
Major Surface Glycoprotein G Polyclonal Antibody

(Catalog # A50853)

Background

Attaches the virion to the host cell membrane by interacting with heparan sulfate, initiating the infection. Interacts with host CX3CR1, the receptor for the CX3C chemokine fractalkine, to modulate the immune response and facilitate infection. Unlike the other paramyxovirus attachment proteins, lacks both neuraminidase and hemagglutinating activities. Secreted glycoprotein G helps RSV escape antibody-dependent restriction of replication by acting as an antigen decoy and by modulating the activity of leukocytes bearing Fcγ receptors.

Description

Major Surface Glycoprotein G Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human respiratory syncytial virus A

Isotype

IgG

Uniprot ID

P03423

Purification

>95%, Protein G purified

Immunogen

Recombinant Human respiratory syncytial virus A Major surface glycoprotein G protein (67-298AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Attachment glycoprotein G, Membrane-bound glycoprotein, mG, G

Application

ELISA; Recommended dilution: ELISA 1:1000-5000
