

Purified Mouse Anti-human Ig Kappa Light Chain Antibody *A8B5, monoclonal, Cross Adsorbed*

Catalog number: V1031305

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 | A8B5 |

Biological Properties

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