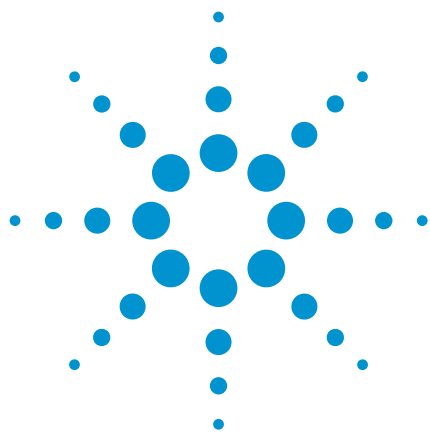




Fatty acids

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

HPLC with an Agilent Polaris C18-A column separates three 18-carbon fatty acids in under 52 seconds using an isocratic eluent.



Agilent Technologies

Conditions

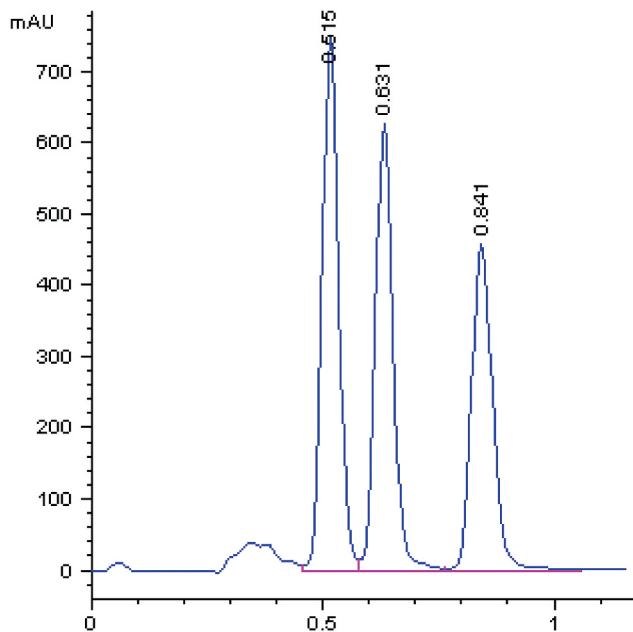
Column : Agilent Polaris C18-A, 4.6 x 30 mm, 3 μ m
Part no. : A2001030X046
Eluent : A: H₂O + 0.1% HCOOH B: MeCN + 0.1% HCOOH A:B = 5:95
Flow : 1.5 mL/min
Detection : UV @ 205 nm
Sample : 1. linolenic acid, 2. linoleic acid, 3. oleic acid
Injection Volume : 2 μ L

Three 18-carbon fatty acids are separated in under 52 seconds using an isocratic eluent.

oleic = CH₃(CH₂)₇CH=CH(CH₂)₇COOH

linoleic = CH₃(CH₂)₄CH=CHCH₂CH=CH(CH₂)₇COOH

linolenic = CH₃CH₂CH=CHCH₂CH=CHCH₂CH=CH(CH₂)₇COOH



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

A01936



Agilent Technologies