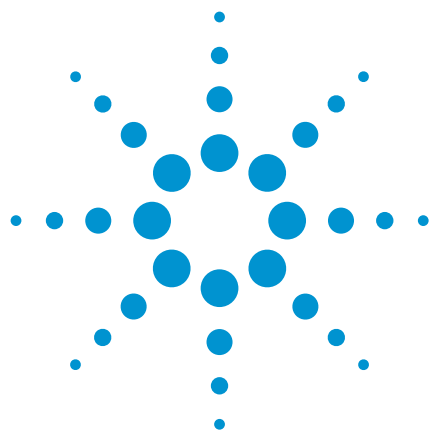




Phenols

Analysis of ethyl derivatives of alkylphenols

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

GC/MS of ten ethyl derivatives of alkylphenols takes less than 18 minutes with an Agilent CP-Sil 8 CB Low Bleed/MS column.



Agilent Technologies

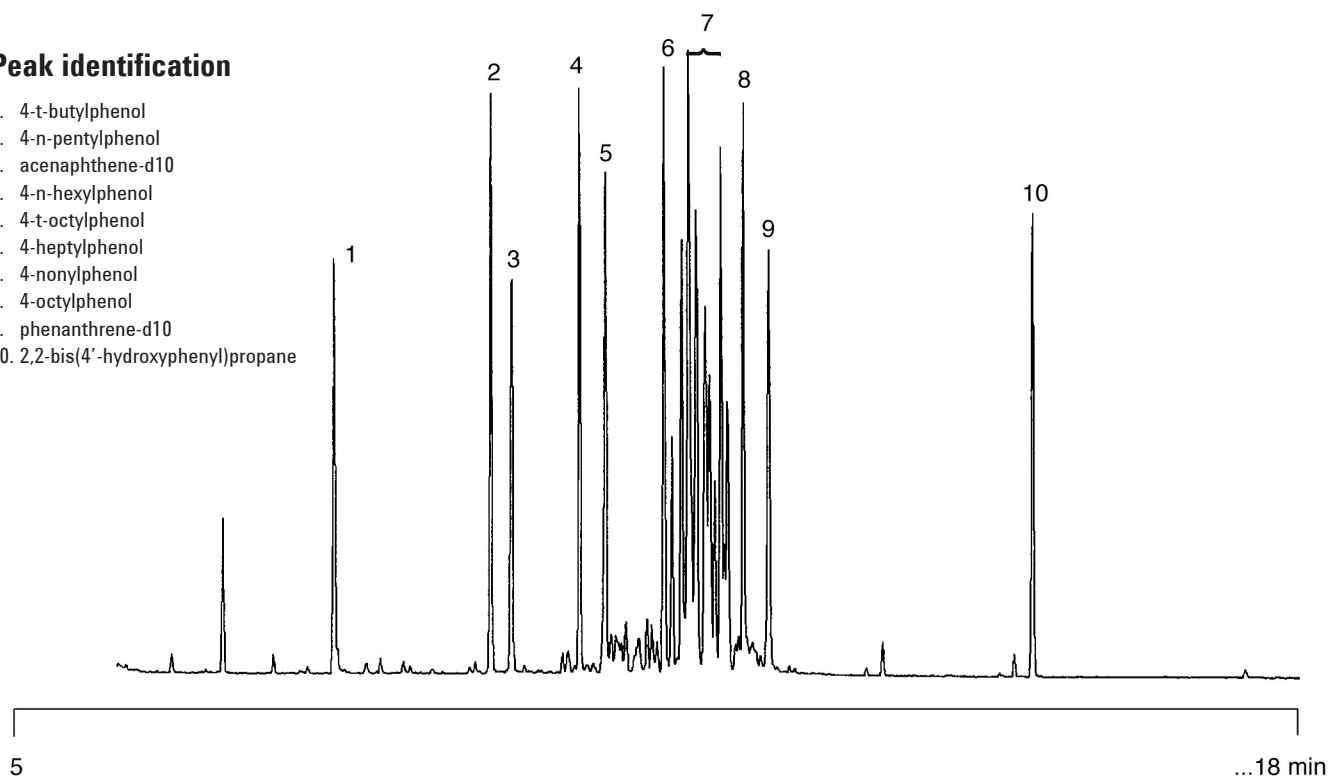
Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB, 0.25 mm x 30 m fused silica
WCOT (df = 0.25 μ m) (Part no. CP5860)
Temperature : 50 °C (1 min) $\rightarrow$ 300 °C, 20 °C/min
Carrier Gas : He, 75 kPa (0.75 bar, 11 psi)
Injector : Splitless, 1min
T = 250 °C
Detector : MSD, TIC
T = 230 °C
Sample Size : 1 μ L
Solvent Sample : hexane

Courtesy : S. Hishimoto, GL Sciences, Japan

Peak identification

1. 4-t-butylphenol
2. 4-n-pentylphenol
3. acenaphthene-d10
4. 4-n-hexylphenol
5. 4-t-octylphenol
6. 4-heptylphenol
7. 4-nonylphenol
8. 4-octylphenol
9. phenanthrene-d10
10. 2,2-bis(4'-hydroxyphenyl)propane



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

A01654



Agilent Technologies