
cagA Polyclonal Antibody

(Catalog # A64204)

Background

May be necessary for the transcription, folding, export, or function of the cytotoxin.

Description

cagA 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

P80200

Purification

>95%, Protein G purified

Immunogen

Recombinant Helicobacter pylori Cytotoxicity-associated immunodominant antigen protein (918-1147AA)

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

Cytotoxicity-associated immunodominant antigen, 120 kDa protein, cagA, cai

Application

ELISA; Recommended dilution: ELISA 1:1000
