

Purified Mouse Anti-human/non-human primates LAR Antibody *W7C6, monoclonal*Catalog number: V1031630
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, non-human primates
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	LAR
Clone	W7C6

Biological Properties

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

Applications

Protein-tyrosine-phosphatase is a transmembrane protein with a molecular weight of 167 kDa, located in the integral component of membrane of cells. In humans, protein-tyrosine-phosphatase has been associated with essential functions like protein tyrosine phosphatase activity. Sequencing of protein-tyrosine-phosphatase has demonstrated it contains 10 types of conserved structural units: Ig-like, fibronectin type-III, TYR_phosphatase_2 and tyrosine-protein phosphatase domain.