



Halogenated hydrocarbons

Analysis of O₂ and N₂ in CF₄

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Molsieve 5Å column module and Agilent 490 Micro GC separates oxygen and nitrogen in carbon tetrafluoride in 70 seconds.



Agilent Technologies

Conditions

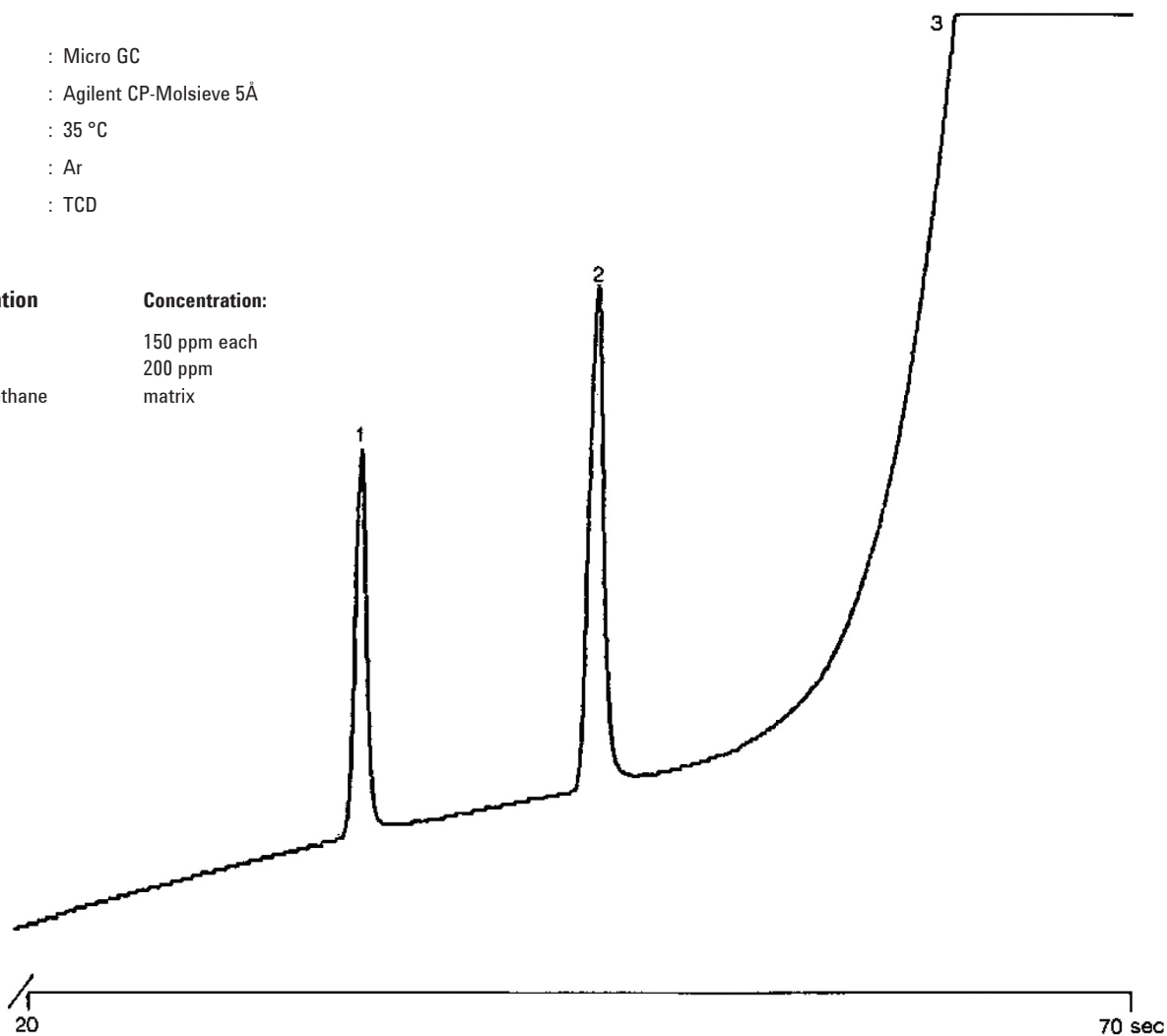
Technique : Micro GC
Column : Agilent CP-Molsieve 5Å
Temperature : 35 °C
Carrier Gas : Ar
Detector : TCD

Peak identification

1. oxygen
2. nitrogen
3. tetrafluoromethane

Concentration:

150 ppm each
200 ppm
matrix



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

A00863



Agilent Technologies