



Anesthetics

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column module and Agilent 490 Micro GC determines impurities in nitrogen in 150 seconds.



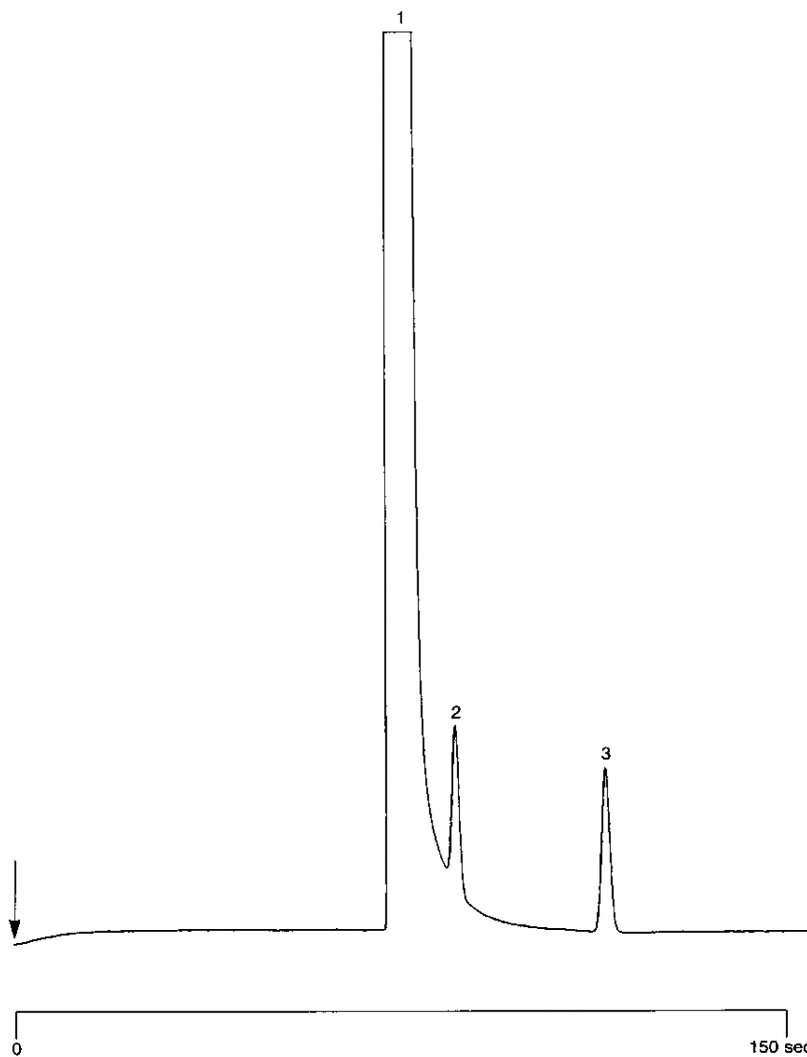
Agilent Technologies

Conditions

Technique : Micro-GC
Column : Agilent PoraPLOT Q
Temperature : 30 °C
Carrier Gas : H₂
Detector : TCD

Peak identification

1.	nitrogen	matrix
2.	methane	35ppm
3.	carbon dioxide	39ppm



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

A01243



Agilent Technologies