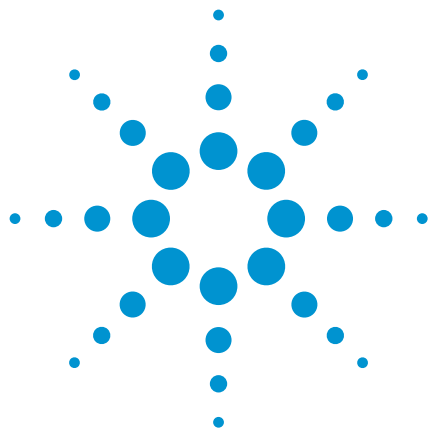




Fatty acids

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

HPLC with Agilent Pursuit C18 5 μm material in a ChromSep cartridge separates three fatty acids in four minutes.



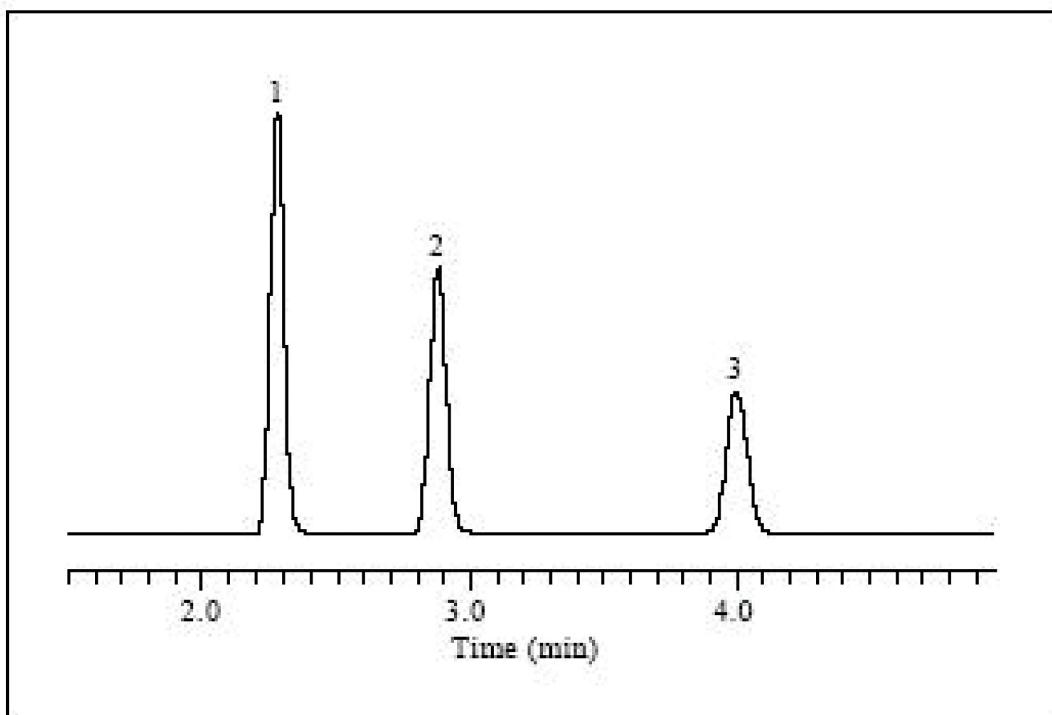
Agilent Technologies

Conditions

Technique : HPLC
Column : Agilent Pursuit C18, 4.6 x 150 mm, 5 μ m
Part no. : Conventional A3000150X046
ChromSep Cartridge System A3000150C046
Temperature : Ambient
Eluent : A: Water + 0.1% HCOOH
B: MeCN + 0.1% HCOOH
A:B = 5:95
Flow rate : 1.5 mL/min
Detection : UV @ 205 nm

Peak identification

1. linolenic acid
2. linoleic acid
3. oleic 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

A02317



Agilent Technologies