

H. Pylori Stain Kit (Microorganism Stain) ab150673

1 References 2 Images

Overview

Product name	H. Pylori Stain Kit (Microorganism Stain)
Product overview	The <i>Helicobacter pylori</i> Rapid Stain (Microorganism Stain) is designed for demonstrating <i>Helicobacter pylori</i> infected tissue. Kit may be used on formalin fixed, paraffin-embedded tissue as well as frozen sections.

Other products for staining tissue sections

Find more kits and reagents in the [special stains guide](#), or products for antigen retrieval, blocking, signal amplification, visualization, counterstaining, and mounting in the [IHC kits and reagents guide](#).

Notes

Staining Interpretation

<i>Helicobacter pylori</i>	Blue
Mucin	Yellow
Background	Light Blue

Control Tissue: *Helicobacter pylori* infected stomach

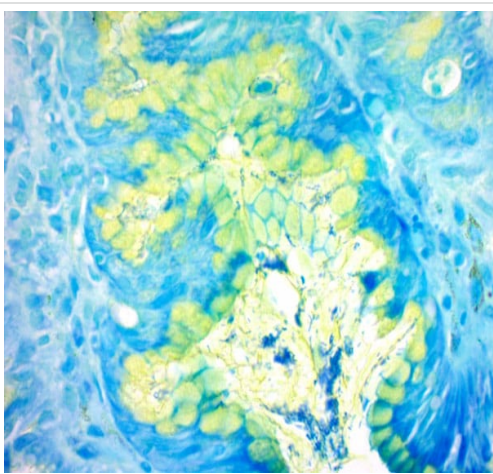
Properties

Components	100 tests
Alcian Yellow solution	1 x 125ml
Periodic acid solution	1 x 500ml
Sodium Hydroxide Solution (3%)	1 x 8ml
Sodium Metabisulfite Solution	1 x 125ml
Toluidine Blue Solution	2 x 6ml

## Relevance

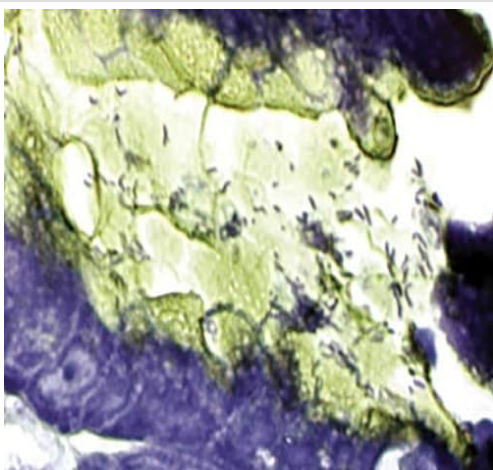
The spiral shaped bacterium *Helicobacter pylori* is strongly associated with inflammation of the stomach and is also implicated in the development of gastric malignancy. *H. pylori* is known to cause peptic ulcers and chronic gastritis in human. It is associated with duodenal ulcers and may be involved in development of adenocarcinoma and low-grade lymphoma of mucosa associated lymphoid tissue in the stomach. More recently this bacterium has also been implicated with a number of vascular disorders including heart disease. It is not clear how *H. pylori* is transmitted or why some patients become symptomatic while others do not. The bacteria are most likely spread from person to person through fecal-to-oral or oral-to-oral routes. Possible environmental reservoirs include contaminated water sources. Serological tests that measure specific *H. pylori* IgG antibodies can determine if a person has been infected although these methods do have drawbacks and limitations.

## Images



ab150673 H. Pylori Stain Kit (Microorganism Stain) staining formalin-fixed-paraffin embedded human stomach.

ab150673 - H. Pylori Stain Kit (Microorganism Stain)



Staining using ab150673 - H. pylori Rapid Stain Kit.

ab150673 - H. Pylori Stain Kit (Microorganism Stain)

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## **Our Abpromise to you: Quality guaranteed and expert technical support**

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
  
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

## **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors