

Application Report 137

Analysis of Acebutolol and Pindolol Antihypertensives Using Ascentis™ RP-Amide

This application demonstrates the suitability of Ascentis RP-Amide for the efficient separation of the antihypertensives acebutolol and pindolol by HPLC.

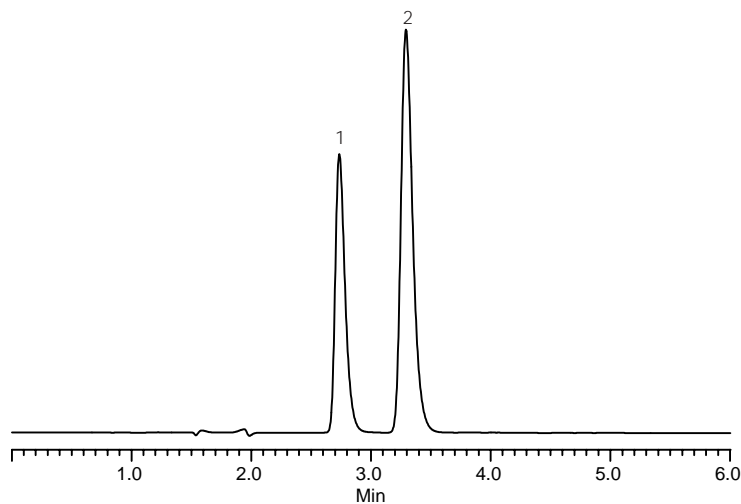
Key Words

Ascentis RP-Amide, acebutolol, 34381-68-5, A3669, pindolol, 13523-86-9, 28410-6, 565324-U, antihypertensive

Author: Hugh Cramer

Acquisition System: Shimadzu
Mongaup

Notebook Reference: 1551



G002600

Conditions

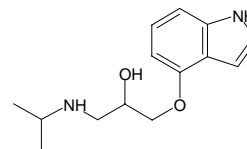
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 80:20, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):acetonitrile
flow rate: 1 mL/min.
temp.: 35 °C
det.: UV, 254 nm
injection: 10 µL
sample: 50 µg/mL each in mobile phase

Peak IDs

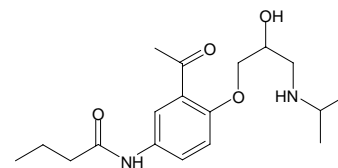
1. Pindolol
2. Acebutolol

Structures

Pindolol - G002379



Acebutolol - G002380



Ascentis is a trademark of Sigma-Aldrich Co.