

**Purified Mouse Anti-human LST1 Antibody**  
**\*LST1/02, monoclonal\***

Catalog number: V1031695

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          | LST1       |
| Clone              | LST1/02    |

**Biological Properties**

---

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

**Applications**

---

Leukocyte-specific transcript 1 protein is a transmembrane protein with a molecular weight of 11 kDa, expressed in the integral component of membrane, cytoplasm and Golgi apparatus of cells. In humans, leukocyte-specific transcript 1 protein is an integral part of organismal processes, for example, cell morphogenesis, immune response and dendrite development. Leukocyte-specific transcript 1 protein is thought to be essential to cell shape, and additionally, is an inhibitor of lymphocyte proliferation.