



Gases

Analysis of gases and water

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q UltiMetal column separates gases in a water sample in ten minutes.



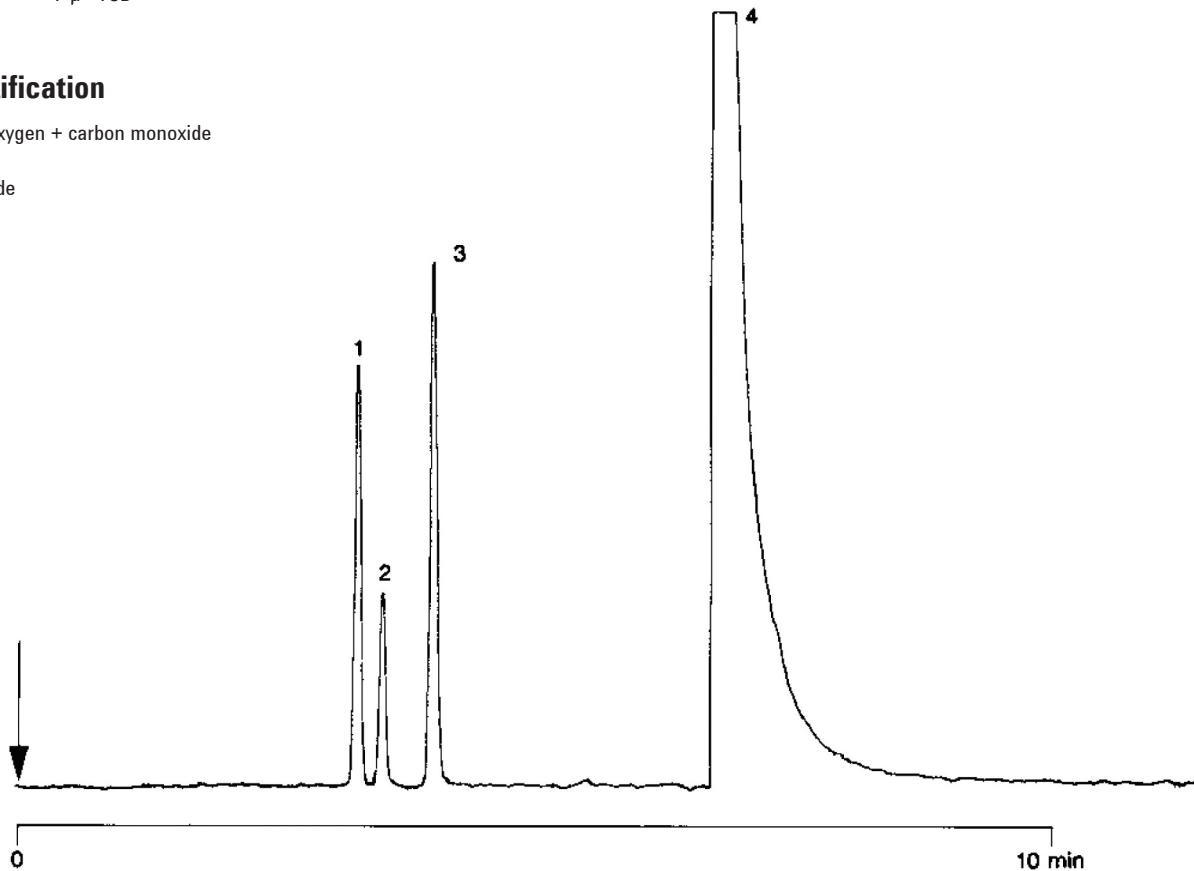
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q Ultimetel, 0.53 mm x 25 m PLOT
PoraPLOT (df = 20 µm) (Part no. CP6954)
Temperature : 75 °C
Carrier Gas : H₂
Injector : Split
Detector : µ - TCD

Peak identification

1. nitrogen + oxygen + carbon monoxide
2. methane
3. carbon dioxide
4. 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

A00992



Agilent Technologies