



# *Helicobacter pylori* IgG ELISA Kit

Catalog Number KA0220

96 assays

Version: 03

Intended for research use only

[www.abnova.com](http://www.abnova.com)

## Table of Contents

<b>Introduction .....</b>	<b>3</b>
Intended Use .....	3
Background .....	3
Principle of the Assay .....	3
<b>General Information .....</b>	<b>4</b>
Materials Supplied .....	4
Storage Instruction .....	4
Materials Required but Not Supplied .....	4
Precautions for Use .....	4
<b>Assay Protocol .....</b>	<b>6</b>
Reagent Preparation .....	6
Sample Preparation .....	6
Assay Procedure .....	6
<b>Data Analysis .....</b>	<b>7</b>
Calculation of Results .....	7
Performance Characteristics .....	8
<b>Resources .....</b>	<b>9</b>
References .....	9
Plate Layout .....	10

## **Introduction**

### **Intended Use**

The *Helicobacter pylori* IgG ELISA Kit is intended for the quantitative determination of IgG antibodies to *Helicobacter pylori* in human serum.

### **Background**

*Helicobacter pylori* is a spiral bacterium cultured from human gastric mucosa discovered by B.J. Marshall in 1982. Studies have indicated that the presence of *H. pylori* is associated with a variety of gastrointestinal diseases including gastritis, duodenal and gastric ulcers, non-ulcer dyspepsia, and gastric adenocarcinoma and lymphoma. The organism is present in 95-98% of patients with duodenal ulcers and 60-90% of patients with gastric ulcers. The studies have also demonstrated that removal of the organism by anti-microbial therapy is correlated with the resolution of symptoms and cure of diseases.

Patients who present clinical symptoms relating to the gastrointestinal tract can be diagnosed for *H. pylori* infection by two methods:

- (1) Invasive techniques – include biopsy followed by culture or histologic examination of biopsy specimen or direct detection of urease activity.
- (2) Non-invasive techniques – include urea breath tests and serological methods.

All of the testing performed on biopsy samples is subject to errors related to sampling and interference of contaminated bacteria. An ELISA test of the presence of *H. pylori* specific IgG antibody is the technique of choice for serologic tests because of its accuracy and simplicity.

### **Principle of the Assay**

Purified *H. pylori* antigen is coated on the surface of microwells. Diluted serum is added to the wells, and the *H. pylori* IgG-specific antibody, if present, binds to the antigen. All unbound materials are washed away. Enzyme conjugate is added, which binds to the antibody-antigen complex. Excess enzyme conjugate is washed off and a solution of TMB Reagent is added. The enzyme conjugate catalytic reaction is stopped at a specific time. The intensity of the color generated is proportional to the amount of IgG-specific antibody in the sample. The results are read by a microwell reader compared in a parallel manner with calibrator and controls.

## General Information

### Materials Supplied

List of component

Component	Amount
Purified <i>H. pylori</i> antigen coated microtiter plate	96 wells
Enzyme Conjugate Reagent (red color)	13 mL
Sample Diluent (green color)	22 mL
<i>H. pylori</i> Negative Control, < 6.25 U/mL	100 µL
<i>H. pylori</i> Standards, 0, 6.25, 12.5, 25, 50, and 100 U/mL	100 µL each
<i>H. pylori</i> Positive Control, > 100 U/mL	100 µL
Wash Buffer (20×)	50 mL
TMB Reagent (One-Step)	11 mL
Stop Solution (1 N HCl)	11 mL

### Storage Instruction

Unopened test kits should be stored at 2-8°C upon receipt and the microtiter plate should be kept in a sealed bag with desiccants to minimize exposure to damp air. Opened test kits will remain stable until the expiration date shown, provided it is stored as described above. A microtiter plate reader with a bandwidth of 10 nm or less and an optical density range of 0-2 OD or greater at 450 nm wavelength is acceptable for use in absorbance measurement.

### Materials Required but Not Supplied

- ✓ Distilled water.
- ✓ Precision pipettes: 5 µL, 100 µL and 200 µL.
- ✓ Disposable pipette tips.
- ✓ Vortex mixer or equivalent.
- ✓ Absorbent paper or paper towel.

### Precautions for Use

For research use only. Not for use in diagnostic procedures.

- Limitations of the procedure
- ✓ Reliable and reproducible results will be obtained when the assay procedure is carried out with a complete understanding of the package insert instructions and with adherence to good laboratory practice.

- ✓ The wash procedure is critical. Insufficient washing will result in poor precision and falsely elevated absorbance readings.
- ✓ Serum samples demonstrating gross lipemia, gross hemolysis, or turbidity should not be used with this test.
- ✓ For professional use only.

## Assay Protocol

### Reagent Preparation

- ✓ All reagents should be allowed to reach room temperature (18-25°C) before use.
- ✓ Dilute 1 volume of Wash Buffer (20×) with 19 volumes of distilled water. For example, dilute 50 mL of Wash Buffer (20×) into distilled water to prepare 1000 mL of Wash Buffer (1×). Wash Buffer is stable for 1 month at 2-8°C. Mix well before use.

### Sample Preparation

- ✓ Serum should be prepared from a whole blood specimen obtained by acceptable techniques. This kit is for use with serum samples without additives only.
- ✓ Specimens may be refrigerated at 2-8°C for up to 7 days or frozen for up to 6 months. Avoid repetitive freezing and thawing of serum sample.

