



Gases

Analysis of impurities in ethylene

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q UltiMetal column separates gaseous impurities in ethylene in less than three minutes.



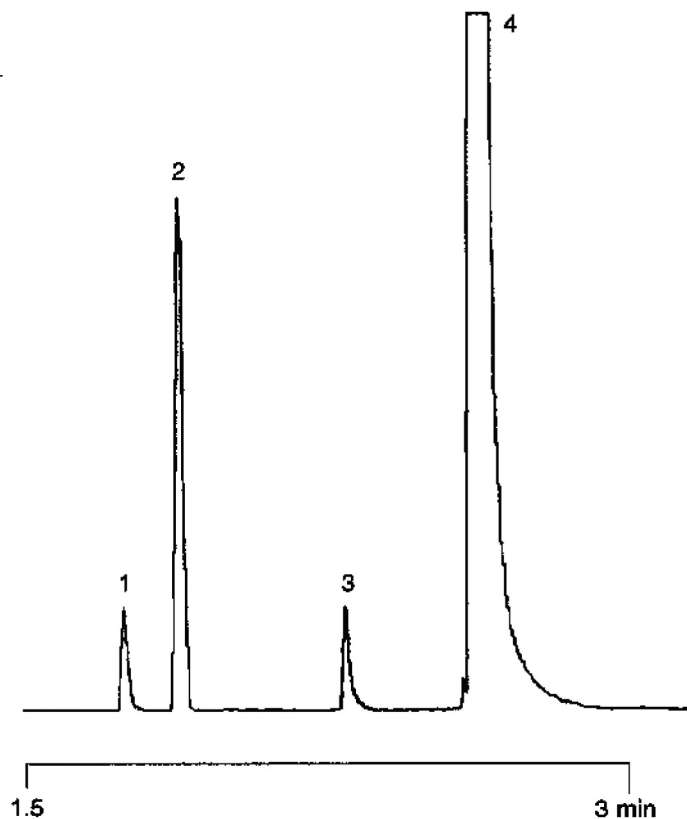
Agilent Technologies

Conditions

Technique : GC-wide-bore
Column : Agilent PoraPLOT Q Ultimet, 0.53 mm x 25 m PLOT
(df = 20 μ m) (Part no. CP6954)
Temperature : 35 $^{\circ}$ C
Carrier Gas : He, 20 kPa (0.2 bar, 2.9 psi)
Injector : Valco valve, 250 μ L
T = 50 $^{\circ}$ C
Detector : FID with a Ni-catalyst methanizer
T = 200 $^{\circ}$ C
Sample Size : 250 μ L

Peak identification

1. carbon monoxide	3 ppm
2. methane	23 ppm
3. carbon dioxide	3 ppm
4. ethylene	99%



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

A01273



Agilent Technologies