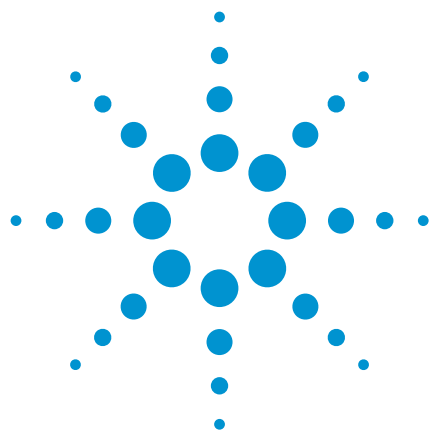




Residual solvents

Analysis of residual solvents on video tape (Brand 2)

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 5 CB column separates eight residual solvents in a sample of video tape in 70 minutes.



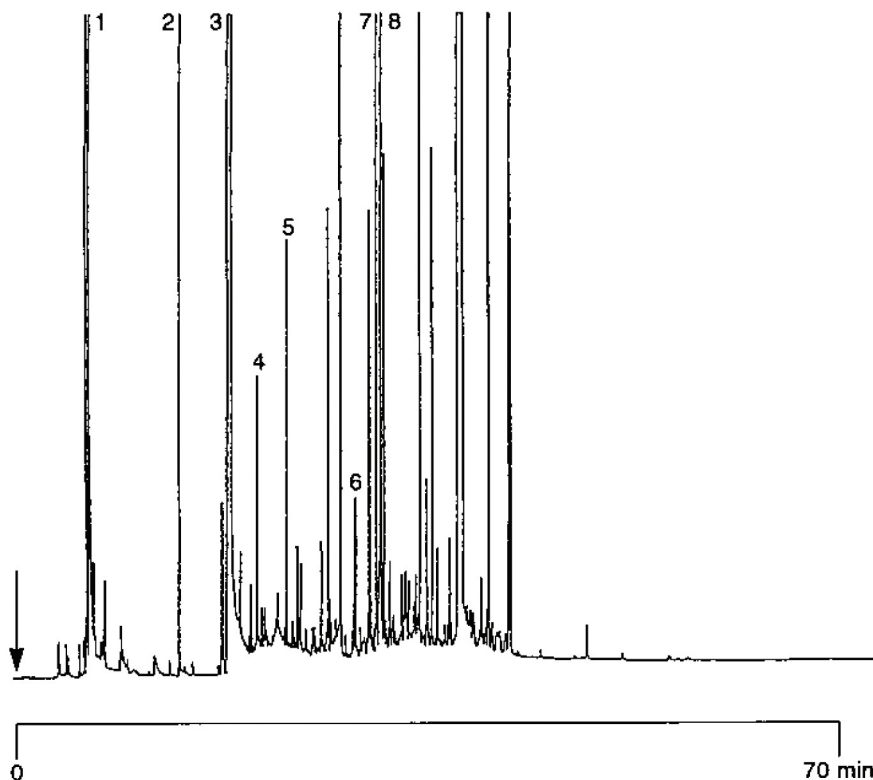
Agilent Technologies

Conditions

Technique : GC-TCT / Dynamic Headspace
Column : Agilent CP-Sil 5 CB, 0.25 mm x 50 m fused silica
WCOT CP-Sil 5 CB (df = 0.4 μ m) (Part no. CP7719)
Temperature : 40°C (10 min) $\rightarrow$ 280°C, 10°C/min
Injector : TCT
Cold trap : CP-Sil 8 CB
0.53 mm; df = 5 μ m
Cryofocussing : -130 °C
Precool time : 3 min
Desorb time : 5 min
Desorb flow : 9.5 mL/min
Desorption oven temp. : 150 °C
Injection temp. : 210 °C
Injection time : 1 min
GC-injection block temp. : 250 °C
Detector : FID
T = 250 °C
Sample Size : 2 mg (2 mm x 40 mm)
Sample Temperature : 150 °C
Concentration Range : ppb/ppm
Courtesy : GL Sciences, Japan

Peak identification

1. 2-butanone
2. toluene
3. cyclohexanone
4. styrene
5. nitrobenzene
6. 1-chloronaphthalene
7. butylated hydroxytoluene (BHT)
8. 1,6-dioxacyclododecane-7,12-dione



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

A01048



Agilent Technologies