



EAGLE
BIOSCIENCES

Pepsinogen I ELISA Assay Kit

Catalog Number:

PP131-K01 (1 x 96 wells)

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

v. 12 (22 APR 24)

EAGLE BIOSCIENCES, INC.

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

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

WWW.EAGLEBIO.COM



INTENDED USE

The Eagle Biosciences Human Pepsinogen I ELISA Assay Kit (enzyme-linked immunoassay kit is intended for the quantitative determination of human pepsinogen I levels in serum. The Eagle Biosciences Human Pepsinogen I ELISA Assay Kit is for research use only and not to be used in diagnostic procedures.

INTRODUCTION

Pepsinogen consists of a single polypeptide chain of 375 amino acids with an average molecular weight of 42 kDa. Pepsinogen I is synthesized at gastric chief cells and mucous neck cells, while pepsinogen II is produced not only by gastric chief cells, mucous neck cells, but also by clear mucous cells of antrum, etc. The clinical applications of measuring pepsinogen I and pepsinogen II are of useful aid in diagnosing severe atrophic gastritis and stomach cancer. It was suggested that the measurement of serum pepsinogens served as a "serological biopsy" for predicting the presence of atrophic gastritis or superficial gastritis.

Atrophic Gastritis: It was found that a serum pepsinogen I level falling to less than 20 ng/ml was highly specific for severe atrophic gastritis. It is also observed that serum pepsinogen I levels fell with increasing severity of mucosal damage in atrophic gastritis. The diagnostic sensitivity and specificity of serum pepsinogen I level for advanced atrophic corpus gastritis are about 92% and 90% respectively. On the other hand, the decrease in serum pepsinogen I levels in patients with pernicious anemia and atrophic gastritis was found to be associated with normal or raised pepsinogen II levels. Therefore, a pepsinogen I/pepsinogen II ratio is significantly lower than those with superficial gastritis or normal remnant mucosa.

Stomach Cancer: Low serum pepsinogen I levels were found in patients with gastric cancer, with a threefold higher incidence. Other studies have concluded that low serum pepsinogen I levels may identify persons at increased risk for intestinal types of stomach cancer.

Duodenal Ulcer: A low serum pepsinogen I level can exclude a diagnosis of duodenal ulcer. Although a high pepsinogen I level has less clinical useful for establishing the diagnosis of a duodenal ulcer, the combination of hypergastrinemia and a highly elevated serum pepsinogen I strongly suggests the possibility of the Zollinger-Ellison syndrome.

PRINCIPLE OF THE ASSAY

The Eagle Biosciences Human Pepsinogen I ELISA Assay Kit is designed, developed and produced for the quantitative measurement of human pepsinogen I level in serum sample. The assay utilizes the two-site "sandwich" technique with two selected monoclonal antibodies that bind to different epitopes of human pepsinogen I without any cross-reaction to human pepsinogen II.

Assay standards, controls and patient serum samples containing human pepsinogen I is added directly to microtiter wells of microplate that was coated with a streptavidin. Simultaneously, a biotinylated antibody and a horseradish peroxidase conjugated antibody are added to each well. After the first incubation period, the wall of microtiter well captures the biotinylated antibody as well as an immunocomplex in the form of "streptavidin – biotin-antibody – pepsinogen I– HRP-antibody". Unbound proteins as well as unbound HRP conjugated antibody in each microtiter well are removed in the subsequent washing step. The well is 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 the pepsinogen I on the wall of the microtiter well is directly proportional to the amount of pepsinogen I in the sample. A calibration curve is generated by plotting the absorbance versus the respective human pepsinogen I concentration for each standard on Point-to-Point, CubicSpline or 4-Parameter plot. The concentration of human pepsinogen I in test samples is determined directly from this standard curve.

REAGENTS: Preparation and Storage

This test 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 the expiration date.

Allow all reagents to come to room temperature prior to use. Reagents from different kit lot numbers should not be combined or interchanged.

1. Streptavidin Coated Microplate Microplate coated with streptavidin.

Qty: 1 x 96 well microplate

Storage: 2 – 8°C

Preparation: Ready to Use

2. Pepsinogen I Tracer Antibody

HRP-conjugated anti-human tracer antibody in a stabilized protein matrix.

