
Scytovirin Polyclonal Antibody

(Catalog # A64238)

Background

Has strong anti-HIV activity against T-tropic strains of HIV-1 and weaker activity against M-tropic strains of HIV-1. Inhibits HIV-1 fusion and infection of CD4 LTR beta-gal cells in vitro. Inhibits fusion of HIV infected CEM-SS cells with uninfected CEM-SS cells, and fusion of HIV-1 Env expressing HL2/3 cells with CD4 LTR beta-gal cells. Binds to HIV gp120, HIV gp160 and to a lesser extent HIV gp41. Binding to HIV gp120 is glycosylation dependent. Binds with high specificity to the tetrasaccharide Man-alpha-1,2-Man-alpha-1,6-Man-alpha-1,6-Man and also binds the higher-order oligosaccharides oligomannose 8 and oligomannose 9. Does not bind to monosaccharides, complex or hybrid N-linked oligosaccharides or chitin.

Description

Scytovirin Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Scytonema varium

Isotype

IgG

Uniprot ID

P86041

Purification

>95%, Protein G purified

Immunogen

Recombinant Scytonema varium Scytovirin protein (1-95AA)

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

Scytovirin1

Application

ELISA
