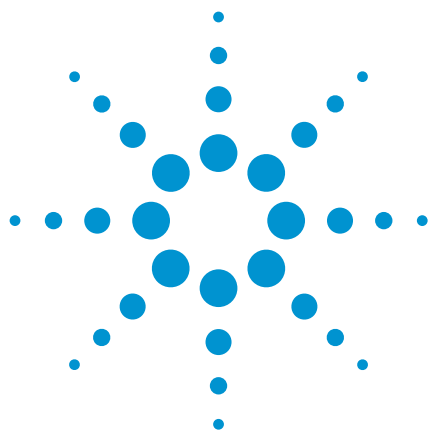




Oxygenates

Analysis of impurities in propylene oxide

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Propox column separates eight impurities in propylene oxide in 11 minutes



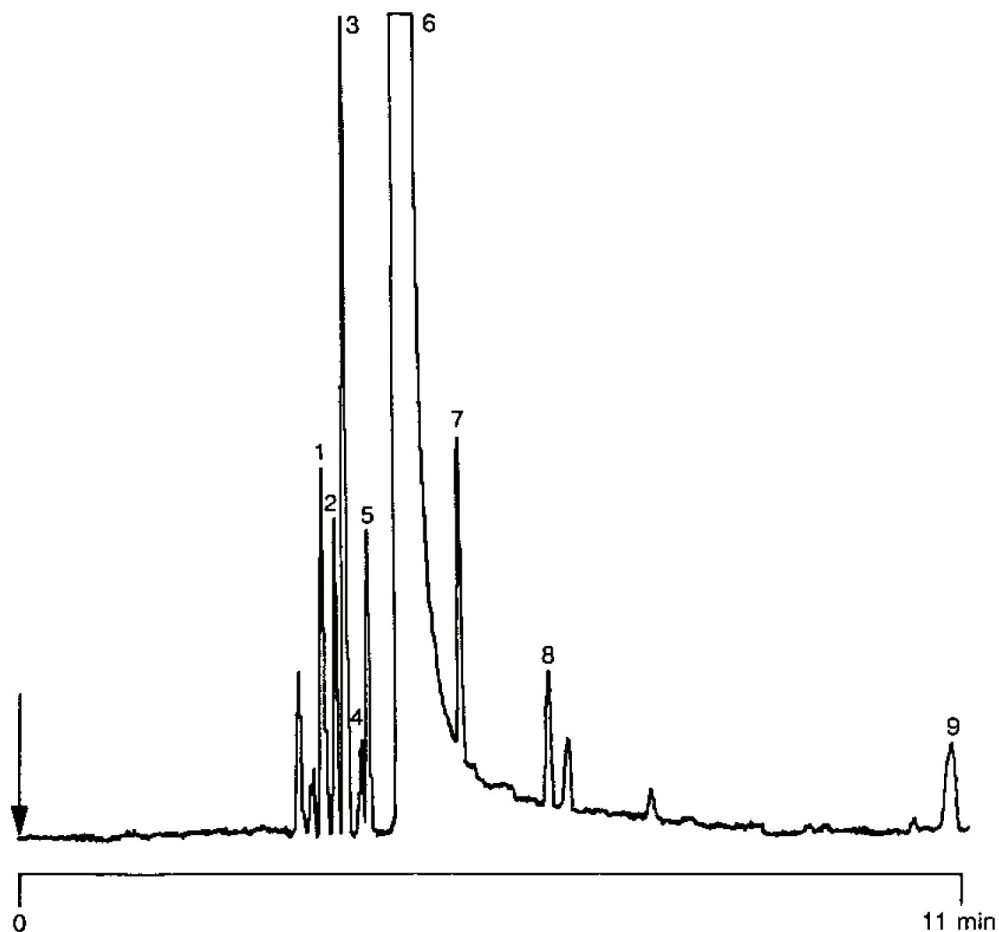
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Propox, 0.53 mm x 50 m fused silica
WCOT CP-Propox
with 0.53 mm x 2.5 m ID retention gap
(Part no. CP8100)
Temperature : 30 °C
Carrier Gas : H₂
Injector : Split, 40 mL/min
T = 100 °C
Detector : FID
T = 275 °C
Sample Size : 0.4 µL

Peak identification

1. 2-chloropropene
2. cis-1-chloropropene
3. trans-1-chloropropene
4. ethylene oxide
5. acetaldehyde
6. propylene oxide
7. propionaldehyde
8. acetone
9. 1,2-dichloropropane



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

A00966



Agilent Technologies