

**PE Mouse Anti-human IgD Antibody *IA6-2,
monoclonal***Catalog number: V1031425
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	IgD
Clone	IA6-2
Conjugate	PE

Biological Properties

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

Applications

Immunoglobulin δ heavy chain is a 65 kDa protein that can be located in the plasma membrane, extracellular space and immunoglobulin complex of cells. Sequencing of immunoglobulin δ heavy chain has exemplified it contains 4 conserved structural units: Ig-like 1, Ig-like 2, Ig-like 3 and Ig-like 4 domain. Immunoglobulin δ heavy chain is thought to be essential to organismal processes, for instance, adaptive immune response and monocyte activation. It is involved in the positive regulation of interleukin-1 β secretion, interleukin-1 secretion and tumor necrosis factor secretion.