



Phenols

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-FFAP column separates eight phenols in six minutes.



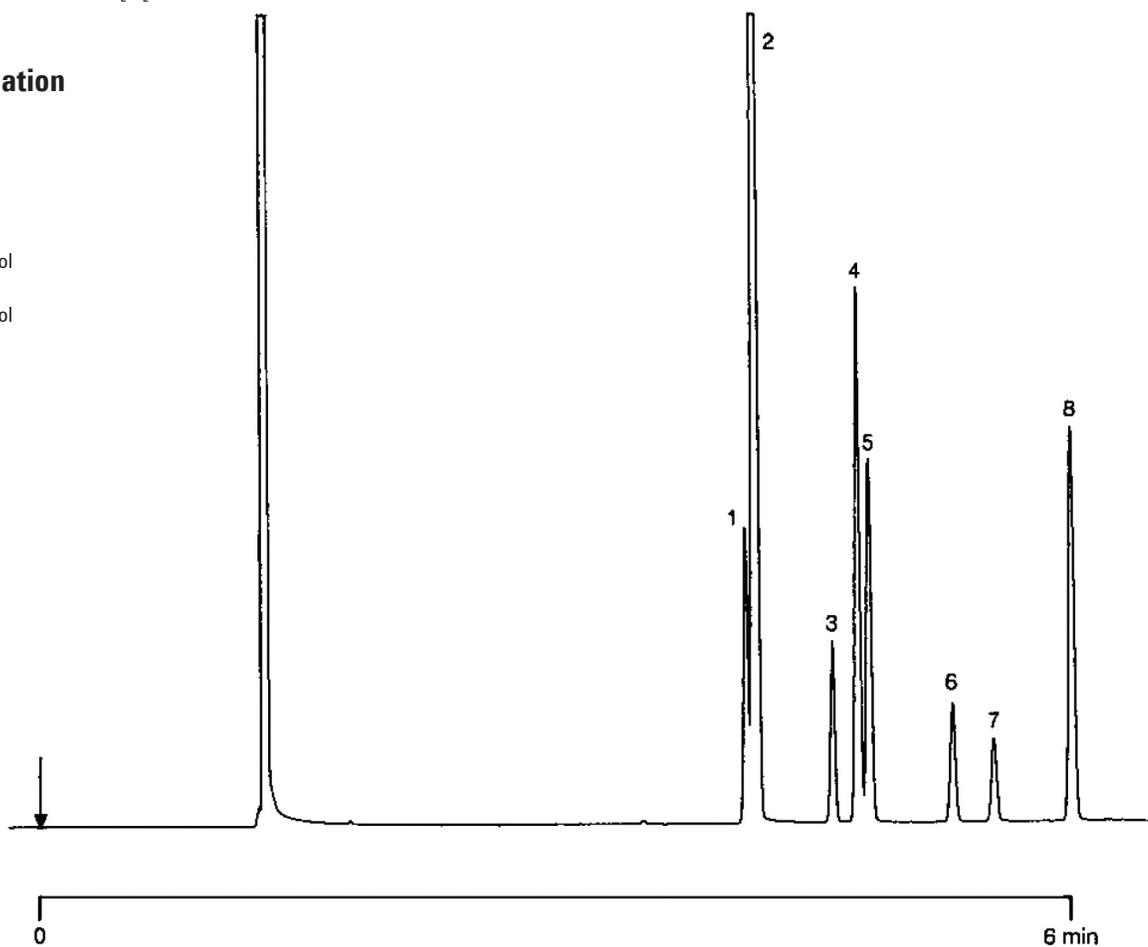
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-FFAP, 0.32 mm x 25 m fused silica WCOT
FFAP-CB (df = 0.3 μ m) (Part no. CP7485)
Temperature : 160 °C + 7½ °C/min → 220 °C
Carrier Gas : H₂, 30 kPa (0.3 bar, 4.5 psi)
Injector : Split
Detector : FID
Sample Size : 0.1% in CH₂Cl₂

Peak identification

1. o-cresol
2. phenol
3. o-ethylphenol
4. p-cresol
5. m-cresol
6. 2,3-dimethylphenol
7. p-ethylphenol
8. 2,4-dimethylphenol



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

A00361



Agilent Technologies