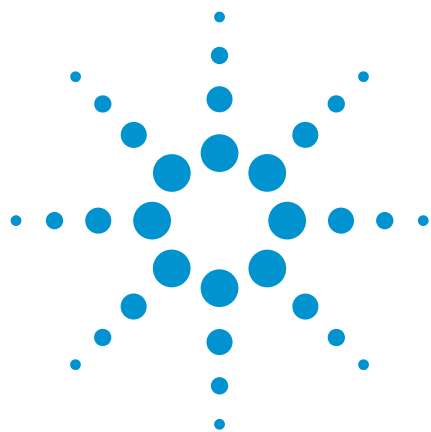




Analysis of nitrofuran metabolites by positive ion electrospray LC/MS/MS

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Nitrofuran antibiotics are veterinary drugs. The use of these in food production is already banned in the United States, Australia, Canada, Japan, Singapore, Bangladesh, and the European Union because of a possible increased cancer-risk through long-term consumption.

Trace amounts of nitrofurans contamination has been found in warm-water prawns, shrimp, and chicken. The contamination may occur from deliberate, direct misuse of the drugs as well as from contaminated feed.

The analyses of nitrofurans are based on the detection of the metabolites of the parent drugs. Since the parent drugs are very rapidly metabolized and the tissue-bound metabolites are detectable for several weeks after administration, the metabolites are a much better marker for the detection of the abuse of these antibiotics.

The metabolites of furazolidone, furaltadone, nitrofurantoin and nitrofurazone are 3-amino-2-oxazolidinone (AOZ), 5-methylmorpholino-3-amino-2-oxazolidinone (AMOZ), 1-aminohydantoin (AH) and semicarbazide (SC), respectively. Nitrobenzaldehyde, the most commonly used derivatizing reagent for these metabolites, forms a nitrophenyl derivative for LC/MS analyses.



Agilent Technologies

Instrumentation

- Agilent ProStar 430 AutoSampler
- Agilent ProStar 210 Isocratic Solvent Delivery Modules
- Agilent 1200L LC/MS equipped with ESI source

Materials and Reagents

- A 5 pg/ μ L standard mixture of 2-nitrobenzylaldehyde derivative of AOZ, AMOZ, AH, and SC was kindly provided by Thai Unique, Thailand.
- All other chemicals are reagent grade or HPLC grade.

Conditions

Column	: Agilent Polaris C18-A, 2 x 50 mm, 3 μ m (Agilent Part Number: A2001050X020)			
Mixer	: 250 μ L static mixer			
Solvent A	: water			
Solvent B	: methanol			
LC Program	: Time (min:sec)	% A	% B	Flow (mL/min)
	0:00	90	10	0.2
	1:00	5	95	
	2:00	5	95	
	2:21	90	10	
	6:00	90	10	0.2
Injection Volume	: 40 μ L			
Sample Solvent	: 10 mM ammonium acetate			

MS Parameters

Collision Gas	: 2.1 mTorr Argon
API Drying Gas	: 20 psi at 300 °C
API Nebulizing Gas	: 50 psi
Scan Time	: 1 s
SIM Width	: 0.7 amu total
Needle	: -4400V
Shield	: -300V
Capillary	: -40V
Detector	: 2000V

Results

The LC method used a 10 minute run cycle. It is clear from the Multiple Reaction Monitoring (MRM) data produced that the 1200L LC/MS system can qualitatively and quantitatively detect nitrofurans to the ultra trace levels required.

10 mM ammonium acetate was used in the sample solution to reduce formation of sodium adducts and to enhance the detection limit for the product ions.

The Limit of Detection (LOD) for SC, AOZ, AH, and AMOZ is about 60 fg/ μ L*, 10 fg/ μ L, 20 fg/ μ L, and 10 fg/ μ L in solution, respectively. The linearity of the detector response was excellent (Figure 1). The R^2 for 209>166, 236>134, 249>134 and 335>291 are 0.999, 0.998, 0.999 and 0.999, respectively. Typical chromatograms obtained are given in Figure 2.

