



## **Gases**

# Analysis of a breath sample

## Application Note

BioPharma

### **Authors**

Agilent Technologies, Inc.

### **Introduction**

Gas chromatography with an Agilent PoraPLOT Q column separates carbon dioxide and water in breath in two minutes.



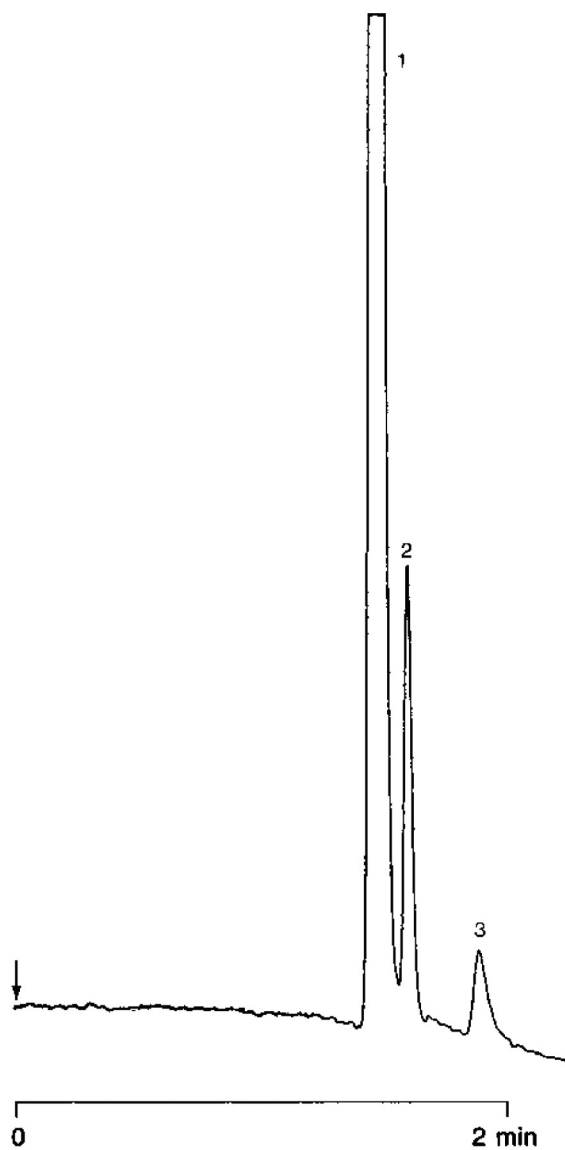
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent PoraPLOT Q, 0.32 mm x 10 m fused silica  
PoraPLOT Q (10  $\mu$ m) (Part no. CP7550)  
Temperature : 150 °C  
Carrier Gas : H<sub>2</sub>, 40 kPa (0.4 bar, 5.8 psi)  
Injector : Splitter  
Detector :  $\mu$  TCD

## Peak identification

1. air
2. carbon dioxide
3. water



[www.agilent.com/chem](http://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

A00279



**Agilent Technologies**