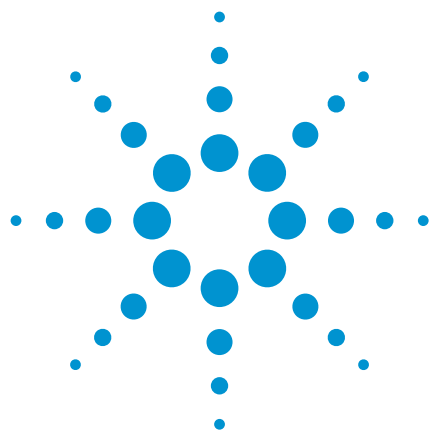




## **Sulfur compounds**

# **Analysis of low-boiling sulfur-containing compounds**

## **Application Note**

Energy & Fuels

### **Authors**

Agilent Technologies, Inc.

### **Introduction**

Gas chromatography with an Agilent CP-Sil 8 CB column separates 12 low-boiling, sulfur-containing compounds in 17 minutes.



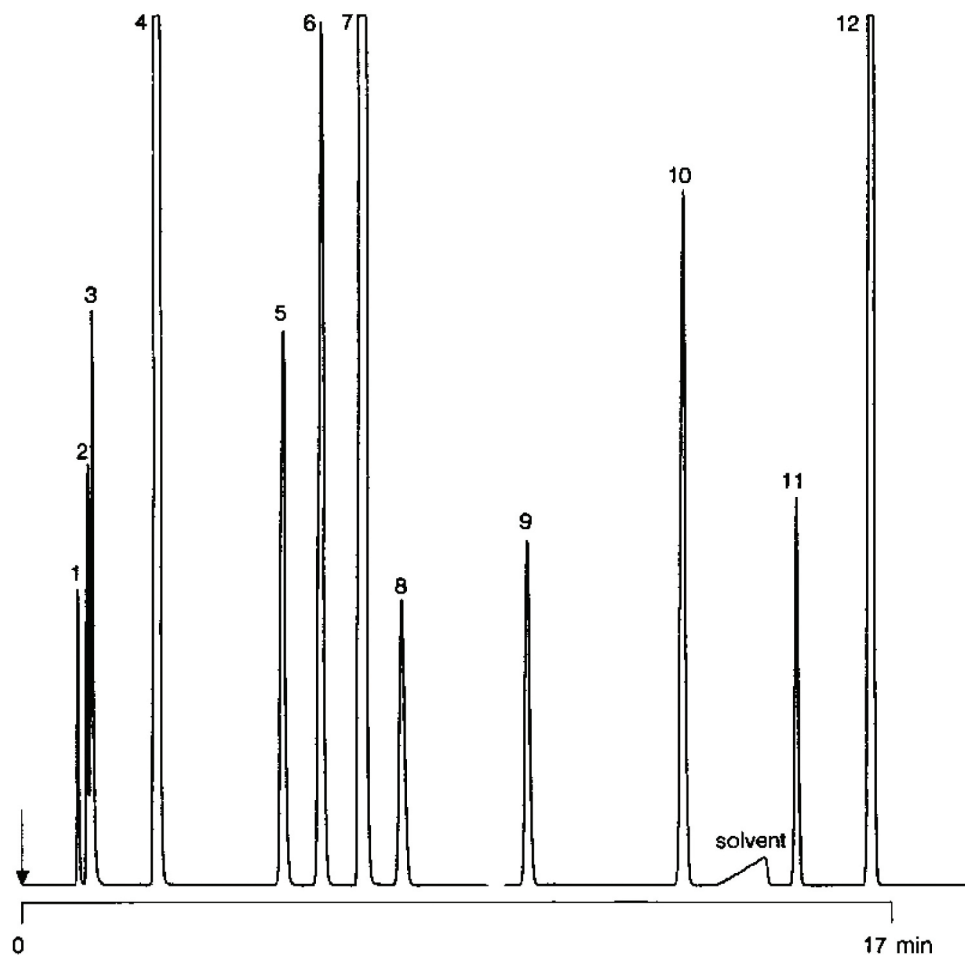
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-Sil 8 CB, 0.53 mm x 25 m fused silica  
WCOT CP-Sil 8 CB (df = 5.0  $\mu$ m) (Part no. CP7656)  
Temperature : 30 °C  $\rightarrow$  110 °C (5 °C/min)  
Carrier Gas : H<sub>2</sub> (3.8 mL/min)  
Injector : T = 200 °C  
Detector : FPD  
T = 200 °C

## Peak identification

1. hydrogen sulfide
2. carbonyl sulfide
3. sulfur dioxide
4. methanethiol
5. methyl sulfide
6. ethanethiol
7. carbon disulfide
8. 2-propanethiol
9. 1-propanethiol
10. thiophene
11. 1-butanethiol
12. methyl disulfide



[www.agilent.com/chem](http://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

A00419



**Agilent Technologies**