



Triglycerides

Analysis of butter fat

Application Note

Food Testing & Agriculture

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-TAP CB for Glycerides column separates 11 components in butter fat in 29 minutes.



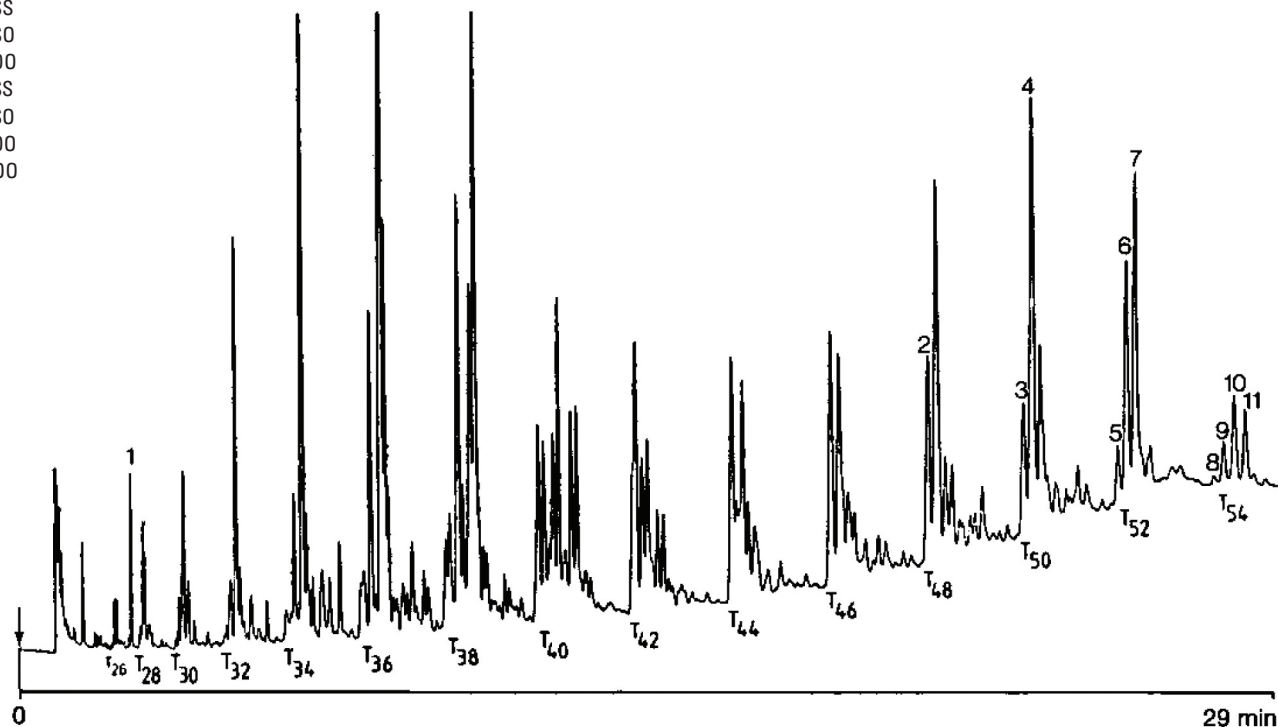
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-TAP CB, 0.25 mm x 25 m WCOT fused silica coated with TAP (0.10 μ m) (Part no. CP7483)
Temperature : 280 °C (1 min) $\rightarrow$ 355 °C, 3 °C/min
Carrier Gas : H₂, 100 kPa (1 bar, 15 psi)
Injector : On-column
Injection : 0.2 μ L of 0.05% butter fat in hexane
Detector : FID

Peak identification

1. cholesterol	P : palmitic acid, hexadecanoic acid	C16:0
2. PPP	S : stearic acid, octadecanoic acid	C18:0
3. PPS	O : oleic acid, cis-9-octadecenoic acid	C18:1
4. PPO		
5. PSS		
6. PSO		
7. POO		
8. SSS		
9. SSO		
10. SOO		
11. OOO		



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

A00223



Agilent Technologies