



Permanent gases

Analysis of gases in water

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Fast gas chromatography with an Agilent PoraPLOT U column separates six permanent gases in water in four minutes.



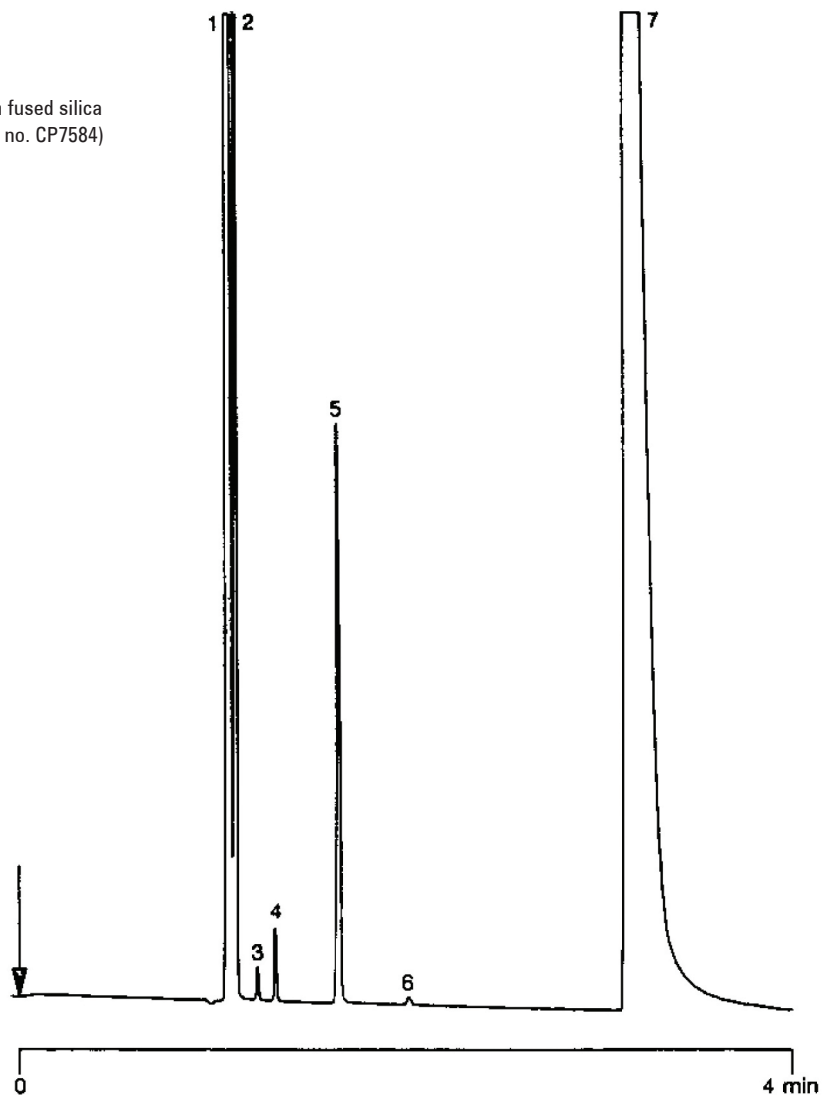
Agilent Technologies

Conditions

Technique : GC-Wide-bore
Column : Agilent PoraPLOT U, 0.53 mm x 25 m fused silica
PLOT PoraPLOT U (df - 20 μ m) (Part no. CP7584)
Temperature : 100 °C
Carrier Gas : H₂, 30 kPa (0.3 bar, 4.3 psi)
Injector : Split, 100 mL/min
Detector : TCD
Sample Size : Headspace 0.1 - 10%

Peak identification

1. air
2. methane
3. carbon dioxide
4. ethane
5. hydrogen sulfide
6. propane
7. water



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

A00684



Agilent Technologies