



Permanent gases and hydrocarbons

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates ten permanent gases and hydrocarbons in 15 minutes.



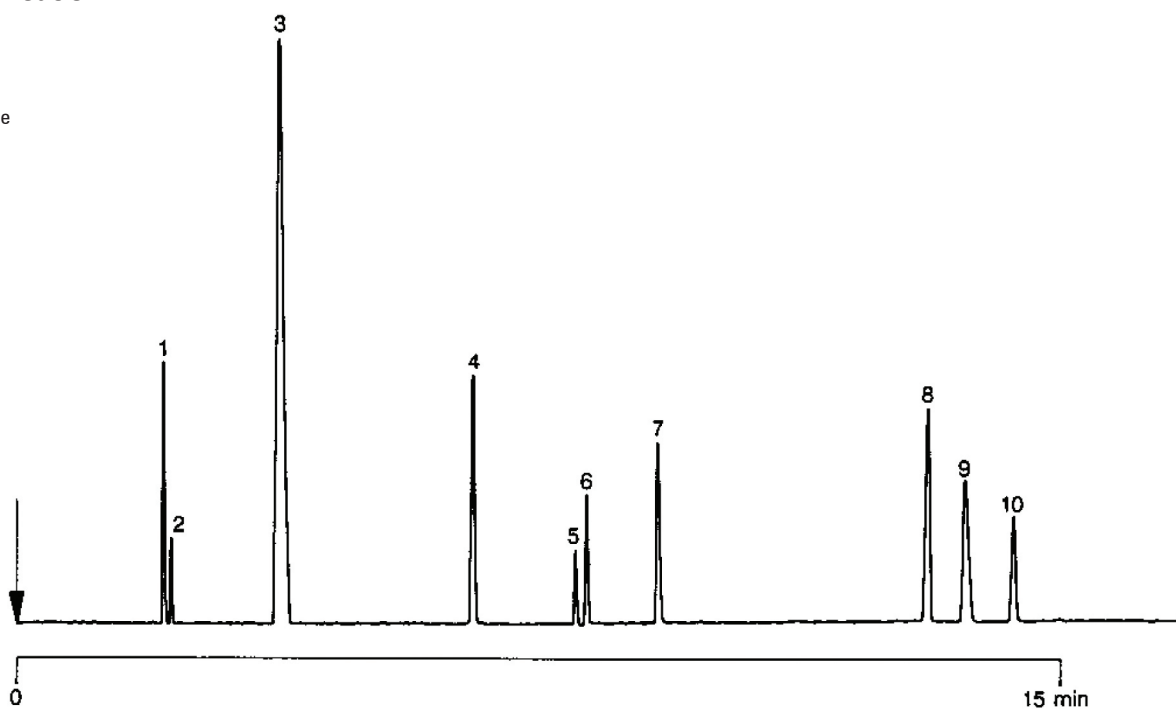
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q, 0.32 mm x 25 m fused silica
PLOT PoraPLOT (df = 10 μ m) (Part no. CP7551)
Temperature : -50 °C (1.5 min) \rightarrow 150 °C, 20 °C/min
Carrier Gas : H₂, 50 kPa (0.5 bar, 7.1 psi)
Injector : Split
Detector : μ -TCD
Sample : 2 to 5% in CH₄

Peak identification

1. nitrogen
2. oxygen
3. methane
4. carbon dioxide
5. ethylene
6. acetylene
7. ethane
8. propylene
9. propane
10. propyne



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

A00680



Agilent Technologies