

Application Report 94

Analysis of Antihypertensive Propranolol and Hydroxy Metabolite Using Ascentis™ C18

This application demonstrates the suitability of Ascentis C18 for the efficient separation of the antihypertensives 4-hydroxypropranolol and propranolol by HPLC.

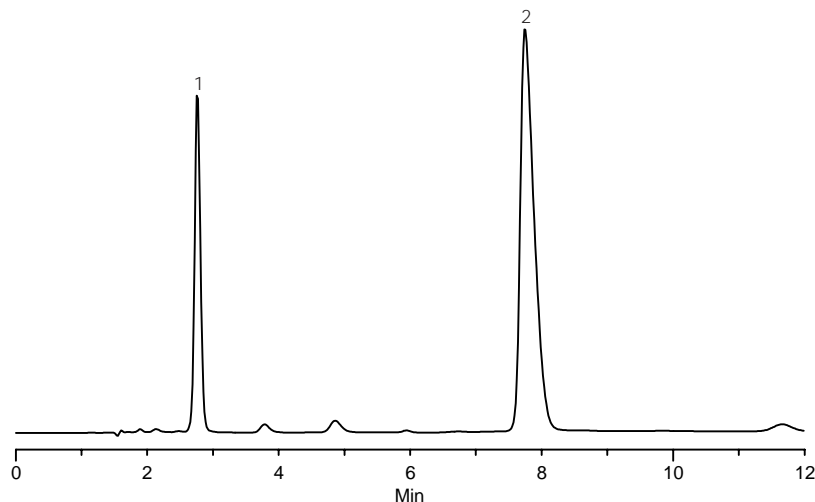
Key Words

Ascentis C18, 4-hydroxypropranolol, propranolol, 318-98-9, 22,298-4, 581324-U, antihypertensive

Author: Hugh Cramer

Acquisition System: Waters Alliance
2690 ID 9371

Notebook Reference: 1486-8?



G002372

Conditions

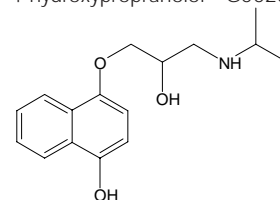
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)
mobile phase: 55:45, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):methanol
flow rate: 1 mL/min.
temp.: 35 °C
det.: UV at 220 nm
injection: 10 µL
sample: 50 µg/mL each in mobile phase

Peak IDs

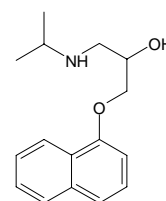
1. 4-hydroxypropranolol
2. Propranolol

Structures

4-hydroxypropranolol - G002374



Propranolol - G002373



Ascentis is a registered trademark of Sigma-Aldrich Co.