



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357) Email: [btinfo@biotium.com](mailto:btinfo@biotium.com)

# Prostate Specific Antigen / PSA Monoclonal Mouse Antibody (KLK3/1248)

## 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.

### Product attributes

Antibody number	#1248
Antibody reactivity (target)	Prostate Specific Antigen
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	KLK3/1248
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	Recombinant full-length human KLK3 protein
Antibody target cellular	Secreted (extracellular)
Verified antibody applications	IHC (FFPE) (verified)
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
Cell/tissue expression	Prostate
Antibody research areas	Cancer, Endocrinology

This datasheet was generated on September 1, 2024 at 04:24:59 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/prostate-specific-antigen-psa-monoclonal-mouse-antibody-klk3-1248/>