



IMMUNOASSAYS AND SERVICES
BIOGENIC AMINES & NEUROSCIENCE | ENDOCRINOLOGY | FOOD SAFETY

LABOR DIAGNOSTIKA NORD GmbH & Co.KG | Am Eichenhain 1 | 48531 Nordhorn | Germany | Tel. +49 5921 8197-0 | Fax +49 5921 8197-222 | info@ldn.de | www.ldn.de

Instructions for use
HistaSure™ ELISA Fast Track



Distributed in the US and Canada by:

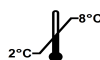
EAGLE BIOSCIENCES, INC.
20A NW BLVD, SUITE 112 NASHUA, NH 03063
P: 617-419-2019 F: 617-419-1110
WWW.EAGLEBIO.COM — INFO@EAGLEBIO.COM



EAGLE
BIOSCIENCES

REF

FC E-3600



1. Intended use and principle of the test

The **HistaSure™ ELISA** Fast Track is intended for the rapid semi-quantitative or quantitative determination of histamine in different scombroid fish types such as tuna, mahi mahi, sardines and for the determination of histamine in fishmeal.

Related Products:

HistaSure™ XL ELISA FC E-3900 (480 determinations)
--

The assay kit provides materials for the determination of derivatized histamine in food extracts. The derivatization is part of the preparation of the samples. By use of the acylation reagent, histamine is quantitatively derivatized into N-acylhistamine. The competitive Histamine ELISA kit uses the microtiter plate format. Histamine is bound to the solid phase of the microtiter plate. Acylated histamine and solid phase bound histamine compete for a fixed number of antiserum binding sites. When the system is in equilibrium, free antigen and free antigen-antiserum-peroxidase complexes are removed by washing. The substrate TMB/peroxidase reaction is monitored at 450 nm. The amount of antibody bound to the solid phase histamine is inversely proportional to the histamine concentration of the sample.

2. Introduction

Histamine testing in fresh fish is a possible control strategy that can be used by seafood processors in their HACCP program to address the hazard of scombrototoxin formation. Histamine is a product of decomposition of histidine caused by the growth of certain bacteria in seafood. The amount of the amine that forms is a function of bacterial species, the temperature and time of exposure, and may exceed 1,000 ppm (mg/kg). Fish containing high levels of histamine has been associated with many examples of poisoning commonly referred to as "scombroid poisoning," a major health problem for consumers. Scombrototoxic fish usually contains levels of histamine in excess of 200 ppm but such fish may be randomly dispersed within a lot. For large fish, histamine is found at variable levels even within individual fish. Quality control measures designed to minimize the occurrence of scombrototoxic fish require the determination of histamine levels in the range of approximately 10 to 200 ppm. Good quality fish contain less than 10 ppm histamine, a level of 30 ppm indicates significant deterioration, and 50 ppm is considered to be evidence of definite decomposition. The defect action level (DAL), the level at which regulatory actions are taken for histamine is 50 ppm (P. L. Rogers, W. F. Staruszkiewicz, Journal of Aquatic Food Product Technology, Vol. 9 (2) 2000 p. 5 - 17).

