



Alcohols

Analysis of butanols in wine

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 butanols in wine in seven minutes.



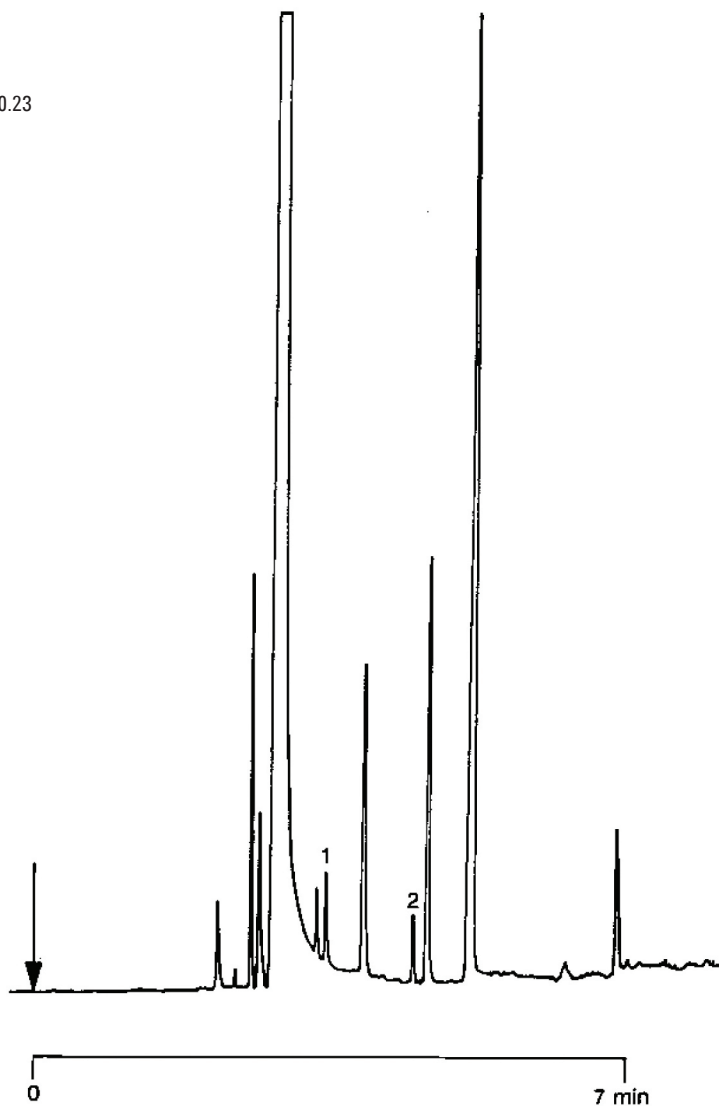
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Wax 57 CB for glycols and alcohols, 0.23 mm x 25 m fused silica WCOT CP-Wax 57 CB (df = 0.2 µm) (Part no. CP7615)
Temperature : 60 °C → 200 °C, 10 °C/min
Carrier Gas : He, 100 kPa (1.0 bar, 14 psi)
Injector : Splitter, 100 mL/min
T = 250 °C
Detector : FID, 16 x 10⁻¹² Afs
T = 250 °C
Sample Size : 1.0 µL

Peak identification

1. 2-butanol
2. 1-butanol



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

A00693



Agilent Technologies