



EAGLE
BIOSCIENCES

Fecal *H. Pylori* Antigen Quantitative ELISA Assay Kit

Catalog Number:

HPL35-K01 (1 x 96 wells)

For Research Use Only. Not for use in diagnostic procedures.

v. 3.0 11.2019

EAGLE BIOSCIENCES, INC.

20A Northwest Blvd., Suite 112, Nashua, NH 03063

Phone: 617-419-2019 Fax: 617-419-1110

WWW.EAGLEBIO.COM



INTENDED USE

Eagle Biosciences *Helicobacter pylori* Antigen ELISA Assay Kit is intended for the quantitative and qualitative detection of *Helicobacter pylori* antigen in feces. This kit is for research use only. The *H. pylori* ELISA Assay Kit is for research use only and not intended for diagnostic procedures.

BACKGROUND

H. pylori (previously known as *Campylobacter pyloridis*) is a type of bacteria that infects the stomach and is a common cause of peptic ulcers. *H. pylori* bacteria can be passed from person to person through direct contact with saliva, vomit or fecal matter. *H. pylori* can also be spread through contaminated food or water.

The infection is normally acquired during childhood. *H. pylori* usually goes undiagnosed until symptoms of a peptic ulcer occur. *H. pylori* infection is quite common and is present in about half the people in the world.

PRINCIPLE OF THE ASSAY

This *Helicobacter pylori* Antigen ELISA Assay Kit "sandwich" ELISA is designed, developed and produced for the quantitative and qualitative measurement of *H. pylori* antigen in stool specimen. The assay utilizes the microplate-based enzyme immunoassay technique by coating highly purified antibody onto the wall of microtiter wells.

Assay calibrators and extracted fecal specimen are added to microtiter wells of microplate that was coated with a highly purified monoclonal *H. pylori* antibody on its wall. During the assay, the *H. pylori* antigen will be bound to the antibody coated plate after an incubation period. The unbound material is washed away and another HRP-conjugated monoclonal antibody which specifically recognizes the protein of *H. pylori* is added for further immunoreactions. After an incubation period, the immunocomplex of "H. pylori Antibody – *H. pylori* Antigen – HRP-conjugated Anti-*H. pylori* Tracer Antibody" is formed if *H. pylori* antigen is present in the test sample. The unbound tracer antibody and other proteins in buffer matrix are removed in the subsequent washing step. HRP conjugated tracer antibody bound to the well is then incubated with a substrate solution in a timed reaction and then measured in a spectrophotometric microplate reader. The enzymatic activity of the tracer antibody bound to *H. pylori* proteins captured on the wall of each microtiter well is directly proportional to the amount of *H. pylori* antigen level in each test specimen.

REAGENTS: Preparation and Storage

The *H. pylori* ELISA Assay Kit must be stored at 2 – 8 °C upon receipt. For the expiration date of the kit refer to the label on the kit box. All components are stable until this expiration date.

Prior to use, allow all reagents to equalize to room temperature. Reagents from different kit lot numbers should not be combined or interchanged.

1. *H.pylori* Antibody Coated Microplate

Microplate coated with H.pylori antibody.

Qty: 1 x 96 well microplate

Storage: 2 – 8°C

Preparation: Ready to Use.



2. **Anti-H.pylori Tracer Antibody**
HRP-conjugated monoclonal H.pylori antibody in a stabilized protein matrix.
Qty: 1 x 12 mL
Storage: 2 – 8°C
Preparation: Ready to Use.
3. **ELISA HRP Substrate**
Tetramethylbenzidine (TMB) with stabilized hydrogen peroxide.
Qty: 1 x 12 mL
Storage: 2 – 8°C
Preparation: Ready to Use.
4. **ELISA Stop Solution**
0.5 M sulfuric acid.
Qty: 1 x 12 mL
Storage: 2 – 25°C
Preparation: Ready to Use.
5. **H.pylori Antigen Calibrators Levels 1 to 6**
Calibrator in bovine serum albumin-based matrix with a proclin preservative.
Qty: 6 x Vials
Storage: 2 – 8°C , <-20°C for long term storage
Do not exceed 3 freeze-thaw cycles.
Preparation: Ready to Use.
6. **H.pylori Antigen Controls**
Calibrator in bovine serum albumin-based matrix with a proclin preservative.
Qty: 2 x Vials
Storage: 2 – 8°C , <-20°C for long term storage
Do not exceed 3 freeze-thaw cycles.
Preparation: Ready to Use.
7. **ELISA Wash Concentrate**
Surfactant in a phosphate buffered saline with non-azide preservative.
Qty: 1 x 30 mL
Storage: 2 – 25°C
Preparation: 30X Concentrate. The contents must be diluted with 870 mL distilled water and mixed well before use.
8. **H.pylori Concentrated Assay Buffer**
Qty: 1 x 30 mL
Storage: 2 – 8°C
Preparation: 4X Concentrate. The contents must be diluted with 90 mL distilled water and mixed well before use.

