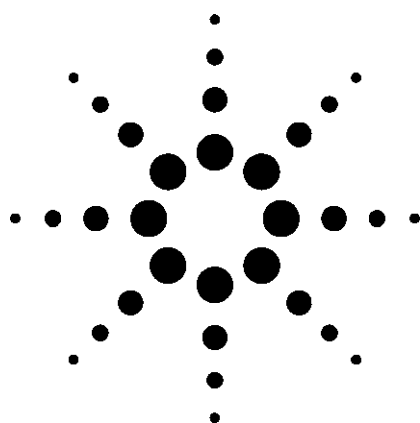


Separation of Aflatoxins by HPLC

Application



Environmental, Food Safety

Authors

Coral Barbas
Facultad de CC Experimentales y de la Salud
Universidad San Pablo-CEU
Urbanización Montepríncipe
Boadilla del Monte, 28668 Madrid
Spain

Andre Dams*
Agilent Technologies, Inc.
Amstelveen
The Netherlands

Ronald E. Majors
Agilent Technologies, Inc.
2850 Centerville Road
Wilmington, DE 19808-1610
USA

*Present address:
Andre Dams
Analytical Consultancy, Amstelveen
The Netherlands

Abstract

Four target aflatoxins (B₁, B₂, G₁, and G₂) were separated by HPLC using an isocratic ternary mixture of water, methanol and acetonitrile, and detected using UV 365 nm. A baseline separation was achieved in less than 5.5 min.

Introduction

Aflatoxins are mycotoxins that are produced by various *Aspergillus flavus* molds. Not only are these compounds extremely toxic, but they are also mutagenic, teratogenic (causing fetal

abnormalities), and carcinogenic. Unfortunately, *A. flavus* is a common mold found in tropical and subtropical countries and has been found to cause aflatoxin contamination. This contamination is a result of inadequate storage conditions for certain agricultural commodities such as peanuts, cereal seeds, dried fruit, and a wide range of tree nuts such as pistachio, pecans, walnuts, almonds, and herbal seeds such as red and black pepper, cloves, and cinnamon. Because of their toxic and carcinogenic nature, there is a very low minimum detectable quantity (MDQ) for aflatoxin contamination in food.

Chemical Nature

Although 18 different aflatoxins have been identified, the four most prevalent aflatoxins are B₁, B₂, G₁, and G₂, whose chemical structures are shown in Figure 1. Aflatoxin B₁ is one of the most potent and abundant environmental mutagens and carcinogens known. Aflatoxins are quite stable in many foods and are fairly resistant to degradation. Collectively, the aflatoxins are chemical derivatives of difurancoumarin. Pure aflatoxin B₁ is a pale-white to yellow crystalline, odorless solid. Aflatoxins are freely soluble in moderately polar solvents such as chloroform, methanol, and dimethyl sulfoxide, and dissolve in water to the extent of 10–20 mg/L. In methanol, they have fairly strong extinction coefficients (around 10,000) at 265 nm and 360–362 nm. They fluoresce under UV radiation, and fluorescence detection is often used for trace analysis in HPLC. Since there are differences in fluorescence yields between B₁ and B₂, and between G₁ and G₂, it can be useful to run both UV and fluorescence detectors in series [1]. Aflatoxins have no polar functional groups, and can be separated by virtue of their hydrophobicity.



Agilent Technologies

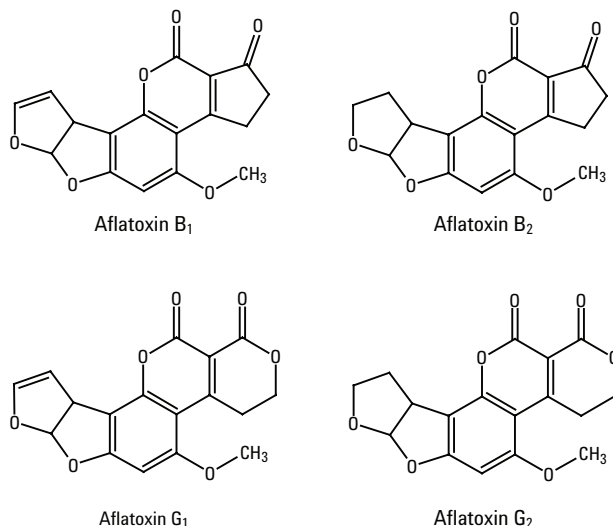


Figure 1. Chemical structures of target aflatoxins.

