

PE Mouse Anti-human PCLO Antibody
PCLO-01, monoclonal

Catalog number: V1031980

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	PCLO
Clone	PCLO-01
Conjugate	PE

Biological Properties

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

Applications

Protein piccolo is a 561 kDa protein that can be expressed in the extracellular exosome, cytoskeleton and GABA-ergic synapse of cells. In Homo sapiens, protein piccolo aids in organismal processes, for example, cAMP-mediated signaling, synaptic vesicle clustering and cytoskeleton organization. Sequencing of protein piccolo has supported it contains 3 types of conserved structural units: PDZ, C2 1 and C2 2 domain. It is the subject of comprehensive study in part because of the fact that it is a component of the presynapse to nucleus signaling pathway. It is an integral part of exocytosis. Protein piccolo binds with profilin, calcium-dependent phospholipid and calcium ion. It is clinically significant because abnormalities in its function have been thought to be involved with diseases like pontocerebellar hypoplasia type 3 (PCH3). Pontocerebellar hypoplasia type 3, an autosomal recessive inherited disorder characterized by cerebral atrophy, macrotia and proptosis, has in particular been of interest to investigators.