



α -Hydroxyacids

Enantiomer separation of hydroxyacids as dioxolanone derivatives

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Enantiomeric separation of eight α -hydroxyacids (dioxolanone derivatives) by gas chromatography with an Agilent CP-Cyclodextrin-B-2,3,6-M-19 column is accomplished in 30 minutes.



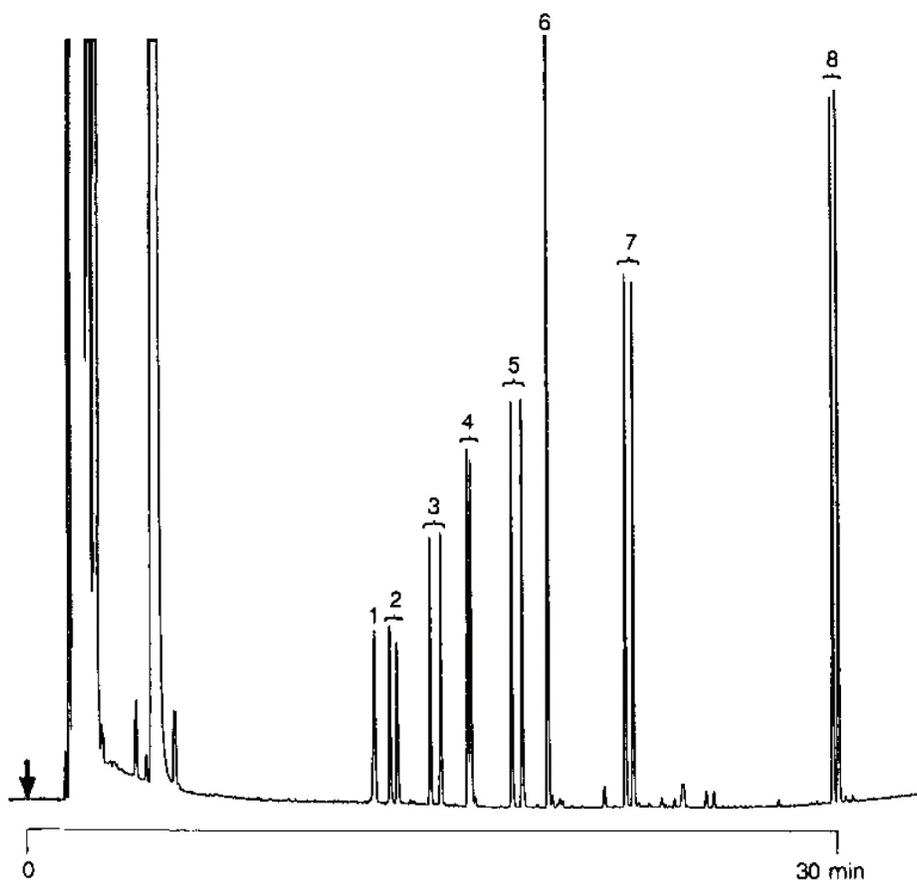
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Cyclodextrin-B-2,3,6-M-19, 0.25 mm x
25 m fused silica WCOT Cyclodextrin-B-2,3,6-M-19
(df = 25 μ m) (Part no. CP7500)
Temperature : 70 °C, 5 min + 5 °C/min \rightarrow 200 °C
Carrier Gas : H₂, 60 kPa (0.6 bar, 8.6 psi)
Injector : Split, 100 mL/min
T = 250 °C
Detector : FID, 4 x 10⁻¹² Afs
T = 275 °C
Concentration Range : 0.01% in hexane

Peak identification

1. glycolic acid
2. lactic acid
3. 2-hydroxy butanoic acid
4. 2-hydroxy-3-methyl butanoic acid
5. 2-hydroxy pentanoic acid
6. 2-hydroxy isohexanoic acid
7. 2-hydroxy hexanoic acid
8. 2-hydroxy octanoic acid



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

A00499



Agilent Technologies