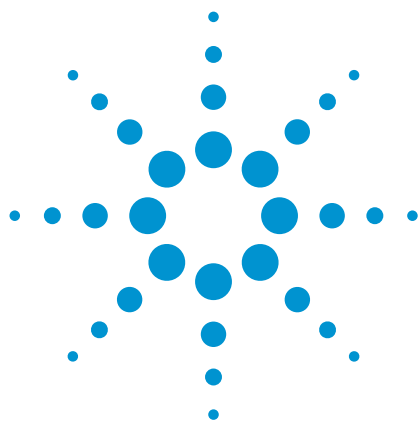




Separation of Mycotoxins with HPLC/UV

Application Note

Pharmaceutical

Author

Institute of Bioanalytical Sciences
Centre of Life Sciences
Anhalt University of Applied Sciences
Bernburg
Germany

Introduction

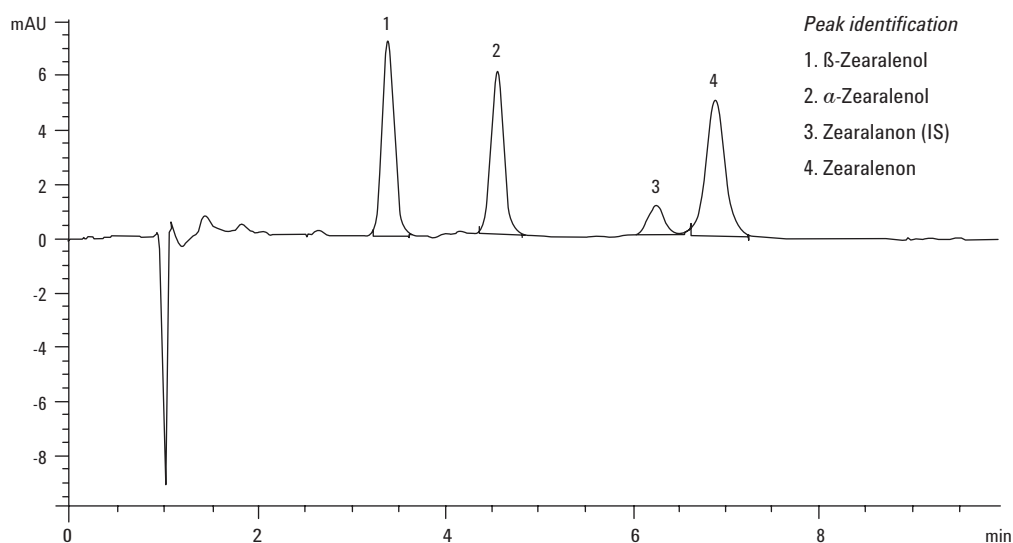
Using HPLC, the Agilent Pursuit XRs Ultra 2.8 Diphenyl column provides good separation of mycotoxins (β -Zearalenol, α -Zearalenol, Zearalanon [IS], and Zearalenon) with isocratic elution and a fast run time.



Agilent Technologies

Conditions

Column	Agilent Pursuit XRs Ultra 2.8 Diphenyl, 2.0 × 100 mm, 2.8 µm (p/n A7521100X020)
Sample size	5 µL
Sample concentration	1 µg/mL
Sample solvent	Acetonitrile
Mobile phases	A: 0.005 M ammonium acetate, 0.1% methanol, 0.02% formic acid, pH 3.6 B: methanol
Gradient	Isocratic A:B 40:60
Flow rate	0.4 mL/min
Temperature	40 °C
Pressure	180 bar
Detector	UV, 236 nm



Conclusion

The Agilent Pursuit XRs Ultra 2.8 Diphenyl column provided good separation with isocratic elution and fast analysis.

For More Information

These data represent typical results. For more information on our products and services, visit our Web site at www.agilent.com/chem.

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. This publication was originally published in 2010.

© Agilent Technologies, Inc., 2011
Published in USA, July 14, 2011
SI-01994



Agilent Technologies