

Application Report 107

Analysis of Cephalosporins Cefoxitin and Cephalothin Using Ascentis™ C18

This application demonstrates the suitability of Ascentis C18 for the efficient separation of cefoxitin and cephalothin by HPLC.

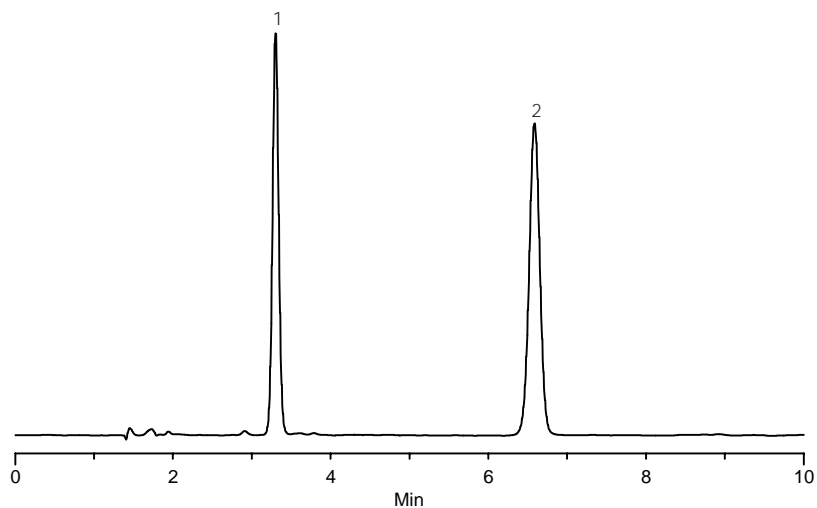
Key Words

Ascentis C18, cefoxitin, 33564-30-6, C4786, cephalothin, 58-71-9, C4520, 581324-U, cephalosporins

Author: Hugh Cramer

Acquisition System: Waters Alliance
2690 ID 9371

Notebook Reference: 1495-18



G002411

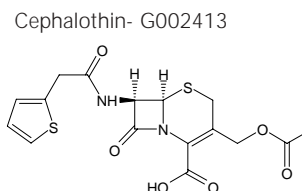
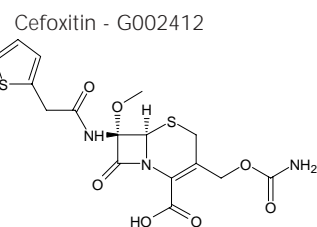
Conditions

column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)
mobile phase: 75:25, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):acetonitrile
flow rate: 1 mL/min.
temp.: 35 °C
det.: UV at 254 nm
injection: 10 µL
sample: 50 µg/mL each in mobile phase

Peak IDs

1. Cefoxitin
2. Cephalothin

Structures



Ascentis is a trademark of Sigma-Aldrich Co.