
VP40 Polyclonal Antibody

(Catalog # A66074)

Background

Promotes virus assembly and budding by interacting with host proteins of the multivesicular body pathway. May facilitate virus budding by interacting with the nucleocapsid and the plasma membrane. Specific interactions with membrane-associated GP and VP24 during the budding process may also occur. The hexamer form seems to be involved in budding. The octamer form binds RNA, and may play a role in genome replication.

Description

VP40 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Reston ebolavirus

Isotype

IgG

Uniprot ID

Q8JPX9

Purification

>95%, Protein G purified

Immunogen

Recombinant Reston ebolavirus Matrix protein VP40 protein (1-331AA)

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

Matrix protein VP40, Membrane-associated protein VP40, VP40

Application

ELISA
