

**Purified Mouse Anti-human CPNE7
Antibody *CPNE7-01, monoclonal***

Catalog number: V103375

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	CPNE7
Clone	CPNE7-01

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	WB

Applications

Copine-7 is a protein with a molecular weight of 70 kDa, found in the nucleus, plasma membrane and cytoplasm of cells. In humans, copine-7 has been associated with key functions such as transporter activity. Sequencing of copine-7 has shown it contains 3 types of conserved structural units: C2 1, C2 2 and VWFA domain. Copine-7 binds to calcium-dependent phospholipid and metal ion, and also, is an integral part of organismal processes, for example, glycerophospholipid biosynthetic process, cellular response to calcium ion and lipid metabolic process.