



Gases

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

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



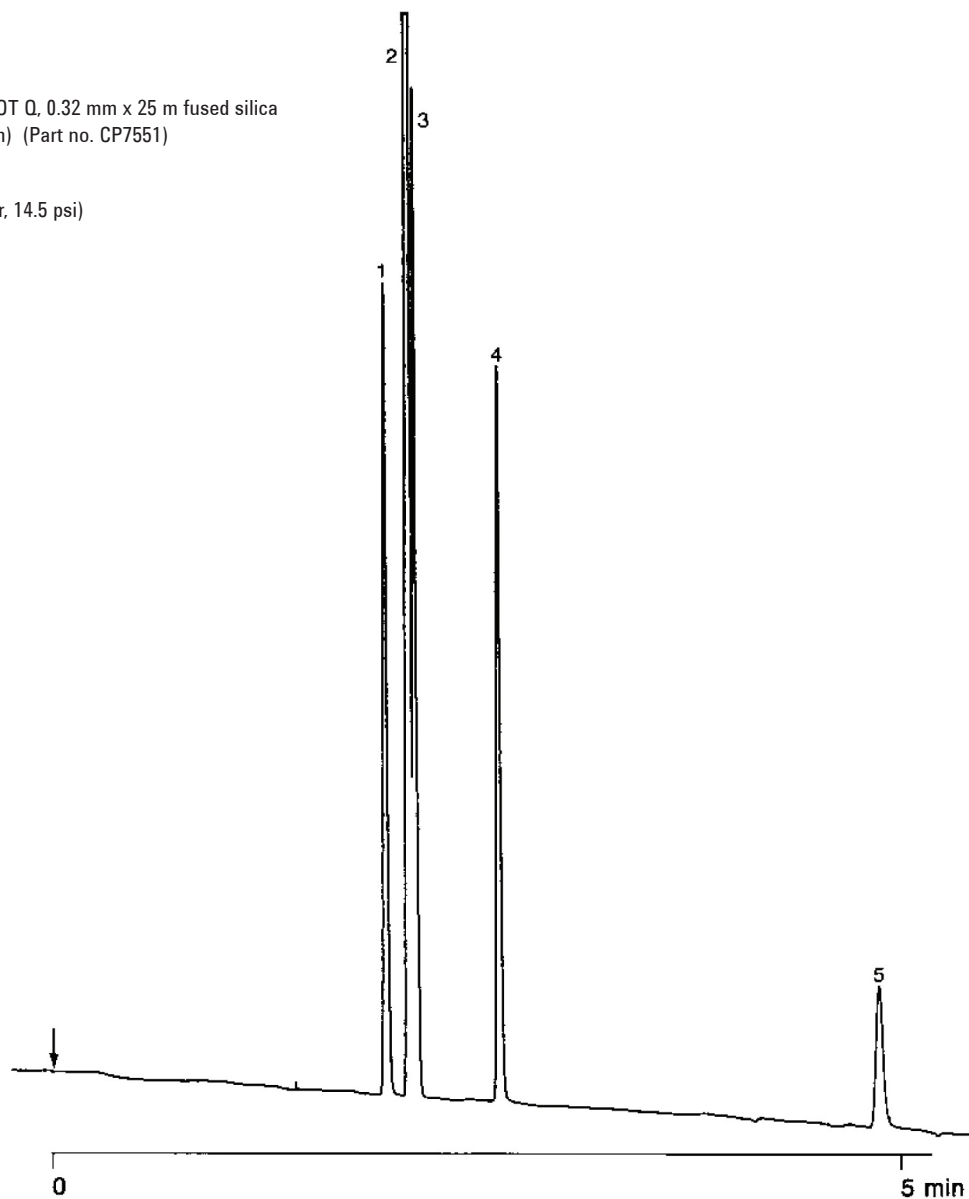
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-PoraPLOT Q, 0.32 mm x 25 m fused silica
PoraPLOT Q (10 μ m) (Part no. CP7551)
Temperature : 0 °C
Carrier Gas : H₂, 100 kPa (1.0 bar, 14.5 psi)
Injector : Splitter
Detector : μ TCD

Peak identification

1. helium
2. nitrogen
3. oxygen
4. methane
5. carbon dioxide



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

A00307



Agilent Technologies