

**PE-Cy5 Mouse Anti-human IgE
Antibody *BE5, monoclonal***Catalog number: V1031470
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 | IgE |
| Clone | BE5 |
| Conjugate | PE-Cy5 |

Biological Properties

| | |
|-------------|---|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE-Cy5 under optimal conditions |
| Application | FC (QC TESTED) |

Applications

Immunoglobulin ϵ heavy chain is a 70 kDa protein that can be expressed in the immunoglobulin complex, plasma membrane and extracellular region of cells. In humans, immunoglobulin ϵ heavy chain plays an important role in organismal processes, namely, adaptive immune response. Sequencing of immunoglobulin ϵ heavy chain has exemplified it contains 5 conserved structural units: Ig-like 1, Ig-like 2, Ig-like 3, Ig-like 4 and Ig-like 5 domain.