



Sulfur compounds

Analysis of sulfur gases

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

GC/MS with an Agilent CP-SilicaPLOT column separates seven sulfur compounds in 25 minutes.



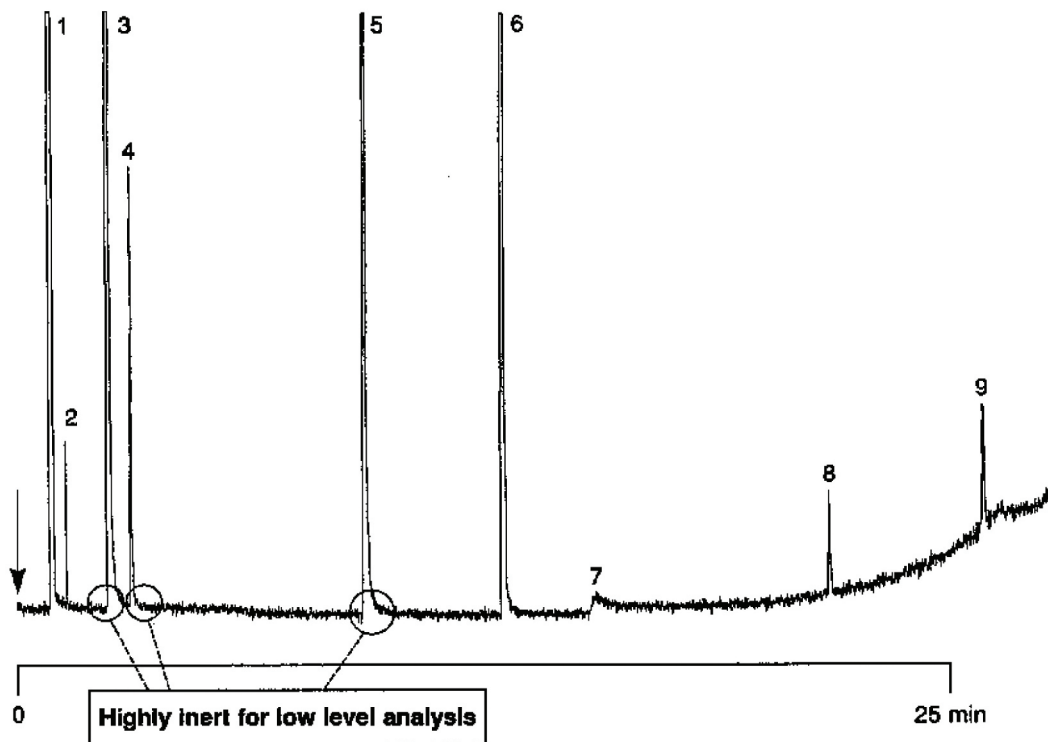
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-SilicaPLOT, 0.32 mm x 30 m fused silica
PLOT CP-SilicaPLOT (df = 4 μ m) (Part no. CP8567)
Temperature : 40 °C (5 min) $\rightarrow$ 200 °C, 10 °C/min
Carrier Gas : He, 30 kPa (0.3 bar, 4.3 psi)
Injector : T = 200 °C
Detector : MS-TIC
Sample Size : 100 μ L
Concentration Range : ca. 1000 ppm in air

Peak identification

1. air
2. carbon dioxide
3. carbonyl sulfide
4. hydrogen sulfide
5. sulfur dioxide
6. methyl sulfide
7. water
8. dimethyl disulfide
9. dimethyl trisulfide



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

A01257



Agilent Technologies