

Application Report 303

Analysis of Anticonvulsants Using Ascentis™ C8

This application demonstrates the suitability of the Ascentis C8 for the analysis of the anticonvulsants felbamate and lamotrigine. Structures along with the optimized chromatogram obtained on the Ascentis C8 are presented.

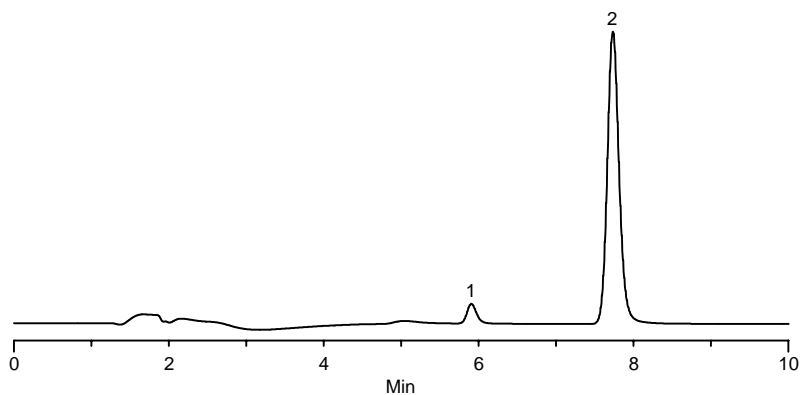
Key Words

felbamate, 25451-15-4, F0778, lamotrigine, 84057-84-1, L3791, Ascentis C8, 581402-U

Author: Hugh M. Cramer

Acquisition System: Shimadzu
Barkaboom Mountain

Notebook Reference: 1558



G003163

Conditions

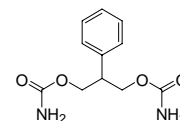
column: Ascentis C8, 15 cm x 2.1 mm I.D., 3 µm particles (581402-U)
mobile phase: 60:40, water with 0.1% ammonium acetate (34674):acetonitrile with 0.1% ammonium acetate (34669)
flow rate: 0.2 mL/min.
temp.: 30 °C
det.: UV at 220 nm
injection: 2 µL
sample: as indicated in mobile phase

Peak IDs

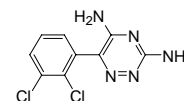
1. Felbamate (50 µg/mL)
2. Lamotrigine (50 µg/mL)

Structures

Felbamate - G003140



Lamotrigine - G003139



Ascentis is a trademark of Sigma-Aldrich Co.