

MaxSignal® IAC for T-2/HT-2 Toxin

Catalog #FOOD-1512-01

ISO
9001
QUALITY
ASSURANCE

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

©PerkinElmer, Inc. • 1512-01 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 T-2/HT-2 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 © 2020-2021 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.
940 Winter Street
Waltham, MA 02451 USA

P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2020-2021, 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-1505-01	MaxSignal® IAC 6-in-1 Combo Mycotoxin Immunoaffinity Column	10 Tests

GENERAL INFORMATION

Purpose

The immunoaffinity column can selectively adsorb T-2 and HT-2 toxins from the sample solution, thereby purifying the sample. The purified sample solution can be used for LC-MS/MS analysis after concentrated with nitrogen gas and resuspended or for HPLC analysis after derivatization.

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

Overview

T-2 toxin (T-2) is a mycotoxin produced by a variety of *Fusaria*.

It mainly pollutes food crops such as wheat, barley, and corn, and their products, and is relatively harmful to human health and cattle industry. HT-2 toxin is the main metabolite of T-2 toxin. These two toxins are strong inhibitors of protein synthesis in the body and mainly affect the functions of blood, liver, kidney, pancreatic muscle and lymphocytes. The general clinical symptoms after poisoning are anorexia, vomiting, diarrhea, growth arrest, as well as reproduction and neurological dysfunction, etc.

Principle

This IAC functions by an antigen-antibody interaction. All four main toxin antibodies are embedded throughout the column. After a sample is extracted and filtered, it is slowly passed through the IAC. The toxoids bind to their corresponding antibody in the column. The IAC is then washed to remove unbound substances. The toxins are then eluted with the eluent, concentrated using nitrogen gas, then injected into an analytical instrument for detection.

KIT CONTENTS, STORAGE, & SHELF LIFE

Each kit contains T-2/HT-2 toxin 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 Certificate of Analysis.

Required Materials Not Provided with the Kit

- HPLC
- 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 (CH₃OH): Chromatography Grade
- Acetonitrile (CH₃CN): Analytical Grade
- Disodium hydrogen phosphate dodecahydrate (Na₂HPO₄ · 12H₂O): Analytical Grade
- Acetic acid (CH₃COOH): Chromatography Grade
- Potassium dihydrogen phosphate (KH₂PO₄): Analytical Grade
- Potassium chloride (KCl): Analytical Grade
- Sodium chloride (NaCl): Analytical Grade
- Tween-20® (C₅₈H₁₁₄O₂₆): Analytical Grade
- Polyethylene glycol: Analytical Grade; polyethylene glycol is used to improve the filtration speed, if the supernatant is obtained by centrifugation, this reagent can be omitted
- 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: the total capacity of the T-2/HT-2 column is 2000 ng, 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 retest.
- **WARNING:** T-2/HT-2 toxin is cytotoxic; 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.

REAGENT PREPARATION

- 1. Preparation of Extraction Solution: 80% v/v Methanol-water**
Combine 800 mL of methanol and 200 mL of distilled/deionized water. Bring to 1 L final volume with distilled/deionized water. Mix well.
- 2. Preparation of Diluent Solution: 0.05M PBS, pH 7.3**
Weigh out 8 g of NaCl, 0.2 g of KCl, 0.2 g of KH_2PO_4 and 1.16 g of $\text{Na}_2\text{HPO}_4 \cdot 12\text{H}_2\text{O}$ into a large, graduated bottle. Dissolve with 800 mL of distilled/deionized water, then bring to a final volume of 1 L. Mix well.
- 3. Preparation of Eluent Solution: 2% v/v acetic acid-methanol**
Combine 2 mL of acetic acid and 98 mL of methanol. Mix well.

SAMPLE PREPARATION

1. Weigh out $25\text{ g} \pm 0.01\text{ g}$ of sample in a bottle. Add 5 g of sodium chloride (NaCl) and 125 mL of Extraction Solution. Solid samples should be homogenized to pass through a 2-mm sieve before use.
2. Homogenize, such as vortex, at high speed ($\geq 10,000\text{ RPM}$) for 1 minute, or shake vigorously on a shaker (200-300 RPM) for 20 minutes.
3. Combine 10 mL of the filtrate with add 40 mL of Diluent, then filter with microfiber filter paper, and collect the filtrate.
4. Use 25 mL of the sample with immunoaffinity column for purification.

Dilution factor = 1

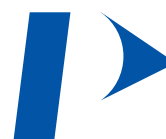
OPERATING PROCEDURE

1. 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.
2. Transfer the appropriate amount of the solution processed in Sample Preparation to fill the syringe.
3. 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.
4. After the liquid has completely flowed through add 10 mL of distilled/deionized water at a rate of 2-3 drops per second. Repeat this step one more time.
5. After the liquid has completely flowed through add 1 mL of 2% acetic acid-methanol solution, collect the eluent with a sample bottle at a flow rate of 1 drop per second, concentrate by passing a slow stream of nitrogen at 50°C to evaporate any residual solvent.
6. LC-MS/MS testing can be conducted, or HPLC testing can be conducted after derivatization in accordance with the GB 5009.118—2016, GB/T 28718-2012 PLC standards.

INTERPRETATION OF RESULTS

T-2/HT-2 Toxin Concentration = Detected Concentration of T-2/HT-2 Toxin
x Dilution Factor

NOTES



PerkinElmer®
For the Better