3. Procedural Cautions, Guidelines and Warnings

- (1) This kit is for professional use only. Users should have a thorough understanding of this protocol for the successful use of this kit. Only the test instruction provided with the kit is valid and has to be used to run the assay. Reliable performance will only be attained by strict and careful adherence to the instructions provided.
- (2) In order to reduce exposure to potentially harmful substances, wear lab coats, disposable latex gloves and protective glasses where necessary.
- (3) All kit reagents and specimens should be brought to room temperature (20 - 25 °C) and mixed gently but thoroughly before use.
- (4) When the use of water is specified for dilution or reconstitution, use deionized, distilled, or ultra-pure water.
- (5) The microplate contains snap-off strips. Unused wells must be stored at 2 °C to 8 °C in the sealed foil pouch and used in the frame provided. Wells are for single use only.
- (6) Once the test has been started, all steps should be completed without interruption. Make sure that required reagents, materials and devices are prepared ready at the appropriate time.
- (7) Incubation times do influence the results. All wells should be handled in the same order and time sequences.
- (8) To prevent contamination of reagents, use a new disposable pipette tip for dispensing each reagent, sample, standard and control.
- (9) Do not mix various lot numbers of kit components within a test and do not use reagents beyond expiry date as shown on the kit labels.
- (10) Avoid contact with Stop Solution containing 0.25 M H₂SO₄. It may cause skin irritation and burns. In case of contact with eyes or skin, flush immediately with water.
- (11) Some reagents contain sodium azide (NaN₃) as preservatives. In case of contact with eyes or skin, flush immediately with water. NaN₃ may react with lead and copper plumbing to form explosive metal azides. When disposing reagents, flush with a large volume of water to avoid azide build-up.
- (12) TMB substrate has an irritant effect on skin and mucosa. In case of possible contact, wash eyes with an abundant volume of water and skin with soap and abundant water. Wash contaminated objects before reusing them.

(13) For information on hazardous substances included in the kit please refer to Safety Data Sheets (SDS). The Safety Data Sheet for this product is available directly on the website of the manufacturer or upon request.

(14) Kit reagents must be regarded as hazardous waste and disposed of according to national regulations.

4. **Storage and stability**

Store the reagents at 2 - 8 °C until expiration date. Do not use components beyond the expiration date indicated on the kit labels. Do not mix various lots of any kit component within an individual assay.

5. **Materials**

5.1 Contents of the kit

BA D-0035 MB 48 **Master Block** - Ready to use

Content: 1 x 48 wells plate in a resealable pouch

BA E-0030 WASH-CONC 50x **Wash Buffer Concentrate** - Concentrated 50x

Content: Buffer with a non-ionic detergent and physiological pH

Volume: 1 x 20 ml/vial, light purple cap

BA E-0055 SUBSTRATE **Substrate** - Ready to use

Content: Chromogenic substrate containing tetramethylbenzidine, substrate buffer and hydrogen peroxide

Volume: 1 x 12 ml/vial, black cap

BA E-0080 STOP-SOLN **Stop Solution** - Ready to use

Content: 0.25 M sulfuric acid

Volume: 1 x 12 ml/vial, light grey cap

Hazards

identification:



H290 May be corrosive to metals.

H314 Causes severe skin burns and eye damage.

Controls - Ready to use

Cat. no.	Component	Colour/Cap	Concentration	Volume / Vial
FC E-3601	C⇒0 PPM	white	0 ppm	4 ml
FC E-3602	C⇒3 PPM	light yellow	3 ppm	4 ml
FC E-3603	C⇒10 PPM	orange	10 ppm	4 ml
FC E-3604	C⇒20 PPM	light green	20 ppm	4 ml
FC E-3605	C⇒30 PPM	light purple	30 ppm	4 ml
FC E-3606	C⇒50 PPM	dark blue	50 ppm	4 ml
FC E-3607	C⇒100 PPM	light grey	100 ppm	4 ml
FC E-3608	C⇒300 PPM	black	300 ppm	4 ml

Content: Ultrapure water with non-mercury stabilizer, spiked with a defined quantity of Histamine

FC E-3611 ACYL-BUFF **Acylation Buffer** - Ready to use

Content: TRIS buffer with non-mercury preservative

Volume: 2 x 50 ml/vial, white cap

FC E-3612 ACYL-REAG **Acylation Reagent** - Ready to use

Content: Acylation Reagent in DMSO, yellow coloured

Volume: 1 x 3 ml/vial, brown cap

FC E-3631 HIS **Histamine Microtiter Strips**- Ready to use

Content: 1 x 48 well (6x8) microwell plate precoated with antigen in resealable pouch with desiccant

