

Purified Mouse Anti-human IgE Antibody
4H10, monoclonalCatalog number: V1031445
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	4H10

Biological Properties

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

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.