



Gases

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

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



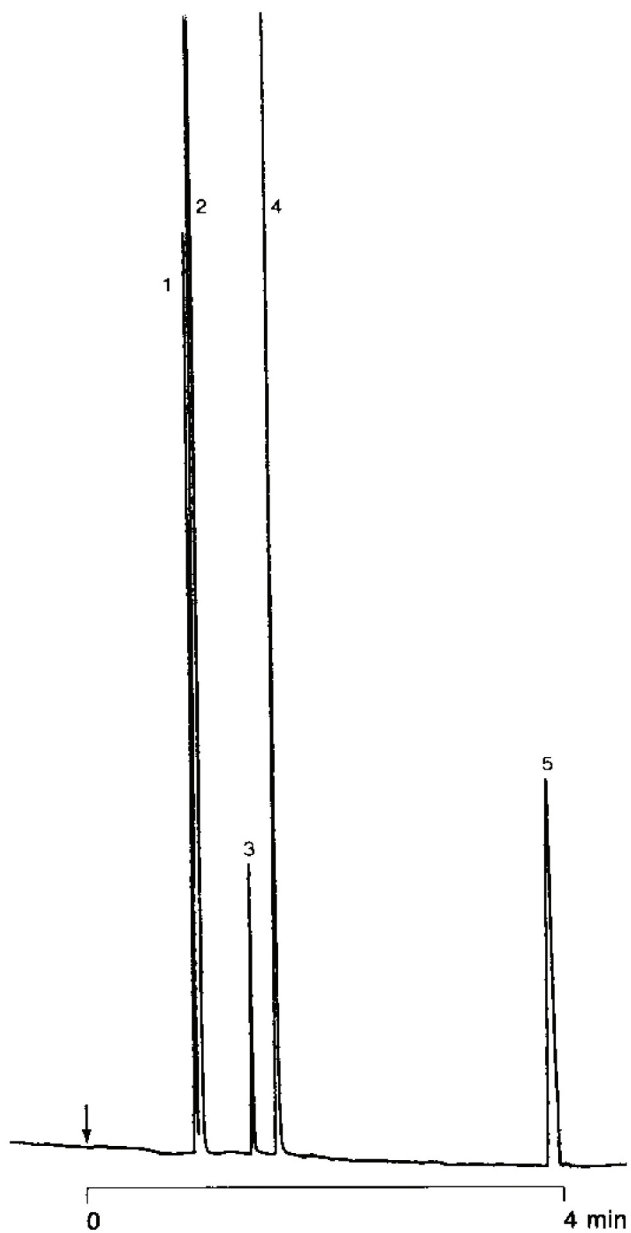
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q, 0.32 mm x 25 m fused silica
PoraPLOT Q
(10 μm) (Part no.CP7551)
Temperature : 25 °C
Carrier Gas : H_2 , 100 kPa (1.0 bar, 14.5 psi)
Injector : Splitter
Detector : μ TCD

Peak identification

1. helium
2. carbon monoxide
3. carbon dioxide
4. nitrous oxide (N_2O)
5. hydrogen sulfide



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

A00294



Agilent Technologies