

**Purified Mouse Anti-human/rabbit/rat
STRO-1 Antibody *STRO-1, monoclonal***Catalog number: V1032165
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	Tris buffered saline (TBS) solution, 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human, rabbit, rat
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	STRO-1
Clone	STRO-1

Biological Properties

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

Applications

Hematopoietic progenitor cell antigen CD34 is a 48 kDa transmembrane protein that can be found in the intercellular bridge, integral component of plasma membrane and cytoplasm of cells. In Homo sapiens, hematopoietic progenitor cell antigen CD34 has been found to be involved in immune response and blood pressure, and in addition, takes part in processes such as organismal processes, namely, vascular wound healing, tissue homeostasis and mesangial cell-matrix adhesion. It plays a role in the downregulation of blood coagulation, gene expression and neuron death but also is a positive regulator of interleukin-10 production, granulocyte colony-stimulating factor production and gene expression. Sequencing of hematopoietic progenitor cell antigen CD34 has shown it contains 2 conserved structural units: cytoplasmic and extracellular domain. Hematopoietic progenitor cell antigen CD34 binds with carbohydrate, transcription factor and sulfate.