

Application Report 218

Analysis of the Anticonvulsants Using Ascentis™ C18

This application demonstrates the suitability of the Ascentis C18 for the analysis of the anticonvulsants clonazepam, clorazepate, and diazepam.

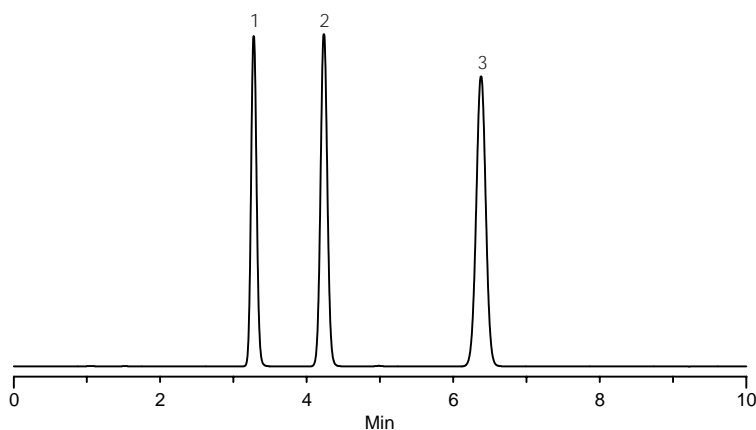
Key Words

clonazepam, 1622-61-3, C4404, clorazepate, 57109-90-7, diazepam, 493-14-5, D9900, Ascentis C18, 581324-U

Author: Hugh M. Cramer

Acquisition System: Mt McKinley
Thermo

Notebook Reference: 1551



G002882

Conditions

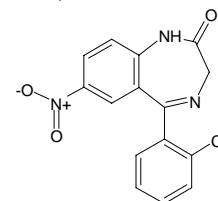
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)
mobile phase: 50:50, water:acetonitrile
flow rate: 1.0 mL/min.
temp.: 35 °C
det.: UV at 254 nm
injection: 10 µL
sample: as indicated in mobile phase

Peak IDs

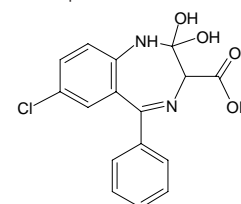
1. Clonazepam (50 µg/mL)
2. Clorazepate (50 µg/mL)
3. Diazepam (50 µg/mL)

Structures

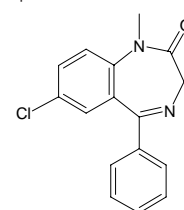
Clonazepam - G002879



Clorazepate - G002880



Diazepam - G002881



Ascentis is a trademark of Sigma-Aldrich Co.