

**FITC Mouse Anti-human Ig Lambda Light Chain Antibody \*4C2, monoclonal, Cross Adsorbed\***Catalog number: V1031375  
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 Lambda Light Chain
Clone	4C2
Conjugate	FITC

**Biological Properties**

---

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

**Applications**

---

Immunoglobulin  $\lambda$ -1 light chain is a 27 kDa protein that can be found in the immunoglobulin complex, plasma membrane and extracellular region of cells. Sequencing of Immunoglobulin  $\lambda$ -1 light chain has shown it contains 2 conserved structural units: Ig-like 1 and Ig-like 2 domain. Immunoglobulin  $\lambda$ -1 light chain is thought to be essential to organismal processes, for instance, adaptive immune response.