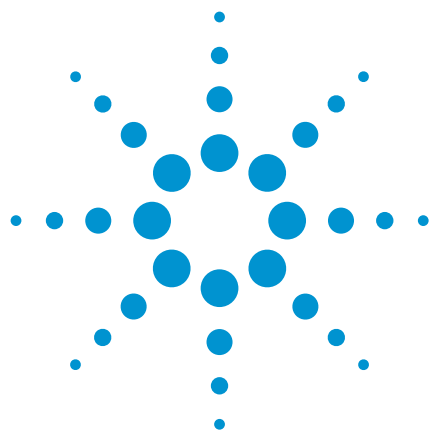




Halohydrins

Separation of positional isomers of chloropropanol

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 13 CB column separates positional isomers of chloropropanol in seven minutes.



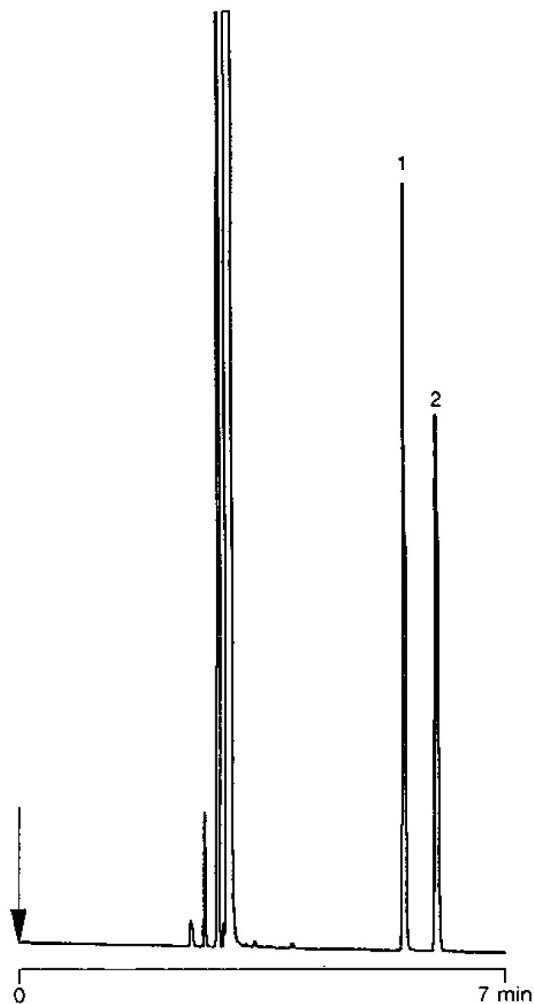
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 13 CB, 0.32 mm x 50 m fused silica
WCOT CP-Sil 13 CB (df = 1.2 µm) (Part no. CP7947)
Temperature : 80 °C
Carrier Gas : H₂, 60 kPa (0.6 bar, 8.6 psi)
Injector : Splitter, 100 mL/min
T = 275 °C
Detector : FID
T = 300 °C
Sample Size : 0.3 µL
Concentration Range : 0.1%
Solvent Sample : diethyl ether

Peak identification

1. 1-chloro-2-propanol
2. 2-chloro-1-propanol



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

A00771



Agilent Technologies