FC E-3640 **HIS** **AB** **CONJ** **Histamine Antiserum Conjugate**- Ready to use
Content: Goat-anti Histamine IgG conjugated with peroxidase
Volume: 1 x 6 ml/vial, red cap

5.2 Additional materials and equipment required but not provided with the kit

- Precision pipette (50 µl)
- Pipette tips (50 µl)
- Manual repetitive pipette (e.g. the Brand HandyStep® S)
- Precision Dispenser Tips (5 ml, 25 ml; e.g. the PLASTIBRAND® PD-Tips)
- Grinder (mill) or house hold blender
- Graduated plastic cylinder (250 ml)
- Water (deionized, distilled, or ultra-pure)
- Scale (capable of weighing 5 – 50 grams, precision 0.1 gram)
- Funnel and filter paper (or alternatively a centrifuge)
- Timer
- Waterproof marker
- Absorbent material (paper towel)
- Microplate Vibration Shaker (shaking amplitude 2 mm; approx. 600 rpm, (e.g. PSU-2T Minishaker *))
- ELISA reader capable of reading absorbance at 450 nm (required for semi-quantitative and quantitative determination)
- Washing device (plate washer or manually)

**Available upon request!*

6. Test procedure

6.1 Preparation of reagents

Wash Buffer

Dilute the 20 ml Wash Buffer Concentrate with water (deionized, distilled, or ultra-pure) to a final volume of 1000 ml. Storage: up to 6 months at 2 - 8 °C.

Acylation Reagent

The Acylation Reagent has a freezing point of 18.5 °C. To ensure that the Acylation Reagent is liquid when being used, it must be ensured that the Acylation Reagent has reached room temperature and forms a homogeneous, crystal-free solution before being used.

Histamine Microtiter Strips

In rare cases residues of the blocking and stabilizing reagent can be seen in the wells as small, white dots or lines. These residues do not influence the quality of the product.

6.2 Sample preparation

The following protocols for the sample preparations are based on the AOAC Official Method 937.07
Sampling should be performed according to national regulation.

A. FRESH FISH • FROZEN FISH

- Keep (fresh) fish frozen prior to analysis.
- Thaw samples under refrigeration or in cold water. Do not thaw the samples in a heated water bath. Discard draining.
- Once thawed, store the samples refrigerated (2 - 8 °C) prior to testing.

whole fish:

Clean, scale and eviscerate fish. In case of small fish 6 in. (≤ 15 cm), use 5 – 10 fish. In case of large fish, from each of ≥ 3 fish cut 3 cross-sectional slices 1 in. (2.5 cm) thick, 1 slice from just back of pectoral fins, 1 slice halfway between first slice and vent, and 1 slice just back of vent. Remove bone. Blend combined samples until homogenous.

fish filet:

Use entire piece. Blend until homogenous.

B. CANNED FISH and other CANNED MARINE PRODUCTS

Place entire content of the can (meat and liquid) in a blender and blend until homogenous.

C. CANNED MARINE PRODUCTS PACKED in OIL, SAUCE, BRINE or BROTH

Drain for 2 minutes on number 8 sieve or dab away the fluid with a paper towel. Place the meat in a blender and blend until homogenous.

D. FISHMEAL

Mix sample until homogenous.

6.3 Extraction

- **Weigh 10 g** of prepared fish sample / fish meal, **add 240 ml water** (deionized, distilled, or ultra-pure) and **homogenize*** for 1-2 minutes in a grinder or blender.
*): Instead of homogenization **fish meal samples** are stirred for 10 minutes at room temperature.
- **Filter** the homogenate through folded filter paper (alternatively an aliquot of the homogenate can be centrifuged for 5 minutes at maximum speed). *If a **lipid layer** forms remove it by suction!*
- Use 50 µl of the **sample extract** for the acylation.

6.4 Histamine ELISA

For the subsequent steps (Acylation and ELISA) allow all reagents and samples to reach room temperature.

