



LC/MS performance test mix

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

An LC/MS test mix is used to demonstrate the HPLC performance of Agilent Pursuit C8 5 μm material in a ChromSep cartridge.



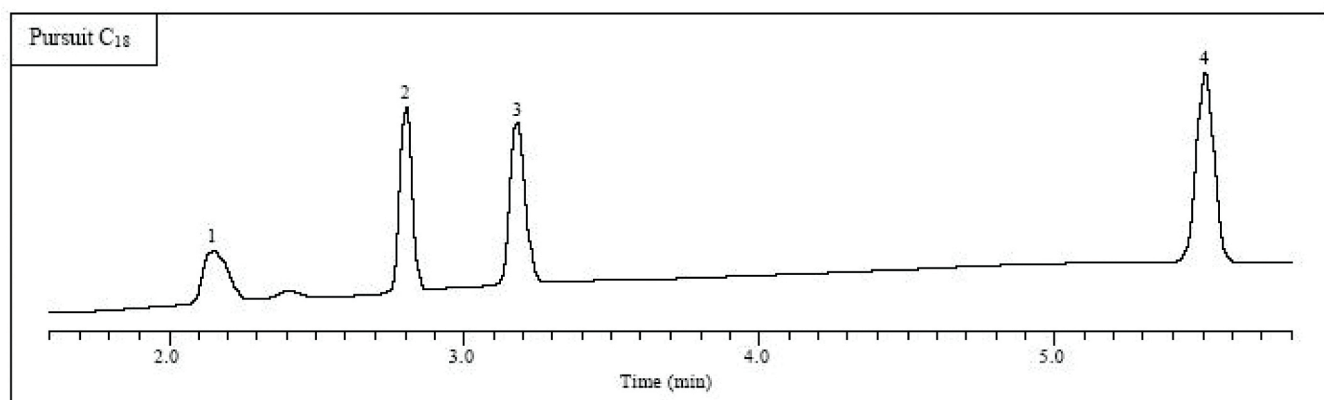
Agilent Technologies

Conditions

Technique : HPLC
Column : Agilent Pursuit C8, 4.6 x 150 mm, 5 µm
Part no. : Conventional A3000030X020
ChromSep Cartridge System A30000200C020
Temperature : Ambient
Eluent : A: 0.05% Formic acid in water
B: 0.05% Formic acid in MeCN
Gradient : 5 - 90% B in 3 minutes
Flow rate : 0.6 mL/min
Detection : UV @ 220 nm

Peak identification

1. aspartame
2. cortisone
3. reserpine
4. dioctyl phthalate



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

A02318



Agilent Technologies