



Trihalomethanes

Application Note

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 13 CB for Halocarbons column separates four trihalomethanes according to EPA 501 in four minutes.



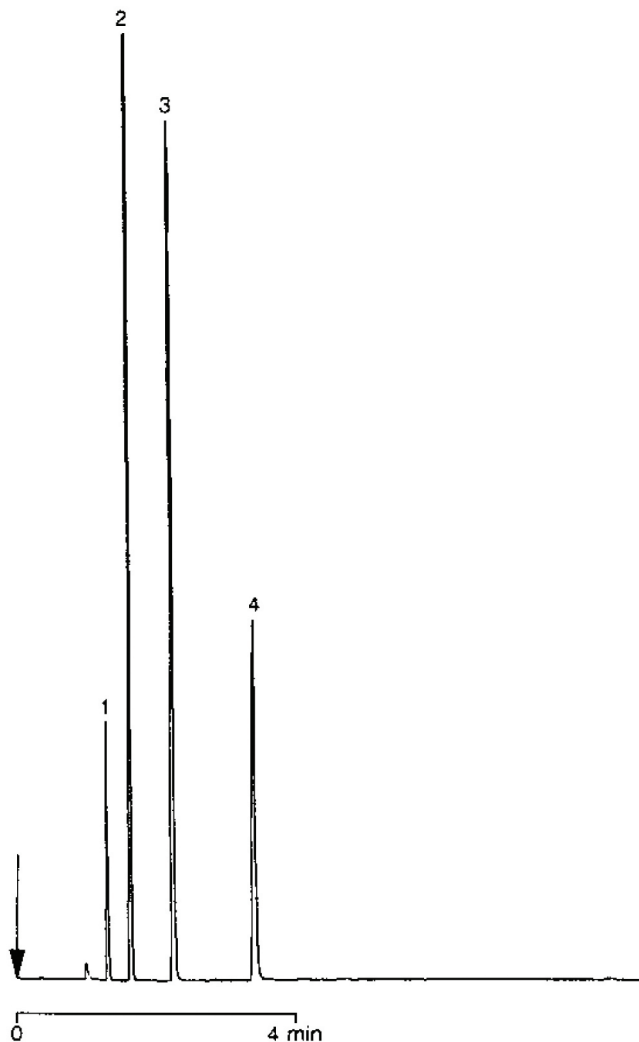
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 13 CB for Halocarbons, 0.32 mm x 25 m fused silica WCOT CP-Sil 13 CB (df = 1.2 μ m)
(Part no. CP7506)
Temperature : 110 °C (isothermal)
Carrier Gas : He, 100 kPa (1 bar, 14 psi)
Injector : Split, 100 mL/min
T = 265 °C
Detector : ECD
T = 300 °C

Peak identification

1. chloroform
2. bromodichloromethane
3. dibromochloromethane
4. bromoform



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

A00776



Agilent Technologies