
VP3 Polyclonal Antibody, HRP Conjugated

(Catalog # A55763)

Background

May act as transcriptional regulator. Induces apoptosis in infected cells. Element of infectious replication cycle.

Description

VP3 Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

Formulation

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

Specificity

Chicken anemia virus

Isotype

IgG

Uniprot ID

Q99152

Purification

>95%, Protein G purified

Immunogen

Recombinant Chicken anemia virus Apoptin protein (1-121AA)

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

VP3

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
