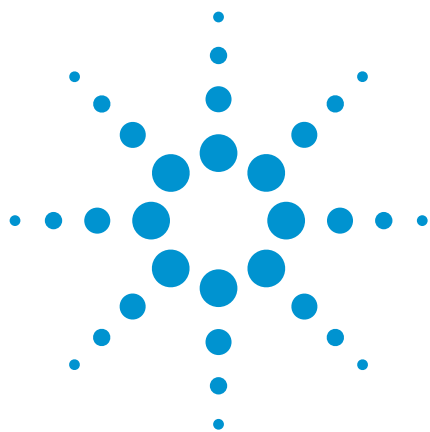




Sulfur compounds

Analysis of aromatic sulfur compounds

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 5 CB-MS column separates three aromatic sulfur compounds in 17 minutes.



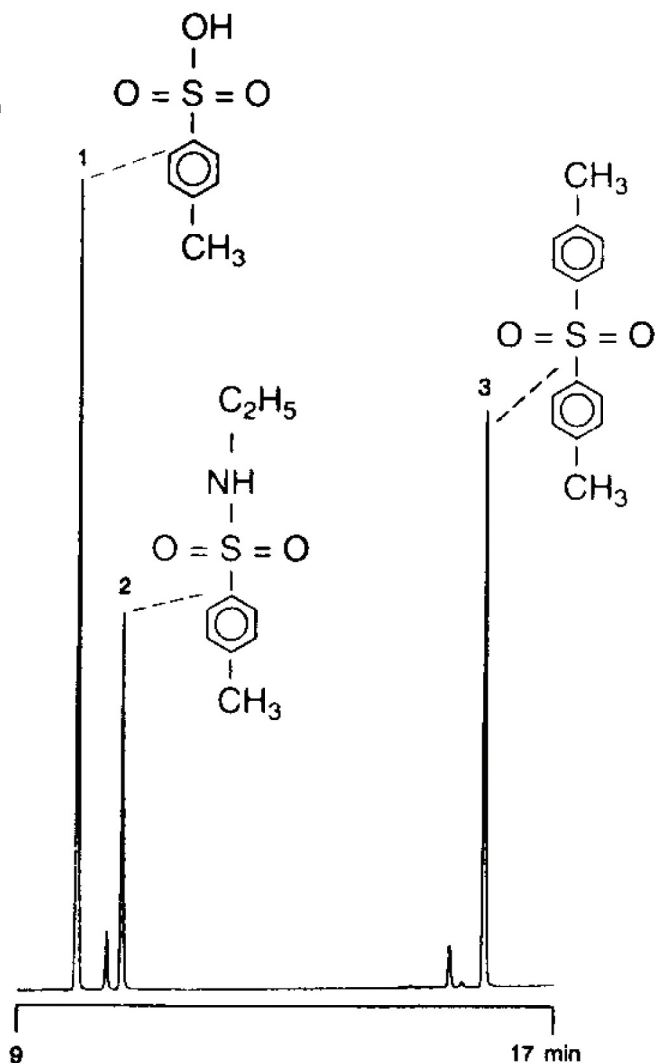
Agilent Technologies

Conditions

Technique	: GC-capillary
Column	: Agilent CP-Sil 5 CB-MS, 0.32 mm x 25 m fused silica WCOT CP-Sil 5 CB-MS (df = 1.2 µm) (Part no. CP7852)
Temperature	: 100°C → 320°C, 10°C/min; 320°C (5 min)
Carrier Gas	: He, 70 cm/s
Injector	: Splitter, 100 mL/min T = 250 °C
Detector	: FID T = 330 °C
Sample Size	: 0.5 µL
Concentration Range	: 1 mg/mL
Solvent Sample	: toluene
Courtesy	: K. Schutjes, Akzo Nobel, Central Research B.V., Deventer, The Netherlands

Peak identification

1. p-toluenesulfonic acid
2. N-ethyl-p-toluenesulfonamide
3. p,p'-ditolylsulfone



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

A00986



Agilent Technologies