

# Application Report 95

## Analysis of Acebutolol and Pindolol Antihypertensives Using Ascentis™ C18

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

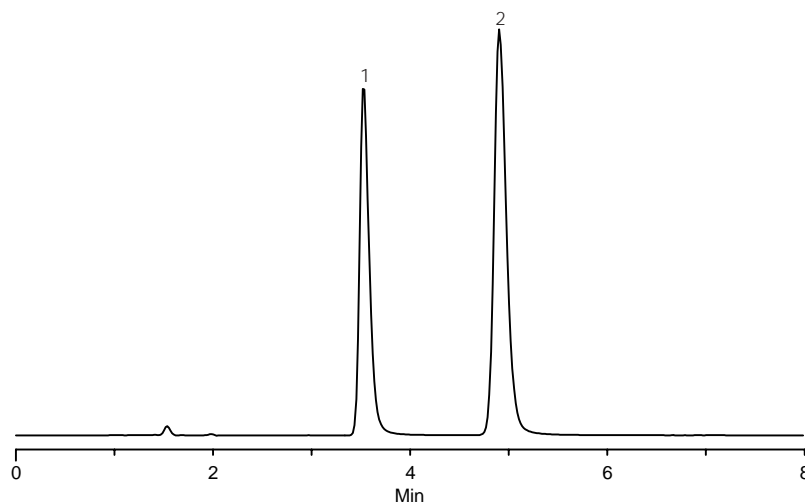
### Key Words

Ascentis C18, acebutolol, 34381-68-5, A3669, pindolol, 13523-86-9, 28410-6, 581324-U, antihypertensive

Author: Hugh Cramer

Acquisition System: Waters Alliance  
2690 ID 9371

Notebook Reference: 1486-74



G002378

### Conditions

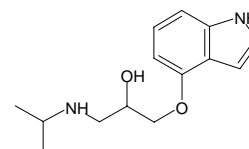
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)  
mobile phase: 80:20, 10 mM dibasic potassium phosphate (pH 7.0 with potassium hydroxide):acetonitrile  
flow rate: 1 mL/min.  
temp.: 35 °C  
det.: UV at 254 nm  
injection: 10 µL  
sample: 50 µg/mL each in mobile phase

### Peak IDs

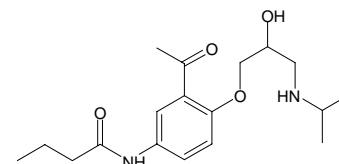
1. Pindolol
2. Acebutolol

### Structures

Pindolol - G002379



Acebutolol - G002380



Ascentis is a registered trademark of Sigma-Aldrich Co.