
vacA Polyclonal Antibody

(Catalog # A64398)

Background

Induces vacuolation of eukaryotic cells. Causes ulceration and gastric lesions.

Description

vacA Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Helicobacter pylori

Isotype

IgG

Uniprot ID

P55981

Purification

>95%, Protein G purified

Immunogen

Recombinant Helicobacter pylori Vacuolating cytotoxin autotransporter protein (37-245AA)

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

Vacuolating cytotoxin autotransporterCleaved into the following 2 chains:Vacuolating cytotoxinVacuolating cytotoxin translocator, vacA, HP_0887

Application

ELISA
