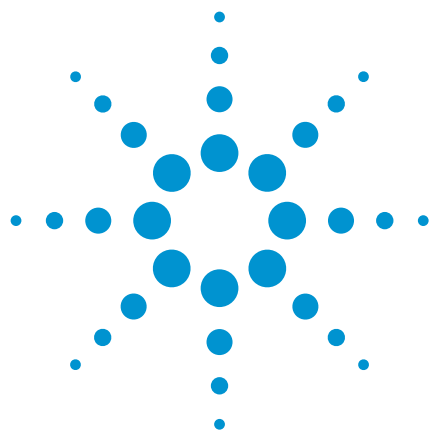




Fatty acid methyl esters, $C_{12} - C_{24}$

Analysis of fish oil on a wide-bore fused silica capillary column

Application Note

Food Testing & Agriculture

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 52 CB column separates 21 C_{12} to C_{24} fatty acid methyl esters in fish oil in 20 minutes.



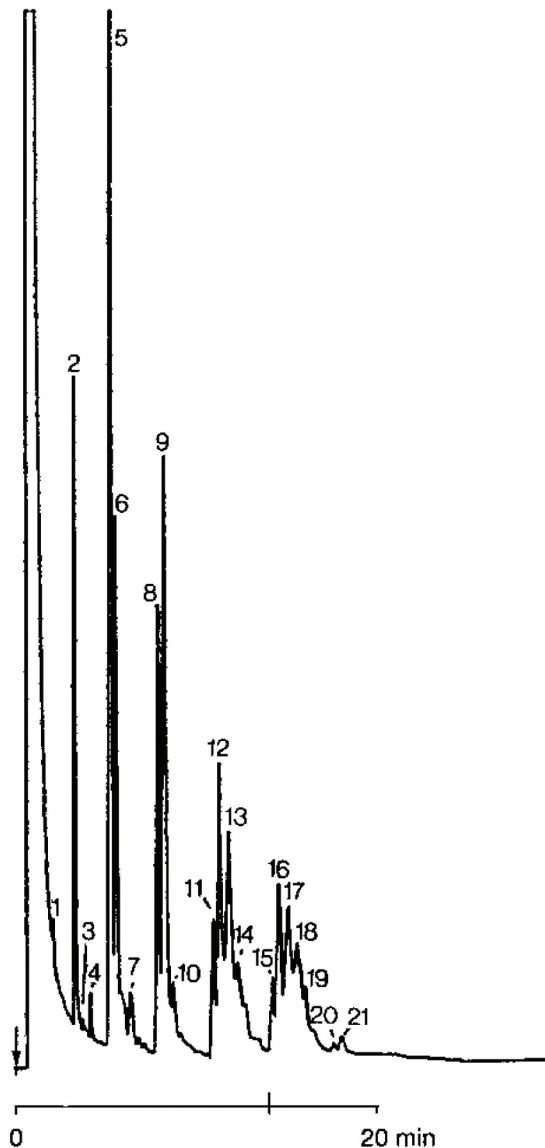
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 52 CB, 0.53 mm x 10 m fused silica
WCOT CP-Sil 52 CB (2.0 μ m) (Part no. CP7648)
Temperature : 180 °C $\rightarrow$ 245 °C, 4 °C/min
Carrier Gas : N₂, 25 cm/s (3 mL/min)
Injector : Direct
T = 260 °C
Detector : FID, 50 x 10⁻¹² Afs
T = 280 °C
Sample Size : 0.1 μ L

Peak identification

1. C12:0
2. C14:0
3. C14:1
4. C15:0
5. C16:0
6. C16:1
7. C17:0
8. C18:0
9. C18:1
10. C18:2
11. C20:0
12. C20:1
13. C20:2
14. C20:3
15. C22:0
16. C22:1
17. C22:2
18. C22:3
19. C22:4
20. C24:0
21. C24:1



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

A00169



Agilent Technologies