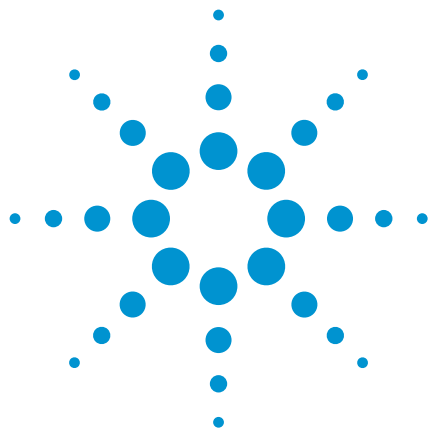




Triglycerides

Analysis of triglycerides in fat and oils

Application Note

Food Testing & Agriculture

Authors

Agilent Technologies, Inc.

Introduction

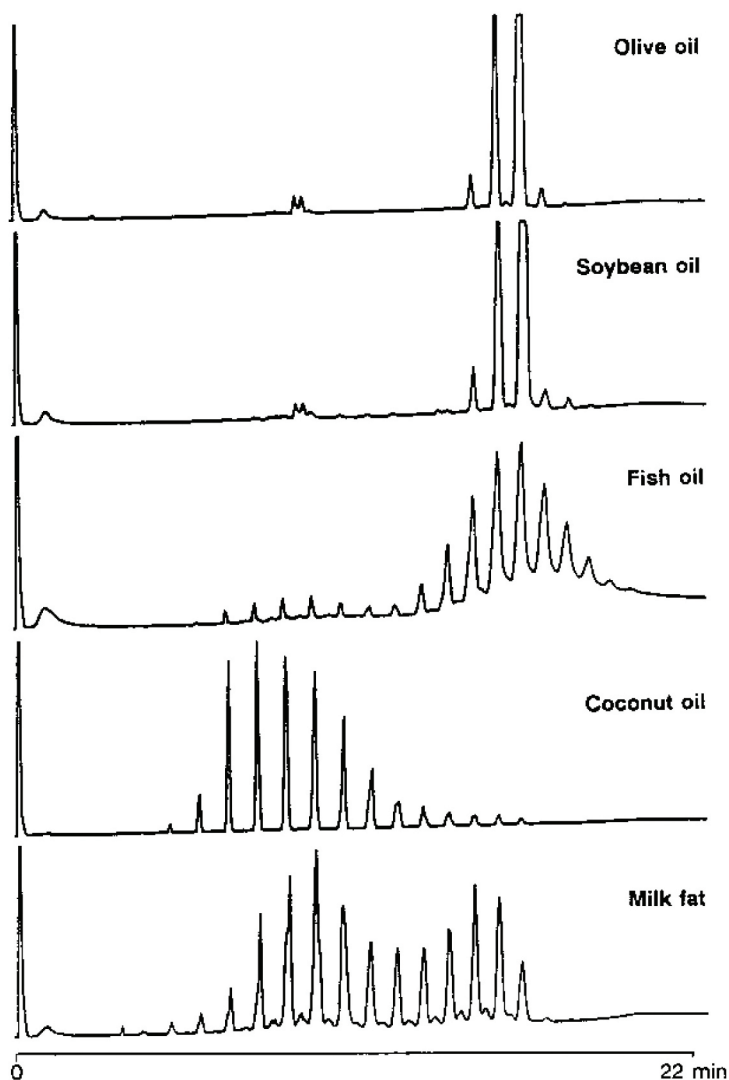
Gas chromatography with an Agilent CP-SimDist column separates triglycerides in olive, soyabean, fish, and coconut oils, and milk fat, in 22 minutes.



Agilent Technologies

Conditions

Technique	: GC-capillary
Column	: Agilent CP-SimDist CB, 0.53 mm x 5 m fused silica WCOT CP-SimDist CB (df = 0.10 µm) (Part no. CP7522)
Temperature	: 200 °C (1 min) → 270 °C, 10 °C/min; 270 °C → 370 °C, 8 °C/min
Carrier Gas	: He
Injector	: PTV, 80 mL/min splitflow T = 390 °C
Detector	: FID T = 380 °C
Sample Size	: 0.6 µL
Concentration range	: 2.5% oil/fat in heptane
Courtesy	: Norweigan Dairies Association, Research Center, Voll, Norwar



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

A00756



Agilent Technologies