SAFETY PRECAUTIONS

The reagents of the *H. pylori* ELISA Assay Kit must be used in a laboratory and are for professional use only. Materials sourced for reagents containing bovine serum albumin were derived in the contiguous 48 United States and obtained only from healthy donor animals



maintained under veterinary supervision and found free of contagious diseases. Wear gloves while performing this assay and handle these reagents as if they are potentially infectious. Avoid contact with reagents containing TMB, hydrogen peroxide, or sulfuric acid. TMB may cause irritation to skin and mucous membranes and cause an allergic skin reaction. TMB is a suspected carcinogen. Sulfuric acid may cause severe irritation on contact with skin. Do not get in eyes, on skin, or on clothing. Do not ingest or inhale fumes. On contact, flush with copious amounts of water for at least 15 minutes. Use Good Laboratory Practices.

MATERIALS REQUIRED BUT NOT PROVIDED

- Precision single channel pipettes capable of delivering 10 μ L, 50 μ L, 100 μ L, and 1000 μ L.
- Repeating dispenser suitable for delivering 100 μ L.
- Disposable pipette tips suitable for above volume dispensing.
- Disposable 12 x 75 mm or 13 x 100 glass or plastic tubes.
- Disposable plastic 1000 mL bottle with cap.
- Aluminum foil.
- Deionized or distilled water.
- Plastic microtiter well cover or polyethylene film.
- ELISA multichannel wash bottle or automatic (semi-automatic) washing system.
- Spectrophotometric microplate reader capable of reading absorbance at 450 nm.

SPECIMEN COLLECTION AND STORAGE

Fresh fecal sample should be collected into a stool sample collection container. It is required to collect a minimum of 1-2 mL liquid stool sample or 1-2 g solid sample. The collected fecal sample must be transported to the lab in a frozen condition (-20°C). If the stool sample is collected and tested the same day, it is allowed to be stored at 2-8°C.

ASSAY PROCEDURE

Reagent Preparation

1. Prior to use allow all reagents to come to room temperature (20-25 °C). Reagents from different kit lot numbers should not be combined or interchanged.
2. ELISA Wash Concentrate must be diluted to working solution prior use. Please see REAGENTS section for details.

Quantitative Assay Procedure

1. Place a sufficient number of microwell strips in a holder to run calibrators controls, and samples in duplicate.



2. Test Configuration

Row	Strip 1	Strip 2	Strip 3
A	Calibrator Level 1	Calibrator Level 5	SAMPLE 1
B	Calibrator Level 1	Calibrator Level 5	SAMPLE 1
C	Calibrator Level 2	Calibrator Level 6	SAMPLE 2
D	Calibrator Level 2	Calibrator Level 6	SAMPLE 2
E	Calibrator Level 3	Control 1	SAMPLE 3
F	Calibrator Level 3	Control 1	SAMPLE 3
G	Calibrator Level 4	Control 2	SAMPLE 4
H	Calibrator Level 4	Control 2	SAMPLE 4

3. Add **100 µL** of calibrators, controls and samples into the designated microwells.

Note: if the collection tubes from CAL35-C50 is used, add two drops of extracted fecal sample into each well.

4. Cover the plate with one plate sealer and aluminum foil. Incubate at **room temperature (20-25 °C)** for **60 minutes**.
5. Remove the plate sealer. Aspirate the contents of each well. Wash each well **5 times** by dispensing **350 µL** of diluted wash solution into each well, and then completely aspirate the contents. Alternatively, an automated microplate washer can be used.
6. Add **100 µL** of Anti-H.pylori Tracer Antibody to each well. Mix by gently tapping the plate.
7. Cover the plate with one plate sealer and aluminum foil. Incubate at **room temperature (20-25 °C)** for **30 minutes**.
8. Remove the plate sealer. Aspirate the contents of each well. Wash each well **5 times** by dispensing **350 µL** of diluted wash solution into each well, and then completely aspirate the contents. Alternatively, an automated microplate washer can be used.
9. Add **100 µL** of ELISA HRP Substrate into each of the wells. Mix by gently tapping the plate.
10. Cover the plate with one plate sealer and aluminum foil. Incubate at **room temperature (20-25 °C)** for **20 minutes**.
11. Remove the aluminum foil and plate sealer. Add **100 µL** of ELISA Stop Solution into each of the wells. Mix by gently tapping the plate.
12. Read the absorbance at **450/620 nm** within **10 minutes** with a microplate reader.

