

Purified Mouse Anti-human IgM Antibody
MA2, monoclonalCatalog number: V1031515
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	Tris buffered saline (TBS) solution, 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Secondary
Clonality	Monoclonal
Host	Mouse
Immunogen	IgM
Clone	MA2

Biological Properties

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

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.