

Purified Mouse Anti-human CLIC5 Antibody
CLIC5-02, monoclonalCatalog number: V103330
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	CLIC5
Clone	CLIC5-02

Biological Properties

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

Applications

Chloride intracellular channel protein is a protein with a molecular weight of 28 kDa, located in the chloride channel complex and cytoplasm of cells. In humans, chloride intracellular channel protein is thought to be essential to organismal processes, for instance, sensory perception of sound. Chloride intracellular channel protein has been thought to be involved with important functions like chloride channel and voltage-gated ion channel activity. Sequencing of Chloride intracellular channel protein has shown evidence of it contains a primary structural unit, the GST C-terminal domain. Chloride intracellular channel protein is an integral part of ion transmembrane transport.