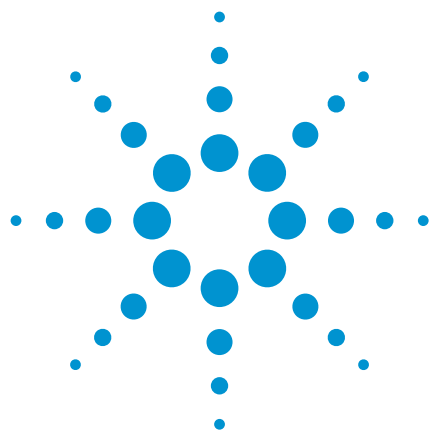




Fatty acid methyl esters, $C_{14} - C_{22}$

Analysis of rapeseed oil

Application Note

Food Testing & Agriculture

Authors

Agilent Technologies, Inc.

Introduction

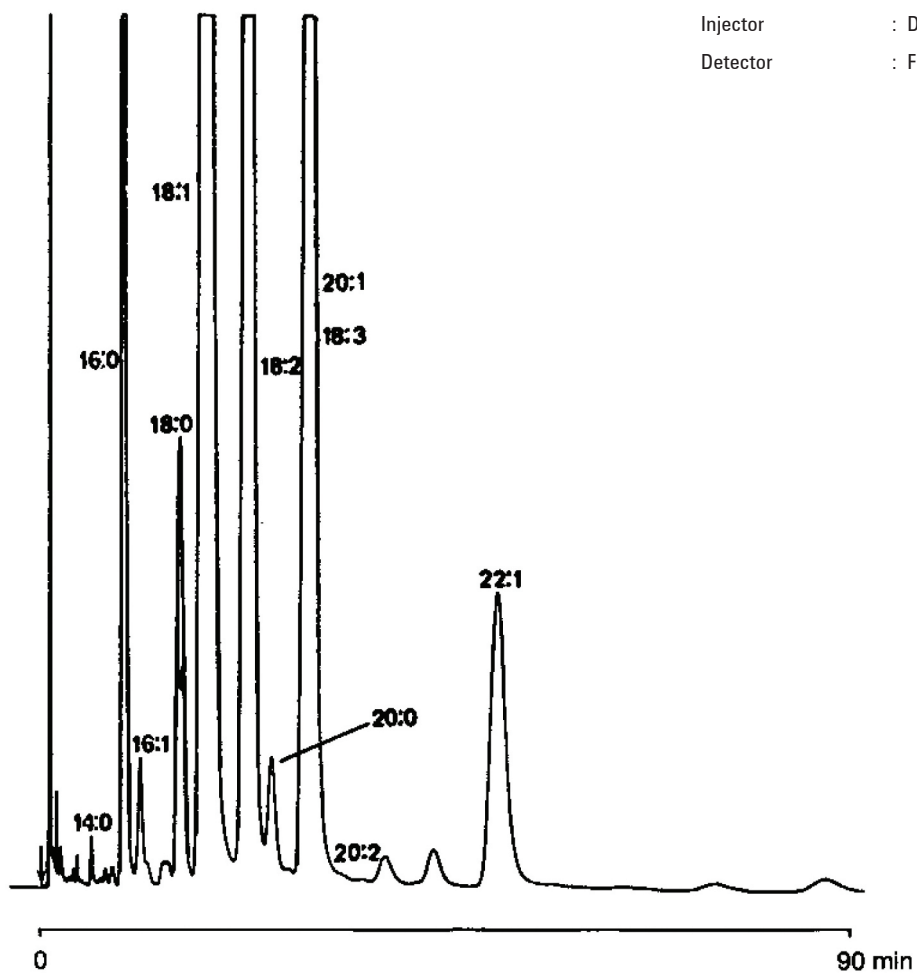
Gas chromatography with packed, custom-made Agilent CP-Sil 84 separates C_{14} to C_{22} fatty acid methyl esters in rapeseed oil in 90 minutes.



Agilent Technologies

Conditions

Technique : GC-packed
Column : Agilent CP-Sil 84, 1/8 in x 2 m SS 15% CP-Sil 84
Chromosorb WHP 100/120 (Part no. CP3601)
Temperature : 185 °C
Carrier Gas : N₂, 12 mL/min
Injector : Direct
Detector : FID, 16 x 10 Afs



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

A00202



Agilent Technologies