

Application Report 96

Analysis of Antihypertensives Using Ascentis™ C18

This application demonstrates the suitability of Ascentis C18 for the efficient separation of the antihypertensives acebutolol, alprenolol, labetalol, and pindolol by HPLC.

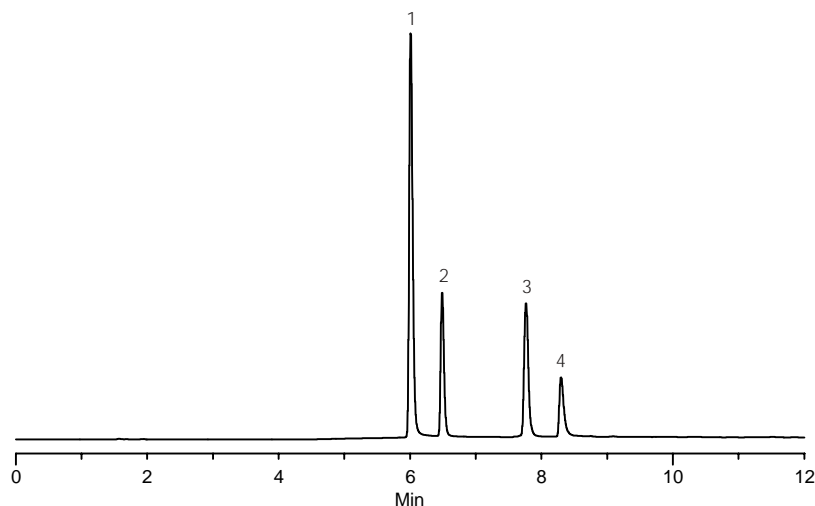
Key Words

Ascentis C18, acebutolol, 34381-68-5, A3669, alprenolol, 13707-88-5, A8676, labetalol, 32780-64-6, L1011, pindolol, 13523-86-9, 28410-6, 581324-U, antihypertensive

Author: Hugh Cramer

Acquisition System: Waters Alliance
2690 ID 9371

Notebook Reference: 1486-74



G002383

Conditions

column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)
mobile phase A: 10 mM monobasic potassium phosphate (pH 7.0 with potassium hydroxide)
mobile phase B: 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
gradient:

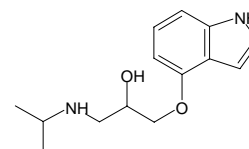
Min.	%A	%B
0	90	10
2	90	10
8	40	60
12	40	60

Peak IDs

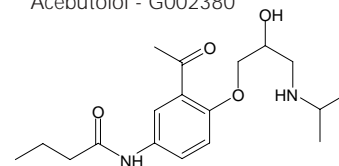
1. Pindolol
2. Acebutolol
3. Labetalol
4. Alprenolol

Structures

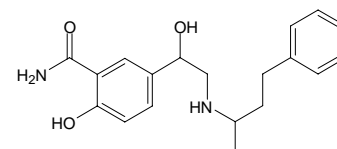
Pindolol - G002379



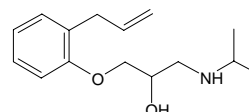
Acebutolol - G002380



Labetalol - G002384



Alprenolol - G002382



Ascentis is a registered trademark of Sigma-Aldrich Co.