



Gases

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Separation of hydrogen and helium by GC on an Agilent CP-Molsieve column in 7.5 minutes.



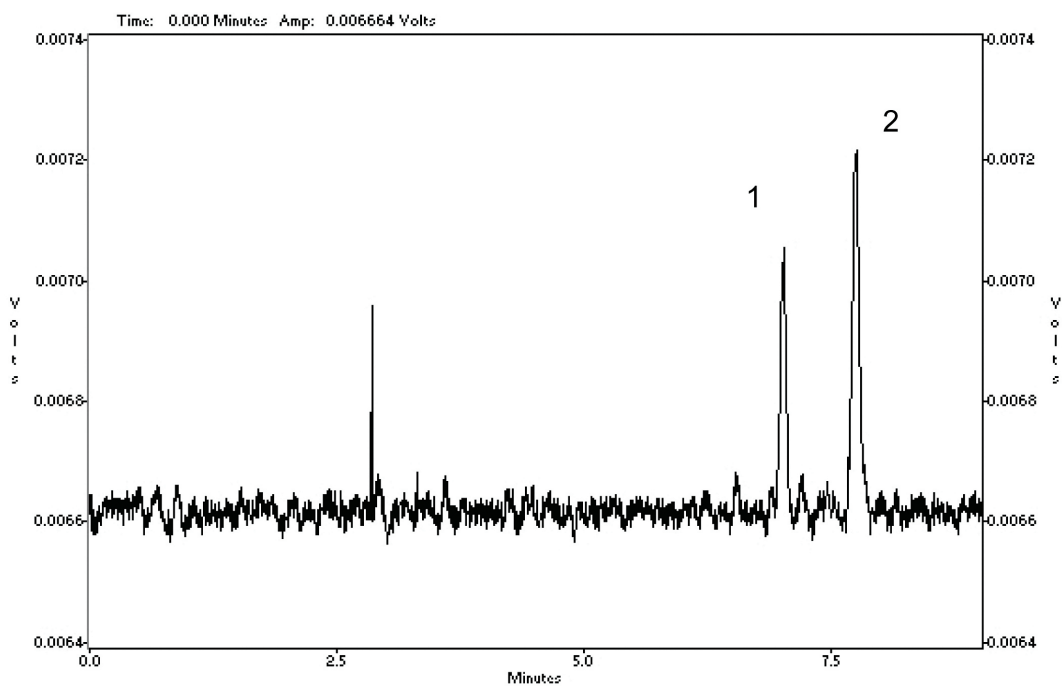
Agilent Technologies

Conditions

Technique : GC
Column : Agilent CP-Molsieve 5Å, 0.53 mm x 50 m fused silica
(df = 50 µm) (Part no. CP7539)
Temperature : 40 °C
Carrier Gas : Nitrogen, 50kPa (7.2 psi)
Injector : Splitter, 40 mL/min
Detector : µ-TCD, 200 °C
Sample Size : 40 µL
Concentration Range : 1% in nitrogen

Peak identification

1. helium
2. hydrogen



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

A02077



Agilent Technologies