

**Purified Mouse Anti-human/mouse/rat SYK  
Antibody \*SYK-01, monoclonal\***

Catalog number: V1032195

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, mouse, rat
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	SYK
Clone	SYK-01

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	IP, WB, IHC(P), ICC

**Applications**

---

Tyrosine-protein kinase SYK (also known as Spleen tyrosine kinase or p72-Syk) is a protein with a molecular weight of 72 kDa, located in the plasma membrane, cytoplasm and nucleus of cells. In humans, tyrosine-protein kinase SYK is an integral part of organismal processes, in particular, viral process, adaptive immune response and lymph vessel development. It is the subject of comprehensive application because it plays a role in the B cell receptor signaling pathway, interleukin-3-mediated signaling pathway and stimulatory C-type lectin receptor signaling pathway. Sequencing of tyrosine-protein kinase SYK has exemplified it contains 3 types of conserved structural units: SH2 2, SH2 1 and protein kinase domain. Tyrosine-protein kinase SYK reacts with phosphatase, Toll-like receptor and protein kinase, and moreover, plays an important role in platelet activation, phagocytosis and arachidonic acid secretion. It has been thought to be involved with essential functions such as protein tyrosine kinase, protein serine/threonine kinase and non-membrane spanning protein tyrosine kinase activity.