

**APC-Cy λ 153;7 Mouse Anti-human Ig
Lambda Light Chain Antibody *1-155-2,
monoclonal***Catalog number: V1031395
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	1-155-2
Conjugate	APC-Cy λ 153;7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC-Cy λ 153;7 under optimal conditions
Application	FC (QC TESTED)

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.