

Purified Rat Anti-mouse LPAM-1 Antibody
DATK32, monoclonal

Catalog number: V1031680

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	Mouse
Class	Primary
Clonality	Monoclonal
Host	Rat
Immunogen	LPAM-1
Clone	DATK32

Biological Properties

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

Applications

Integrin β -7 is a transmembrane protein with a molecular weight of 101 kDa, expressed in the focal adhesion, integrin complex and plasma membrane of cells. In humans, integrin β -7 is thought to be essential to immune response. It reacts with integrin, metal ion and cell adhesion molecule. It is the subject of extensive research due to the fact that it is a member of the integrin-mediated signaling pathway. It has been closely linked to essential functions such as virus receptor and signaling receptor activity. Sequencing of Integrin β -7 has exemplified it contains 3 conserved structural units: VWFA, extracellular and cytoplasmic domain. It aids in organismal processes, namely, cell migration, T cell migration and extracellular matrix organization.