A. SEMI-QUANTITATIVE DETERMINATION

For the semi-quantitative determination select the desired cut-off you need from the controls provided with the kit. The kit controls have the following concentrations:

control 3, 10, 20, 30, 50, 100 or 300 ppm.

B. QUANTITATIVE DETERMINATION

For the quantitative determination use the following controls provided with the kit:

control 0, 3, 10, 30, 100 and 300 ppm.

These 6 controls are used to establish the standard curve (please refer to section 7.)

6.4.1 Acylation

1. Pipette 50 µl of control(s) and sample extracts into the respective wells of the Master Block .
2. Add 1.5 ml of Acylation Buffer in 1 (!) pipetting step to all wells. <i>The use of a repetitive pipette together with a new Precision Dispenser Tip (25 ml, please refer to chapter 5.), is recommended.</i>
3. Add 50 µl of Acylation Reagent to all wells. (Colour change from yellow to pink!) Continue without any delay with step 4.! ⚠ <i>The use of a repetitive pipette together with a new Precision Dispenser Tip (2.5 ml, please refer to chapter 5.), is recommended.</i>
4. Incubate 5 min at RT (20 - 25 °C) on a shaker (approx. 600 rpm). ⚠ Make sure that mixing is complete (slight pink colour).
5. Take 50 µl for the ELISA

6.4.2 Histamine ELISA

1. Pipette 50 µl of the acylated control(s) and samples into the wells of the Histamine Microtiter Strips .
2. Pipette 100 µl of the Histamine Antiserum Conjugate into all wells. <i>The use of a repetitive pipette together with a new Precision Dispenser Tip (5 ml, please refer to chapter 5.), is recommended.</i>
3. Incubate 10 min at RT (20 - 25 °C) on a shaker (approx. 600 rpm).
4. Discard or aspirate the contents of the wells. Wash the plate 3 x by adding 300 µl of Wash Buffer , discarding the content and blotting dry each time by tapping the inverted plate on absorbent material.
5. Pipette 100 µl of the Substrate into all wells. <i>The use of a repetitive pipette together with a new Precision Dispenser Tip (5 ml, please refer to chapter 5.), is recommended.</i>
6. Incubate for 10 min at RT (20 - 25 °C) on a shaker (approx. 600 rpm). ⚠ Avoid exposure to direct sunlight!
7. Add 100 µl of the Stop Solution to each well and shake the microtiter plate to ensure a homogeneous distribution of the solution. <i>The use of a repetitive pipette together with a new Precision Dispenser Tip (5 ml, please refer to chapter 5.), is recommended.</i>
8. Read the absorbencies of the solutions in the wells within 10 minutes using a microplate reader set to 450 nm .

7. Calculation of results

A. SEMI-QUANTITATIVE RESULTS:

If the absorbance of the sample is **higher** than that of the selected **Cut-Off Control**, the sample has **passed**.

If the absorbance of the sample is **lower** than that of the selected **Cut-Off Control**, the sample has **failed**.

B. QUANTITATIVE RESULTS:

The absorbance readings of the 6 controls (0, 3, 10, 30, 100 and 300 ppm) are used to establish a standard curve.

Plot the absorbance readings of the controls (y-axis, linear) against the corresponding control concentrations (x-axis, log) using a concentration of 0.001 ppm for the 0-control (*this alignment is mandatory because of the logarithmic presentation of the data*). For the curve fitting a non-linear regression has to be applied.

The concentrations of the samples can be read **directly** from this standard curve.

⚠ *If a sample is off-curve it has to be diluted with water 1:10 and re-assayed. The result obtained has to be multiplied by the dilution factor of 10.*

8. Warranty

This test kit was produced according to the latest developments in technology and subjected to stringent internal and external quality control checks. Any alteration of the test kit or the test procedure as well as the usage of reagents from different charges may have a negative influence on the test results and are therefore not covered by warranty. The manufacturer is not liable for damages occurring during transit.