1. Qualitative Assay Procedure

1. Place a sufficient number of microwell strips in a holder to run positive control [H.pylori Antigen Calibrator Level 6] negative control [diluted 1X Assay Buffer], and samples in duplicate.
2. Test Configuration



Row	Strip 1	Strip 2	Strip 3
A	Negative Control	SAMPLE 3	SAMPLE 7
B	Negative Control	SAMPLE 3	SAMPLE 7
C	Positive Control	SAMPLE 4	SAMPLE 8
D	Positive Control	SAMPLE 4	SAMPLE 8
E	SAMPLE 1	SAMPLE 5	SAMPLE 9
F	SAMPLE 1	SAMPLE 5	SAMPLE 9
G	SAMPLE 2	SAMPLE 6	SAMPLE 10
H	SAMPLE 2	SAMPLE 6	SAMPLE 10

1. Add **100 µL** of positive control [H.pylori Antigen Calibrator Level 6], negative control [diluted 1X Assay Buffer], and samples into the designated microwells.

Note: if the collection tubes from CAL35-C50 is used, add two drops of extracted fecal sample into each well.

2. Cover the plate with one plate sealer and aluminum foil. Incubate at **room temperature (20-25 °C)** for **60 minutes**.
3. Remove the plate sealer. Aspirate the contents of each well. Wash each well **5 times** by dispensing **350 µL** of diluted wash solution into each well, and then completely aspirate the contents. Alternatively, an automated microplate washer can be used.
4. Add **100 µL** of Anti-H.pylori Tracer Antibody to each well. Mix by gently tapping the plate.
5. Cover the plate with one plate sealer and aluminum foil. Incubate at **room temperature (20-25 °C)** for **30 minutes**.
6. Remove the plate sealer. Aspirate the contents of each well. Wash each well **5 times** by dispensing **350 µL** of diluted wash solution into each well, and then completely aspirate the contents. Alternatively, an automated microplate washer can be used.
7. Add **100 µL** of ELISA HRP Substrate into each of the wells. Mix by gently tapping the plate.
8. Cover the plate with one plate sealer and aluminum foil. Incubate at **room temperature (20-25 °C)** for **20 minutes**.
9. Remove the aluminum foil and plate sealer. Add **100 µL** of ELISA Stop Solution into each of the wells. Mix by gently tapping the plate.
10. Read the absorbance at **450nm** within **10 minutes** with a microplate reader.

PROCEDURAL NOTES

1. It is recommended that all calibrators and unknown samples be assayed in duplicate in the Eagle Biosciences *H. pylori* ELISA Assay Kit. The average absorbance reading of each duplicate should be used for data reduction and the calculation of results.
2. Keep light-sensitive reagents in the original amber bottles.
3. Store any unused antibody-coated strips in the foil zipper bag with desiccant to protect from moisture.
4. Careful technique and use of properly calibrated pipetting devices are necessary to ensure reproducibility of the *H. pylori* ELISA Assay test.
5. Incubation times or temperatures other than those stated in this insert may affect the results.
6. All reagents of the *H. pylori* ELISA Assay Kit should be mixed gently and thoroughly prior to use. Avoid foaming.



INTERPRETION OF RESULTS

1. Quantitative Measurement

1. Calculate the average absorbance for each pair of duplicate test results.
2. Subtract the average absorbance of the calibrator 1 (0 ng/mL) from the average absorbance of all other readings to obtain corrected absorbance.
3. The calibration curve is generated by the corrected absorbance of all calibrator levels on the ordinate against the calibrator concentration on the abscissa using point-to-point or log-log paper. Appropriate computer assisted data reduction programs may also be used for the calculation of results.
4. The *H. Pylori* concentrations for the unknown samples are read directly from the calibration curve using their respective corrected absorbance.

