

**Purified Mouse Anti-human/mouse NCK1
Antibody *EM-06, monoclonal***

Catalog number: V1031820

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, mouse
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	NCK1
Clone	EM-06

Biological Properties

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

Applications

Cytoplasmic protein NCK1 is a 43 kDa protein that can be found in the endoplasmic reticulum, cytoplasm and ribosome of cells. It is also known as NCK adaptor protein 1 and SH2/SH3 adaptor protein NCK- α . In humans, NCK adaptor protein 1 takes part in processes such as cell migration. It is thought to be essential to organismal processes, namely, signal complex assembly, lamellipodium assembly and response to other organism. It has been thought to be involved with vital functions like signaling receptor complex adaptor, cytoskeletal anchor and protein-macromolecule adaptor activity. Sequencing of NCK adaptor protein 1 has exemplified it contains 4 types of conserved structural units: SH3 1, SH3 2, SH3 3, and SH2 domain. NCK adaptor protein 1 binds with cadherin, ephrin receptor and eukaryotic initiation factor eIF2.