

PRODUCT INFORMATION

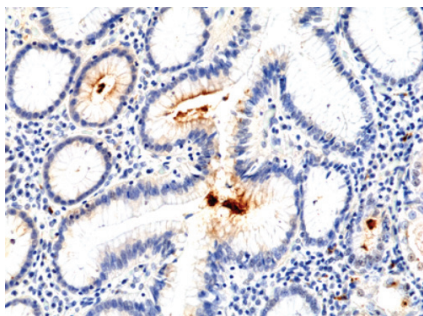


Helicobacter pylori Rabbit Monoclonal Antibody (RM400) Item No. 32325

Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	<i>H. pylori</i>
Immunogen:	Proteins purified from <i>H. pylori</i>
Cross Reactivity:	(+) <i>H. pylori</i>
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:	RM400
Host:	Rabbit
Isotype:	IgG
Application:	Immunohistochemistry; the recommended starting dilution is 1:100-1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Immunohistochemical staining of formalin-fixed and paraffin-embedded human stomach tissue using *Helicobacter pylori* Rabbit Monoclonal Antibody (RM400) at a dilution of 1:100.

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, 02/19/2024

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.¹ Transmission of *H. pylori* can occur directly through person-to-person contact, as well as through contaminated drinking water, raw vegetables, or animal milk.² *H. pylori* infection can lead to chronic gastritis, stomach inflammation and oxidative stress, peptic ulcer disease, and gastric cancer.¹ It is considered an oncogenic agent and, although a low percentage of patients infected with *H. pylori* develop stomach cancer, it is responsible for the majority of stomach cancers. In mouse models of *H. pylori* infection, antibiotics can reduce and prevent gastritis and pre-malignant changes. Cayman's *Helicobacter pylori* Rabbit Monoclonal Antibody (RM400) can be used for immunohistochemistry (IHC).

References

1. Chang, A.H. and Parsonnet, J. Role of bacteria in oncogenesis. *Clin. Microbiol. Rev.* **23(4)**, 837-857 (2010).
2. Mezmale, L., Coelho, L.G., Bordin, D., *et al.* Review: Epidemiology of *Helicobacter pylori*. *Helicobacter* **25(Suppl 1)**, e12734 (2020).

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