

Purified Mouse Anti-human/mouse/rat Clathrin Heavy Chain Antibody *BF-06, monoclonal*

Catalog number: V103325

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	Stabilizing Tris buffered saline (TBS) solution, 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human, mouse, rat, cow, pig
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Clathrin Heavy Chain
Clone	BF-06

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	FC (QC TESTED), IP, WB, ICC, ELISA

Applications

Clathrin heavy chain is a protein with a molecular weight of 187 kDa, expressed in the cytosol, actin cortical patch and clathrin complex of cells. In *Saccharomyces cerevisiae* (strain ATCC 204508 / S288c), clathrin heavy chain has been closely linked to essential functions such as structural molecule activity. Clathrin heavy chain is an integral part of organismal processes, for example, receptor-mediated endocytosis, intracellular protein transport and Golgi to endosome transport. It binds with clathrin light chain. Clathrin heavy chain has been closely linked to critical functions such as structural molecule activity.