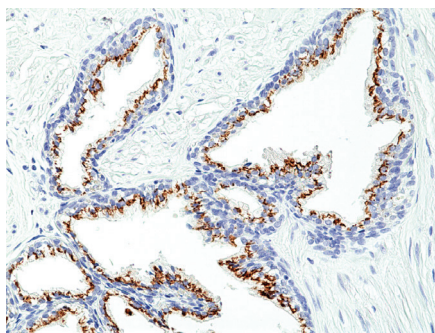
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	p501s, PCANAP6, Prostate Cancer-associated Protein 6
Immunogen:	Peptide corresponding to the C-terminus of human Prostein
Cross Reactivity:	(+) Prostein
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:	RM426
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC); the recommended starting dilution is 1:100-1:400. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Immunohistochemical staining of formalin-fixed and paraffin-embedded human prostate cancer tissue using Prostein (C-Term) Rabbit Monoclonal Antibody (Clone RM426) at a 1:100 dilution.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human prostate tissue using Prostein (C-Term) Rabbit Monoclonal Antibody (Clone RM426) at a 1:100 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, 01/19/2024

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

Prostein, also known as prostate cancer-associated protein 6 (PCANAP6), is a prostate-specific protein encoded by *SLC45A3* that is expressed in normal and malignant prostate tissues.^{1,2} It is composed of 11 transmembrane domains and localizes to the Golgi apparatus.^{3,4} Prostein mRNA and protein expression is increased by androgen stimulation in LNCaP prostate cancer cells, and prostein translation is reduced by miRNA-126* in the same model.^{1,5} It has been used as a biomarker for prostate cancer and to determine the origin of metastatic tumors.⁶ Reduced prostein expression is associated with *SLC45A3-ERG* gene fusions and reduced recurrence-free survival time in patients with prostate cancer.⁷ Cayman's Prostein (C-Term) Rabbit Monoclonal Antibody (Clone RM426) can be used for immunohistochemistry (IHC).

References

1. Xu, J., Kalos, M., Stolk, J.A., et al. Identification and characterization of prostein, a novel prostate-specific protein. *Cancer Res.* **61**(4), 1563-1568 (2001).
2. Kalos, M., Askaa, J., Hylander, B.L., et al. Prostein expression is highly restricted to normal and malignant prostate tissues. *Prostate* **60**(3), 246-256 (2004).
3. Pin, E., Henjes, F., Hong, M.-G., et al. Identification of a novel autoimmune peptide epitope of prostein in prostate cancer. *J. Proteome Res.* **16**(1), 204-216 (2017).
4. Kristiansen, I., Stephan, C., Jung, K., et al. Sensitivity of HOXB13 as a diagnostic immunohistochemical marker of prostatic origin in prostate cancer metastases: Comparison to PSA, prostein, androgen receptor, *ERG*, *NKX3.1*, *PSAP*, and *PSMA*. *Int. J. Mol. Sci.* **18**(6), 1151 (2017).
5. Musiyenko, A., Bitko, V., and Barik, S. Ectopic expression of miR-126*, an intronic product of the vascular endothelial EGF-like 7 gene, regulates prostein translation and invasiveness of prostate cancer LNCaP cells. *J. Mol. Med. (Berl.)* **86**(3), 313-322 (2008).
6. Queisser, A., Hagedorn, S.A., Braun, M., et al. Comparison of different prostatic markers in lymph node and distant metastases of prostate cancer. *Mod. Pathol.* **28**(1), 138-145 (2015).
7. Perner, S., Rupp, N.J., Braun, M., et al. Loss of *SLC45A3* protein (prostein) expression in prostate cancer is associated with *SLC45A3-ERG* gene rearrangement and an unfavorable clinical course. *Int. J. Cancer* **132**(4), 807-812 (2013).

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