

Sulfur compounds

Analysis of sulfur compounds in wine

Application Note

Food Testing & Agriculture

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates 13 sulfur compounds in a wine in 33 minutes.



Agilent Technologies

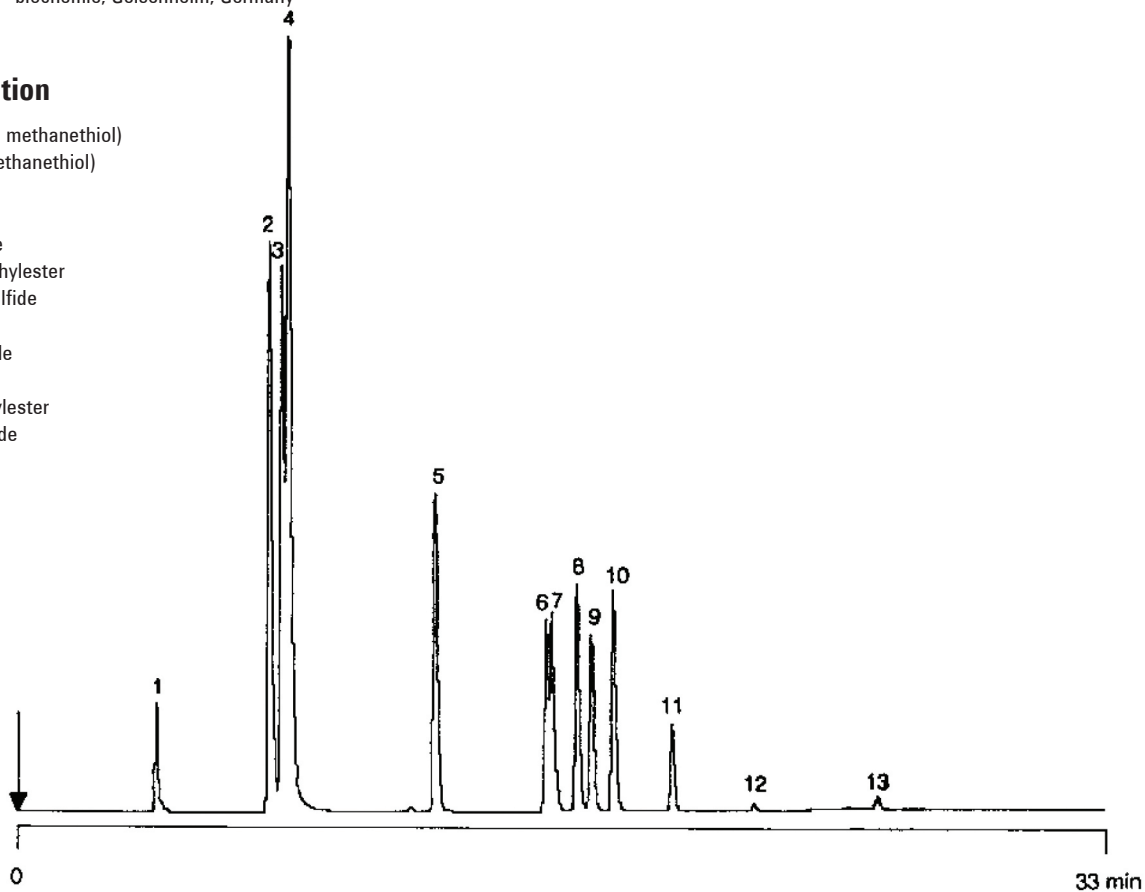
Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q, 0.32 mm x 10 m fused silica
PoraPLOT Q (df = 10 µm) (Part no. CP7550)
Temperature : 120 °C (2 min) 5 °C/min → 250 °C (15 min)
Carrier Gas : N₂, 30 kPa (0.3 bar, 4.2 psi), 8 cm/s
Injector Method : Purge and Trap
Detector : MS

Courtesy : Mrs D. Rauhut, Institut für mikrobiologie und
biochemie, Geisenheim, Germany

Peak identification

1. methyl mercaptan (methanethiol)
2. ethyl mercaptane (ethanethiol)
3. dimethyl sulfide
4. carbon disulfide
5. ethyl methyl sulfide
6. thioacetic acid methylester
7. isopropyl methyl sulfide
8. diethyl sulfide
9. methyl propyl sulfide
10. dimethyldisulfide
11. thioacetic acid ethylester
12. ethyl methyl disulfide
13. diethyl disulfide



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

A00444



Agilent Technologies