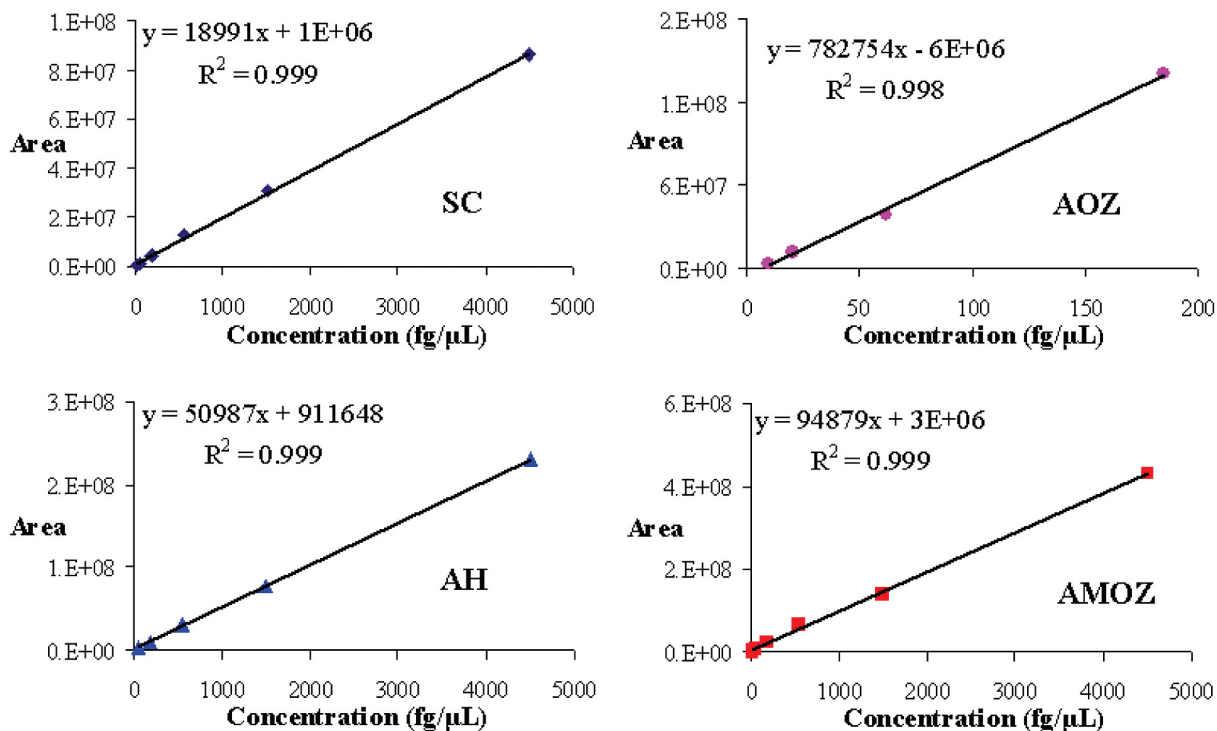


Figure 1. Standard Calibration Curves

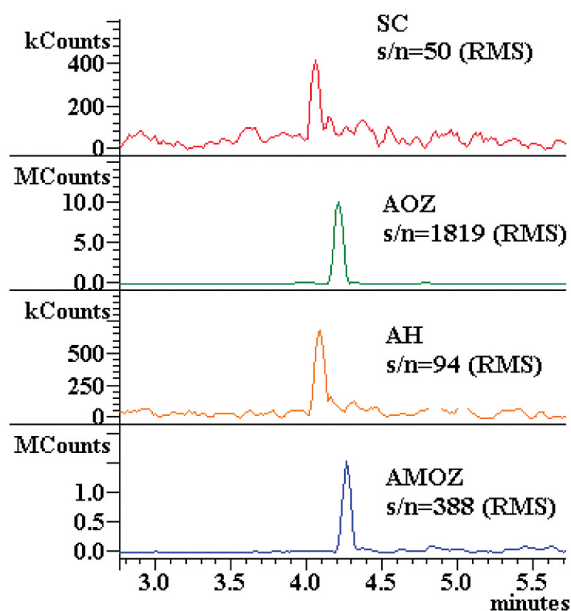


Figure 2. MRM Chromatograms

Conclusion

The Agilent 1200L LC/MS/MS system demonstrates excellent performance for the analyses of nitrofurans metabolites and is able to assist in the detection of potentially harmful substances in food imports/exports to improve food safety throughout the world.

* pg/μL=ppb and fg/μL=ppt

www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Printed in the USA

31 October, 2011

First published prior to 11 May, 2010

A02052



Agilent Technologies