Qty: 1 x 0.6 mL

Storage: 2 – 8°C

Preparation: 21X Concentrate. The contents must be diluted with tracer antibody diluent and mixed well before use.

3. Tracer Antibody Diluent

Buffer for antibody dilution according to the assay procedures.

Qty: 1 x 12 mL

Storage: 2 – 8°C

Preparation: Ready to use

4. Pepsinogen I Capture Antibody

Biotinylated anti-human pepsinogen I capture antibody in a stabilized protein matrix.

Qty: 1 x 0.6 mL

Storage: 2 – 8°C

Preparation: 21X Concentrate. The contents must be diluted with tracer antibody diluent and mixed well before use.

5. 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.

6. ELISA HRP Substrate

Tetramethylbenzidine (TMB) with stabilized hydrogen peroxide.

Qty: 1 x 12 mL



Storage: 2 – 8°C

Preparation: Ready to Use

7. ELISA Stop Solution

0.5 M sulfuric acid

Qty: 1 x 12 mL

Storage: 2 – 25°C

Preparation: Ready to Use

8. Pepsinogen I Calibrators Levels 1 to 6

Lyophilized human pepsinogen I in a bovine serum albumin-based matrix with a non-azide preservative. Refer to vials for exact concentration.

Qty: 6 x vials

Storage: 2 – 8°C , <-20°C for long term storage Do not exceed 3 freeze-thaw cycles.

Preparation: Must be reconstituted with 0.5 mL of demineralized water, allowed to sit for 10 minutes, and then mixed by inversions or gentle vortexing. Make sure that all solids are dissolved completely prior to use.

9. Pepsinogen I Controls

Lyophilized human pepsinogen I in a bovine serum albumin-based matrix with a non-azide preservative. Refer to vials for exact concentration.

Qty: 2 x vials

Storage: 2 – 8°C , <-20°C for long term storage Do not exceed 3 freeze-thaw cycles.

Preparation: Must be reconstituted with 0.5 mL of demineralized water, allowed to sit for 10 minutes, and then mixed by inversions or gentle vortexing. Make sure that all solids are dissolved completely prior to use.

SAFETY PRECAUTIONS

The Human Pepsinogen I ELISA Assay Kit reagents are for Research Use Only. Source material which contains reagents of bovine serum albumin was derived in the contiguous 48 United States. It was 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 were 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. 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 20 µL, 25 µL, 100 µL, and 1000 µL etc.
- Repeating dispenser suitable for delivering 100 µL.
- 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 caps.
- 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 450nm.

SPECIMEN COLLECTION & STORAGE

Only 50 µL of human serum is required for human pepsinogen I measurement in duplicate. No special preparation of individual is necessary prior to specimen collection. However, a 10 hour fasting serum sample is recommended for the test. Samples should not be taken from patients taking biotin-containing multivitamins or dietary supplements at least 48 hours prior to specimen collection. Whole blood should be collected and must be allowed to clot for minimum 30 minutes at room temperature before the serum is separated by centrifugation (850 – 1500xg for 10 minutes). The serum should be separated from the clot within three hours of blood collection and transferred to a clean test tube. Serum samples should be stored at – 20°C or below until measurement. Avoid more than three freeze-thaw cycles of specimen.

ASSAY PROCEDURE

Reagent Preparation

1. Prior to use allow all reagents to come to room temperature. Reagents from different kit lot numbers should not be combined or interchanged.
2. ELISA Wash Concentrate must be diluted to working solution prior to use. Please see REAGENTS section for details.
3. Reconstitute all assay calibrators level 1 to level 6 and controls by adding 0.5 mL of demineralized water to each vial. Allow the calibrators and controls to sit undisturbed for 10 minutes, and then mix well by inversions or gentle vortexing. One must make sure that all solid is dissolved completely prior to use. These reconstituted calibrators and controls may be stored at 2 – 8°C for up to 3 days or at –20°C or below for long-term storage. Do not exceed 3 freeze-thaw cycles.

