

**Purified Mouse Anti-herpes simplex virus
Glycoprotein C (HSV-1) Antibody *T96,
monoclonal***Catalog number: V103750
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	Herpes simplex virus
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Glycoprotein C (HSV-1)
Clone	T96

Biological Properties

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

Applications

Envelope glycoprotein C is a 55 kDa transmembrane protein that can be located in the virion membrane and integral component of membrane of cells. Sequencing of envelope glycoprotein C has demonstrated it contains 3 types of conserved structural units: virion surface, cytoplasmic and Ig-like domain. Envelope glycoprotein C aids in organismal processes, for instance, suppression by virus of host complement activation, viral entry into host cell and adhesion receptor-mediated virion attachment to host cell, and in addition, recognizes complement.