9. Application lists for different kind of fish samples

All fish samples tested so far are suitable for the **HistaSure™ ELISA^{Fast Track}**. The lists below depict some major applications in different matrices.

	Fish species	Presentation
Species validated through AOAC Certification	Tuna	canned chunk light
		fresh/frozen yellow fin
	Mahi Mahi	fresh/frozen
	Sardines	canned in oil
	Fishmeal	

	Fish species	Presentation
Species validated through in-house testing	Mackerel	smoked
	Anchovy	fresh
		brined
		in sauce
	Shad	dry salted
		fermented
	Herring	smoked
	Salmon	smoked
	Bonito	lakerda
	Swordfish	fresh
Marlin	fresh	

Fish products validated through in-house testing	Maldive fish
	Fish sauce (sardines/anchovy)
	Oyster sauce (oyster/sardines)

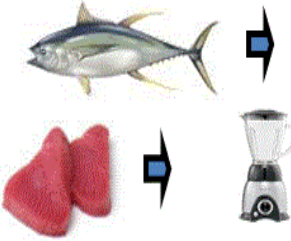
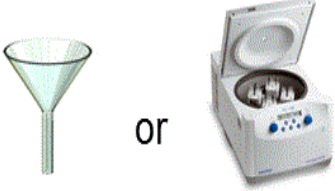
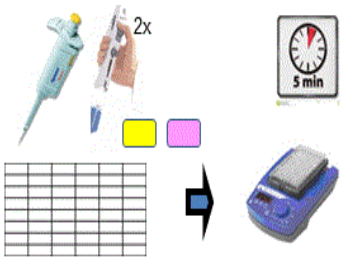
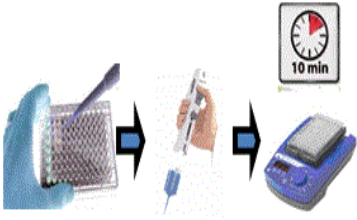
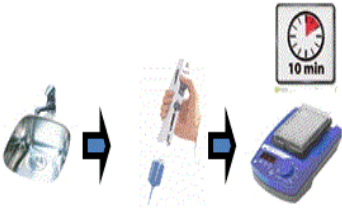
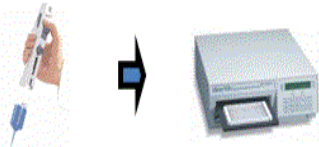
10. Assay characteristics

AOAC performance tested method for fresh/frozen yellowfin tuna, canned tuna-chunk light in water, frozen mahi mahi, canned sardines in oil and fish meal.

Analytical Specificity (Cross Reactivity)	Substance			Cross Reactivity (%) Histamine	
	Histamine			100	
	L-Tryptophan			nd	
	Tryptamine			nd	
	3-Methylhistamine			0.44	
	L-Histidine			nd	
	L-Tyrosine			nd	
	L-Phenylalanine			nd	
	Tyramine			0.69	
	Cadaverine			0.40	
	Spermine			nd	
	Putrescine			nd	
	Trimethylamine			nd	
	nd = not detectable				
Accuracy and Precision					
Recovery				Intra CV	
Sample	Fortification range (ppm)	Mean recovery (%)	Recovery range (%)	Mean CV (%) (n=7)	Range CV (%)
Fresh/Frozen	5.25 – 218.6	91.8	85.2 – 99.6	7.58	4.13 – 10.8
Canned Tuna	5.49 – 267.3	99.8	94.7 – 106.5	8.74	3.17 – 13.9
Frozen Mahi	6.38 – 199.0	87.3	79.1 – 103.1	6.19	2.97 – 9.56
Canned	5.27 – 249.9	87.6	76.8 – 99.7	5.65	1.80 – 9.22
Fish meal	9.4 – 244.2	86.0	79.0 – 93.9	4.89	2.09 – 7.92
Mean (ppm)		n		Inter CV %	
Lot to Lot		25.8		3 lots	
				4.16	
Detection limits					
LOD (Limit of Detection) 0.44 ppm			LOQ (Limit of Quantification) 1.31 ppm		
Method Comparison					
LDN HistaSure Elisa vs AOAC 977.13 fluorometric method: Fresh/Frozen Tuna, Canned Tuna, Frozen Mahi Mahi and Canned Sardines					

