

PRODUCT INFORMATION

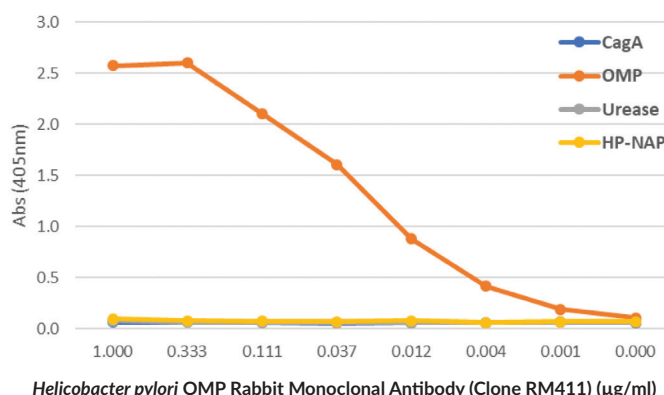


Helicobacter pylori OMP Rabbit Monoclonal Antibody (Clone RM411) Item No. 32342

Overview and Properties

Contents:	This vial contains 100 µg of protein A-affinity purified monoclonal antibody.
Synonyms:	<i>H. pylori</i> , Outer Membrane Protein
Immunogen:	Proteins purified from <i>H. pylori</i>
Cross Reactivity:	(+) OMP; (-) CagA, HP-NAP, Urease
Species Reactivity:	(+) <i>H. pylori</i>
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM411
Host:	Rabbit
Isotype:	IgG
Application:	ELISA; the recommended starting concentration is 0.05-0.5 µg/ml. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Helicobacter pylori OMP Rabbit Monoclonal Antibody (Clone RM411) (µg/ml)

ELISA of *Helicobacter Pylori* protein using *Helicobacter pylori* OMP Rabbit Monoclonal Antibody (Clone RM411). The plate was coated with 1 µg/ml of CagA, OMP, Urease, or HP-NAP of *H. Pylori*. *Helicobacter pylori* OMP Rabbit Monoclonal Antibody (Clone RM411) was used as the primary antibody and an alkaline phosphatase-conjugated anti-rabbit IgG was used as the secondary antibody.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/30/2021

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Helicobacter pylori (*H. pylori*) is a Gram-negative bacterium that can infect and colonize the stomach, leading to chronic gastritis, stomach inflammation and oxidative stress, peptic ulcer disease, and gastric cancer.¹ *H. pylori* expresses greater than 30 outer membrane proteins (OMPs) on the surface of its outer membrane encoded by two gene families, helicobacter outer membrane porins (*hop*) and hop-related (*hor*).²⁻⁴ OMPs have roles in *H. pylori* attachment and colonization, the expression of pathogenic factors, antibiotic resistance, and the maintenance of the outer membrane integrity. Cayman's *Helicobacter pylori* OMP Rabbit Monoclonal Antibody (Clone RM411) can be used for ELISA. The antibody recognizes a 23 kDa outer membrane protein.

References

1. Chang, A.H. and Parsonnet, J. Role of bacteria in oncogenesis. *Clin. Microbiol. Rev.* **23**(4), 837-857 (2010).
2. Matsuo, Y., Kido, Y., Yamaoka, Y., et al. *Helicobacter pylori* outer membrane protein-related pathogenesis. *Toxins (Basel)* **9**(3), 101 (2017).
3. Xu, C., Soyfoo, D.M., Wu, Y., et al. Virulence of *Helicobacter pylori* outer membrane proteins: An updated review. *Eur. J. Clin. Microbiol. Infect. Dis.* **39**(10), 1821-1830 (2020).
4. Odenbreit, S., Swoboda, K., Barwig, I., et al. Outer membrane protein expression profile in *Helicobacter pylori* clinical isolates. *Infect. Immun.* **77**(9), 3782-3790 (2009).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM