



Alcohols and esters

Analysis of 'fire-water'

Application Note

Food Testing & Agriculture

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 57 CB for Glycols and Alcohols column separates 21 alcohols and esters in 'fire water' in 21 minutes.



Agilent Technologies

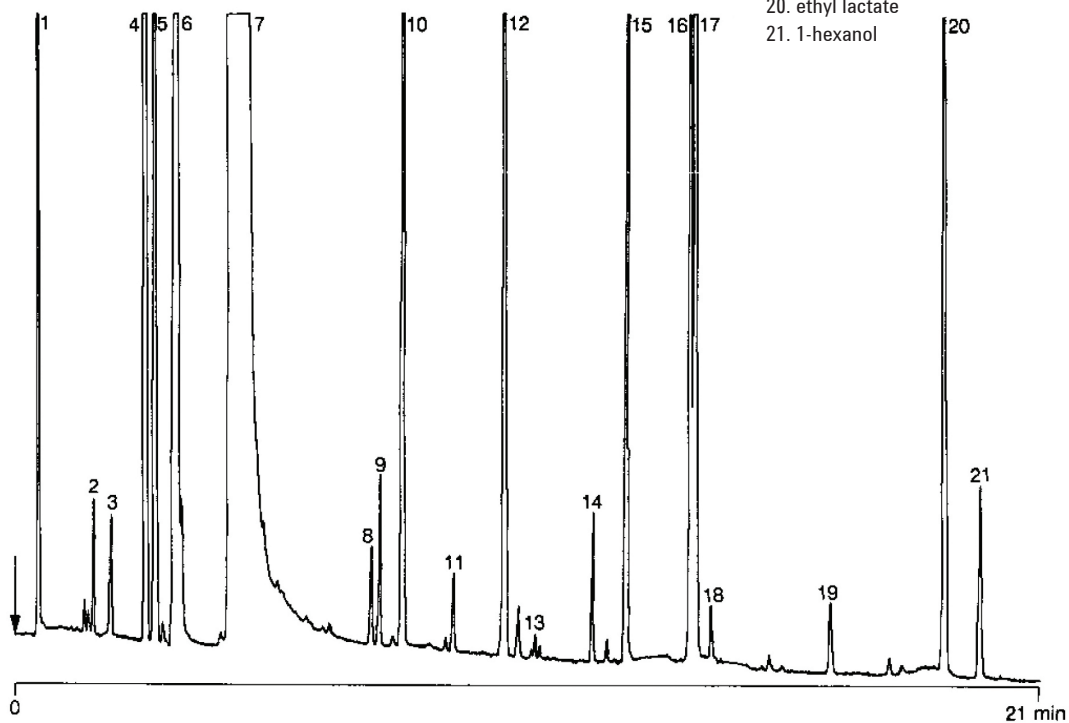
Conditions

Technique : GC-capillary
Column : Agilent CP-Wax 57 CB, 0.25 mm x 50 m fused silica
WCOT CP-Wax 57 CB (df = 0.2 µm)
(Part no. CP97723)
Temperature : 40 °C → 100 °C, 7 °C/min; 100 °C (4 min)
Carrier Gas : He, 140 kPa (1.4 bar, 20 psi)
Injector : Split, 1:50
T = 210 °C
Detector : FID
T = 210 °C
Sample Size : 1 µL

Courtesy : Maria de Jesus Caetano Tavares,
Wine and Vineyard Institute,
Sacavém, Portugal

Peak identification

1. acetaldehyde
2. methyl acetate
3. acrolein (2-propenal)
4. ethyl acetate
5. acetal
6. methanol
7. ethanol
8. ethyl butyrate
9. 2-butanol
10. propanol
11. hexanal
12. isobutanol
13. allyl alcohol
14. n-butanol
15. 4-methyl-2-pentanol (I.S.)
16. 2-methyl-1-butanol
17. 3-methyl-1-butanol
18. 2-hexanol
19. acetoin (3-hydroxy-2-butanone)
20. ethyl lactate
21. 1-hexanol



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

A00619



Agilent Technologies