



Gases

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column module and Agilent 490 Micro GC separates nitrous oxide in air in 75 seconds.



Agilent Technologies

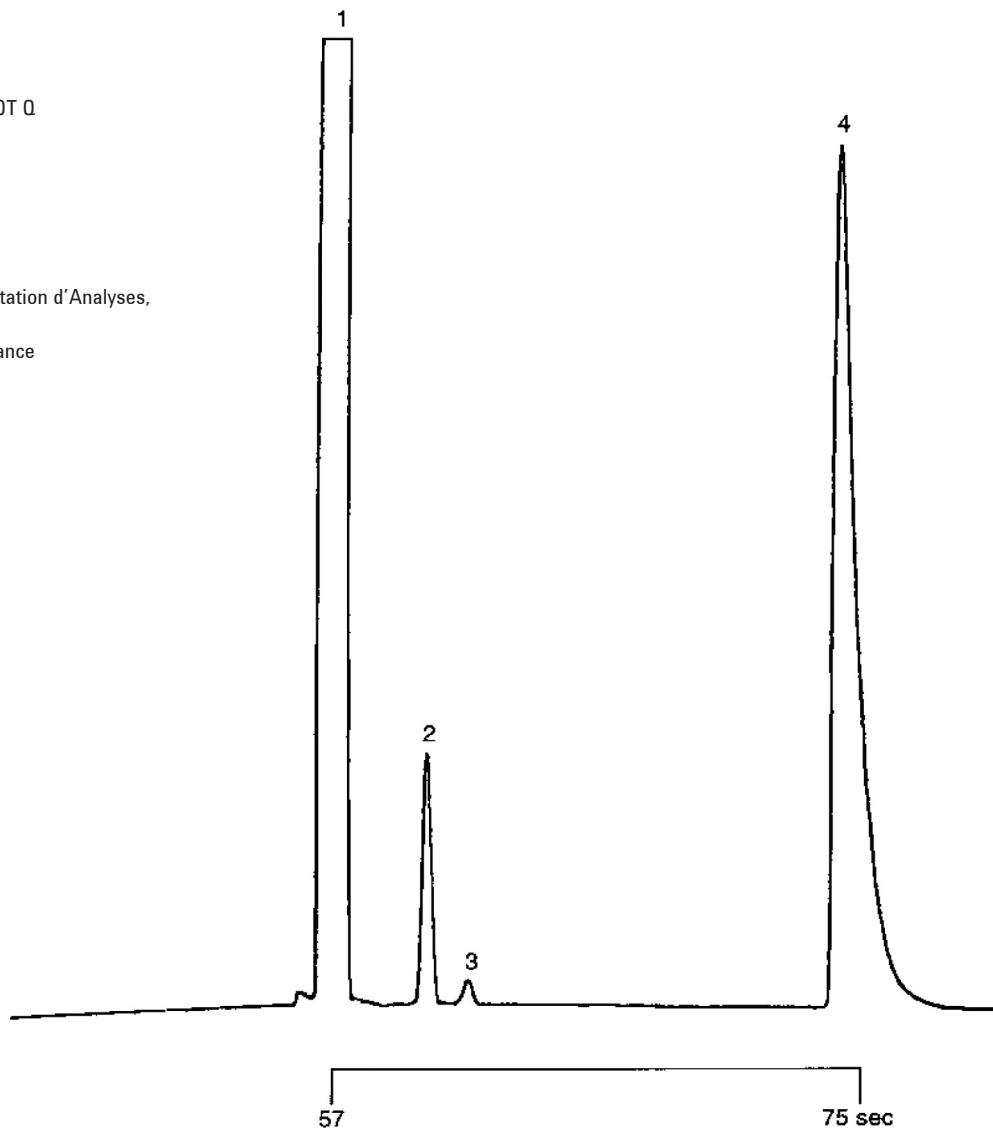
Conditions

Technique : Micro-GC
Column : Agilent PoraPLOT Q
Temperature : 60 °C
Carrier Gas : He
Detector : TCD

Courtesy : Société de Prestation d'Analyses,
Quad-lab,
Vert-le-Petit, France

Peak identification

1. air
2. carbon dioxide
3. nitrous oxide 40 ppm
4. water



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

A01240



Agilent Technologies