

Application Report 277

Analysis of Analgesics Acetaminophen and Gentisic Acid Using Ascentis™ RP-Amide

This application demonstrates the suitability of Ascentis RP-Amide for the efficient separation of the analgesics acetaminophen and gentisic acid by HPLC.

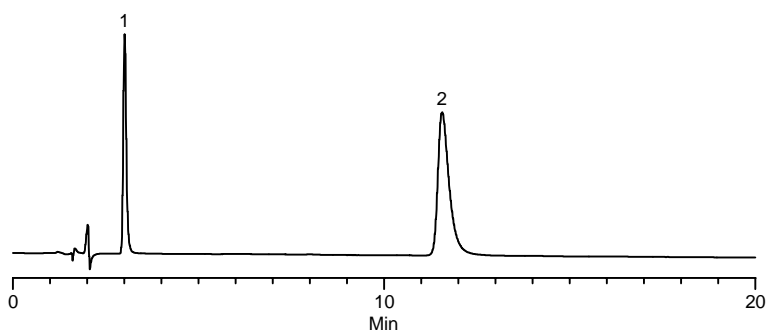
Key Words

Ascentis RP-Amide, 565324-U, acetaminophen, A7085, 103-90-2, gentisic acid, G5254, 490-79-9

Author: Jacinth A.M. McKenzie

Acquisition System: Hitachi LC

Notebook Reference: 1550-07



G003035

Conditions

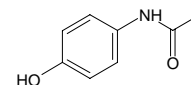
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 80:20, 20 mM phosphoric acid (pH 2.0 unadjusted):acetonitrile
flow rate: 1.0 mL/min.
temp.: 35°C
det.: UV at 220 nm
injection: 10 µL
sample: as indicated (in 20 mM phosphoric acid, pH 2.0)

Peak IDs

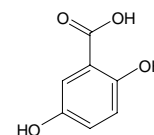
1. Acetaminophen (50 µg/mL)
2. Gentisic acid (50 µg/mL)

Structures

Acetaminophen - G002257



Gentisic acid - G002341



Ascentis is a trademark of Sigma-Aldrich Co.