
ipaH9.8 Polyclonal Antibody

(Catalog # A55475)

Background

Effector proteins function to alter host cell physiology and promote bacterial survival in host tissues. This protein is an E3 ubiquitin ligase that interferes with host's ubiquitination pathway and modulates the acute inflammatory responses, thus facilitating bacterial colonization within the host cell.

Description

ipaH9.8 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Shigella flexneri serotype X

Isotype

IgG

Uniprot ID

D2AJU0

Purification

>95%, Protein G purified

Immunogen

Recombinant Shigella flexneri serotype X E3 ubiquitin-protein ligase ipaH9.8 protein (1-545AA)

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

Invasion plasmid antigen ipaH9.8, ipaH9.8, SFxv_5076

Application

ELISA; Recommended dilution: 1:1000 – 1:5000
