

# Application Report 93

## Analysis of Antihypertensive Metoprolol and Hydroxy Metabolite Using Ascentis™ C18

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

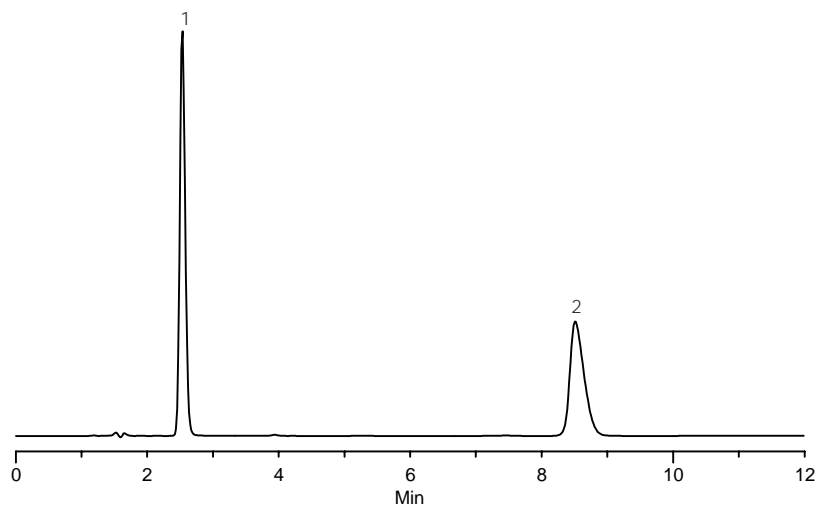
### Key Words

Ascentis C18, hydroxymetoprolol, metoprolol, 56392-17-7, M5391, 581324-U, antihypertensive

Author: Hugh Cramer

Acquisition System: Waters Alliance  
2690 ID 9371

Notebook Reference: 1486-82



G002369

### Conditions

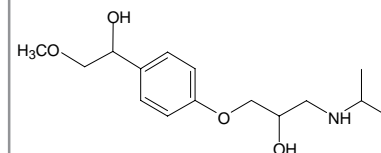
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)  
mobile phase: 70:30, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):methanol  
flow rate: 1 mL/min.  
temp.: 35 °C  
det.: UV at 220 nm  
injection: 10 µL  
sample: 50 µg/mL each in mobile phase

### Peak IDs

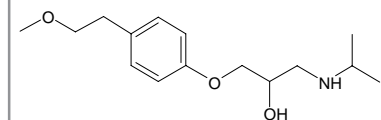
1. Hydroxymetoprolol
2. Metoprolol

### Structures

Hydroxymetoprolol - G002371



Metoprolol - G002370



Ascentis is a registered trademark of Sigma-Aldrich Co.