

Purified Mouse Anti-human IgA Antibody
AD3, monoclonalCatalog number: V1031405
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	IgA
Clone	AD3

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	IHC(P), ELISA

Applications

Immunoglobulin heavy constant α 1 (also known as Ig α -1 chain C region TRO, Ig α -1 chain C region BUR or Ig α -1 chain C region) is a protein with a molecular weight of 44 kDa, located in the external side of plasma membrane, immunoglobulin complex, circulating and extracellular exosome of cells. Sequencing of Ig α -1 chain C region BUR has shown it contains 3 conserved structural units: Ig-like 1, Ig-like 2 and Ig-like 3 domain. It recognizes antigen. Ig α -1 chain C region BUR is the subject of comprehensive study stemming from the fact that it plays a role in the B cell receptor signaling pathway and complement activation, classical pathway. It is an integral part of organismal processes, for instance, retina homeostasis, leukocyte migration and antibacterial humoral response. It is an enhancer of respiratory burst and B cell activation.