2. Qualitative Measurement

1. Visual

1. Positive or reactive: Any sample well that is obviously more yellow than the negative control well.
2. Negative or non-reactive: Any sample well that is not obviously more yellow than the negative control well.

Note: The negative control, as well as some patient samples, may show some slight yellow color. A sample well must be obviously darker or more yellow than the negative control well, when it is interpreted as a positive result.

2. ELISA Reader

1. Calculate the average absorbance for each pair of duplicate test results.
2. Calculate the cut-offs:
 - Positive cut-off: $1.1 \times (\text{mean extinction of the negative control} + 0.10)$.
 - Negative cut-off: $0.9 \times (\text{mean extinction of the negative control} + 0.10)$.
3. Interpret test results
 - Positive: patient sample extinction is greater than the positive cut-off.
 - Negative: patient sample extinction is less than the negative cut-off.
 - Equivocal: sample extinction is between the positive and negative cut-off.
4. Assay Quality Control
 - Positive control must show an average OD reading greater than 0.8.
 - Negative control must show an average OD reading less than 0.18.

LIMITATIONS OF THE PROCEDURE

1. The results obtained with this *H. pylori* Antigen Test Kit serve only as an aid to diagnosis and should not be interpreted as diagnostic in themselves without taking other clinical findings such as stomach endoscope and biopsy, etc.
2. Single *H. pylori* negative results in untreated patients do not rule out *H. pylori* infection.
3. For unknown sample value read directly from the assay that is greater than the highest calibrator, it is recommended to measure a further diluted sample for more accurate measurement.
4. Bacterial or fungal contamination of serum specimens or reagents, or cross-contamination between reagents may cause erroneous results.

QUALITY CONTROL

To assure the validity of the results each assay should include adequate controls with known *H. pylori* antigen levels. We recommend that all assays include the laboratory's own controls.



EXPECTED VALUES

1. Quantitative Measurement

Stool from 25 normal adults were measured with this ELISA. We found that normal people show undetectable H. pylori antigen in the extracted stool sample according to the sample collection, extraction and assay procedures described in this insert. The suggested positive cut-off for fecal H. pylori antigen is 3 ng/mL.

2. Qualitative Measurement

3. Stool samples from 29 negative specimens and 17 positive specimens were tested with this ELISA.

Samples EagleBio's ELISA	True Positive	True Negative	Total
Positive	17	0	17
Negative	0	29	29
Total	17	29	46

Sensitivity: 100% (17/17)

Specificity: 100% (29/29)

Accuracy: 100% (46/46)

EXAMPLE DATA

1. Quantitative Measurement

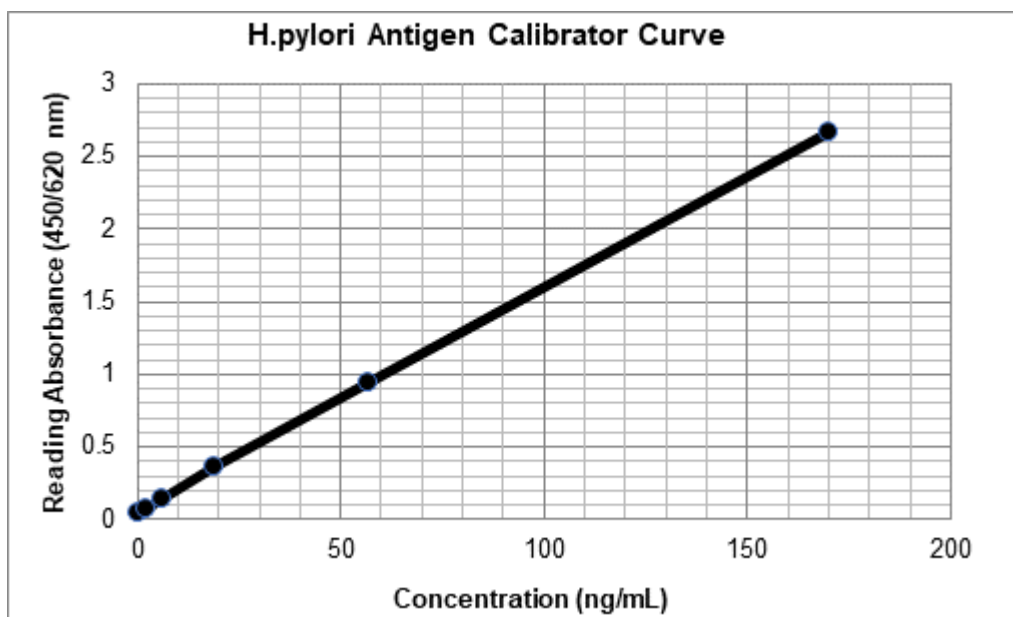
A typical absorbance data and the resulting calibration curve from Fecal H. Pylori antigen ELISA are represented.

