

Purified Mouse Anti-herpes simplex virus Glycoprotein G (HSV-2) Antibody *T303, monoclonal*

Catalog number: V103755

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 G (HSV-2)
Clone	T303

Biological Properties

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

Applications

Major surface glycoprotein G (sometimes called attachment glycoprotein G, mG or membrane-bound glycoprotein) is a transmembrane protein with a molecular weight of 38 kDa, expressed in the extracellular region, host cell surface and integral component of membrane of cells. Sequencing of major surface glycoprotein G has demonstrated it contains 2 conserved structural units: intravirion and virion surface domain. Major surface glycoprotein G is thought to be essential to organismal processes, for instance, virion attachment to host cell, mitigation of host immune response by virus and viral entry into host cell.