

Application Report 302

Analysis of Organic Acids Using Ascentis™ C8

This application demonstrates the suitability of the Ascentis C8 for the analysis of the organic acids syringic and homoveratric. Structures along with the optimized chromatogram obtained on the Ascentis C8 are presented.

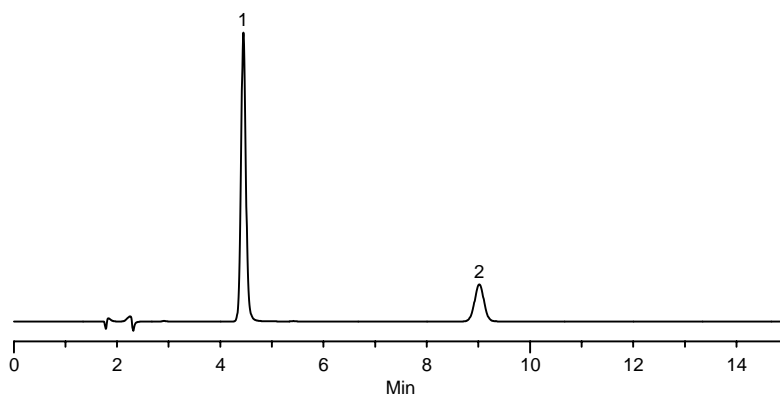
Key Words

syringic acid, 530-57-4, S6881, homoveratric acid, 93-40-3, 53650, Ascentis C8, 581424-U

Author: Hugh M. Cramer

Acquisition System: Hitachi Mt Fuji

Notebook Reference: 1558



G003160

Conditions

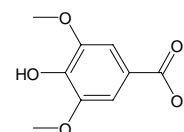
column: Ascentis C8, 15 cm x 4.6 mm I.D., 5 µm particles (581424-U)
mobile phase: 80:20, 10 mM ammonium phosphate monobasic (pH 2.50 with phosphoric acid):acetonitrile
flow rate: 1.0 mL/min.
temp.: 35 °C
det.: UV at 220 nm
injection: 10 µL
sample: as indicated in mobile phase

Peak IDs

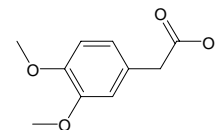
1. Syringic acid (50 µg/mL)
2. Homoveratric acid (50 µg/mL)

Structures

Syringic acid - G003017



Homoveratric acid - G003161



Ascentis is a trademark of Sigma-Aldrich Co.