
III Polyclonal Antibody

(Catalog # A64208)

Background

Plays essential roles both in the penetration of the viral genome into the bacterial host via pilus retraction and in the extrusion process. During the initial step of infection, G3P mediates adsorption of the phage to its primary receptor, the tip of host F-pilus. Subsequent interaction with the host entry receptor tolA induces penetration of the viral DNA into the host cytoplasm. In the extrusion process, G3P mediates the release of the membrane-anchored virion from the cell via its C-terminal domain.

Description

III Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Enterobacteria phage M13

Isotype

IgG

Uniprot ID

P69168

Purification

>95%, Protein G purified

Immunogen

Recombinant Enterobacteria phage M13 Attachment protein G3P protein (19-424AA)

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 protein G3P, Gene 3 protein, G3P, Minor coat protein, III

Application

ELISA
