



Halogenated compounds

Analysis of 1,2-dibromomethane and
1,2-dibromo-3-chloropropane to EPA 504

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 13 CB for Halocarbons column separates 1,2-dibromomethane and 1,2-dibromo-3-chloropropane according to EPA in seven minutes.



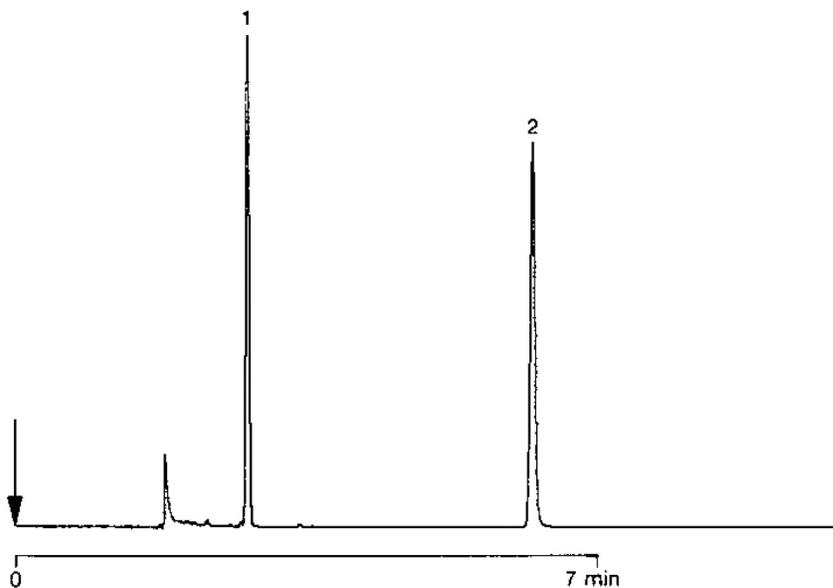
Agilent Technologies

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 : 150 °C (isothermal)
Carrier Gas : He, 50 kPa (0.5 bar, 7.2 psi)
Injector : Split, 100 mL/min
T = 265 °C
Detector : ECD
T = 300 °C

Peak identification

- 1,2-dibromoethane
- 1,2-dibromo-3-chloropropane



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

A00777



Agilent Technologies