

**Biotin Mouse Anti-human Albumin  
Antibody \*AL-01, monoclonal, Cross  
Adsorbed\***Catalog number: V103075  
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	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Albumin
Clone	AL-01
Conjugate	Biotin

**Biological Properties**

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against non-human primates serum and then conjugated with Biotin under optimal conditions
Application	IHC(P), ELISA, RIA, WB

**Applications**

Albumin is a protein with a molecular weight of 80 kDa ( 65-67 kDa ), found in the extracellular space, cytoplasm and nucleus of cells. In humans, albumin has been closely linked to important functions such as antioxidant activity. Sequencing of albumin has shown it contains 3 conserved structural units: albumin 1, albumin 2 and albumin 3 domain. It binds with zinc ion, fatty acid and copper ion. Albumin negatively regulates programmed cell death and apoptotic process. It has been found to be involved in organismal processes, namely, cellular response to starvation, receptor-mediated endocytosis and post-translational protein modification. Mutations and abnormalities in albumin have been associated with a number of diseases, for example, analbuminemia (ANALBA) and familial dysalbuminemic hyperthyroxinemia,(FDH). Analbuminemia, an autosomal recessive inheritance disorder characterized by hypoalbuminemia, hypotension and lipodystrophy, has in particular been of interest to scientists.