

# Application Report 91

## Analysis of the Antiemetic Meclizine Using Ascentis™ C18

This application demonstrates the suitability of Ascentis C18 for the efficient elution of the antiemetic meclizine by HPLC.

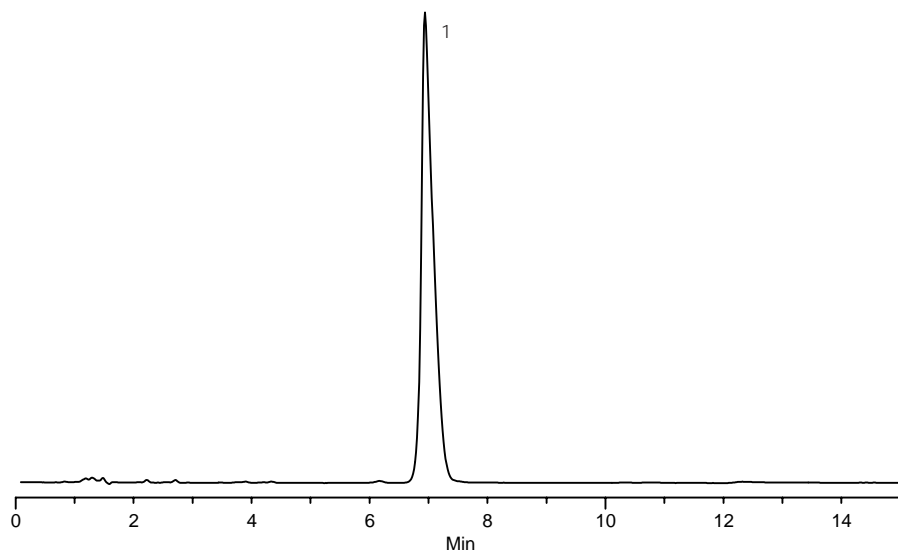
### Key Words

Ascentis C18, meclizine, 1104-22-9, M1637, 581324-U, antiemetic

Author: Hugh Cramer

Acquisition System: Waters Alliance  
2690 ID 9371

Notebook Reference: 1486-61



G002358

### Conditions

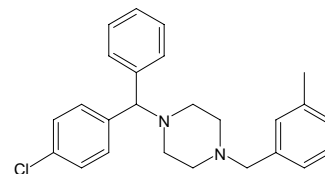
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)  
mobile phase: 50:50, 10 mM monobasic potassium phosphate  
(pH 3.0 with phosphoric acid):acetonitrile  
flow rate: 1 mL/min.  
temp.: 35 °C  
det.: UV at 230 nm  
injection: 10 µL  
sample: 50 µg/mL in mobile phase

### Peak IDs

1. Meclizine

### Structures

Meclizine - G002359



Ascentis is a registered trademark of Sigma-Aldrich Co.