



Gases

Analysis of oxygen and argon content in high purity nitrogen

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Molsieve column identifies the oxygen and argon content in a sample of high purity nitrogen in less than two minutes.



Agilent Technologies

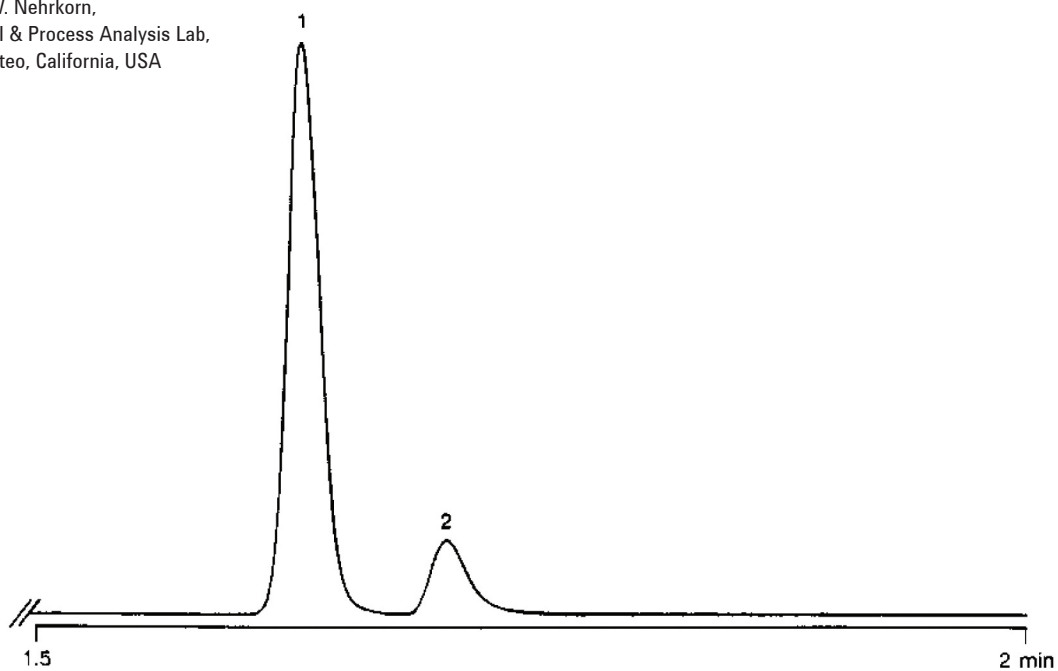
Conditions

Technique : GC-wide-bore
Column : Agilent CP-Molsieve 5Å, 0.53 mm x 25 m fused silica
PLOT CP-Molsieve 5Å (df = 50 µm)
(Part no. CP7538)
Temperature : 35 °C
Carrier Gas : He, 55 kPa (0.55 bar, 8 psi)
Injector : Splitter: 1:10
Detector : TCD
Sample Size : 2 mL gas sampling loop

Courtesy : David W. Nehrkorn,
Material & Process Analysis Lab,
San Mateo, California, USA

Peak identification

1. argon	600 ppm
2. oxygen	67 ppm



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

A00767



Agilent Technologies