



Haloethers

Analysis of haloethers according to EPA 611

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB column separates five haloethers according to EPA 611 in 25 minutes.



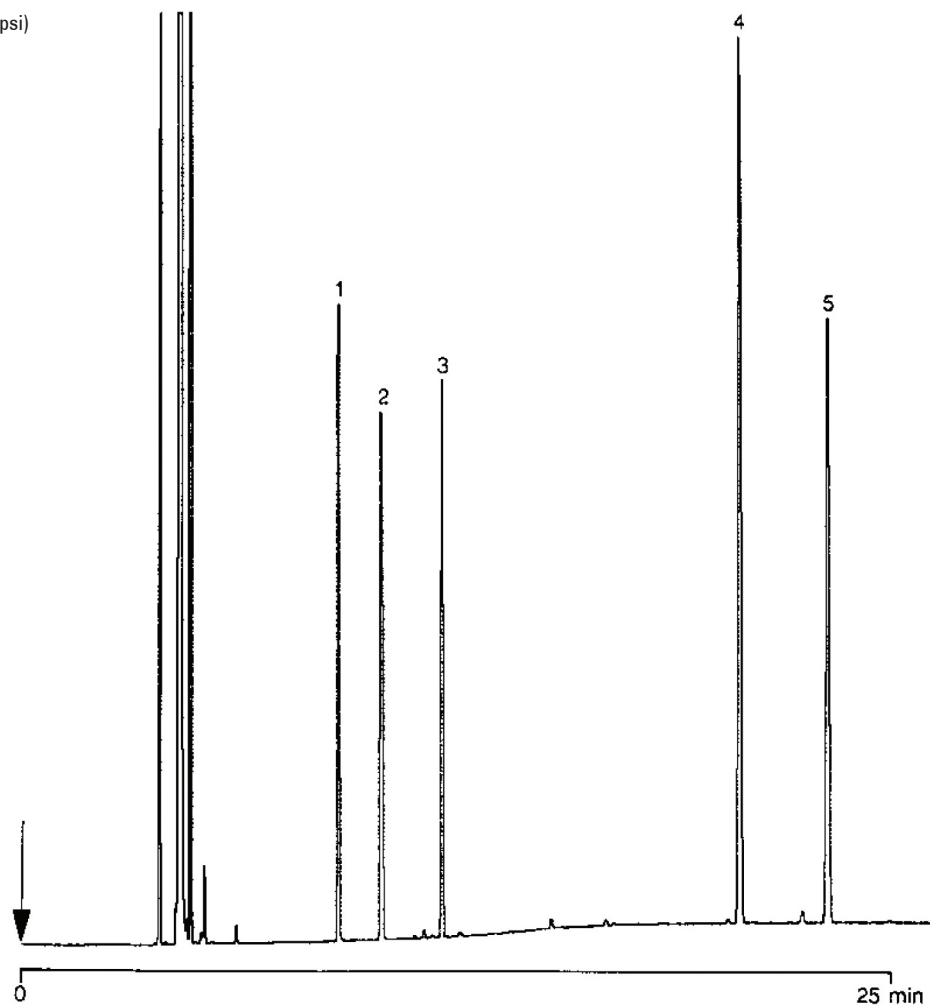
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB, 0.32 mm x 50 m fused silica
WCOT CP-Sil 8 CB (df = 1.2 μ m)
(Part no. CP7771)
Temperature : 100 °C $\rightarrow$ 250 °C, 10 °C/min
Carrier Gas : He, 90 kPa (0.9 bar, 13 psi)
Injector : Split, 100 mL/min
T = 265 °C
Detector : FID
T = 300 °C

Peak identification

1. bis(2-chloroethyl)ether
2. bis(2-chloroisopropyl)ether
3. bis(2-chloroethoxy)methane
4. 4-chlorophenyl phenylether
5. 4-bromophenyl phenylether



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

A00787



Agilent Technologies