



# *Helicobacter pylori* IgA ELISA Kit

Catalog Number KA0964

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 .....	7
<b>Resources .....</b>	<b>9</b>
References .....	9
Plate Layout .....	10

## **Introduction**

### **Intended Use**

The *Helicobacter pylori* IgA ELISA Kit is intended for the detection of IgA antibody to *H. pylori* in human serum and plasma.

### **Background**

*H. pylori* is detectable in nearly 100% of adult patients with duodenal ulcer and about 80% of patients with gastric ulcer. An association between *H. pylori* and gastric cancer is confirmed. In developing countries, where most children become infected by the age of 10, gastric cancer rates are very high. In the USA and other developed countries, standards of hygiene and the increasing socioeconomic status of the population have reduced the incidence of infection, and in parallel, the rates of peptic ulcers and gastric cancer have declined. There is excellent correlation between the clinical presentation of gastritis, the presence of *H. pylori* in the stomach and elevated serum *H. pylori* IgG and IgA antibodies. ELISA sensitivity and specificity are 90%, and the predictive value of a negative result for is very high. *H. pylori* IgG and/or IgA antibodies falls significantly after successful antibacterial therapy. Eradication of *H. pylori* is associated with a significant reduction in duodenal ulcer recurrence. *pylori* strains are classified into two broad groups - those that express both VacA and CagA (type I) and those that produce neither (type II). Type I strains are predominate in patients with ulcers and cancer. Up to 50% of adults is infected with *H. pylori*, but most of them are asymptomatic and will not develop ulcer. The reason is they are infected with type II. 80-100% of patients with duodenal ulcer disease produce CagA antibodies against a 128 kd antigen compared with 60-63% of *H. pylori*-infected persons with gastritis only, indicating that serologic responses to the 128 kd protein are more prevalent among *H. pylori*-infected persons with duodenal ulcers than infected persons without peptic ulceration. In *H. pylori*-infected patients who develop gastric cancer, serum IgG against CagA 94% sensitive and 93% specific, indicating that detection of antibodies to CagA is useful marker for diagnosis of duodenal ulcer and gastric cancer.

### **Principle of the Assay**

Diluted patient serum is added to wells coated with purified antigen. IgA specific antibody, if present, binds to the antigen. All unbound materials are washed away and the enzyme conjugate is added to bind to the antibody-antigen complex, if present. Excess enzyme conjugate is washed off and substrate is added. The plate is incubated to allow the hydrolysis of the substrate by the enzyme. The intensity of the color generated is proportional to the amount of IgA specific antibody in the sample.

## General Information

### Materials Supplied

List of component

MATERIALS PROVIDED	96 Tests
Microwell coated with <i>H. pylori</i> antigen	12 x 8 x 1
Sample Diluent: 1 bottle (ready to use)	22 ml
Calibrator: 1 vial (ready to use)	1 ml
Positive Control: 1 vial (ready to use)	1 ml
Negative Control: 1 vial (ready to use)	1 ml
Enzyme conjugate: 1 bottle (ready to use)	12 ml
TMB Substrate: 1 bottle (ready to use)	12 ml
Stop Solution: 1 bottle (ready to use)	12 ml
Wash concentrate 20X: 1 bottle	25 ml

### Storage Instruction

- ✓ Store the kit at 2-8°C.
- ✓ Keep microwells sealed in a dry bag with desiccants.
- ✓ The reagents are stable until expiration of the kit.
- ✓ Do not expose test reagents to heat, sun, or strong light.

### Materials Required but Not Supplied

- ✓ Distilled or deionized water
- ✓ Precision pipettes
- ✓ Disposable pipette tips
- ✓ ELISA reader capable of reading absorbance at 450nm
- ✓ Absorbance paper or paper towel
- ✓ Graph paper

### Precautions for Use

- ✓ Precautions
  - Potential biohazardous materials:
 

The calibrator and controls contain human source components which have been tested and found non-reactive for hepatitis B surface antigen as well as HIV antibody with FDA licensed reagents. However, there is no test method that can offer complete assurance that HIV, Hepatitis B virus or other

infectious agents are absent. These reagents should be handled at the Biosafety Level 2, as recommended in the Centers for Disease Control/National Institutes of Health manual, "Biosafety in Microbiological and Biomedical Laboratories" 1984.

- This kit is designed for research use only.
- Optimal results will be obtained by strict adherence to the test protocol. Precise pipetting as well as following the exact time and temperature requirements is essential.
- Do not pipette by mouth. Do not smoke, eat, or drink in the areas in which specimens or kit reagents are handled.
- The components in this kit are intended for use as an integral unit. The components of different lots should not be mixed.
- Control sera and sample diluent contain preserved with sodium azide. Sodium azide may react with lead and copper plumbing to form explosive metal azide. On disposal, flush with a large volume of water.

✓ Limitations of the Test

Lipemic or hemolyzed samples may cause erroneous results.

## **Assay Protocol**

### **Reagent Preparation**

Prepare 1X Wash buffer by adding the contents of the bottle (25 ml, 20X) to 475 ml of distilled or deionized water. Store at room temperature (18-26°C).

