

Application Report 344

LC-MS Analysis of Benzodiazepines Using Ascentis™ C8

Two benzodiazepines are separated on Ascentis C8 and detected by mass spectrometry.

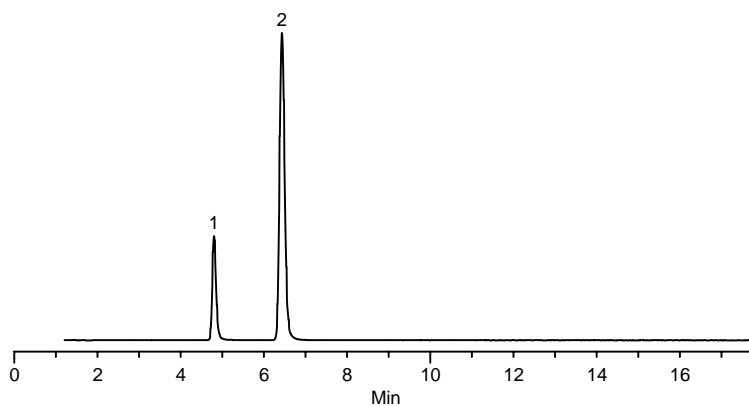
Key Words

LC-MS, Ascentis C8, drugs of abuse, 581402-U, lorazepam, flunitrazepam, 34967, benzodiazepines, F9261, 1622-62-4, L1764, 846-49-1, 581402-U

Author: Carmen T. Santasania

Acquisition System: W2795

Notebook Reference: 1570-13



G003313

Conditions

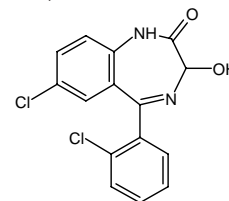
column: Ascentis C8, 15 cm x 2.1 mm I.D., 3 μm particles (581402-U)
mobile phase: 55:45, 10 mM ammonium acetate (pH 6.98 unadjusted):acetonitrile
flow rate: 0.2 mL/min.
temp.: 35 °C
det.: MS, ESI (+), Selected Ion Recording Mode
injection: 5 μL
sample: 1 μg/mL each in 10 mM ammonium acetate (pH unadjusted)

Peak IDs

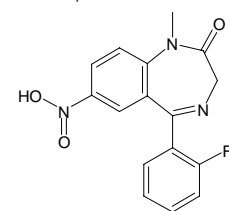
1. Lorazepam (M+H)⁺= 321.01
2. Flunitrazepam (M+H)⁺= 314.08

Structures

Lorazepam - G003312



Flunitrazepam - G003311



Ascentis is a trademark of Sigma-Aldrich Co.