



# Gases

## Application Note

Environmental

### Authors

Agilent Technologies, Inc.

### Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates nitrogen and oxygen in two minutes.



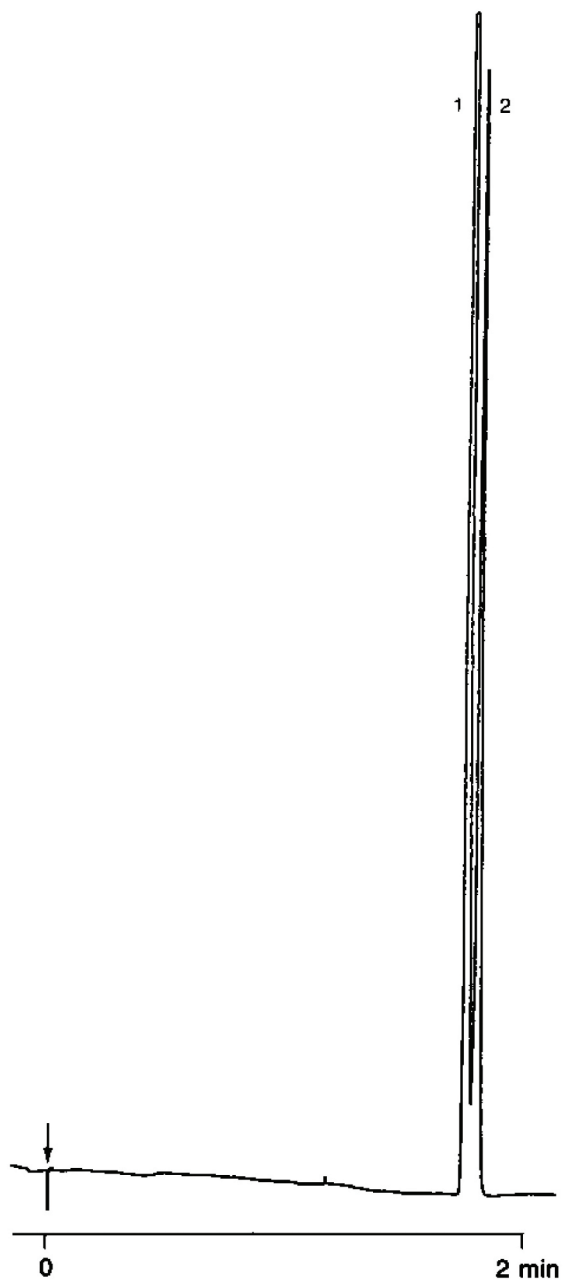
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent PoraPLOT Q, 0.32 mm x 25 m fused silica  
PoraPLOT Q (10  $\mu\text{m}$ ) (Part no. CP7551)  
Temperature : 0  $^{\circ}\text{C}$   
Carrier Gas :  $\text{H}_2$ , 100 kPa (1.0 bar, 14.5 psi)  
Injector : Splitter  
Detector :  $\mu$  TCD

## Peak identification

1. nitrogen
2. oxygen



[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

A000292



**Agilent Technologies**