

**PE Mouse Anti-human Ig Lambda Light Chain Antibody \*1-155-2, monoclonal\***Catalog number: V1031390  
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          | PE                    |

**Biological Properties**

|             |                                                                                                   |
|-------------|---------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE 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.