

**Purified Mouse Anti-human/mouse/pig
Kinesin Heavy Chain Antibody *KN-03,
monoclonal***Catalog number: V1031605
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	Tris buffered saline (TBS) solution, 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human, mouse, pig, rat
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Kinesin Heavy Chain
Clone	KN-03

Biological Properties

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

Applications

Kinesin-1 heavy chain is a 127 kDa protein that can be expressed in the vesicle, centriolar satellite and microtubule of cells. It is alternatively called ubiquitous kinesin heavy chain and conventional kinesin heavy chain. In humans, kinesin-1 heavy chain has been associated with key functions such as ATPase and microtubule motor activity. It binds to ATP, JUN kinase and microtubule. Sequencing of Kinesin-1 heavy chain has exemplified it contains a primary structural unit, the kinesin motor domain. It takes part in processes such as organismal processes. Kinesin-1 heavy chain aids in membrane potential, and in addition, positively regulates potassium ion transport, protein localization to plasma membrane and vesicle fusion.