

Purified Mouse Anti-human LCK Antibody
LCK-01, monoclonal, Cross Adsorbed

Catalog number: V1031650

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	LCK
Clone	LCK-01

Biological Properties

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against mouse serum and then conjugated with under optimal conditions
Application	FC, IP, WB, ICC

Applications

Tyrosine-protein kinase Lck is a 58 kDa protein that can be located in the pericentriolar material, extracellular exosome and cytosol of cells. It is alternatively called Lymphocyte cell-specific protein-tyrosine kinase, Protein YT16 and T cell-specific protein-tyrosine kinase. In Homo sapiens, tyrosine-protein kinase Lck has been thought to be involved with key functions such as protein tyrosine kinase, protein serine/threonine phosphatase and non-membrane spanning protein tyrosine kinase activity. It is an integral part of lymphocyte activation, cell population proliferation and defense response to virus. Tyrosine-protein kinase Lck is the subject of intensive examination in part because of the fact that it is a component of the positive regulation of T cell receptor signaling pathway, T cell receptor signaling pathway and positive regulation of intrinsic apoptotic signaling pathway. It is involved in the positive regulation of T cell activation, protein kinase B signaling and leukocyte cell-cell adhesion, and in addition, plays an important role in organismal processes, namely, cell differentiation, protein phosphorylation and T cell costimulation. Sequencing of tyrosine-protein kinase Lck has exemplified it contains 3 types of conserved structural units: protein kinase, SH2 and SH3 domain. Tyrosine-protein kinase Lck reacts with ATP, T cell receptor and protein phosphatase. Mutations and abnormalities in tyrosine-protein kinase Lck have been associated with a number of diseases, for example, immunodeficiency 22 (IMD22). Immunodeficiency 22, an autosomal recessive inheritance disorder characterized by failure to thrive, autoimmunity and immunodeficiency, has in specific been of interest to scientists.