



Hydrocarbons, $C_1 - C_3$, oxygenated hydrocarbons, C_2

Analysis of ethylene oxide

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT U column analyzes ethylene oxide spiked with light hydrocarbons, vinyl chloride and acetaldehyde in 15 minutes.



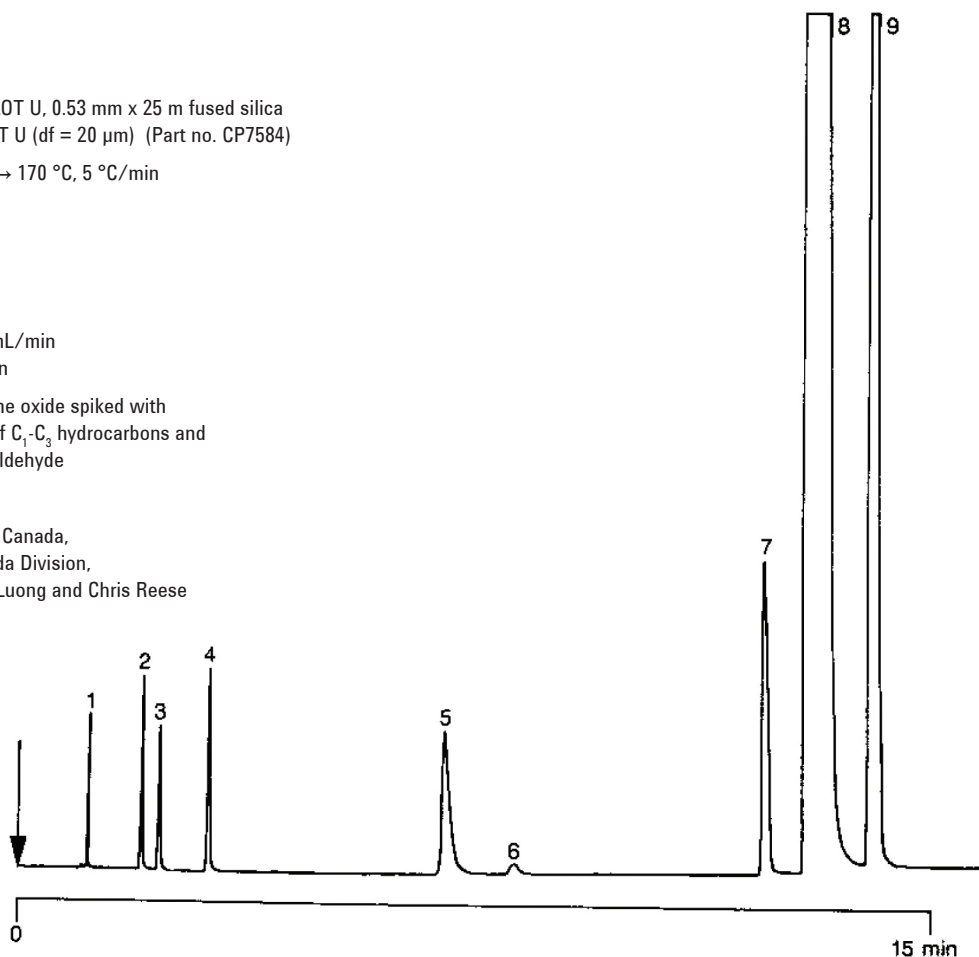
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT U, 0.53 mm x 25 m fused silica
PLOT PoraPLOT U (df = 20 μ m) (Part no. CP7584)
Temperature : 40 °C (2 min) \rightarrow 170 °C, 5 °C/min
Carrier Gas : He, 25 mL/min
Injector : Splitless
T = 200 °C
Detector : FID
hydrogen, 30 mL/min
air, 400 mL/min
Sample : 1 mL of ethylene oxide spiked with
50 ppm (v/v) of C₁-C₃ hydrocarbons and
1% (v/v) acetaldehyde
Courtesy : Dow Chemical Canada,
Western Canada Division,
R&D Lab, Jim Luong and Chris Reese

Peak identification

1. methane
2. ethylene
3. ethane
4. acetylene
5. propylene/propane
6. cyclopropane
7. vinylchloride
8. ethylene oxide
9. acetaldehyde



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

A00674



Agilent Technologies