

Application Report 228

LC-MS Analysis of Fatty Acids Using Ascentis™ RP-Amide

This application demonstrates the suitability of the Ascentis RP-Amide for the analysis of cholesterol using an MS detector.

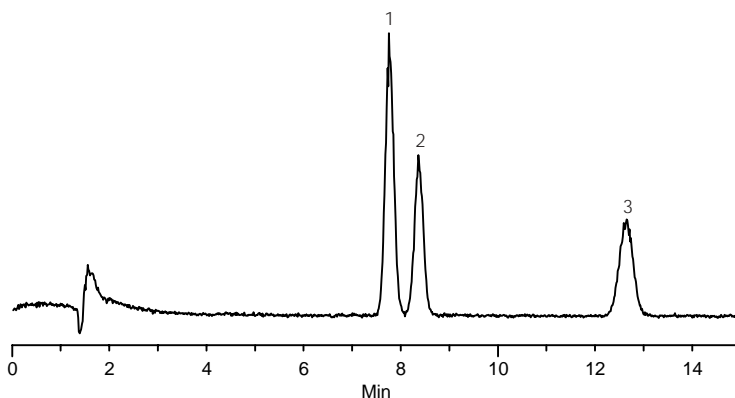
Key Words

Ascentis RP-Amide, 565324-U, palmitic acid, 57-10-3, 258725, stearic acid, 57-11-4, S4751, oleic acid, 112-80-1, O1008

Author: Hugh M. Cramer

Acquisition System: Waters Micro Mass ZQ

Notebook Reference: 1551



G002905

Conditions

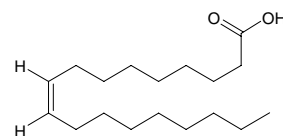
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 5:95, 0.1% formic acid in water:acetonitrile
flow rate: 1.0 mL/min., split to detector
temp.: 35 °C
det.: ESI(-)
injection: 5 µL
sample: 10 µg/mL in acetonitrile

Peak IDs

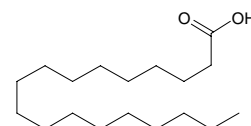
1. Oleic acid (m/z 281)
2. Palmitic acid (m/z 255)
3. Stearic acid (m/z 283)

Structures

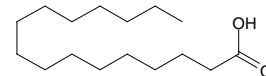
Oleic acid - G002902



Palmitic acid - G002903



Stearic acid - G002904



Ascentis is a trademark of Sigma-Aldrich Co.