

Antibiotics

Potential LC/MS method for analysis of antibiotics

Application Note

BioPharma

Authors

Agilent Technologies, Inc.

Introduction

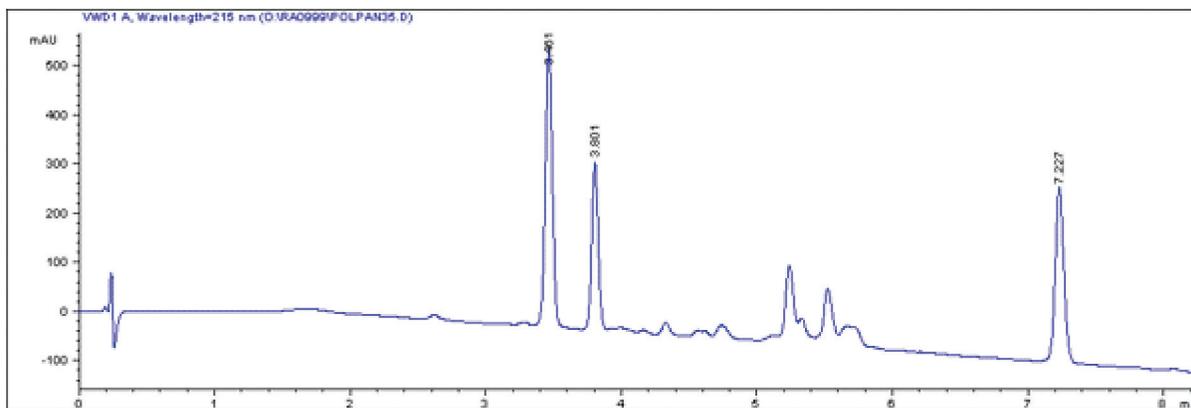
HPLC with an Agilent Polaris C18-A column separates three antibiotics in seven minutes.



Agilent Technologies

Conditions

Column : Agilent Polaris C18-A, 2.0 x 50 mm, 5 µm
Part no. : A2000050X020
Eluent : A: 0.1% formic acid
 B: MeCN+0.1% formic acid
Gradient : 0 to 50%B in 10 min
Flow : 0.8 mL/min
Detection : UV @ 215nm
Sample : 1. cephaloridine
 2. penicillin G
 3. penicillin V



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

A01930



Agilent Technologies