Manual Assay Procedure

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

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



2. Prepare the antibody working solution by 1:21 fold dilution of the tracer antibody and capture antibody with the diluent. For each strip, it is required to mix **1 mL** of the diluent with **50 µL** of the tracer antibody and **50 µL** capture antibody in a clean test tube. *Note: This antibody working solution should be freshly prepared.*
3. Add **25 µL** of calibrators, controls, and samples into the designated microwells. Mix by gently tapping the plate.
4. Add **100 µL** of antibody working solution to each microwell.
5. Cover the plate with one plate sealer and aluminum foil. Incubate at **room temperature (20-25 °C)** for **60 minutes**.
6. Remove the plate sealer. Aspirate the contents of each microwell. Wash each microwell **5 times** by dispensing **350 µL** of diluted wash solution into each microwell, and then completely aspirate the contents. Alternatively, an automated microplate washer can be used.
7. Add **100 µL** of substrate into each microwell. 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 and add **100 µL** of Stop Solution into each of the microwells. Mix by gently tapping the plate.
10. Read the absorbance at **450 nm** within **10 minutes** with a microplate reader.

Automated Assay Procedure

1. Prepare the antibody working solution by 1:21 fold dilution of the tracer antibody and capture antibody with the diluent. For each strip, it is required to mix **1 mL** of the diluent with **50 µL** of the tracer antibody and **50 µL** capture antibody) in a clean test tube. *Note: This antibody working solution should be freshly prepared.*
2. Add **25 µL** of calibrators, controls, and samples into the designated microwells.
3. Add **100 µL** of antibody working solution to each microwell.
4. Incubate plate with initial shaking for 1 minutes and further incubation at **37°C** for **45 minutes**.
5. Aspirate the contents of each microwell. Wash each microwell **5 times** by dispensing **350 µL** of diluted wash solution into each microwell, and then completely aspirate the contents.
6. Add **100 µL** of substrate into each microwell.
7. Incubate plate at **37°C** for **15 minutes**.
8. Add **100 µL** of Stop Solution into each of the microwells.
9. Read the absorbance at **450 nm**.

Note: The above automated ELISA procedure has been performed on DS2 system. A satisfactory patient sample correlation was observed between the manual and automated assay procedures ($r = 0.943$, slope = 1.0958). One may adjust the procedure according to different automated ELISA system used in each laboratory.

PROCEDURAL NOTES

1. It is recommended that all standards, controls and unknown samples be assayed in duplicate. 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 bottles and avoid unnecessary exposure to the light.



3. Store any unused antibody coated strips in the foil Ziploc bag with desiccant to protect from moisture.
4. Careful technique and use of properly calibrated pipetting devices are necessary to ensure reproducibility of the test.
5. Incubation times or temperatures other than those stated in this insert may affect the results.
6. Avoid air bubbles in the microwell as this could result in lower binding efficiency and higher CV% of duplicate reading
7. All reagents should be mix gently and thoroughly prior use. Avoid foaming.

INTERPRETATION OF RESULTS

1. Calculate the average absorbance for each pair of duplicate test results.
2. Subtract the average absorbance of the calibrator level 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. It is recommended to use following curve fits: (1) Point-to-Point, or (2) 4-Parameter or (3) CubicSpline.
5. The human pepsinogen I concentrations for the controls and patient samples are read directly from the calibration curve using their respective corrected absorbance.

LIMITATIONS OF THE PROCEDURE

1. Since there is no Gold Standard concentration available for human pepsinogen I measurement, the values of assay calibrators were established by diluting a highly purified human pepsinogen I in a protein matrix.
2. For unknown sample value read directly from the assay that is greater than 300 ng/mL, it is recommended to measure a further diluted sample for more accurate measurement.
3. If there is not a microplate reader in your laboratory able to read beyond 2.0 at OD 450 nm, adjust the computer program for an assay without the calibrator level 6 from the calibrator set.
4. Bacterial or fungal contamination of serum specimens or reagents, or cross-contamination between reagents may cause erroneous results.
5. Water deionized with polyester resins may inactive the horseradish peroxidase enzyme.

QUALITY CONTROL

To assure the validity of the results each assay should include adequate controls with known pepsinogen I levels. We recommend that all assays include the laboratory's own human serum based pepsinogen I controls in addition to those provided with this kit.

EXPECTED VALUES

Seventy-three normal adult sera were measured with this human pepsinogen I ELISA. The expected normal range is listed in the following table with different percentile cut-off and the median level of this group of population is 62.8 ng/mL.



