



Aldehydes, C₁ Formalin headspace gas analysis

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Sil 5 CB column separates components in formalin headspace gas in eight minutes.



Agilent Technologies

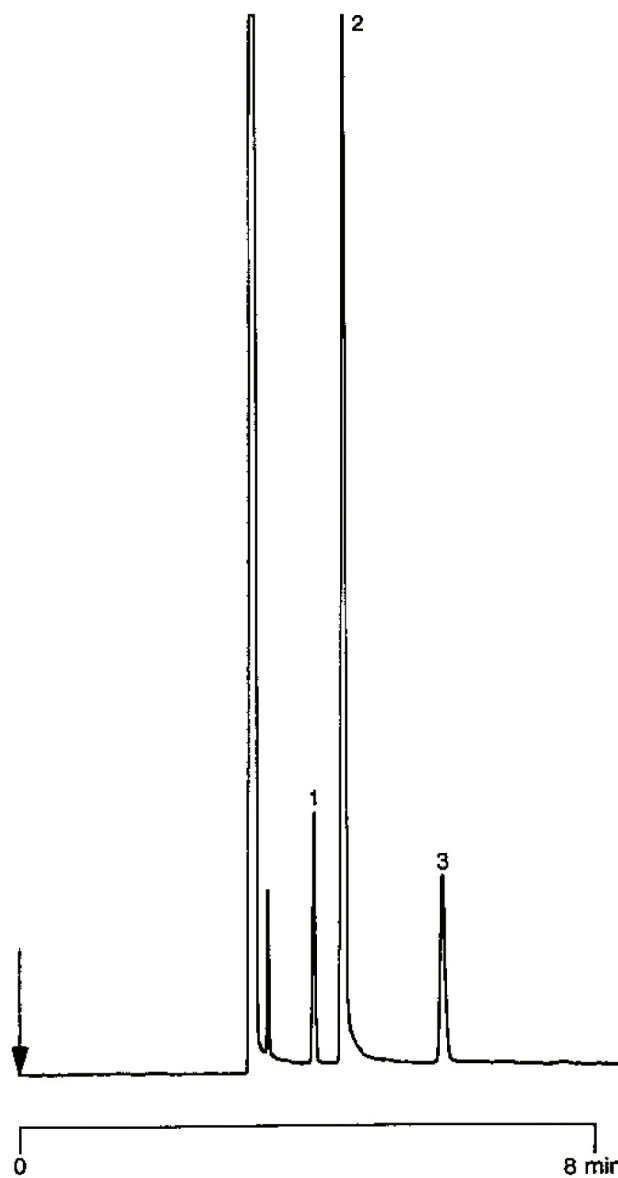
Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5 CB fused silica WCOT CP-Sil 5 CB
(df = 5.0 μ m) (Part no. CP7690)
Temperature : 40 °C
Carrier Gas : He, 120 kPa (1.2 bar, 17 psi), 2 mL/min
Injector : Splitter, 1:20
T = 200 °C
Detector : μ -TCD, 8 x 10 Amp
T = 220 °C, filaments T = 300 °C
Sample Size : 0.4 mL

Courtesy : GL Sciences, Japan

Peak identification

1. formaldehyde
2. water
3. methanol



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

A00624



Agilent Technologies