

**Purified Mouse Anti-herpes simplex virus
Glycoprotein B (HSV-1/HSV-2) Antibody
*T111, monoclonal***

Catalog number: V103745
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 B (HSV-1/HSV-2)
Clone	T111

Biological Properties

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

Applications

Envelope glycoprotein B (gB) is a 100 kDa transmembrane protein that can be found in the viral envelope, virion membrane and host cell plasma membrane of cells. Sequencing of envelope glycoprotein B has shown evidence of it contains 2 conserved structural units: intravirion and virion surface domain. Envelope glycoprotein B takes part in organismal processes, namely, virion attachment to host cell and viral entry into host cell.