Percentile Cut-off	Normal Range (ng/mL)
95%	25 – 200
90%	30 – 150
85%	40 – 120
80%	40 – 100

It is highly recommended that each laboratory should establish their own normal range for pepsinogen I based on local populations.

Patients with atrophic gastritis, as well as patients with stomach cancer would have a pepsinogen I level below 20 ng/mL. However, gastroendoscope and tissue biopsy should be used as final and confirmative diagnostic method.

EXAMPLE DATA

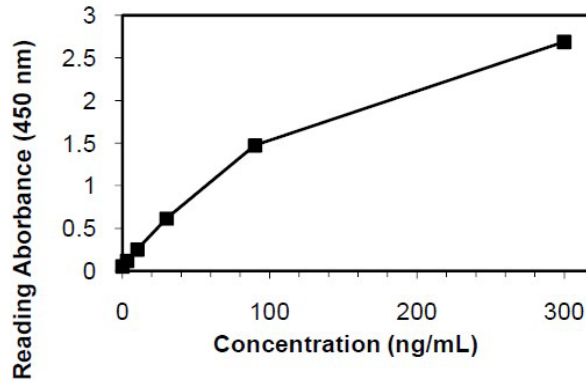
A typical absorbance data and the resulting calibration curves from are represented.

Microwell ID	Reading Absorbance (450 nm)			Concentration (ng/mL)
	Readings	Average	Corrected	
Calibrator Level 1: 0 ng/mL	0.053	0.052	0.000	
	0.050			
Calibrator Level 2: 3 ng/mL	0.119	0.119	0.067	
	0.118			
Calibrator Level 3: 10 ng/mL	0.262	0.254	0.202	
	0.246			
Calibrator Level 4: 30 ng/mL	0.616	0.619	0.567	
	0.622			
Calibrator Level 5: 90 ng/mL	1.565	1.476	1.424	
	1.387			
Calibrator Level 6: 300 ng/mL	2.766	2.685	2.633	
	2.604			
Control 1	0.373	0.368	0.316	16.2
	0.363			
Control 2	1.692	1.640	1.588	118
	1.587			

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



Example Data



PERFORMANCE CHARACTERISTICS

Sensitivity

The sensitivity of this human pepsinogen I ELISA is 0.1 ng/mL as determined by measuring zero calibrator 16 times in the same assay and calculating the detection limit at 3 standard deviations above the pepsinogen I zero calibrator. The assay analytical sensitivity is approximately 0.5 ng/mL.

Hook Effect

It was determined that this pepsinogen I ELISA did not show any high dose "hook" effect up to 10,000 ng/mL of pepsinogen I

Reproducibility and Precision

The intra-assay precision is validated by measuring two samples in a single assay with 20 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)	18.2	121.1	17.5	123.7
CV (%)	5.3	4.8	6.9	5.7

Linearity

Two human serum samples were diluted with assay buffer and assayed. The results are as follows:



Sample	Observed (ng/mL)	Expected (ng/mL)	Recovery (%)
Sample 1	31.90	-	-
50%	16.21	15.95	102
25%	7.95	7.78	102
12.5%	3.73	3.99	93
6.25%	2.11	1.99	106
Sample 2	252.00	-	-
50%	125.27	126.00	99
25%	64.12	63.00	102
12.5%	31.25	31.50	99
6.25%	16.92	15.75	107

Spike Recovery

Two patient samples were spiked with various amounts of human pepsinogen I and assayed. The results are as follows:

Samples	Observed (ng/mL)	Expected (ng/mL)	Recovery (%)
Sample 1	18.6	-	-
10	12.6	14.3	88
30	25.1	24.3	103
90	56.2	54.3	103
Sample 2	121.1	-	-
10	61.3	65.6	93
30	70.9	75.6	94
90	104.7	105.6	99

Specificity

This assay measures human pepsinogen I without any cross-reaction to human pepsinogen II.

