

Purified Mouse Anti-human Ig Lambda Light Chain Antibody *Rs4, monoclonal*

Catalog number: V1031350

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	Secondary
Clonality	Monoclonal
Host	Mouse
Immunogen	Ig Lambda Light Chain
Clone	Rs4

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	ICC, ELISA

Applications

Immunoglobulin λ -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 λ -1 light chain has shown it contains 2 conserved structural units: Ig-like 1 and Ig-like 2 domain. Immunoglobulin λ -1 light chain is thought to be essential to organismal processes, for instance, adaptive immune response.