

**Biotin Mouse Anti-human IgE Antibody**  
**\*BE5, monoclonal\***Catalog number: V1031450  
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	Biotin

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with Biotin under optimal conditions
Application	ELISA

**Applications**

Immunoglobulin  $\epsilon$  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  $\epsilon$  heavy chain plays an important role in organismal processes, namely, adaptive immune response. Sequencing of immunoglobulin  $\epsilon$  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.