

## MaxSignal® IAC for Total Aflatoxin Aflatoxin B1/B2/G1/G2/M1/M2 Immunoaffinity Column

Catalog #FOOD-1513-01

**ISO 9001** Quality Assurance  
Manufactured in compliance with our ISO 9001 certified quality management system.

©PerkinElmer, Inc. • 1513-01 Version 2019-1

### NOTICES

Except as specifically set forth in its terms and conditions of sale, PerkinElmer, Inc. and its subsidiaries ("PerkinElmer"), make no warranty of any kind, either express or implied, with regard to this document or the use of the product, including, but not limited to, the implied warranties of merchantability and fitness for a particular purpose. PerkinElmer shall not be liable for any omissions, or errors or inaccuracies contained herein. PerkinElmer shall not be liable for any damages, including special, consequential or incidental damages in connection with furnishing, performance or use of this material or the product.

PerkinElmer has made every effort to ensure that this document is accurate. PerkinElmer disclaims liability for any inaccuracies or omissions that may have occurred. Information in this document is subject to change without notice and does not represent a commitment on the part of PerkinElmer. This document supersedes and replaces all information supplied prior to the publication hereof. The material in this document is for informational purposes only. PerkinElmer makes no commitment to update or keep current the information in this document and reserves the right to make improvements to this document and/or to the products described in this document, at any time without notice. If you find information in this document that is incorrect, misleading, or incomplete, PerkinElmer would appreciate your comments and suggestions.

MaxSignal® IAC for Total Aflatoxin is intended as a screening tool for research use only. This product is not intended for clinical diagnostic use.

### COPYRIGHT INFORMATION

This document, including all photographs and illustrations, contains proprietary information that is protected by copyright. All rights are reserved. No part of this publication may be reproduced in any form whatsoever or translated into any language without the prior, written permission of PerkinElmer. Copyright © 2020 PerkinElmer, Inc. Produced in the U.S.A.

### TRADEMARKS

Registered names, trademarks, etc. used in this document, even when not specifically marked as such, are protected by law. MaxSignal® is a registered trademark of PerkinElmer.

PerkinElmer, Inc. P: (800) 762-4000 or  
940 Winter Street (+1) 203-925-4602  
Waltham, MA 02451 USA [www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

## RELATED PRODUCTS

CATALOG #	PRODUCT	QTY
FOOD-1000-00	AuroFlow™ X	00 strips
FOOD-1000-00	MaxSignal® X	00

## GENERAL INFORMATION

### Purpose

The immunoaffinity column can selectively adsorb aflatoxins from the sample solution, thereby purifying the sample. The purified sample solution can then be directly used for HPLC analysis.

Affinity columns can be used in combination with HPLC to achieve rapid testing, and to increase signal-to-noise ratio and improve the accuracy of the detection method.

This affinity column can be used for the extraction and purification of aflatoxins (B<sub>1</sub>, B<sub>2</sub>, G<sub>1</sub>, G<sub>2</sub>, M<sub>1</sub>, M<sub>2</sub>) in samples.

### Overview

Aflatoxins are toxic metabolites of a class of fungi (such as *Aspergillus flavus* and *Aspergillus parasiticus*). They are highly carcinogenic and are found mainly in grains, peanuts, nuts, cottonseeds, animal feed, vegetable oils, as well as animal tissues and blood. Among them, Aflatoxin B<sub>1</sub> (AFT B<sub>1</sub>) ranks first in terms of toxicity, carcinogenicity and frequency of contamination. Aflatoxin M<sub>1</sub> is a hydroxylated metabolite of Aflatoxin B<sub>1</sub> and a strong carcinogen. Cow milk and its products are prone to Aflatoxin M<sub>1</sub> contamination.

### Principle

The basis of the measurement is the antigen-antibody reaction. Antibodies are connected to the column and the aflatoxin in the sample is extracted, filtered, and diluted, and then passed slowly through the aflatoxin immunoaffinity column. The toxins bind to the antibodies in the column and the immunoaffinity column is then washed to remove other unrelated substances that have not been bound. Aflatoxin is eluted with methanol, and then injected into the analytical instrument for detection.

