

**PE Mouse Anti-human Ig Kappa Light Chain  
Antibody \*A8B5, monoclonal, Cross  
Adsorbed\***Catalog number: V1031335  
Unit size: 100 tests**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	Secondary
Clonality	Monoclonal
Host	Mouse
Immunogen	Ig Kappa Light Chain
Clone	A8B5
Conjugate	PE

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against rabbit, hamster, sheep, guinea pig, goat serum and then conjugated with PE under optimal conditions
Application	FC (QC TESTED)

**Applications**

---

Immunoglobulin  $\kappa$  light chain is a 27 kDa protein that can be located in the plasma membrane, immunoglobulin complex and extracellular region of cells. In Homo sapiens, immunoglobulin  $\kappa$  light chain plays an important role in organismal processes, for example, adaptive immune response. Sequencing of immunoglobulin  $\kappa$  light chain has demonstrated it contains 2 conserved structural units: Ig-like 1 and Ig-like 2 domain.