



Anesthetics

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column and Agilent 490 Micro GC analyzes nitrous oxide in air in 35 seconds.



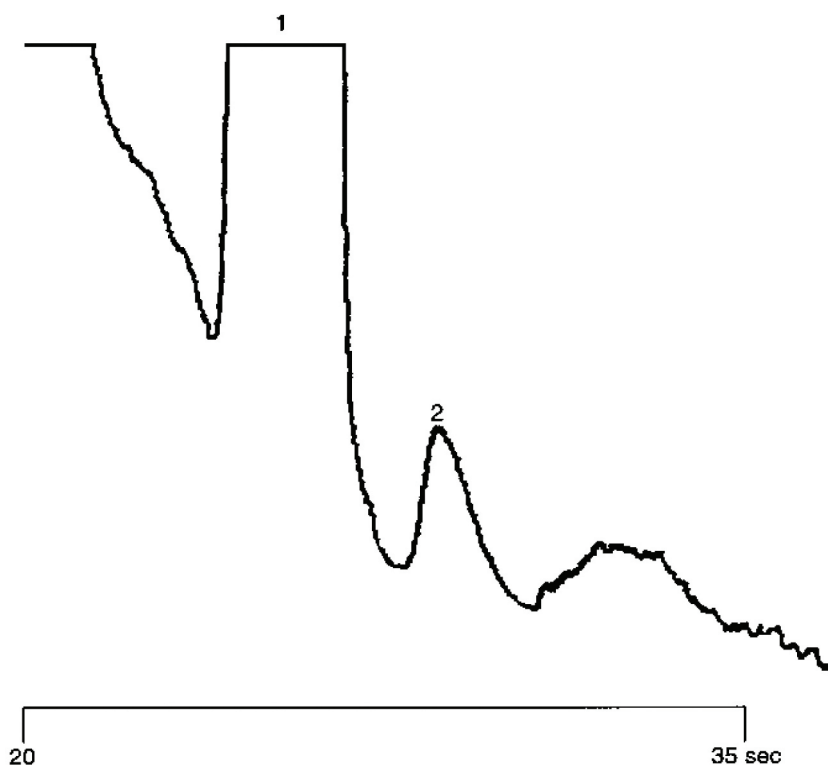
Agilent Technologies

Conditions

Technique : Micro-GC
Column : Agilent PoraPLOT Q
Temperature : 35 °C
Carrier Gas : He
Detector : TCD

Peak identification

1. carbon dioxide
2. nitrous oxide 1ppm



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

A01244



Agilent Technologies