

# Halogenated hydrocarbons

## Separation of CFC (chlorofluorocarbons)

### Application Note

Environmental

#### Authors

Agilent Technologies, Inc.

#### Introduction

Gas chromatography with an Agilent CP-SilicaPLOT column separates three chlorofluorocarbons and methane in six minutes.



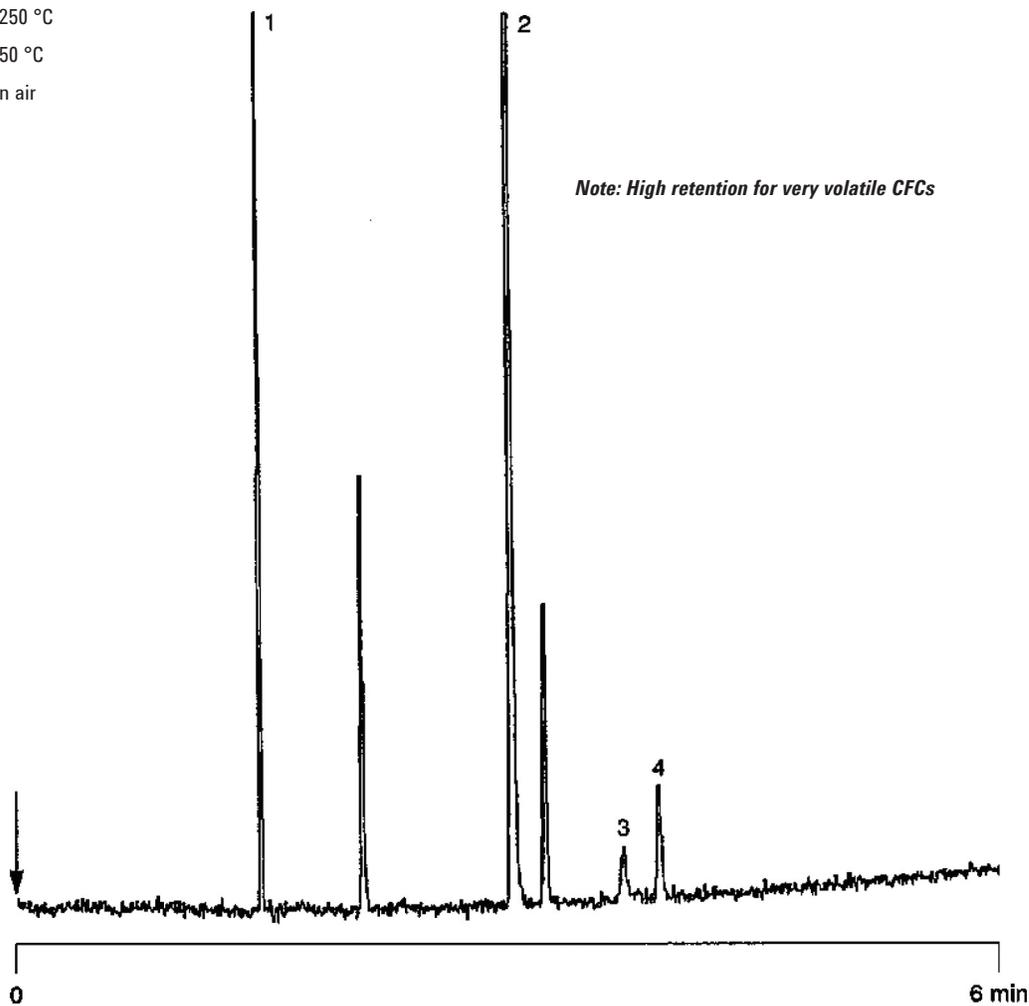
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-SilicaPLOT, 0.32 mm x 30 m fused silica  
PLOT CP-SilicaPLOT (df = 4  $\mu$ m) (Part no. CP8567)  
Temperature : 90 °C (3 min)  $\rightarrow$  200 °C, 10 °C/min  
Carrier Gas : He, 100 kPa (1 bar, 14 psi)  
Injector : Split, T = 250 °C  
Detector : FID, T = 250 °C  
Concentration Range : 100 ppm in air

## Peak identification

1. methane
2. CFC 32
3. CFC 12
4. CFC 22



[www.agilent.com/chem](http://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

A01259



**Agilent Technologies**