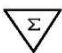

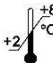







Flow Chart HistaSure™ ELISA^{Fast Track} FC E-3600 / FC E-3900 Quantitative / Semi-Quantitative

- For a quantitative result use the following controls provided in the kit: control 0, 3, 10, 30, 100 and 300 ppm
- For a semi-quantitative result select the desired cut-off you need from the controls provided with the kit

<p>1. Extraction</p> <ul style="list-style-type: none"> Weigh 10 grams of fish or fishmeal Add 240 ml of water Homogenize 1 -2 min. (fishmeal stir 10 min) 		<p>2. Filter or centrifuge</p> <ul style="list-style-type: none"> Filter the homogenate or centrifuge an aliquot of the homogenate (5 min max speed) If a lipid layer forms, remove by suction! 															
<p>3. Acylation</p> <ul style="list-style-type: none"> Pipet 50 µl control(s) / sample extract into the masterblock Add 1.5 ml Acylation Buffer Add 50 µl Acylation Reagent Incubate for 5 min on a shaker 		<p>4. ELISA: step 1</p> <ul style="list-style-type: none"> Pipet 50 µl acylated control(s)/sample into the ELISA wells Add 100 µl Antiserum Conjugate Incubate for 10 min on a shaker 															
<p>5. ELISA: step 2 and 3</p> <ul style="list-style-type: none"> Discard the content of the wells and wash each well 3x Add 100 µl Substrate Incubate for 10 min on a shaker 		<p>6. ELISA: step 4</p> <ul style="list-style-type: none"> Add 100 µl of Stop Solution Shake shortly Read the plate at 450 nm 															
<p>Calculation of Results: Quantitative</p> <ul style="list-style-type: none"> Plot the absorbance readings of the calibrators (y-axis, linear) against the corresponding calibrator concentrations (x-axis, log). <table border="1" style="width: 100%; border-collapse: collapse; margin: 5px 0;"> <thead> <tr> <th style="text-align: left;">Control Histamine (ppm)</th> <th>0.001</th> <th>3.0</th> <th>10</th> <th>30</th> <th>100</th> <th>300</th> </tr> </thead> <tbody> <tr> <td style="height: 20px;"> </td> <td> </td> <td> </td> <td> </td> <td> </td> <td> </td> <td> </td> </tr> </tbody> </table> <ul style="list-style-type: none"> Use non-linear regression for curve fitting Read the concentrations of the samples 		Control Histamine (ppm)	0.001	3.0	10	30	100	300								<p>Calculation of Results: Semi-quantitative</p> <ul style="list-style-type: none"> If the absorbance of the samples > then that of selected cut-off, the sample has passed If the absorbance of the samples < then that of selected cut-off, the sample has failed 	
Control Histamine (ppm)	0.001	3.0	10	30	100	300											

⚠ The liability of the manufacturer shall be limited to the replacement of defective products. The manufacturer takes no liability for any damages or expenses arising directly or indirectly from the use of this product.

Symbols:

	Contains sufficient for <n> tests		Manufacturer		Storage temperature
	Catalogue number		Batch code		Expiry date
	Caution		Content		Consult instructions for use
	For research use only!				

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 insert, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.

distributed in the US/Canada by:



EAGLE
BIOSCIENCES

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH
03063

Phone: 617-419-2019 FAX: 617-419-1110

www.EagleBio.com • info@eaglebio.com