

Application Report 82

Analysis of Antibacterials Using Ascentis™ C18

This application demonstrates the suitability of Ascentis C18 for the analysis of an antibacterial mix consisting of chloramphenicol, nitrofurantoin and furaltadone. Structures along with the optimized chromatogram obtained on the Ascentis C18 are presented.

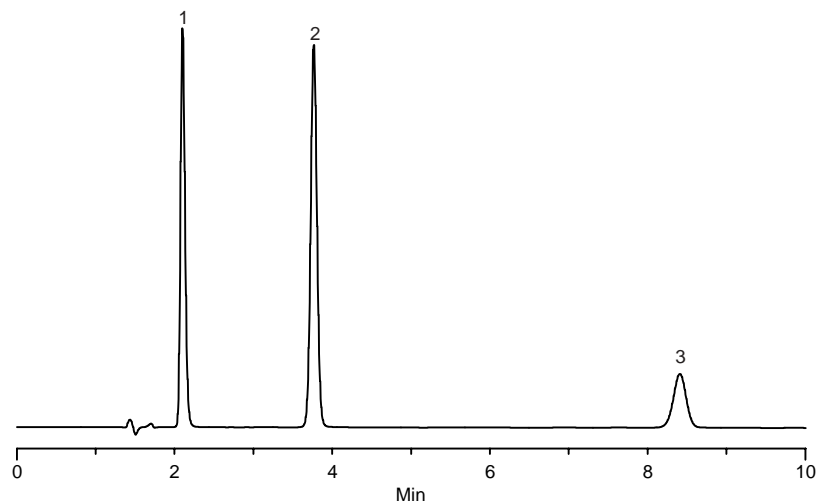
Key Words

chloramphenicol, 56-75-7, C-0378, nitrofurantoin, 67-20-9, N-7878, furaltadone, 139-91-3, F-9130, 581324-U

Author: Shawn R. Wyatt

Acquisition System: Lab 35, Waters Alliance

Notebook Reference: 1445-49



G002325

Conditions

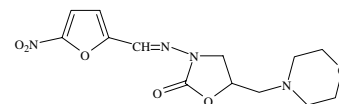
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)
mobile phase: 75:25 water, 10 mM ammonium formate (pH 3.0 with formic acid):acetonitrile
flow rate: 1.0 mL/min.
temp.: 35 °C
det.: UV, 254 nm
injection: 5 µL
sample: 50 µg/mL each in acetonitrile

Peak IDs

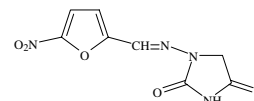
1. Furaltadone
2. Nitrofurantoin
3. Chloramphenicol

Structures

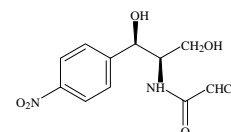
Furaltadone - G002324



Nitrofurantoin - G002223



Chloramphenicol - G002322



Ascentis is a trademark of Sigma-Aldrich Co.