REFERENCES

1. Kuipers EJ. In through the out door: serology for atrophic gastritis. *Eur J Gastroenterol Hepatol.* 2003 Aug;15(8):877-9.
2. Miki K, Morita M, Sasajima M, Hoshina R, Kanda E, Urita Y. Usefulness of gastric cancer screening using the serum pepsinogen test method. *Am J Gastroenterol.* 2003 Apr;98(4):735-9.
3. Miki K. [Serum pepsinogen I/II ratio test] *Nippon Rinsho.* 2003 Jan;61(1):92-5. Japanese.
4. So JB, Yeoh KG, Moochala S, Chachlani N, Ho J, Wong WK, Mack P, Goh PM. Serum pepsinogen levels in gastric cancer patients and their relationship with *Helicobacter pylori* infection: a prospective study. *Gastric Cancer.* 2002;5(4):228-32.
- 5.



- 6.
7. Korstanje A, den Hartog G, Biemond I, Lamers CB. The serological gastric biopsy: a nonendoscopic diagnostic approach in management of the dyspeptic patient: significance for primary care based on a survey of the literature. *Scand J Gastroenterol Suppl.* 2002;(236):22-6. Review.
8. Sipponen P, Harkonen M, Alanko A, Suovaniemi O. Diagnosis of atrophic gastritis from a serum sample. *Clin Lab.* 2002;48(9-10):505-15. Review.
9. Tabata H, Fuchigami T, Kobayashi H, Sakai Y, Nakanishi M, Tomioka K, Nakamura S, Matsumoto T, Fujishima M. Difference in degree of mucosal atrophy between elevated and depressed types of gastric epithelial tumors. *Scand J Gastroenterol.* 2001 Nov;36(11):1134-40.
10. Varis K, Sipponen P, Laxen F, Samloff IM, Huttunen JK, Taylor PR, Heinonen OP, Albanes D, Sande N, Virtamo J, Harkonen M. Implications of serum pepsinogen I in early endoscopic diagnosis of gastric cancer and dysplasia. Helsinki Gastritis Study Group. *Scand J Gastroenterol.* 2000 Sep;35(9):950-6.
11. Fernandez R, Vizoso F, Rodriguez JC, Merino AM, Gonzalez LO, Quintela I, Andicoechea A, Truan N, Diez MC. Expression and prognostic significance of pepsinogen C in gastric carcinoma. *Ann Surg Oncol.* 2000 Aug;7(7):508-14.
12. Kalinovskii VP, Gamaiunova VB, Shumakov AP, Khanson KP. [Radioimmunoassay of serum pepsinogen I in chronic gastritis and stomach cancer] *Vopr Onkol.* 2000;46(2):153-5. Russian.
13. Shumakov AR, Fedorov SN, Kalinovskii VP, Khanson KP. [Evaluation of pepsinogen A expression in stomach cancer] *Vopr Onkol.* 1999;45(3):238-40. Russian.
14. Kitahara F, Kobayashi K, Sato T, Kojima Y, Araki T, Fujino MA. Accuracy of screening for gastric cancer using serum pepsinogen concentrations. *Gut.* 1999 May;44(5):693-7.
15. Samloff IM and Taggart RT. Pepsinogens, pepsins, and peptic ulcer. *Clinical and Investigative Medicine* 1987;10:215-221.
16. Samloff IM. Slow moving protease and the seven pepsinogens. Electrophoretic demonstration of the existence of eight proteolytic fractions in human gastric mucosa. *Gastroenterology* 1969;57:659-669.

SHORT ASSAY PROCEDURE

Manual Assay Procedure	Automated Assay Procedure
<ol style="list-style-type: none">1. Add 25 μL of calibrators, controls, and samples into the designated microwells.2. Add 100 μL of antibody working solution into the designated microwells.3. Mix, cover, and incubate at room temperature (20-25C) for 60 minutes.4. Wash each microwell five times.5. Add 100 μL of substrate to each microwell.6. Cover and incubate at room temperature (20-25 DC) for 20 minutes.7. Add 100 μL of the stop solution to each microwell.8. Read the absorbance at 450 nm.	<ol style="list-style-type: none">1. Add 25 μL of c;11Qt,>rators, controls, and samples into the designated microwell s:2. Add 100 μL of antibody working solution to each microwell.3. Incubate plate with initial shaking for 1 minutes and further incubation at 37°C for 45 minutes.4. Wash each microwell five times.5. Add 100 μL of substrate into each microwell.6. Incubate plate at 37° C for 15 minutes.7. Add 100 μL of Stop Solution into each of the microwells.8. Read the absorbance at 450 nm



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 866-411-8023.