

**Purified Mouse Anti-human SOCS3
Antibody *SO1, monoclonal***

Catalog number: V1032135

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	SOCS3
Clone	SO1

Biological Properties

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

Applications

Suppressor of cytokine signaling 3 (sometimes called SSI-3, Cytokine-inducible SH2 protein 3, STAT-induced STAT inhibitor 3 or CIS-3) is a protein with a molecular weight of 25 kDa, located in the phosphatidylinositol 3-kinase complex and cytosol of cells. In humans, suppressor of cytokine signaling 3 has been associated with key functions such as protein kinase inhibitor and 1-phosphatidylinositol-3-kinase regulator activity. It plays a role in the upregulation of tyrosine phosphorylation of STAT protein and cell differentiation but also is an inhibitor of tyrosine phosphorylation of STAT protein, apoptotic process and inflammatory response. Sequencing of suppressor of cytokine signaling 3 has shown it contains 2 types of conserved structural units: SOCS box and SH2 domain. Suppressor of cytokine signaling 3 binds to phosphotyrosine residue. Suppressor of cytokine signaling 3 is the subject of comprehensive research because it plays a role in the receptor signaling pathway via JAK-STAT, cytokine-mediated signaling pathway and regulation of interferon- γ -mediated signaling pathway. It is an integral part of organismal processes, for example, protein ubiquitination, cellular response to leukemia inhibitory factor and post-translational protein modification.