



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 57 CB for Glycols and Alcohols column separates four impurities in ethanol in 19 minutes.



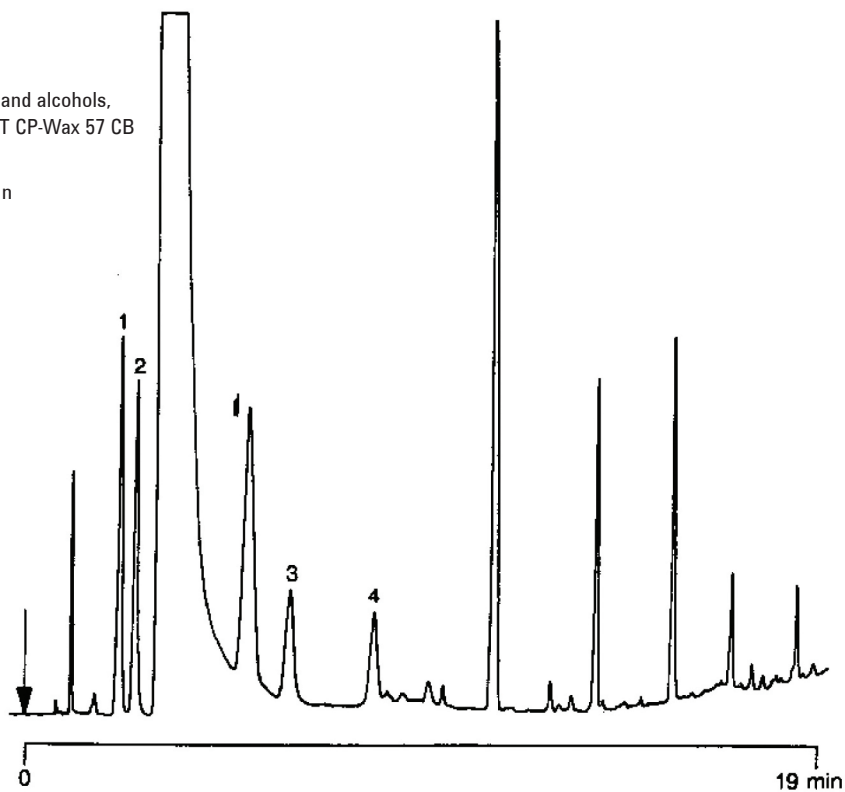
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Wax 57 CB for glycols and alcohols,
0.53 mm x 25 m fused silica WCOT CP-Wax 57 CB
(df = 0.5 μ m) (Part no. CP7617)
Temperature : 50 °C (6 min) $\rightarrow$ 180 °C, 10 °C/min
Carrier Gas : H₂, 10 kPa (0.1 bar, 1 .4 psi)
Injector : Splitter, 10 mL/min
T = 250 °C
Detector : FID, 2 x 10⁻¹¹ A/s
T = 250 °C
Concentration Range : 0.01 - 1.0%

Peak identification

1. ethyl acetate
2. methanol
3. n-propanol
4. isobutanol



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

A00689



Agilent Technologies