



Gases

Analysis of carbon dioxide in air

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates carbon dioxide in air in 20 seconds.



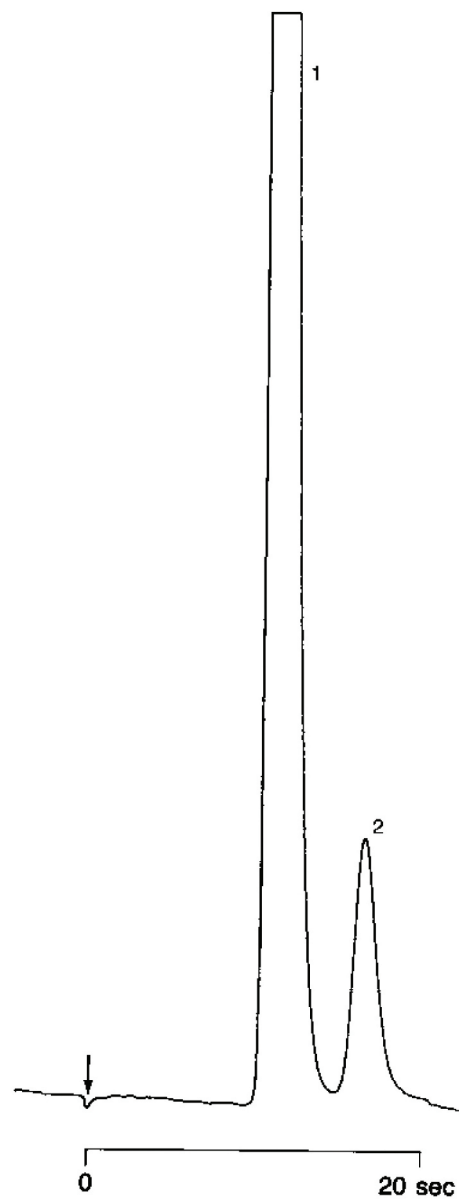
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q, 0.32 mm x 1 m fused silica
PoraPLOT Q (10 μ m) (Custom-made)
Temperature : 30 °C
Carrier Gas : H₂, 10 kPa (0.1 bar, 1.5 psi)
Injector : Splitter
Detector : TCD

Peak identification

1. nitrogen/oxygen
2. 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

A00288



Agilent Technologies