



Aldehydes $C_2 - C_5$

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 5 CB column separates nine C_2 to C_5 aldehydes in nine minutes.



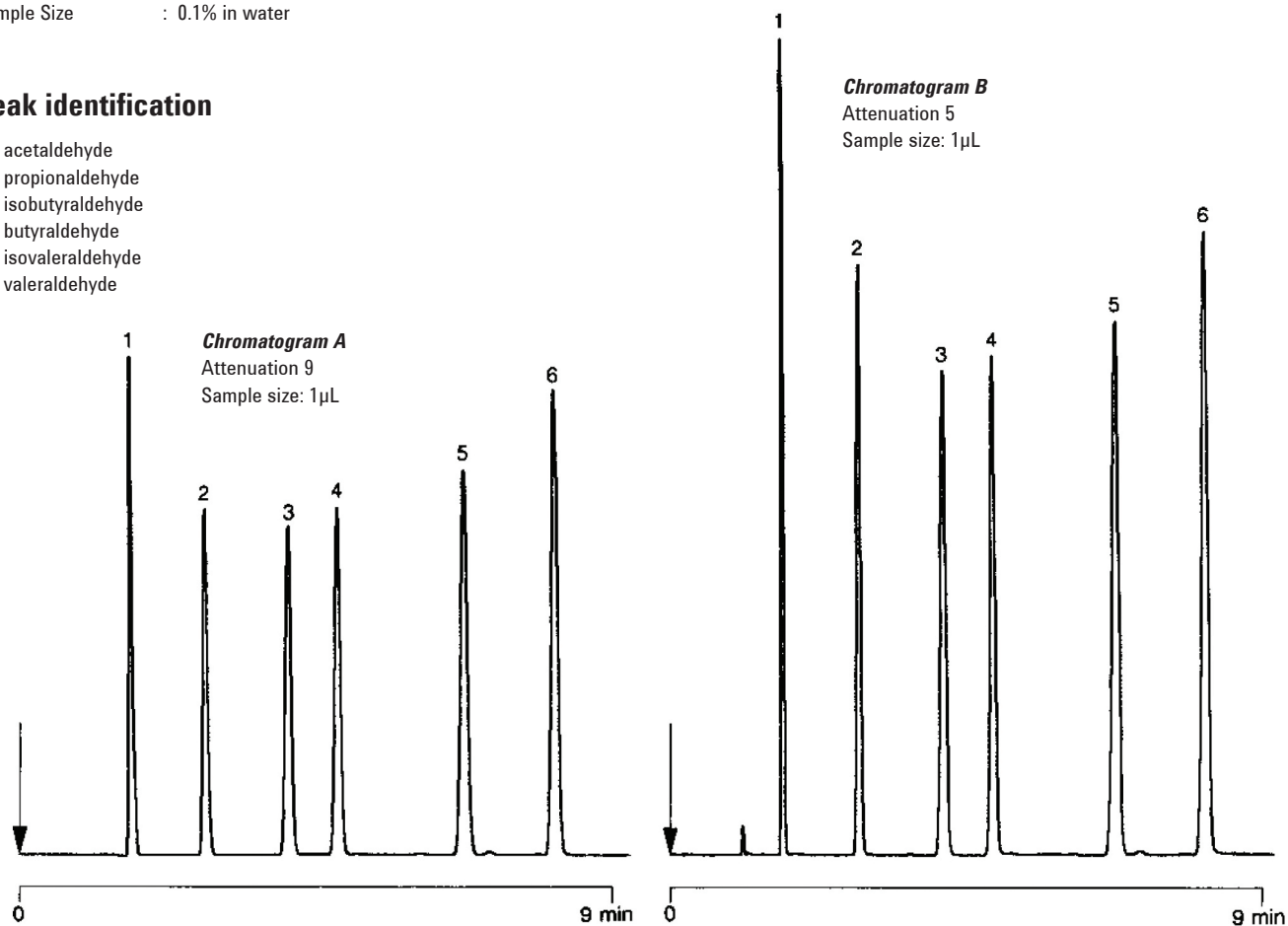
Agilent Technologies

Conditions

Technique : GC-wide-bore
Column : Agilent CP-Sil 5 CB, 0.53 mm x 50 m fused silica
WCOT CP-Sil 5 CB (df = 5 μ m) (Part no. CP7685)
Temperature : 30 °C $\rightarrow$ 80 °C, 5 °C/min
Carrier Gas : He, 100 kPa (1 bar, 14.5 psi), 78 cm/s
Injector : Split, 100 mL/min
T = 150 °C
Detector : FID
T = 200 °C
Sample Size : 0.1% in water

Peak identification

1. acetaldehyde
2. propionaldehyde
3. isobutyraldehyde
4. butyraldehyde
5. isovaleraldehyde
6. valeraldehyde



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

A00844



Agilent Technologies