

**Purified Mouse Anti-hepatitis b virus HBsAg
(HBV) Antibody *HB5, monoclonal***Catalog number: V103785
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	Hepatitis b virus
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	HBsAg (HBV)
Clone	HB5

Biological Properties

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

Applications

Large envelope protein is a 49 kDa transmembrane protein that can be expressed in the virion membrane and integral component of membrane of cells. It is also known as L glycoprotein, major surface antigen and large surface protein. In Hepatitis B virus, major surface antigen is thought to be essential to organismal processes, in particular, caveolin-mediated endocytosis of virus by host cell, virion attachment to host cell and fusion of virus membrane with host endosome membrane. It recognizes N-myristoyl glycine by host. Sequencing of major surface antigen has demonstrated it contains 6 conserved structural units.