## KIT CONTENTS, STORAGE, & SHELF LIFE

Each box contains Aflatoxin immunoaffinity columns of various specifications and 1 instruction manual. Store the entire kit at 2–8°C. Do not use this product past the expiration date indicated on the box label.

### Required Materials Not Provided with the Kit

- HPLC
- Derivatization device: such as optical derivatization device, photochemical derivatization device, iodine derivatization device
- Centrifuge capable of at least 3,000-4,000 x g
- Nitrogen gas evaporator apparatus
- Nitrogen gas tank and pressure regulator
- LC-MS (LC-MS/MS)
- Air-pressure controller
- Air pump
- Balance with 0.01 g readability
- High-speed homogenizer (i.e. rotary shaker, vortexer, stomacher, or equivalent) (maximum speed ≥ 10,000 RPM)
- Grinder
- Sieving screen: 2-mm
- pH meter (or pH test paper)
- Graduated cylinder: 10 mL & 100 mL
- Funnel: 50 mL
- Syringe: 10 mL & 20 mL
- Pipette and pipette tips
- Homogenization flask (or 250-mL conical flask with pestle)
- Sample tubes and bottles
- Qualitative filter paper
- Microfiber filter paper (e.g. Whatman 934-AH)
- Column holder and syringe connector plug (for use with 6-mL immunoaffinity columns)
- Methanol (CH<sub>3</sub>OH): Chromatography Grade
- Acetonitrile (CH<sub>3</sub>CN): Analytical Grade
- Disodium hydrogen phosphate dodecahydrate (Na<sub>2</sub>HPO<sub>4</sub> · 12H<sub>2</sub>O): Analytical Grade
- Acetic acid (CH<sub>3</sub>COOH): Chromatography Grade
- Potassium dihydrogen phosphate (KH<sub>2</sub>PO<sub>4</sub>): Analytical Grade
- Potassium chloride (KCl): Analytical Grade
- Sodium chloride (NaCl): Analytical Grade
- Tween-20® (C<sub>58</sub>H<sub>114</sub>O<sub>26</sub>): Analytical Grade
- Hydrochloric acid (HCl): Analytical Grade
- Sodium hydroxide (NaOH): Analytical Grade
- Distilled/deionized water

## PRECAUTIONS

- Allow the immunoaffinity column to equilibrate to room temperature (20–25°C) before use.
- The immunoaffinity column should be stored at 2–8°C; do not freeze.
- Do not use any expired immunoaffinity column.
- The sample volume can be increased or decreased appropriately as needed. The volume of the extraction solution should be adjusted accordingly.
- The pH of the loading solution onto the immunoaffinity column should be 6–8. If it deviates from this range, the pH should be adjusted with dilute hydrochloric acid or dilute sodium hydroxide.
- Maintaining consistency (such as polarity, pH, and concentration) between the test solvent loaded into any analytical instrument and the mobile phase can help eliminate any adverse solvent effects.

- Column capacity:

Toxin name	Column capacity: ng	Toxin name	Column capacity: ng
AFT B1	300	AFT G2	100
AFT B2		AFT M1	
AFT G1		AFT M2	

When the content of the toxin in the sample divided by the dilution factor is higher than the column capacity, it is necessary to reduce the volume of the sample solution appropriately and test again.

- WARNING:** Aflatoxin is toxic and carcinogenic; protective equipment such as gloves and masks should always be used during handling.
- Vessels and tools used to handle toxin solutions should be completely immersed in a sodium hypochlorite solution (5% v/v) overnight.
- Ensure the LC-MS/MS is clean and the tubing is primed appropriately for each run.
- Follow appropriate instrument precautions if using HPLC.
- If milk powder samples are difficult to dissolve, incubate at 50°C for 30 minutes. If it is still not dissolved, increase the proportion of water appropriately.
- For samples that are particularly turbid after centrifugation, increase the rotation speed or increase the centrifugation time appropriately.

