

**PE Mouse Anti-human LIME1 Antibody**  
**\*LIME-06, monoclonal\***

Catalog number: V1031675

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	LIME1
Clone	LIME-06
Conjugate	PE

**Biological Properties**

---

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

**Applications**

---

Lck-interacting transmembrane adapter 1 is a 30 kDa transmembrane protein that can be expressed in the integral component of membrane and B cell receptor complex of cells. In mice, Lck-interacting transmembrane adapter 1 is the subject of intensive examination because of the fact that it plays a role in the T cell receptor signaling pathway and B cell receptor signaling pathway. It is thought to be essential to organismal processes, for example, adaptive immune response. Lck-interacting transmembrane adapter 1 is an integral part of release of sequestered calcium ion into cytosol, I-kappaB kinase/NF-kappaB signaling and transcription by RNA polymerase II. It recognizes protein kinase and S-palmitoyl cysteine. Sequencing of Lck-interacting transmembrane adapter 1 has supported it contains 2 types of conserved structural units: cytoplasmic and extracellular domain.