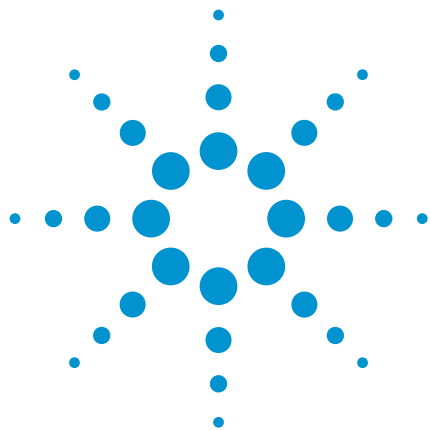




Alcohols, C₁ – C₄

Analysis of impurities in ethanol

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 52 CB column separates three impurities in ethanol in six minutes.



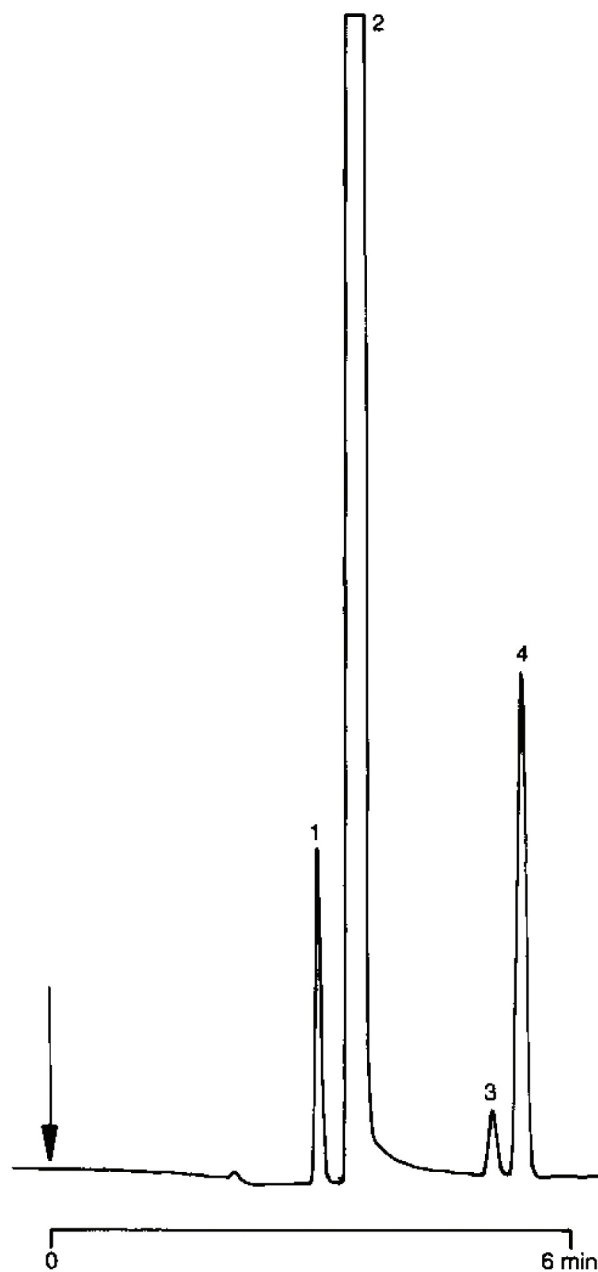
Agilent Technologies

Conditions

Technique : GC-wide-bore
Column : Agilent CP-Wax 52 CB, 0.53 mm x 50 m fused silica
WCOT CP-Wax 52 CB (df = 2 μ m) (Part no. CP7658)
Temperature : 80 °C
Carrier Gas : N₂, 70 kPa (0.7 bar, 10 psi)
Injector : Split, 1:100
T = 150 °C
Detector : FID
T = 250 °C
Sample Size : 0.5 μ L
Concentration Range : 1%
Solvent Sample : heptane

Peak identification

	Concentration range:
1. methanol	$\pm 0.7\%$ v/v
2. ethanol	$\pm 98.0\%$ v/v
3. 2-butanol	$\pm 0.1\%$ v/v
4. 1-propanol	$\pm 1.0\%$ v/v



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

A00900



Agilent Technologies