

Purified Mouse Anti-human SIT Antibody
SIT-01, monoclonal

Catalog number: V1032115

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	SIT
Clone	SIT-01

Biological Properties

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

Applications

Signaling threshold-regulating transmembrane adapter 1 is a 21 kDa transmembrane protein that can be located in the integral component of plasma membrane and extracellular exosome of cells. It is also known as Suppression-inducing transmembrane adapter 1, gp30/40 and SHP2-interacting transmembrane adapter protein. Sequencing of suppression-inducing transmembrane adapter 1 has shown it contains 2 types of conserved structural units: extracellular and cytoplasmic domain. It plays an important role in organismal processes, in particular, T cell homeostasis, adaptive immune response and signal transduction. Signaling threshold-regulating transmembrane adapter 1 aids in T cell activation.