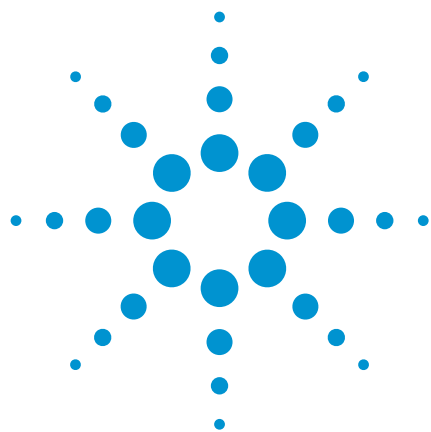




Alcohols and esters

Analysis of volatile compounds in beer

Application Note

Food Testing & Agriculture

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Carbowax 400 for Volatiles in Alcohol column separates 11 alcohols and esters in beer in 12 minutes.



Agilent Technologies

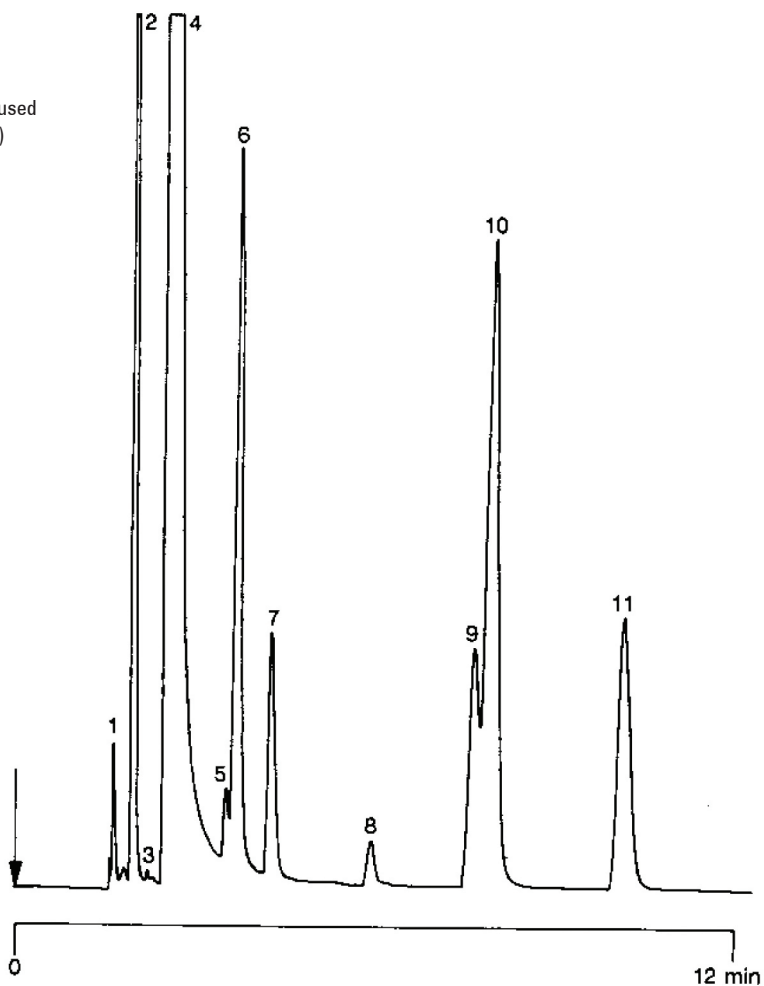
Conditions

Technique : GC-capillary
Column : Agilent CP-Carbowax 400, 0.32mm x 50 m fused silica WCOT CP-Carbowax 400 (df = 0.2 µm) (Part no. CP7527)
Temperature : 60 °C
Carrier Gas : H₂, 70 kPa (0.7 bar, 10 psi)
Injector : Splitter, 40 mL/min (1:20)
T = 125 °C
Detector : FID, 2^d
T = 145 °C
Sample Size : 1 mL headspace

Courtesy : Mr. H. Klein,
Österreichische Brau-AG,
Qualitätssicherung + Entwicklung,
Linz, Österreich

Peak identification

1. acetaldehyde
2. ethyl acetate
3. methanol
4. ethanol
5. n-propanol
6. isoamyl acetate
7. 2-methylpropanol
8. hexyl acetate
9. 2-methyl-1-butanol
10. 3-methyl-1-butanol
11. n-pentanol (I.S.)



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

A00620



Agilent Technologies