

Application Report 305

Analysis of Cardiac Drugs Using Ascentis™ C8

This application demonstrates the suitability of the Ascentis C8 for the analysis of cardiac drugs. Structures along with the optimized chromatogram obtained on the Ascentis C8 are presented.

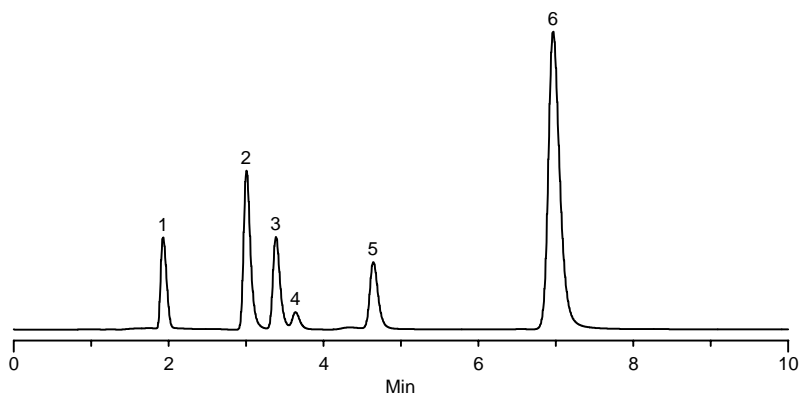
Key Words

acebutolol, 34381-68-5, A3669, metoprolol, 56392-17-7, M5391, timolol, 26921-17-5, T6394, oxprenolol, 6452-73-9, propranolol, 318-98-9, 222984, Ascentis C8, 581402-U

Author: Hugh M. Cramer

Acquisition System: Shimadzu
Barkaboom Mountain

Notebook Reference: 1558



G003172

Conditions

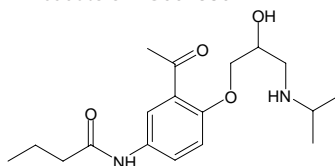
column: Ascentis C8, 15 cm x 2.1 mm I.D., 3 µm particles (581402-U)
mobile phase: 47:53, water with 0.1% ammonium acetate (34674):acetonitrile with 0.1% ammonium acetate (34669)
flow rate: 0.2 mL/min.
temp.: 30 °C
det.: UV at 220 nm
injection: 2 µL
sample: as indicated in mobile phase

Peak IDs

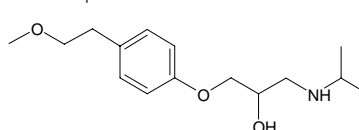
1. Maleic acid (50 µg/mL)
2. Acebutolol (50 µg/mL)
3. Metoprolol (50 µg/mL)
4. Timolol (50 µg/mL)
5. Oxprenolol (50 µg/mL)
6. Propranolol (50 µg/mL)

Structures

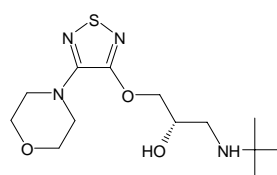
Acebutolol - G002380



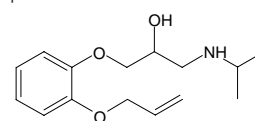
Metoprolol - G002370



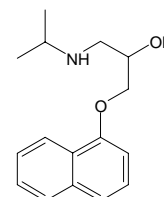
Timolol - G003170



Oxprenolol - G003171



Propranolol - G002373



Ascentis is a trademark of Sigma-Aldrich Co.