

Purified Mouse Anti-human SHIP1 Antibody
SHIP-02, monoclonalCatalog number: V1032105
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	SHIP1
Clone	SHIP-02

Biological Properties

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

Applications

Phosphatidylinositol 3,4,5-trisphosphate 5-phosphatase 1 (sometimes called Phosphatidylinositol 4,5-bisphosphate 5-phosphatase, p150Ship, , hp51CN, Inositol polyphosphate-5-phosphatase of 145 kDa, SH2 domain-containing inositol 5'-phosphatase 1, SIP-145 or Inositol polyphosphate-5-phosphatase D) is a protein with a molecular weight of 133 kDa, found in the plasma membrane, membrane raft and cytosol of cells. In humans, Phosphatidylinositol 3,4,5-trisphosphate 5-phosphatase 1 is thought to be essential to organismal processes, in particular, apoptotic process, immunoglobulin mediated immune response and leukocyte migration. Sequencing of phosphatidylinositol 3,4,5-trisphosphate 5-phosphatase 1 has supported it contains a primary structural unit, the SH2 domain. Phosphatidylinositol 3,4,5-trisphosphate 5-phosphatase 1 has been closely linked to vital functions like inositol-polyphosphate 5-phosphatase, inositol-1,3,4,5-tetrakisphosphate 5-phosphatase and phosphatidylinositol-4,5-bisphosphate 5-phosphatase activity. It is the subject of extensive research in part because of the fact that it is a member of the T cell receptor signaling pathway and cytokine-mediated signaling pathway.