

# PRODUCT INFORMATION

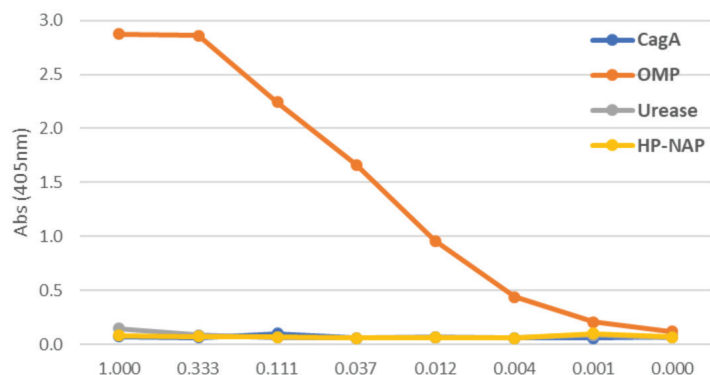


## *Helicobacter pylori* OMP Rabbit Monoclonal Antibody (Clone RM410) Item No. 32341

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µg of protein A-affinity purified monoclonal antibody.
<b>Synonym:</b>	Outer Membrane Protein
<b>Immunogen:</b>	Proteins purified from <i>H. pylori</i>
<b>Cross Reactivity:</b>	(+) OMP; (-) CagA, Urease, HP-NAP
<b>Species Reactivity:</b>	(+) <i>H. pylori</i>
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
<b>Clone:</b>	RM410
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	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 RM410) (µl/ml)

A titer ELISA using *Helicobacter pylori* OMP Rabbit Monoclonal Antibody (Clone RM410). The plate was coated with 1 µg/ml of *H. pylori* CagA, OMP, urease, or NAP. A serial dilution of *Helicobacter pylori* OMP Rabbit Monoclonal Antibody (Clone RM410) 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, 03/09/2021

**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* 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.<sup>1</sup> *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*).<sup>2-4</sup> OMPs have roles in *H. pylori* attachment and colonization, the expression of pathogenic factors, antibiotic resistance, and maintenance of the outer membrane integrity. Cayman's *Helicobacter pylori* OMP Rabbit Monoclonal Antibody (Clone RM410) can be used for ELISA. The antibody recognizes a 23 kDa OMP.

## 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., and Yamaoka, Y. *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