



## **Gases**

# Analysis of formaldehyde in air

## Application Note

Environmental

### **Authors**

Agilent Technologies, Inc.

### **Introduction**

Gas chromatography with an Agilent CP-Sil 5 CB column module and Agilent 490 Micro GC separates formaldehyde in air in 30 seconds.



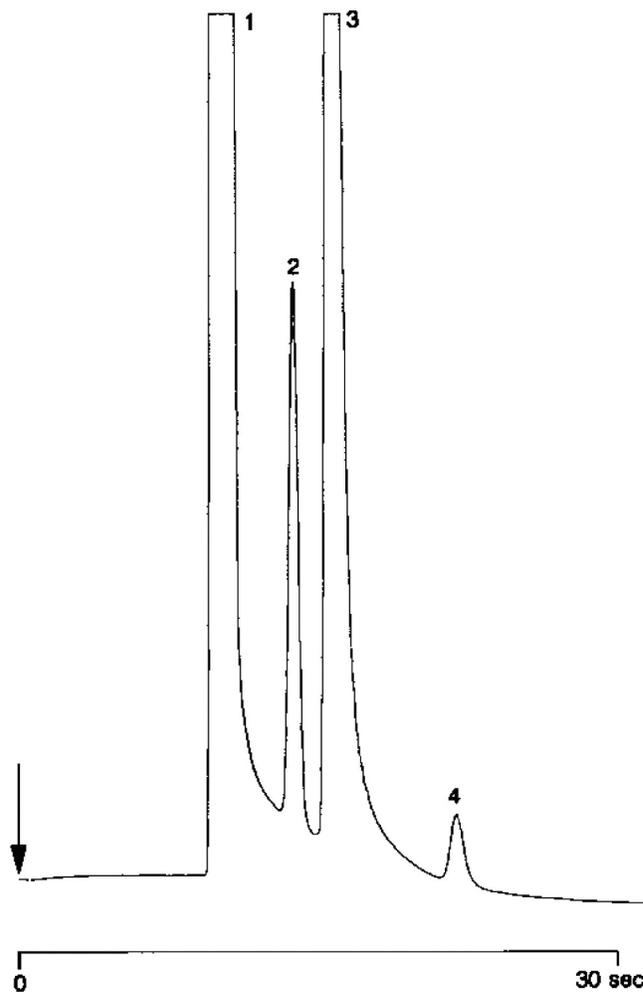
**Agilent Technologies**

## Conditions

Technique : Micro GC  
Column : Agilent CP-Sil 5 CB  
Temperature : 40 °C  
Carrier Gas : He  
Detector : TCD

### Peak identification      Concentration:

- |                 |          |
|-----------------|----------|
| 1. air          |          |
| 2. formaldehyde | 1000 ppm |
| 3. water        |          |
| 4. methanol     |          |



[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

A00989



**Agilent Technologies**