



Ethers, C₄ – C₈

Application Note

Energy & Fuels

Authors

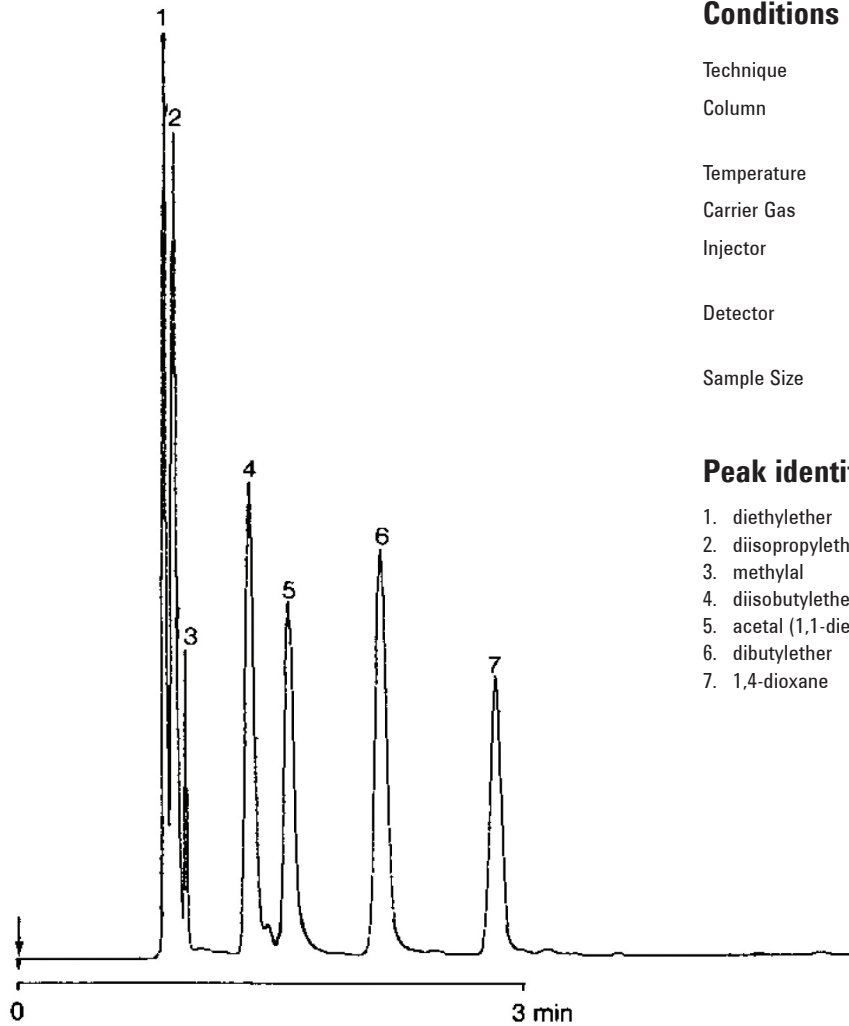
Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 52 CB column separates seven C₄ to C₈ ethers in three minutes.



Agilent Technologies



Conditions

Technique	: GC-capillary
Column	: Agilent CP-Wax 52 CB, 0.53 mm x 25 m fused silica WCOT CP-Wax 52 CB (2.0 μ m) (Part no. CP7658)
Temperature	: 50 °C $\rightarrow$ 200 °C, 10 °C/min
Carrier Gas	: N ₂ , 47 cm/s (10 mL/min)
Injector	: Direct T = 250 °C
Detector	: FID, 100 x 10 ⁻¹² Afs T = 275 °C
Sample Size	: 0.2 μ L

Peak identification

1. diethylether
2. diisopropylether
3. methylal
4. diisobutylether
5. acetal (1,1-diethoxyethane)
6. dibutylether
7. 1,4-dioxane

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

A00180



Agilent Technologies