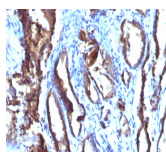


Prostate Specific Antigen / PSA Monoclonal Mouse Antibody (3E6)



Product Description

Recognizes a single protein of 33-34 kDa, identified as the prostate specific antigen (PSA). This MAb is highly specific to PSA and stains prostatic secretory and ductal epithelium in both normal and neoplastic tissues. PSA is a chymotrypsin-like serine protease (kallikrein family) exclusively produced by the prostate epithelium, and abundant in seminal fluid. PSA can be detected in the sera of patients with prostatic carcinoma. It is predominantly complexed to a liver-derived serine protease inhibitor, alpha-1-antichymotrypsin (ACT). A higher proportion of serum PSA is complexed to ACT in prostate cancer than in benign prostate hyperplasia. This MAb makes an excellent pair with MAb A67-B/E3 for PSA tests.

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.

Catalog number key for antibody number 1777, Anti-Prostate Specific Antigen|PSA (#3E6)

Product attributes

Antibody number	#1777
Antibody reactivity (target)	Prostate Specific Antigen, PSA
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	#3E6
Isotype	IgG1, kappa
Molecular weight	33-34 kDa
Synonyms	Antigen; prostate-specific (APS); Gamma-seminoprotein; hK3; Kallikrein related peptidase 3; Kallikrein-3 (KLK3); KLK2A1; P-30 antigen; Semenogelase; Seminin
Human gene symbol	KLK3
Entrez gene ID	354
SwissProt	P07288
Unigene	171995
Immunogen	Human Prostate Specific Antigen (PSA)
Verified antibody applications	IHC (FPPE) (verified)
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Human
Positive control	PC12 cells. Normal prostate or prostate carcinoma.
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	Prostate
Antibody research areas	Cancer, Endocrinology
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
BNC74	CF®740	742/767	633-685	775/50	CF®740 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, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on November 28, 2024 at 06:44:42 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/prostate-specific-antigen-psa-monoclonal-mouse-antibody-3000000/>