

Prostate Specific Antigen Monoclonal Mouse Antibody (1A7)

Product Description

This antibody 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 | #0926 |
| Antibody reactivity (target) | Prostate Specific Antigen |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | 1A7 |
| 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 | PSA (p30) from human sperm plasma |
| Antibody target cellular | Secreted (extracellular) |
| Verified antibody applications | IHC (FFPE) (verified) |
| Species reactivity | Cat, Human |
| Antibody application notes | Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistochemistry (formalin-fixed): 0.5-1.0 ug/mL for 30 minutes at RT, Flow cytometry: 0.5-1 ug/million cells, Immunofluorescence: 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user |
| Positive control | PC-3 or LNCaP cell lysates (WB), Prostate tissue (IHC), LNCaP cells (IF). |
| 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 |
| Tumor expression | Prostate cancer |
| Antibody research areas | Cancer |

This datasheet was generated on September 1, 2024 at 04:25:10 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/prostate-specific-antigen-monoclonal-mouse-antibody-1a7/>