



# Gases

## Application Note

Environmental

### Authors

Agilent Technologies, Inc.

### Introduction

Gas chromatography with an Agilent PoraPLOT U column separates four gases in five minutes.



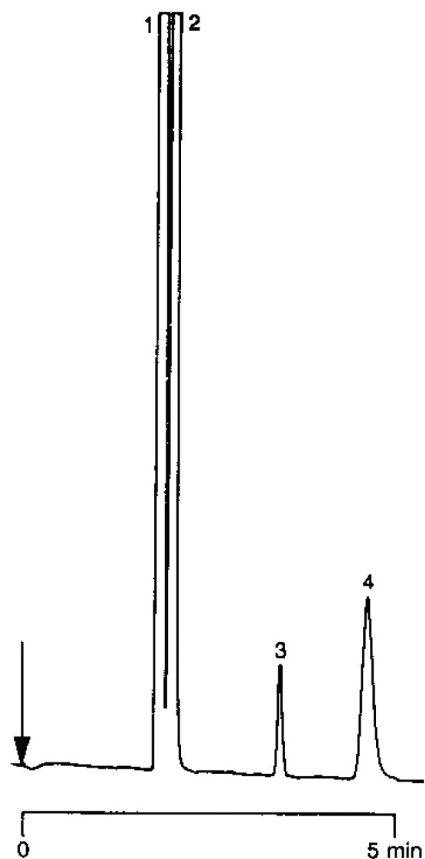
**Agilent Technologies**

## 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 : 26 °C  
Carrier Gas : H<sub>2</sub>  
Injector : Split, 100 mL/min  
Detector : TCD  
Sample Size : 1 - 10%

## Peak identification

1. helium
2. nitrogen + carbon monoxide
3. carbon dioxide
4. ethane



[www.agilent.com/chem](http://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

A00683



**Agilent Technologies**