

Application Report 205

LC-UV Analysis of the Antimicrobial Therapeutic Carbadox Using Ascentis™ RP-Amide Column

Carbadox, an antimicrobial drug and its major metabolite, quinoxaline-2 carboxylic acid are separated on an Ascentis RP-Amide column. Detection was at 320 nm

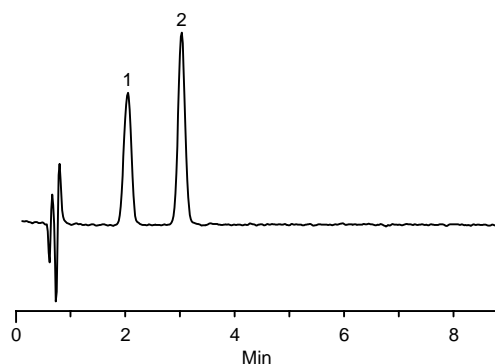
Key Words

carbadox, quinoxaline-2 carboxylic acid, 6804-07-5, 879-65-2, C6770, 29340-7, antimicrobial, metabolites, LC-UV Ascentis RP-Amide, 565323-U

Author: Carmen T. Santasania

Acquisition System: W2790

Notebook Reference: 1548-67



G002847

Conditions

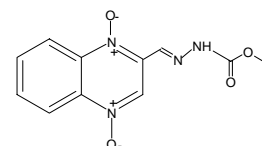
column: Ascentis RP-Amide, 5 cm x 4.6 mm I.D., 5 µm particles (565323-U)
mobile phase: 85:15, 10 mM ammonium formate (pH 2.5 with formic acid):acetonitrile
flow rate: 1.0 mL/min
temp.: 35 °C
det.: UV at 320 nm
injection: 5 µL
sample: carbadox (1 µg/mL) and quinoxaline-2 carboxylic acid (10 µg/mL) in 50:50 10 mM ammonium formate (pH 2.5):acetonitrile

Peak IDs

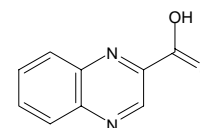
1. Carbadox
2. Quinoxaline-2 carboxylic acid

Structures

Carbadox - G002845



Quinoxaline-2 carboxylic acid - G002846



Ascentis is a trademark of Sigma-Aldrich Co.