## REAGENT PREPARATION

- Preparation of Extraction Solution 3: 70% v/v Methanol-water**  
Combine 700 mL of methanol and 300 mL of distilled/deionized water. Bring to 1 L final volume with distilled/deionized water. Mix well.
- Preparation of Extraction Solution: 80% v/v Acetonitrile-water**  
Combine 800 mL of acetonitrile and 200 mL of distilled/deionized water. Bring to 1 L final volume with distilled/deionized water. Mix well.
- Preparation of 1% Tween-water solution**  
Combine 1 mL of Tween-20, with distilled/deionized water to a final volume of 100 mL.
- Preparation of Phosphate Buffer Solution (PBS)**  
Weigh out 8.00 g of NaCl, 1.20 g of Na<sub>2</sub>HPO<sub>4</sub> (or 2.92 g of Na<sub>2</sub>HPO<sub>4</sub> · 12H<sub>2</sub>O), 0.20 g of KH<sub>2</sub>PO<sub>4</sub> and 0.20 g of KCl to a container and add 900 mL of distilled/deionized water to dissolve the powders, adjust the pH to 7.4 by using HCl, then add distilled/deionized water to a final volume of 1000 mL.

## SAMPLE PREPARATION

### Method 1: Suitable for samples of grain crops such as corn, wheat, nuts, corn oil, flower oil, feed, etc.

- Weigh out 25 g ± 0.01 g of sample in a bottle. Add 5 g of sodium chloride (NaCl) and 125 mL of Extraction Solution 3. Solid samples should be homogenized to pass through a 2-mm sieve before use.
- Homogenize, such as vortex, at high speed (≥ 10,000 RPM) for 1 minute, or shake vigorously on a shaker (200-300 RPM) for 20 minutes. Filter with microfiber filter paper. Collect the filtrate.
- Combine 10 mL of the filtrate with 20 mL of distilled/deionized water to dilute, then filter with microfiber filter paper, and collect the filtrate as sample solution. Use 15 mL of the sample solution with the immunoaffinity column for purification.

Dilution factor = 1

### Method 2: Suitable for samples of flour, wheat flour, chili, pepper, soybean paste, sesame oil and other vegetable oils, straw, peanuts and their products

- Weigh 25 g ± 0.01 g of sample into a bottle. Add 125 mL of Extraction Solution. Solid samples should be homogenized to pass through a 2-mm sieve.
- Homogenize, such as vortex, at high speed (≥ 10,000 RPM) for 1 minute, or shake vigorously on a shaker (200-300 RPM) for 20 minutes. Filter with microfiber filter paper. Collect the filtrate.
- Combine 10 mL of the filtrate with 70 mL of distilled/deionized water to dilute, then filter with microfiber filter paper, and collect the filtrate as sample solution. Use 20 mL of the sample solution with the immunoaffinity column for purification.

Dilution factor = 2

### Method 3: Suitable for Chinese medicine (19 types of Chinese medicine specified in Version 2015 of the Pharmacopoeia)

- Weigh out 15 g ± 0.01 g of sample (solid samples should be ground and passed through a 1-mm sieving screen) and add to 75 mL of Extraction Solution (80% acetonitrile-water solution).
- Homogenize, such as vortex, at high speed (≥ 10,000 RPM) for 1 minute or shake vigorously on a shaker (200-300 RPM) for 20 minutes. Filter with microfiber filter paper. Collect the filtrate.
- Combine 10 mL of the filtrate with 40 mL of 1% Tween-water solution, then filter with microfiber filter paper, and collect the filtrate as sample solution.
- Use 25 mL of the sample solution with the immunoaffinity column for purification.

