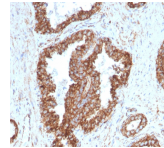


PMEPA1 / TMEPAI Monoclonal Mouse Antibody (PMEPA1/2698)



Product Description

PMEPA1 (prostate transmembrane protein, androgen induced 1) is a 287 amino acid single-pass membrane protein that contains WW-binding motifs and localizes to the cell membrane. Expressed at high levels in prostate, kidney and ovary, PMEPA1 interacts with NEDD4 and may play a role in regulating AR (androgen receptor) levels, specifically in prostate cells. Down regulation of PMEPA1 is observed in prostate tumors, suggesting that PMEPA1 may exhibit activity as a tumor suppressor. Overexpression of this protein may play a role in multiple types of cancer.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Product attributes

Antibody number	#2698
Antibody reactivity (target)	PMEPA1, TMEPAI
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	PMEPA1/2698
Isotype	IgG1, kappa
Molecular weight	32 kDa
Synonyms	PMEPA; PMEPA1; Prostate transmembrane protein; androgen induced 1; Solid tumor-associated 1 protein; STAG1; TMEPAI; Transmembrane prostate androgen-induced protein; Transmembrane, prostate androgen induced RNA
Human gene symbol	PMEPA1
Entrez gene ID	56937
SwissProt	Q96W9
Unigene	517155
Immunogen	Recombinant full-length human PMEPA1 protein
Antibody target cellular localization	Membrane/vesicular
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human
Positive control	Human prostate, breast or renal cell carcinoma tissue (IHC).
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Validated in protein array	Monospecific
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Cell/tissue expression	Kidney, Ovary, Prostate
Tumor expression	Prostate cancer

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, of LI-COR Bioscience.

This datasheet was generated on September 1, 2024 at 12:20:04 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/pmepa1-tmepai-monoclonal-mouse-antibody-pmepa1-2698/>