

Application Report 116

Analysis of Calcium Channel Blocker Verapamil and Methoxy Metabolite using Ascentis™ C18

This application demonstrates the suitability of Ascentis C18 for the efficient separation of (+/-) -verapamil and methoxyverapamil acid by HPLC.

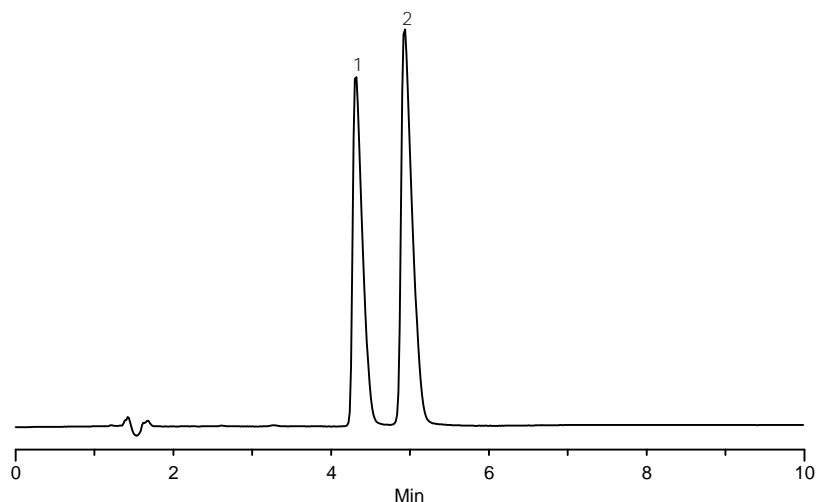
Key Words

Ascentis C18, (+/-)verapamil, 23313-68-0, V4629, methoxyverapamil, 16662-47-8, M5644, 581324-U

Author: Hugh Cramer

Acquisition System: Waters Alliance
2690 ID 9371

Notebook Reference: 1486-92



G002437

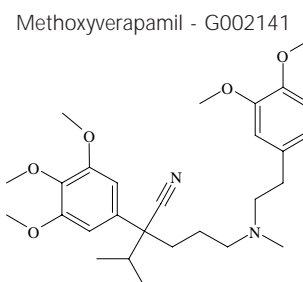
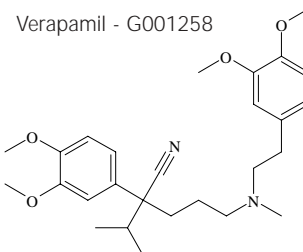
Conditions

column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)
mobile phase: 60:40, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):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. (+/-)verapamil
2. Methoxyverapamil

Structures



Ascentis is a trademark of Sigma-Aldrich Co.