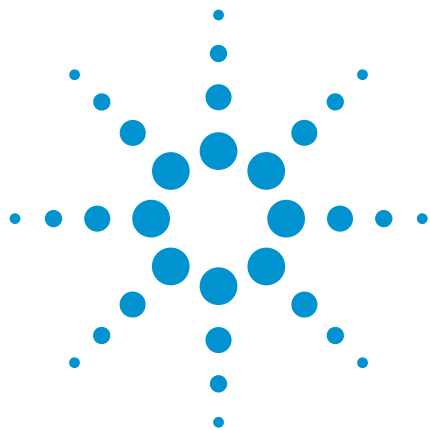




Solvents

Fast analysis of residual solvents via headspace analysis

Application Note

BioPharma

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with a 0.15 mm id Agilent CP-Sil 5 CB column separates residual solvents in less than seven minutes.



Agilent Technologies

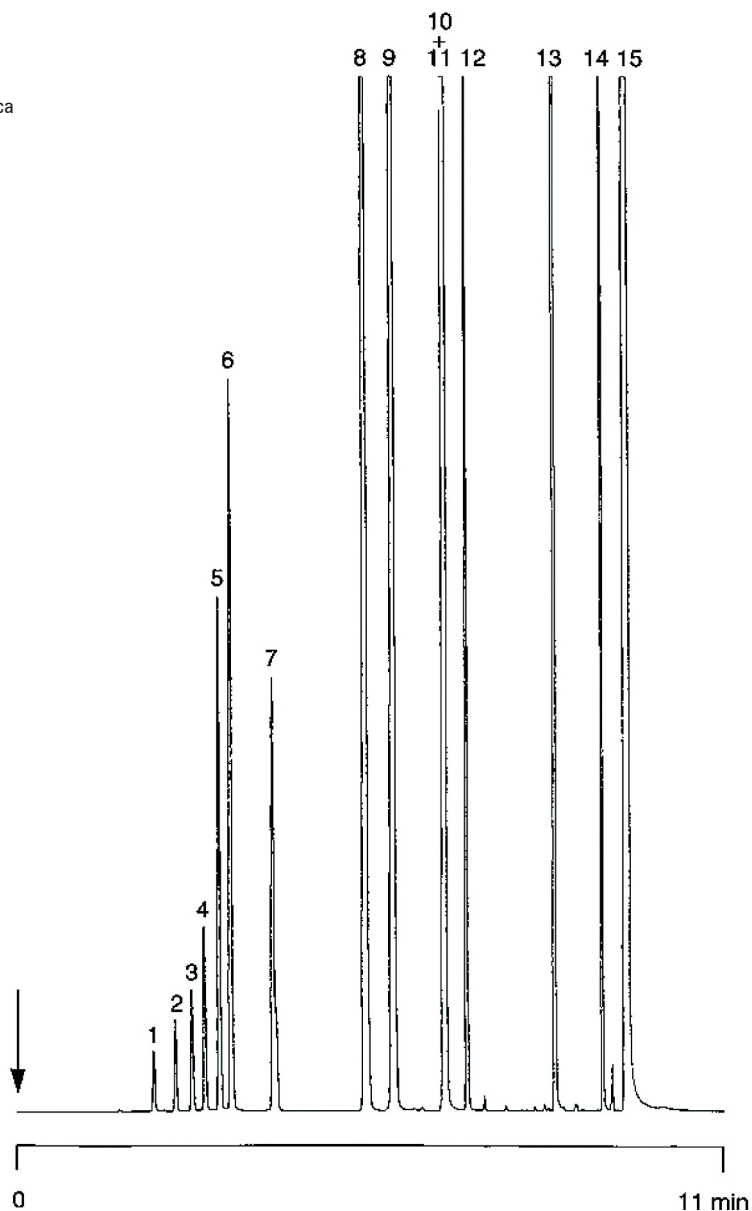
Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5 CB, 0.15 mm x 25 m fused silica
WCOT (df = 1.2 µm) (Part no. CP7693)
Temperature : 40 °C (6 min) → 160 °C, 50 °C/min
Carrier Gas : He, 330 kPa (3.3 bar, 46 psi)
Injector : Split,
T = 250 °C
Detector : FID
T = 250 °C
Sample Size : 1 mL headspace
Concentration Range : 0.01 -1%

Courtesy : I. Dijkstra, Solvay Pharmaceuticals,
Weesp, The Netherlands

Peak identification

1. methanol
2. methyl formate
3. ethanol
4. acetonitrile
5. acetone
6. isopropanol
7. dichloromethane + methylacetate
8. MTBE
9. methyl ethyl ketone
10. ethylacetate
11. hexane
12. tetrahydrofuran
13. heptane
14. toluene
15. DMSO



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

A01502



Agilent Technologies