

Application Report 90

Analysis of Nitrazepam Anticonvulsant Using Ascentis™ C18

This application demonstrates the suitability of Ascentis C18 for the efficient elution of the anticonvulsant nitrazepam by HPLC.

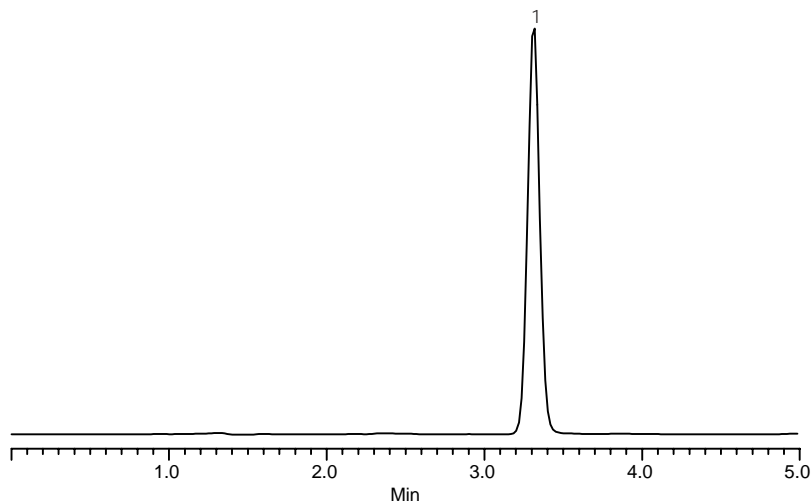
Key Words

Ascentis C18, nitrazepam, 146-22-5, 581324-U, anticonvulsant

Author: Hugh Cramer

Acquisition System: Waters Alliance
2690 ID 9371

Notebook Reference: 1486-61



G002356

Conditions

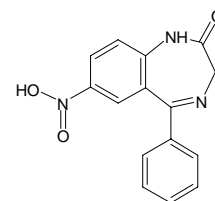
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)
mobile phase: 50:50, 10 mM monobasic potassium phosphate (pH 7.0 with potassium hydroxide):acetonitrile
flow rate: 1 mL/min.
temp.: 35 °C
det.: UV at 220 nm
injection: 10 µL
sample: 50 µg/mL each in mobile phase

Peak IDs

1. Nitrazepam

Structures

Nitrazepam - G002357



Ascentis is a registered trademark of Sigma-Aldrich Co.