

**PE-Cy5 Mouse Anti-human IgM
Antibody *CH2, monoclonal***Catalog number: V1031540
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	IgM
Clone	CH2
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 μ heavy chain is a protein with a molecular weight of 74 kDa, located in the extracellular region, plasma membrane and immunoglobulin complex of cells. In humans, immunoglobulin μ heavy chain is an integral part of organismal processes, for example, adaptive immune response. Sequencing of Immunoglobulin μ heavy chain has shown it contains 5 conserved structural units: Ig-like 1, Ig-like 2, Ig-like 3, Ig-like 4 and Ig-like 5 domain.