

**Purified Mouse Anti-human/mouse/cow
HSP90 β Antibody *MBH90B, monoclonal***Catalog number: V1031235
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, cow
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	HSP90 β
Clone	MBH90B

Biological Properties

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

Applications

Heat shock protein HSP 90- β is a protein with a molecular weight of 96 kDa, found in the protein-containing complex, apical plasma membrane and ficolin-1-rich granule lumen of cells. In humans, heat shock protein HSP 90- β reacts with histone methyltransferase, DNA polymerase and ion channel. It has been closely linked to vital functions like nitric-oxide synthase regulator, protein kinase regulator and protein dimerization activity, and moreover, takes part in processes such as protein ubiquitination, cellular protein localization and cellular response to heat. It is the subject of extensive application because of the fact that it plays a role in the regulation of interferon- γ -mediated signaling pathway, regulation of type I interferon-mediated signaling pathway and Fc- γ receptor signaling pathway involved in phagocytosis, and furthermore, is a negative regulator of protein metabolic process, proteasomal protein catabolic process and transforming growth factor β activation. On the other hand, it also acts to positively regulate protein binding, protein localization to cell surface and nitric oxide biosynthetic process. Heat shock protein HSP 90- β aids in organismal processes, namely, placenta development, response to cocaine and cellular response to drug.