

Gases

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates three gases in two minutes.



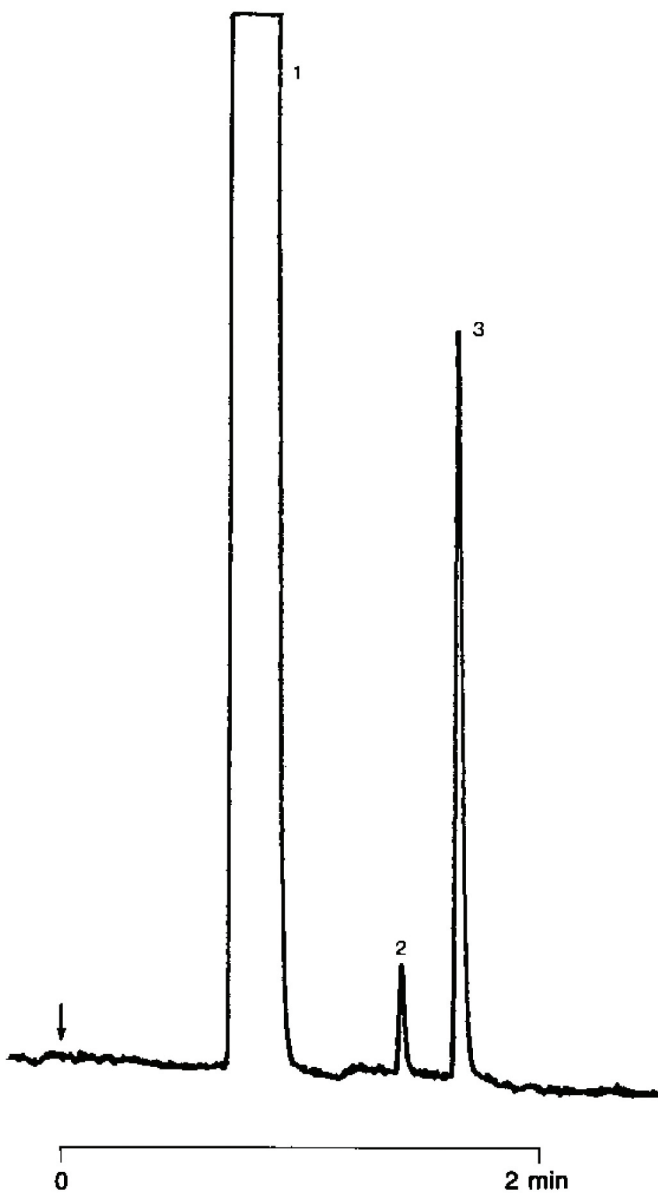
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q, 0.32 mm x 10 m fused silica
PoraPLOT Q (10 μ m) (Part no. CP7550)
Temperature : 25 $^{\circ}$ C
Carrier Gas : H₂, 45 kPa (0.45 bar, 6.5 psi)
Injector : Splitter
Detector : μ TCD

Peak identification

1. carbon monoxide
2. carbon dioxide (100 ppm)
3. nitrous oxide (700 ppm)



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

A00284



Agilent Technologies