

ureA Polyclonal Antibody

(Catalog # A53606)

Background

Ammonia produced by ureolysis increases the gastric pH thereby providing an environment permissive for colonization of the stomach.

Description

ureA 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

P14916

Purification

>95%, Protein G purified

Immunogen

Recombinant Helicobacter pylori Urease subunit alpha protein (1-238AA)

Storage

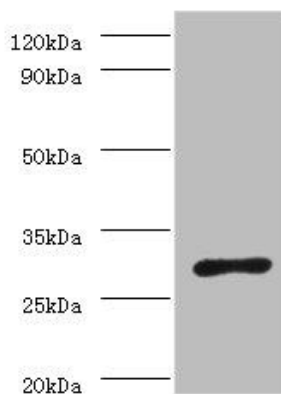
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Tubulin-folding cofactor C TBCC

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: ureA Polyclonal Antibody at 2ug/ml+Helicobacter pylori bacteria liquid
Goat polyclonal to rabbit at 1/10000 dilution
Predicted band size: 27kDa
Observed band size: 27kDa