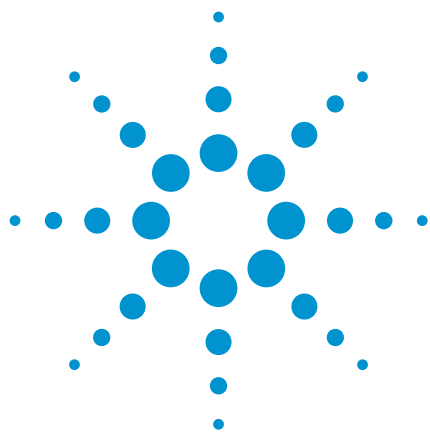




Tocopherols

Application Note

Food Testing & Agriculture

Authors

Agilent Technologies, Inc.

Introduction

HPLC using an Agilent Polaris NH2 column separates three tocopherols in just over four minutes.



Agilent Technologies

Conditions

Column : Agilent Polaris NH2, 4.6 x 150 mm, 5 µm
Part no. A2001030X046

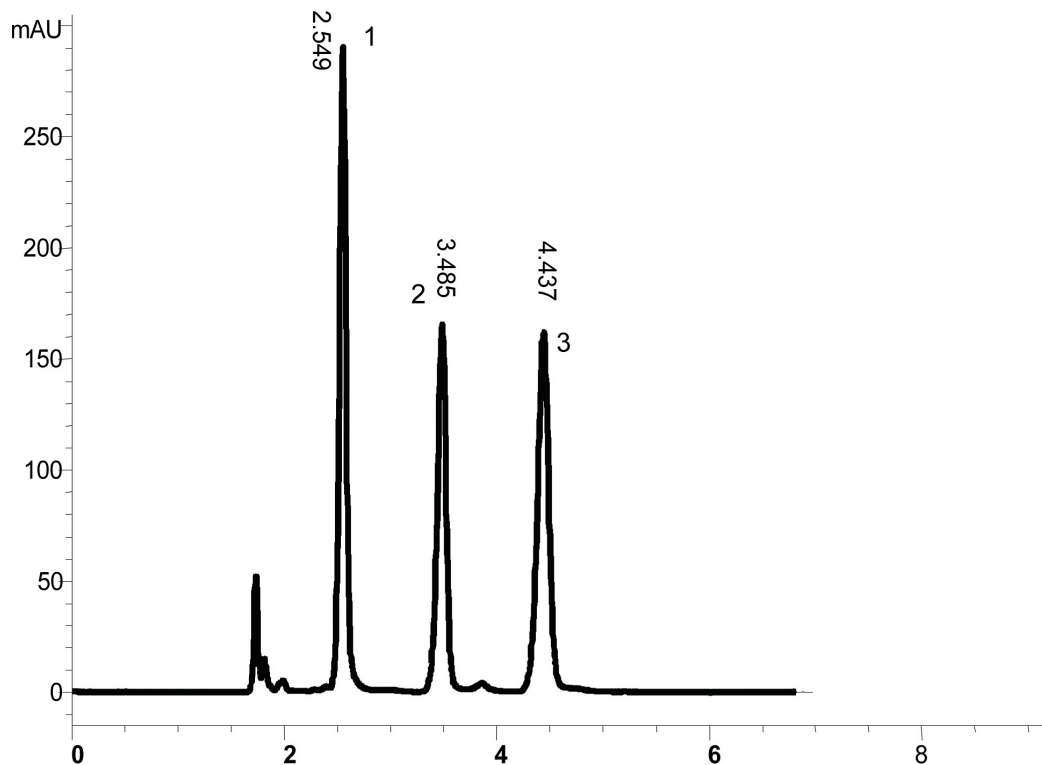
Eluent : A: Heptane
B: CH₂Cl₂ : MeOH 80:20

Flow : 1.0 mL/min

Detection : UV @ 295 nm

Peak identification

1. alpha-tocopherol
2. gamma-tocopherol
3. delta-tocopherol



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

A02000



Agilent Technologies