

Application Report 301

Analysis of Antiasthmatics Using Ascentis™ C8

This application demonstrates the suitability of the Ascentis C8 for the analysis of antiasthmatics. Structures along with the optimized chromatogram obtained on the Ascentis C8 are presented.

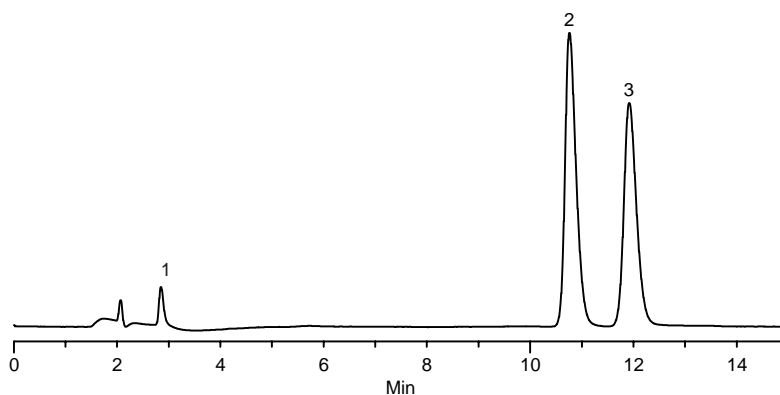
Key Words

formoterol, 73573-87-2, F9552, clenbuterol, 37148-27-9, C5423, Ascentis C8, 581402-U, fumaric acid

Author: Hugh M. Cramer

Acquisition System: Shimadzu
Barkaboom Mountain

Notebook Reference: 1558



G003160

Conditions

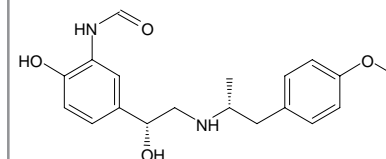
column: Ascentis C8, 15 cm x 2.1 mm I.D., 3 µm particles (581402-U)
mobile phase: 75:25, 10 mM ammonium phosphate (pH 2.5 with phosphoric acid):methanol
flow rate: 0.2 mL/min.
temp.: 35 °C
det.: UV at 220 nm
injection: 2 µL
sample: as indicated in mobile phase

Peak IDs

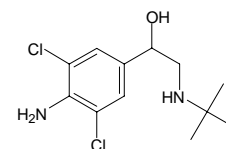
1. Fumaric acid
2. Clenbuterol (50 µg/mL)
3. Formoterol (50 µg/mL)

Structures

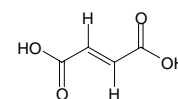
Formoterol - G003159



Clenbuterol - G002641



Fumaric acid - G003318



Ascentis is a trademark of Sigma-Aldrich Co.