

**Biotin Mouse Anti-human PSA Antibody**  
**\*A67-B/E3, monoclonal\***Catalog number: V1032050  
Unit size: 0.1 mg**Product Details**

---

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

**Antibody Properties**

---

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	PSA
Clone	A67-B/E3
Conjugate	Biotin

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with Biotin under optimal conditions
Application	IHC(P), WB, ELISA

**Applications**

---

Prostate-specific antigen (PSA) is a protein with a molecular weight of 29 kDa, expressed in the extracellular exosome, protein-containing complex and extracellular region of cells. In humans, PSA plays an important role in organismal processes, in particular, cellular protein metabolic process, zymogen activation and proteolysis. PSA is a negative regulator of angiogenesis. It has been associated with critical functions like serine-type peptidase, serine-type endopeptidase and endopeptidase activity. Sequencing of PSA has shown it contains a primary structural unit, the peptidase S1 domain. PSA aids in systemic arterial blood pressure.