

Application Report 122

LC-MS Analysis of Rifampicin Using Ascentis™ C18

This application demonstrates efficient analysis of rifampicin on Ascentis C18.

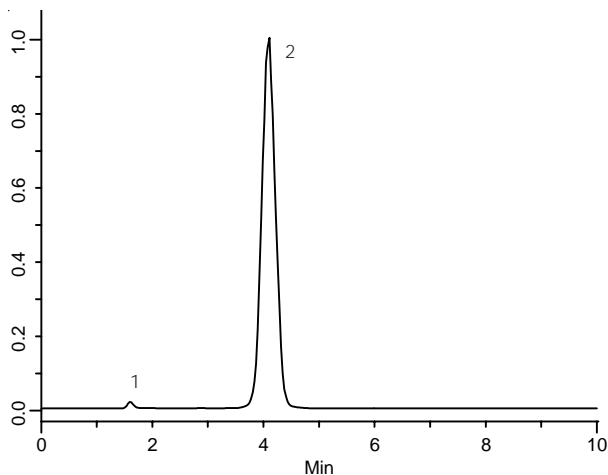
Key Words

Rifamycin AMP; rifampicin, rifampin, R3501; 13292-46-1; antibiotic; antibacterial tuberculostatic; LC-MS; Ascentis C18, 581303-U, mass spectrometry

Author: Hillel Brandes

Acquisition System: Thermo Surveyor & LCQAdvantage

Notebook Reference: 1498



G002447

Conditions

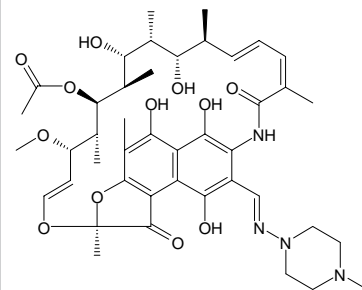
column: Ascentis C18, 5 cm x 2.1 mm I.D., 5 µm particles (581303-U)
mobile phase: 60:40, 33 mM formic acid:acetonitrile
flow rate: 0.2 mL/min.
temp.: 35 °C
det.: ESI (+), full scan
injection: 2 µL
sample: 10 mg/L in 50% methanol

Peak IDs

1. Rifampicin Impurity (m/z 821.1)
2. Rifampicin (m/z 823.1)

Structures

Rifampicin - G002448



Ascentis is a trademark of Sigma-Aldrich Co.