



Oxygenated hydrocarbons, C₃

Determination of ppm levels of water in pure propylene oxide

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT U column determines ppm levels of water in pure propylene oxide in three minutes.



Agilent Technologies

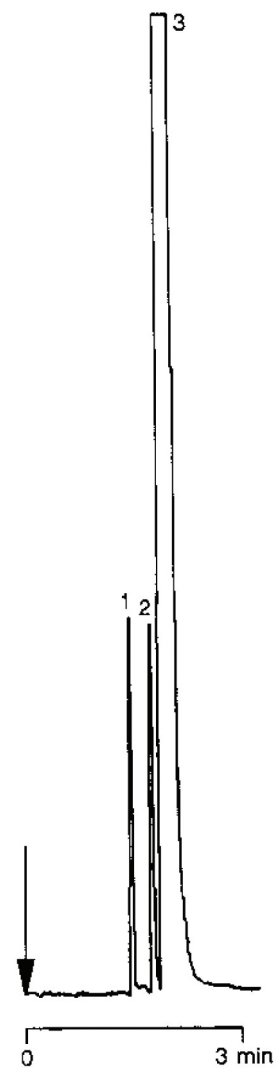
Conditions

Technique : GC-wide-bore
Column : Agilent PoraPLOT U, 0.53 mm x 25 m fused silica
PLOT PoraPLOT U (df = 20 µm) (Part no. CP7584)
Temperature : 190 °C
Carrier Gas : He, 5 mL/min
Injector : Direct
T = 100 °C
Detector : TCD
T = 170 °C
Sample Size : 5 µL

Courtesy : D. Korczewski, Dow Deutschland Inc., Werk Stade,
Germany

Peak identification

1. air
2. water (55 ppm)
3. propylene oxide



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

A00758



Agilent Technologies