

**Purified Mouse Anti-human SCIMP
Antibody *NVL-07, monoclonal***

Catalog number: V1032095

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	SCIMP
Clone	NVL-07

Biological Properties

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

Applications

SLP adapter and CSK-interacting membrane protein is a 17 kDa transmembrane protein that can be found in the filopodium, tetraspanin-enriched microdomain and leading edge membrane of cells. In humans, SLP adapter and CSK-interacting membrane protein is an integral part of organismal processes, for instance, cellular response to molecule of fungal origin and cytokine production involved in immune response. It is the subject of intensive examination stemming from the fact that it plays a role in the toll-like receptor 7 signaling pathway, toll-like receptor 3 signaling pathway and toll-like receptor TLR1:TLR2 signaling pathway. SLP adapter and CSK-interacting membrane protein has been associated with vital functions like molecular adaptor activity. SLP adapter and CSK-interacting membrane protein recognizes microtubule and S-palmitoyl cysteine. It positively regulates interleukin-12 production, apoptotic process and dendritic cell cytokine production.