### **Sample Preparation**

- ✓ Collect blood specimens and separate the serum.
- ✓ Specimens may be refrigerated at 2-8°C for up to seven days or frozen for up to six months. Avoid repetitive freezing and thawing.

### **Assay Procedure**

Bring all specimens and kit reagents to room temperature (18-26°C) and gently mix.

1. Place the desired number of coated strips into the holder.
2. Negative control, positive control, and calibrator are ready to use. Prepare 1:21 dilution of test samples, by adding 10 µl of the sample to 200 µl of sample diluent. Mix well.
3. Dispense 100 µl of diluted sera, calibrator 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. Incubate for 20 minutes at room temperature.
4. Remove liquid from all wells. Wash wells three times with 300 µl of 1X wash buffer. Blot on absorbance paper or paper towel.
5. Dispense 100 µl of enzyme conjugate to each well and incubate for 20 minutes at room temperature.
6. Remove enzyme conjugate from all wells. Wash wells three times with 300 µl of 1X wash buffer. Blot on absorbance paper or paper towel.
7. Dispense 100 µl of TMB substrate and incubate for 10 minutes at room temperature. Add 100 µl of stop solution.
8. Read O.D. at 450 nm using ELISA reader within 15 min. A dual wavelength is recommended with reference filter of 600-650 nm.

## Data Analysis

### Calculation of Results

1. Check Calibrator Factor (CF) value on the calibrator bottle. This value might vary from lot to lot. Make sure you check the value on every kit.
2. Calculate the cut-off value: Calibrator OD x Calibrator Factor (CF).
3. Calculate the Ab (Antibody) Index of each determination by dividing the O.D. value of each sample by cut-off value.

✓ Example of typical results:

Calibrator mean OD = 0.8

Calibrator Factor (CF) = 0.5

Cut-off Value =  $0.8 \times 0.5 = 0.400$

Positive control O.D. = 1.2

Ab Index =  $1.2 / 0.4 = 3$

Patient sample O.D. = 1.6

Ab Index =  $1.6 / 0.4 = 4.0$

✓ Quality Control

The test run may be considered valid provided the following criteria are met:

1. The O.D. of the Calibrator should be greater than 0.250.
2. The Ab index for Negative control should be less than 0.9.
3. The Ab Index for Positive control should be greater than 1.2.

✓ Interpretation

The following is intended as a guide to interpretation of *H. pylori* IgA test results; each laboratory is encouraged to establish its own criteria for test interpretation based on sample populations encountered.

✓ Antibody Index Interpretation

<0.9 No detectable antibody to *H. pylori* IgA by ELISA.

0.9-1.1 Borderline positive. Follow-up testing is recommend if clinically indicated.

>1.1 Detectable antibody to *H. pylori* by ELISA.

### Performance Characteristics

✓ Sensitivity and Specificity

94 sera from patients with suspected *H. pylori* infections were tested by this *H. pylori* IgA ELISA and a reference ELISA method. 14 sera were positive and 73 were negative by both methods (95% agreement).

The results are summarized below:

	<i>H. pylori</i> IgA ELISA		
	+	-	Total
Reference ELISA Kit +	14	3	17
-	4	73	77
Total	18	76	94

✓ Precision

• Intra-Assay Study

Serum	No. of Replicates	Mean	Standard Deviation	Coefficient of Variation %
1	16	1.43	0.096	6.71
2	16	0.98	0.067	6.83
3	16	0.26	0.019	7.30

• Inter-Assay Study

Serum	No. of Replicates	Mean	Standard Deviation	Coefficient of Variation %
1	10	1.29	0.121	9.37
2	10	0.91	0.089	9.78
3	10	0.23	0.025	10.86

## Resources

### References

1. Cutler AF; Prasad VM; Santogade P. Four-year trends in Helicobacter pylori IgG serology following successful eradication. *Am J Med* 1998;105(1):18-20
2. Holtmann G; Talley NJ; Mitchell H; Hazell S. Antibody response to specific H. pylori antigens in functional dyspepsia, duodenal ulcer disease, and health. *Am J Gastroenterol* 1998; 93(8):1222-7.
3. Parsonnet J; Replogle M; Yang S; Hiatt R. Seroprevalence of CagA-positive strains among Helicobacter pylori-infected, healthy young adults. *J Infect Dis* 1997;175(5):1240-2.
4. Klaamas K; Held M; Wadström T; Lipping A; Kurtenkov O. IgG immune response to Helicobacter pylori antigens in patients with gastric cancer as defined by ELISA and immunoblotting. *Int J Cancer* 1996; 67(1):1-5.
5. Matsukura N; Onda M; Tokunaga A; Kato S; Yoshiyuki T; Hasegawa H; Yamashita K; Tomtitchong P; Hayashi A. Role of Helicobacter pylori infection in perforation of peptic ulcer: an age- and gender-matched case-control study. *J Clin Gastroenterol* 1997;10:S235-9.

**Plate Layout**

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