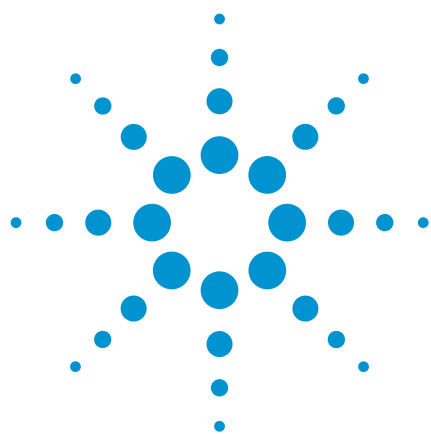




Halogenated hydrocarbons, C₁ – C₃

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates 15 chlorofluorocarbons in 25 minutes.



Agilent Technologies

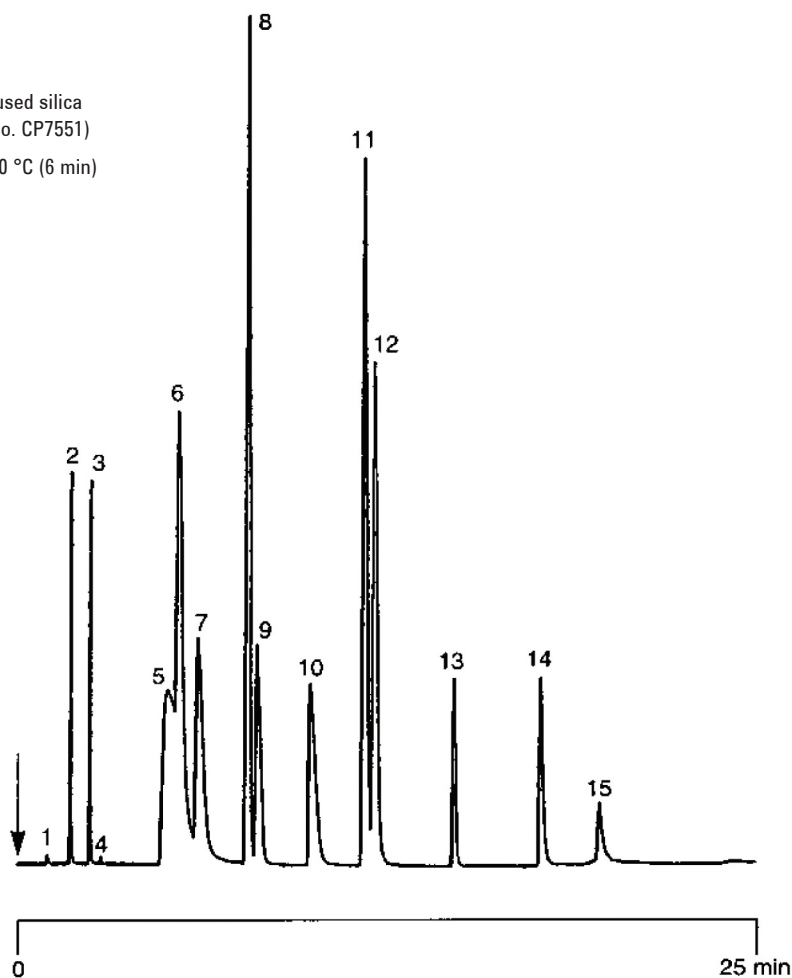
Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q, 0.32 mm x 25 m fused silica
PLOT PoraPLOT Q (df = 10 µm) (Part no. CP7551)
Temperature : 30 °C (5 min) → 200 °C, 12 °C/min; 200 °C (6 min)
Carrier Gas : He, 150 kPa (1.5 bar, 21 psi)
Injector : Split, 100 mL/min
T = 180 °C
Detector : FID
T = 200 °C

Courtesy : Mr Valtz, Cerep,
Ecole des Mines de Paris,
Paris, France

Peak identification

1. CFC 14, tetrafluoromethane
2. CFC 23, trifluoromethane
3. CFC 32, difluoromethane
4. CFC 13, chlorotrifluoromethane
5. CFC 218, octafluoropropane
6. CFC 143a, 1,1,1-trifluoroethane
7. CFC 125, pentafluoroethane
8. CFC 152a, 1,1-difluoroethane
9. CFC 22, chlorodifluoromethane
10. CFC 12, dichlorodifluoromethane
11. CFC 142b, 1-chloro-1,1-difluoroethane
12. CFC 124, 1-chloro-1,2,2,2-tetrafluoroethane
13. trifluoroethanol
14. CFC 113, 1,1,2-trichloro-1,2,2-trifluoroethane
15. FC-75, fluorinert



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

A001110



Agilent Technologies