

Application Report 229

Analysis of the Antihypercholesterolemic Mevinolin (Levostatin) and Degradation Products Using Ascentis™ C18

This application demonstrates the suitability of the Ascentis C18 for the analysis of the antihypercholesterolemic mevinolin and its degradation products using an MS detector.

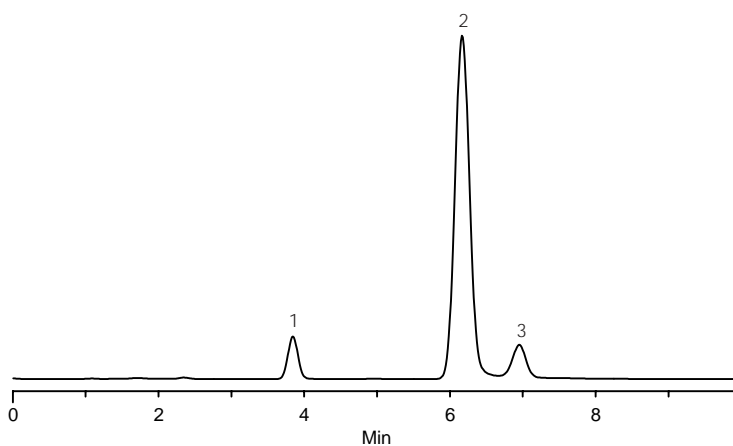
Key Words

Ascentis C18, 581324-U, mevinolin, levostatin, monacolin K, 75330-75-5, M2147, mevinolinic acid, acetonitrile LC-MS, 34967

Author: Hugh M. Cramer

Acquisition System: Waters Micro Mass ZQ

Notebook Reference: 1551



G002911

Conditions

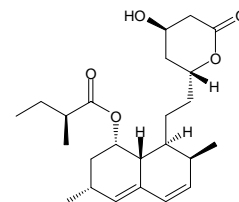
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)
mobile phase: 30:70, 10 mM ammonium formate (pH 3.0 with formic acid):acetonitrile
flow rate: 1.0 mL/min., split to detector
temp.: 35 °C
det.: ESI(+)
injection: 5 µL
sample: 10 µg/mL total in mobile phase

Peak IDs

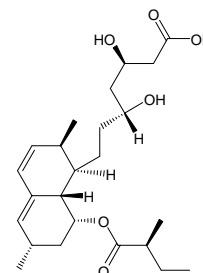
1. Mevinolinic acid (m/z 461, (M+K⁺))
2. Mevinolin (m/z 443, (M+K⁺))
3. Mevinolinic acid methyl ester (m/z 475, (M+K⁺))

Structures

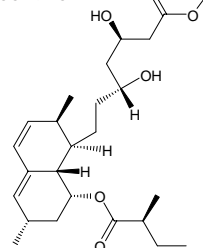
Mevinolinic acid - G002908



Mevinolin - G002909



Mevinolinic acid methyl ester - G002910



Ascentis is a trademark of Sigma-Aldrich Co.