

Application Report 92

Analysis of Antihistamines Using Ascentis™ C18

This application demonstrates the suitability of Ascentis C18 for the efficient separation of (+/-)-chlorpheniramine, (+/-)-brompheniramine, and triprolidine by HPLC.

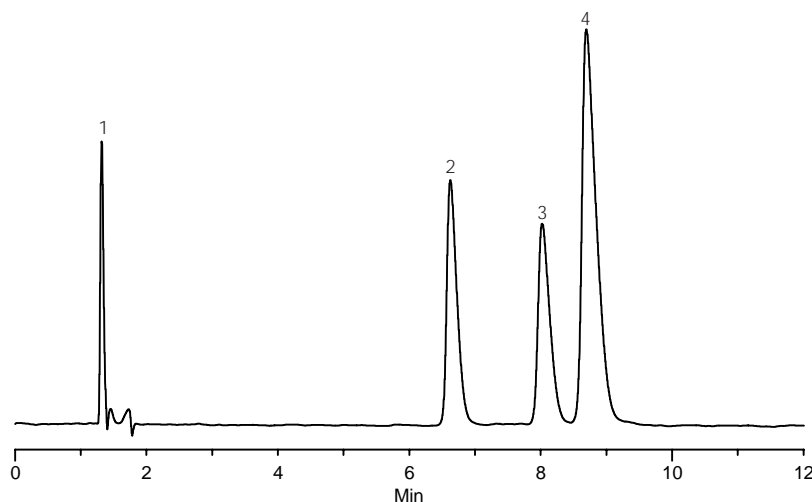
Key Words

Ascentis C18, 581324-U, (+/-)-chlorpheniramine, 7054-11-7, C3025, (+/-)-brompheniramine, 980-71-2, B2390, triprolidine, 6138-79-0, T6764, maleic acid

Author: Hugh Cramer

Acquisition System: Waters Alliance
2690 ID 9371

Notebook Reference: 1486-61



G002365

Conditions

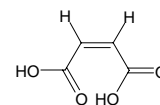
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)
mobile phase: 75:25, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):acetonitrile
flow rate: 1 mL/min.
temp.: 35 °C
det.: UV at 254 nm
injection: 10 µL
sample: 10 µg/mL each in mobile phase

Peak IDs

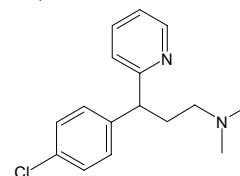
1. Maleic acid
2. (+/-)-chlorpheniramine
3. (+/-)-brompheniramine
4. Triprolidine

Structures

Maleic acid - G002520



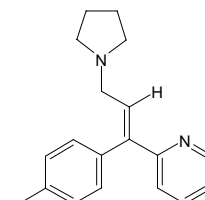
Chlorpheniramine - G002368



Brompheniramine - G002367



Triprolidine - G002366



Ascentis is a registered trademark of Sigma-Aldrich Co.