HPLC Methodology

While thin-layer chromatography was frequently used in the past, HPLC has been used in recent years because of its ease of operation and better quantitation. Most HPLC methods published to date have used reversed-phase HPLC on C18 bonded phases [1–4], where the aflatoxins are separated by their hydrophobicity. Most published separations have been performed on 5- μ m columns of 25-cm in length. The use of smaller particle size packings in shorter columns with faster separation times are now in vogue. These columns show that the same separations can be achieved in less time than on the longer columns with similar resolution. In the present study, we desired to show that using such a column can provide improved results compared to the older methods. See Figure 2.

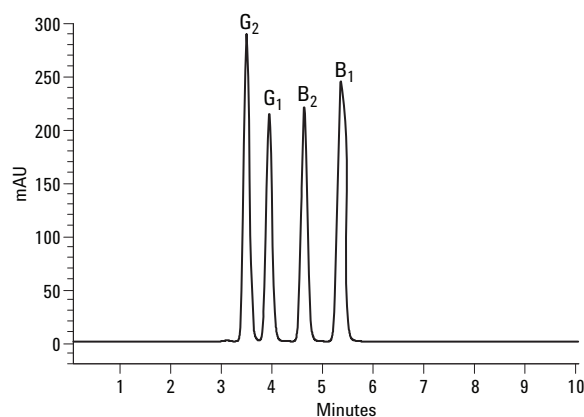


Figure 2. Reversed-phase separation of target aflatoxins using ZORBAX XDB-C18 Rapid Resolution column.

Experimental Conditions

Chemicals: The aflatoxins were purchased from Sigma Aldrich (Madrid, Spain).

HPLC Conditions

Column: ZORBAX Eclipse XDB-C18, 4.6 mm \times 150 mm, 3.5 μ m

Mobile phase: Water/MeOH/ACN; 50/40/10 (V/V/V)

Flow rate: 0.8 mL/min

Temperature: Ambient

Detector: UV 365 nm

Injection volume: 10 μ L (0.044 mg/mL)

Results and Discussion

All four aflatoxins were separated using an isocratic ternary mixture of water, methanol, and acetonitrile. A baseline separation was achieved in less than 5.5 min. Although UV detection was shown here, in some cases, lower levels of detection may be obtained for B₂ and G₂ using fluorescence ($\lambda_{\text{ex}} = 365$ nm, $\lambda_{\text{em}} = 455$ nm) detection. Mass spectroscopic detection has also been used [1].

References

1. R. Schuster, G. Marx, and M. Rothaupt, "Analysis of Mycotoxins by HPLC with Automated Spectroscopic Confirmation by Spectral Library", Agilent Technologies, publication 5091-8692E, www.agilent.com/chem
2. A. Gratzfeld-Heusgen, "HPLC Analysis of Aflatoxins in Pistachio Nuts", Agilent Technologies, publication 5966-0632E, www.agilent.com/chem
3. I. Ferreira, E. Mendes, and M. Oliveira, (2004) *J. Liquid Chromatog.*, **27** (2), 325–334.
4. E. Chiaavaro, C. Dall'Asta, G. Galaverna, A. Biancardi, and E. Gambarelli, (2001) *J. Chromatogr. A*, **937** (1–2), 31–40.

For More Information

For more information on our products and services, visit our Web site at www.agilent.com/chem.

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material. Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc. 2005

Printed in the USA
August 16, 2005
5989-3634EN