Dilution factor = 1

### Method 4: Suitable for tea samples such as black tea, green tea, oolong tea, and Pu'er tea

- Weigh 5 g ± 0.01 g of sample into a bottle. Add 25 mL of Extraction Solution 1. Solid samples should be homogenized to pass through a 1-mm sieve before use.
- Homogenize, such as vortex, at high speed (≥ 10,000 RPM) for 1 minute, or shake vigorously on a shaker (200-300 RPM) for 20 minutes. Filter with microfiber filter paper. Collect the filtrate.
- Combine 10 mL of the filtrate with 70 mL of distilled/deionized water. Mix well. Filter with microfiber filter paper, and collect the filtrate.
- Use 20 mL of the filtrate as the final sample for testing.

Dilution factor = 2

### Method 5: Liquid milk, yogurt

- Weigh out 4 g of sample that has been uniformly mixed (accurate to 0.001 g) and place it into a 50-mL centrifuge tube, add 10 mL of methanol, and vortex for 3 minutes.
- Centrifuge at 3,000-4,000 x g for 10 minutes.
- Transfer 7 mL of the supernatant into a centrifuge tube, add 40 mL of distilled/deionized water. Mix well.

Dilution factor = 0.5

### Method 6: Milk powder, special dietary foods

- Weigh out 1 g of the sample (accurate to 0.001 g) and place it into a 50-mL centrifuge tube, then add 4 mL of hot distilled/deionized water at 50°C and vortex to mix well. If the milk powder cannot be completely dissolved, place the centrifuge tube in a water bath of 50°C to dissolve the milk powder completely. After the sample has cooled to 20°C, add 10 mL of methanol, and vortex for 3 minutes.
- Centrifuge at 3,000-4,000 x g for 10 minutes.
- Transfer 7 mL of the supernatant into a centrifuge tube add 40 mL of distilled/deionized water. Mix well.

Dilution factor = 2

### Method 7: Cheese

- Weigh out 1 g of finely diced sample that has been passed through a round-hole sieve with 1–2 mm diameter (accurate to 0.001 g) and place it into a 50-mL centrifuge tube, add 1 mL of distilled/deionized water and 14 mL of methanol and shake for 30 minutes.
- Centrifuge at 3,000-4,000 x g for 10 minutes.
- Transfer 10 mL of the supernatant to a round bottom flask.
- Subject to rotary evaporation until 2 mL or less remain, dilute with 30 mL of PBS.

Dilution factor = 1.5

## OPERATING PROCEDURE

- Remove the column and place into a column holder. Remove the plunger of a syringe, then attach the syringe through the connector plug above the column to complete the connection. Secure to an air-pressure controller, if available.
- Transfer the appropriate amount of the solution processed in Sample Preparation to fill the syringe.
- Remove the cap under the affinity column (do not discard as this will be used in the next step). Adjust the air-pressure to have a flow rate of 1–2 drops/second.
- After the liquid has flowed through wash with 10 mL of 0.1% Tween-water solution, then wash with 10 mL of distilled/deionized water at a flow rate of 2-3 drops per second;
- After the liquid has flowed through, prepare the sample from Methods 1, 2, 3, or 4, by adding 1 mL of methanol for elution at a flow rate of 1 drop/second, and collect the eluate and adjust the volume to precisely 1 mL. Pass the eluate through a 0.22 µm micropore filter, transfer into a sample bottle, and use it for HPLC analysis.
- Prepare the sample from Methods 5, 6, or 7, after replacing the syringe with a new one, load 2 x 2 mL of acetonitrile for elution and collect the eluate. Place the eluent under a slow stream of nitrogen gas at 50°C to evaporate any residual solvents. Dissolve the dried residue with 1 mL of the appropriate solvent. Filter the solution through a 0.22 µm micropore filter, then transfer to a sample bottle to be used for HPLC analysis.

*Also applicable to GB 5009.22-2016, GB 5009.24-2016 and pharmacopoeia methods.*

## INTERPRETATION OF RESULTS

Aflatoxin Concentration = Detected Concentration x Dilution Factor