Note: This curve should not be used in lieu of calibration curve run with each assay.

Well ID	Reading Absorbance (450/620 nm)			Concentration (ng/mL)
	Readings	Average	Corrected	
Calibrator Level 1: 0 ng/mL	0.047	0.047	0.000	
	0.047			
Calibrator Level 2: 2.1 ng/mL	0.075	0.073	0.026	
	0.071			
Calibrator Level 3: 6.3 ng/mL	0.140	0.142	0.095	
	0.144			
Calibrator Level 4:	0.357	0.363	0.316	
	0.368			



18.9 ng/mL				
Calibrator Level 5: 56.7 ng/mL	0.909	0.939	0.892	
	0.968			
Calibrator Level 6: 170 ng/mL	2.682	2.665	2.618	
	2.647			
Control 1	0.076	0.076	0.029	
	0.074			
Control 2	0.737	0.724	0.677	
	0.711			



2. Qualitative Measurement

	Reading Absorbance (450 nm)	Average
Negative Control	0.049 0.050	0.050
Positive Control	1.332 1.376	1.354

Positive Cut-Off = $1.1 \times (0.050 + 0.10) = 0.165$

Negative Cut-Off = $0.9 \times (0.050 + 0.10) = 0.135$



PERFORMANCE CHARACTERISTICS

Sensitivity

The sensitivity of the Eagle Biosciences Fecal H. pylori Ag ELISA as determined by the 95% confidence limit on 16 duplicate determination of zero calibrator is approximately 0.165 ng/mL.

Specificity

The assay does not cross react to the following organisms: Cryptosporidium parvum, Giardia lamblia, rotavirus and adenovirus.

Reproducibility and Precision

The intra-assay precision is validated by measuring two samples in a single assay with 12 replicate determinations. The inter-assay precision is validated by measuring two samples in duplicate in 12 individual assays. The results are as follows:

Sample	Intra-Assay		Inter-Assay	
	1	2	1	2
Mean (ng/mL)	13.1	1.8	13.9	1.8
CV (%)	5.4	2.8	5.9	5.2

Linearity

Two (2) stool samples were diluted with assay buffer and tested. The results are as follows:

Samples	Observed (ng/mL)	Recovery (%)
Sample A	77.4	-
50%	38.1	98.4
25%	17.5	90.4
Sample B	24.8	-
50%	12.2	98.6
25%	6.3	102.7

Spike Recovery

Two amples were spiked with calibrators each other in equal volume and assayed. The results indicate below:



Samples	Observed (ng/mL)	% Recovery (%)
Sample A	0.3	-
+ Level 2: 1.9 ng/mL	1.1	97.2
+ Level 4: 16.7 ng/mL	7.1	84.2
+ Level 5: 50 ng/mL	20.9	83.2
Sample B	0.2	-
+ Level 2: 1.9 ng/mL	1.0	93.8
+ Level 4: 16.7 ng/mL	7.0	82.0
+ Level 5: 50 ng/mL	21.0	83.1

REFERENCES

1. Janoff EN, Smith PD, Blaser MJ. Acute antibody responses to H. Pylori are depressed in patients with AIDS. J Infect Dis. 1988 Apr;157(4):798-804..
2. Rosenblatt JE, Sloan LM, Schneider SK. Evaluation of an enzyme-linked immunosorbent assay for the detection of H. Pylori in stool specimens. Diagn Microbiol Infect Dis. 1993 May-Jun;16(4):337-41.
3. Stibbs HH, Samadpour M, Manning JF. Enzyme immunoassay for detection of H. Pylori cyst antigens in formalin-fixed and unfixed human stool. J Clin Microbiol. 1988 Sep;26(9):1665-9.
4. Stibbs HH. Monoclonal antibody-based enzyme immunoassay for H. Pylori antigen in human stool. J Clin Microbiol. 1989 Nov;27(11):2582-8.
5. Ungar BL, Yolken RH, Nash TE, Quinn TC. Enzyme-linked immunosorbent assay for the detection of H. Pylori in fecal specimens. J Infect Dis. 1984 Jan;149(1):90-7

Warranty Information

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 617-419-2019.