



# Enantiomers

## Analysis of $\gamma$ -lactones

### Application Note

Materials Testing & Research

#### Authors

Agilent Technologies, Inc.

#### Introduction

Gas chromatography with an Agilent CP-Chirasil-DEX CB column separates enantiomers of eight gamma-lactones in 35 minutes.



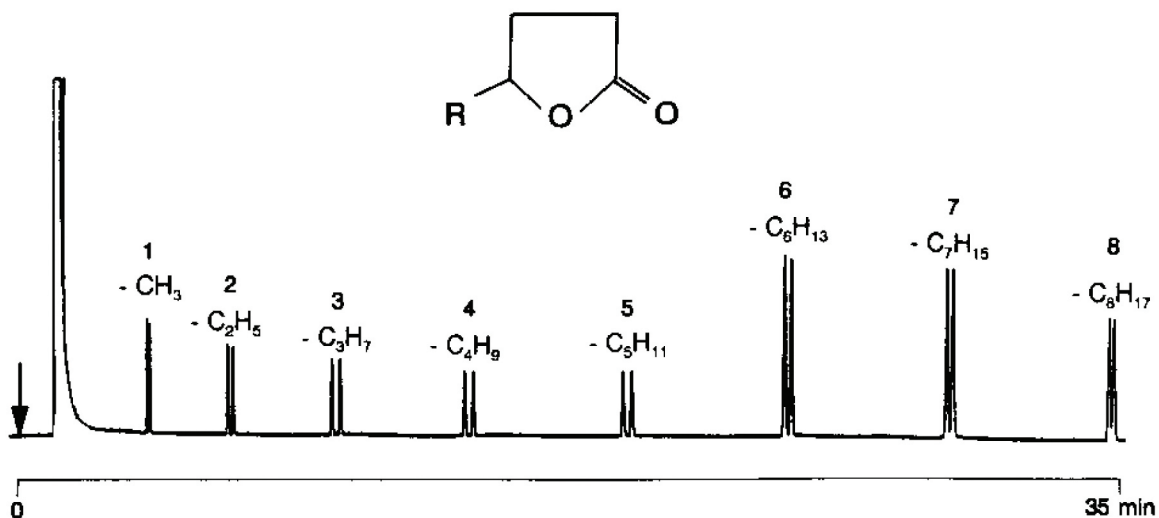
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-Chirasil-DEX CB, 0.25 mm x 25 m fused silica WCOT CP-Chirasil-DEX CB (df = 0.25 µm) (Part no. CP7502)  
Temperature : 100 °C (1 min) → 180 °C, 2 °C/min  
Carrier Gas : H<sub>2</sub>, 65 kPa (0.65 bar, 9.3 psi)  
Injector : Split, 100 mL/min  
T = 250 °C  
Detector : FID  
T = 275 °C

## Peak identification

1. (±)-γ-methyl-γ-butyrolactone
2. (±)-γ-ethyl-γ-butyrolactone
3. (±)-γ-propyl-γ-butyrolactone
4. (±)-γ-butyl-γ-butyrolactone
5. (±)-γ-pentyl-γ-butyrolactone
6. (±)-γ-hexyl-γ-butyrolactone
7. (±)-γ-heptyl-γ-butyrolactone
8. (±)-γ-octyl-γ-butyrolactone



[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

A00969



**Agilent Technologies**