

Application Report 170

Analysis of Calcium Channel Blocker Verapamil and Methoxy Metabolite Using Ascentis™ RP-Amide

This application demonstrates the suitability of Ascentis RP-Amide for the efficient separation of (+/-)verapamil and methoxyverapamil acid by HPLC.

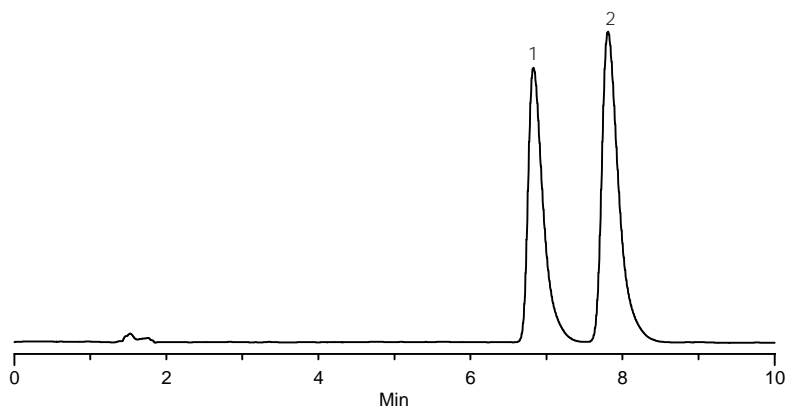
Key Words

Ascentis RP-Amide, (+/-)verapamil, 23313-68-0, V4629, methoxyverapamil, 16662-47-8, M5644, 565324-U

Author: Hugh Cramer

Acquisition System: Shimadzu
Mongaup

Notebook Reference: 1551



G002665

Conditions

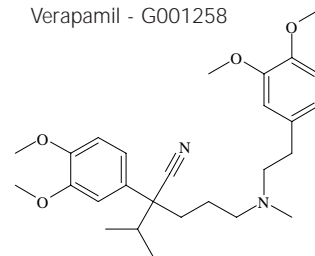
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 70:30, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):acetonitrile
flow rate: 1 mL/min.
temp.: 35 °C
det.: UV, 220 nm
injection: 10 µL
sample: 50 µg/mL each in mobile phase

Peak IDs

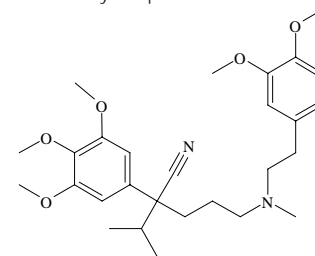
1. (+/-)verapamil
2. Methoxyverapamil

Structures

Verapamil - G001258



Methoxyverapamil - G002141



Ascentis is a trademark of Sigma-Aldrich Co.