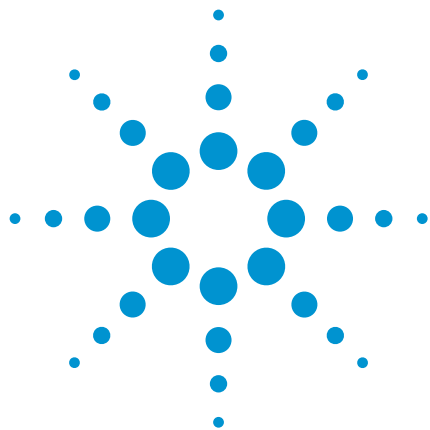




Halogenated hydrocarbons, $C_1 - C_2$

Analysis of CFCs

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates eight chlorofluorocarbons in six minutes.



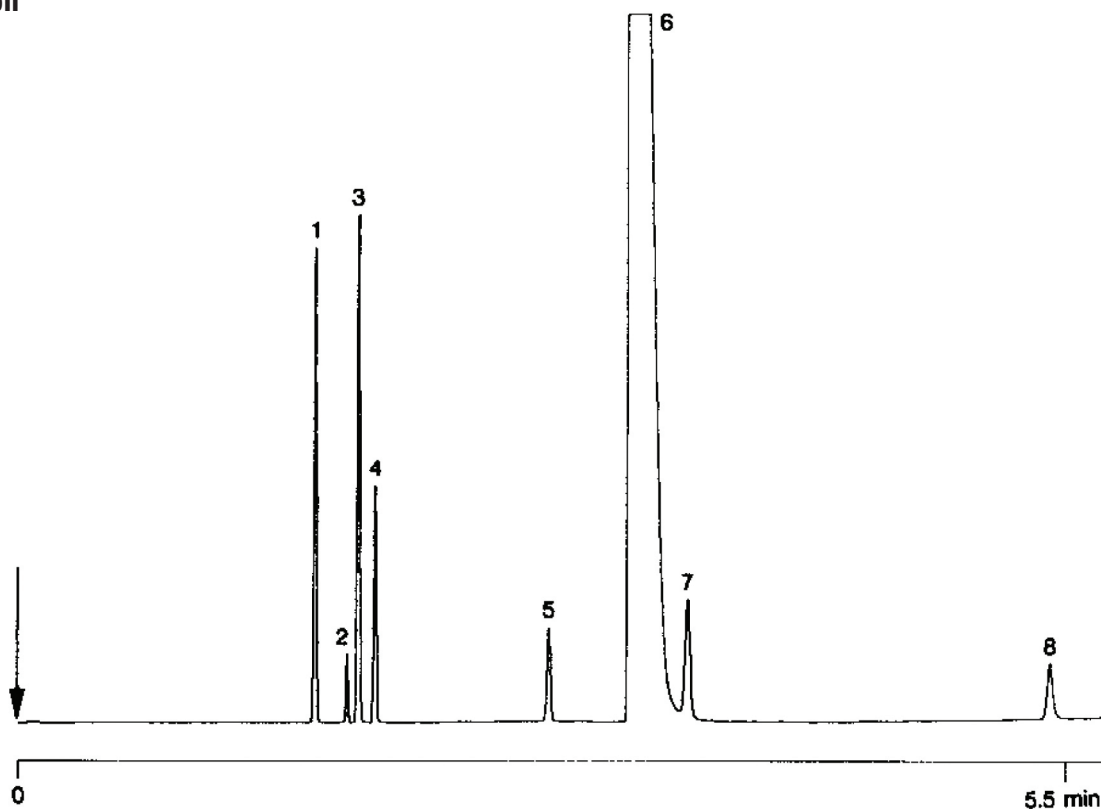
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q, 0.32 mm x 25 m fused silica
PLOT PoraPLOT (df = 20 μ m) (Part no. CP7555)
Temperature : 100 °C $\rightarrow$ 200 °C, 15 °C/min
85 °C (3 min) $\rightarrow$ 140 °C, 10 °C/min
Carrier Gas : He
Injector : Direct
Detector : FID
Concentration Range : ca. 50 ppm in N₂ + ca. 5000 ppm CFC 12

Peak identification

1. methane
2. CFC 23 (CHF₃)
3. CFC 41 (CH₃F)
4. CFC 32 (CH₂F₂)
5. CFC 22 (CHClF₂)
6. CFC 12 (CF₂Cl₂)
7. CFC 31 (CH₂ClF)
8. CFC 21 (CHCl₂F)



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

A00918



Agilent Technologies