### Assay Procedure

1. Secure the desired number of coated wells in the holder.
2. Prepare 1:40 dilution for test samples, all six *H. pylori* standards, negative control, and positive control by adding 5 µL of the sample to 200 µL of sample diluent. Mix well.
3. Dispense 100 µL of diluted sera, six standards, and controls into the appropriate wells. For the reagent blank, dispense 100 µL sample diluent in 1A well position. Tap the holder to remove air bubbles from the liquid and mix well for 10 seconds.
4. Incubate at room temperature for 30 minutes.
5. At the end of the incubation period, remove liquid from all wells. Rinse and flick the microtiter wells 4 times with diluted wash buffer (1×) and then one time with distilled water. (Please do not use tap water.)
6. Dispense 100 µL of enzyme conjugate to each well. Mix gently for 10 seconds.
7. Incubate at room temperature for 30 minutes.
8. Remove enzyme conjugate from all wells. Rinse and flick the microtiter wells 4 times with diluted wash buffer (1×) and then one time with distilled water.
9. Add 100 µL of TMB Reagent to each well. Mix gently for 10 seconds.
10. Incubate at room temperature for 20 minutes.
11. Add 100 µL of Stop Solution to each well including the 2 blanks.
12. Mix gently for 30 seconds. It is important to make sure that all the blue color changes to yellow color completely.
13. Read the optical density at 450 nm within 15 minutes with a microtiter plate reader.

*Important Note: The wash procedure is critical. Insufficient washing will result in improper color development.*

## Data Analysis

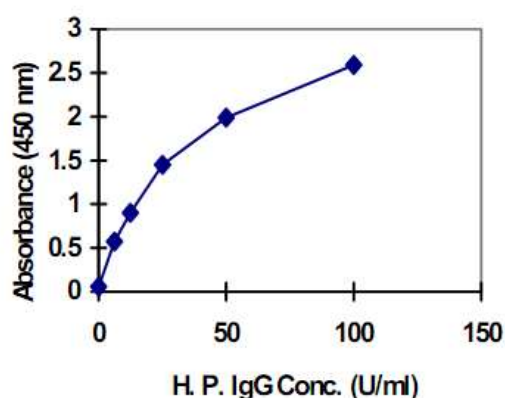
### Calculation of Results

- ✓ Calculate the mean absorbance value (A<sub>450</sub>) for each set of reference standards, controls and samples.
- ✓ Construct a standard curve by plotting the mean absorbance obtained from each reference standard against its concentration in U/mL on graph paper, with absorbance values on the vertical or Y axis, and concentrations on the horizontal or X axis.
- ✓ Use the mean absorbance values for each specimen to determine the corresponding concentration of *H. pylori* IgG in U/mL from the standard curve.

#### ✓ Example of standard curve

Results of a typical standard run with optical density readings at 450 nm shown in the Y axis against *H. pylori* IgG concentrations shown in the X axis. This standard curve is for the purpose of illustration only, and should not be used to calculate unknowns. Each user should obtain his or her own data and standard curve.

<i>H. pylori</i> (U/mL)	Absorbance (450 nm)
0	0.059
6.25	0.573
12.5	0.901
25	1.450
50	1.988
100	2.591



## Performance Characteristics

- Expected values

A cut-off level is set at 20 U/mL for normal samples. Values below 20 U/mL are considered normal. Values above 20 U/mL are regarded as positive. Values above 100 U/mL should be re-assayed at a higher dilution, e.g. 1:802 (first with 1:41, and then 1:20). Results obtained from this 1:802 dilution should be multiplied by 20 to reflect the true *H. pylori* IgG concentration.

The comparison of Abnova ELISA *H. pylori* IgG test to a commercial ELISA kit results are summarized in the following table.

		Reference ELISA			
		N	E	P	Total
Abnova ELISA	N	96(D)	1	4(B)	101
	E	2	2	1	5
	P	3(C)	0	105(A)	108
	Total	101	3	110	214

$$\text{Sensitivity} = A / (A+B) = 107 / 109 = 99\%$$

$$\text{Specificity} = D / (C+D) = 96 / 99 = 97\%$$

$$\text{Accuracy} = (A+D) / (A+B+C+D) = 201 / 208 = 97\%$$

The precision of the assay was evaluated by testing three different sera of 20 replicates over 4 days. The intra-assay and inter-assay C.V. are summarized below.

	7.5 U/mL	22 U/mL	80 U/mL
Intra-assay	9.1%	8.5%	6.4%
Inter-assay	10.5%	8.9%	7.5%



## Resources

### References

- ✓ Marshall, B.J. and J.R. Warren. Unidentified curved bacilli in the stomach of patients with gastritis and Peptic ulceration. *Lancet* 1: 1311-1314, 1984.
- ✓ Ruaws, E.A.J. and G.N.J. Tytgat. Cure of duodenal ulcer associated with eradication of *Helicobacter pylori*, *Lancet* 335: 1233-35, 1990.
- ✓ Perez-Perez, G.I., S.S. Wilkin, M.D. Decker and M.J. Blaswer. Seroprevalence of *Helicobacter pylori* infection in couples. *J. Clin. Microbiol.* 29:642-644, 1991.

**Plate Layout**

12								
11								
10								
9								
8								
7								
6								
5								
4								
3								
2								
1								
	A	B	C	D	E	F	G	H