



Gases

Analysis of laboratory air

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Molsieve 5Å column separates oxygen, nitrogen and argon in laboratory air in 60 seconds.



Agilent Technologies

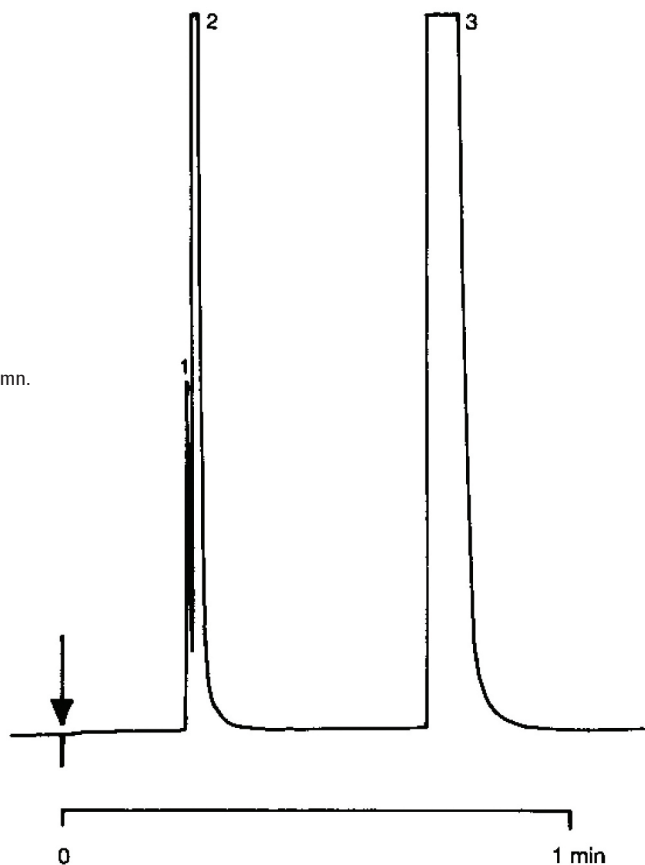
Conditions

Technique : GC-capillary
Column : Agilent CP-Molsieve 5Å, 0.53 mm x 10 m fused silica
PLOT Molsieve 5Å (df = 50 µm) (Part no. CP7537)
Temperature : 25 °C
Carrier Gas : H₂, 45 kPa (0.45 bar, 6.5 psi); 12 mL/min
Injector : Direct
Detector : µ-TCD, Attn. 8 x
Sample Size : 10 µL

Note: compare Application Note 146 for the same analysis on a Narrowbore column.

Peak identification

1. argon
2. oxygen
3. nitrogen



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

A00437



Agilent Technologies