

Purified Mouse Anti-human Ig Kappa Light Chain Antibody *MEM-09, monoclonal*

Catalog number: V1031300

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 Kappa Light Chain
Clone	MEM-09

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	FC (QC TESTED), IP, WB, IHC(P), ICC, ELISA

Applications

Immunoglobulin κ 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 κ light chain plays an important role in organismal processes, for example, adaptive immune response. Sequencing of immunoglobulin κ light chain has demonstrated it contains 2 conserved structural units: Ig-like 1 and Ig-like 2 domain.