

Application Report 146

Author: Jacinth A.M. McKenzie

Acquisition System: Hitachi LC

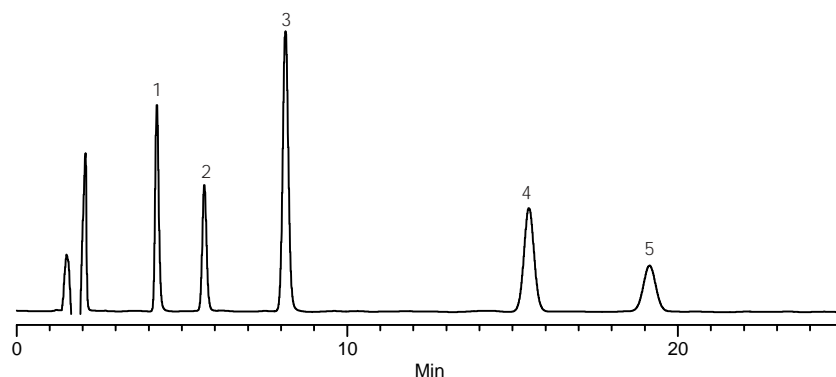
Notebook Reference: 1550-25

Analysis of Benzoic Acid Derivatives Using Ascentis™ RP-Amide

Ascentis RP-Amide is shown to be suitable for the efficient separation of organic acids, benzoic acid, 4-hydroxybenzoic acid, acetylsalicylic acid (aspirin), 2-hydroxybenzoic acid (salicylic acid) and ethyl paraben by HPLC-UV. Structures and a representative chromatogram are presented.

Key Words

Ascentis RP-Amide, 565324-U, organic acids, benzoic acid, 65-85-0, 47849, 4-hydroxybenzoic acid, 99-96-7, 240141, acetylsalicylic acid, 50-78-2, A3160, 2-hydroxybenzoic acid, 69-72-7, ethyl paraben, R472370



G002627

Conditions

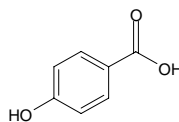
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 70:30, 0.1% TFA in water:0.1% TFA in acetonitrile
flow rate: 1.0 mL/min.
temp.: 30 °C
det.: UV, 220 nm
injection: 10 µL
sample: as indicated in 0.1% TFA in water

Peak IDs

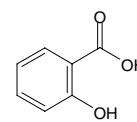
1. 4-hydroxybenzoic acid (50 mg/mL)
2. Acetylsalicylic acid (50 mg/mL)
3. Benzoic acid (50 mg/mL)
4. 2-hydroxybenzoic acid (50 mg/mL)
5. Ethyl paraben (50 mg/mL)

Structures

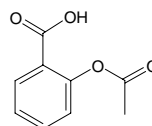
4-hydroxybenzoic acid - G002623



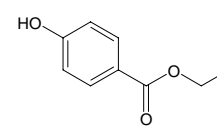
2-hydroxybenzoic acid - G002624



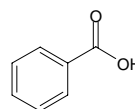
Acetylsalicylic acid - G002625



Ethyl paraben - G002626



Benzoic acid - G002622



Ascentis is a trademark of Sigma-Aldrich Co.