

Application Report 235

Analysis of Cold Remedy Ingredients Using Ascentis™ C18

This application demonstrates the suitability of the Ascentis C18 for the analysis of cold remedy ingredients. Structures along with the optimized chromatogram obtained on the Ascentis C18 are presented.

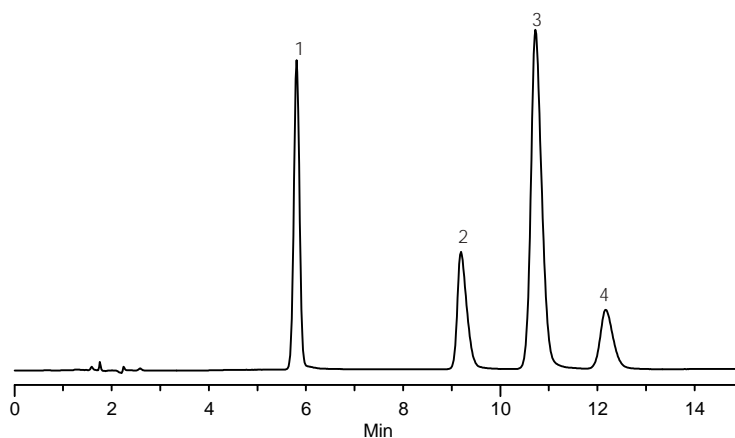
Key Words

acetaminophen, 103-90-2, A7085, pseudoephedrine, 345-78-8, 294616, codeine, 76-57-3, C1653, doxylamine, 562-10-7, D3775, Ascentis C18, 581324-U

Author: Hugh M. Cramer

Acquisition System: Cinnabar
Waters 2690

Notebook Reference: 1551



G002921

Conditions

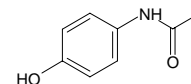
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)
mobile phase: 93:7, 25 mM potassium phosphate (pH 2.5 with phosphoric acid):
acetonitrile
flow rate: 1.0 mL/min.
temp.: 35 °C
det.: UV at 215 nm
injection: 10 µL
sample: as indicated in buffer

Peak IDs

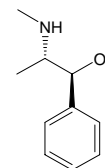
1. Acetaminophen (50 µg/mL)
2. Pseudoephedrine (50 µg/mL)
3. Codeine (50 µg/mL)
4. Doxylamine (50 µg/mL)

Structures

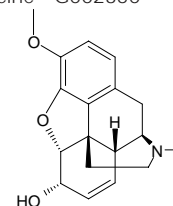
Acetaminophen - G002257



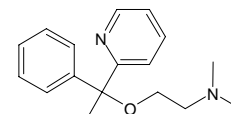
Pseudoephedrine - G002922



Codeine - G002350



Doxylamine - G002923



Ascentis is a trademark of Sigma-Aldrich Co.