

Application Report 236

Analysis of Cold Remedy Ingredients Using Ascentis™ RP-Amide

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

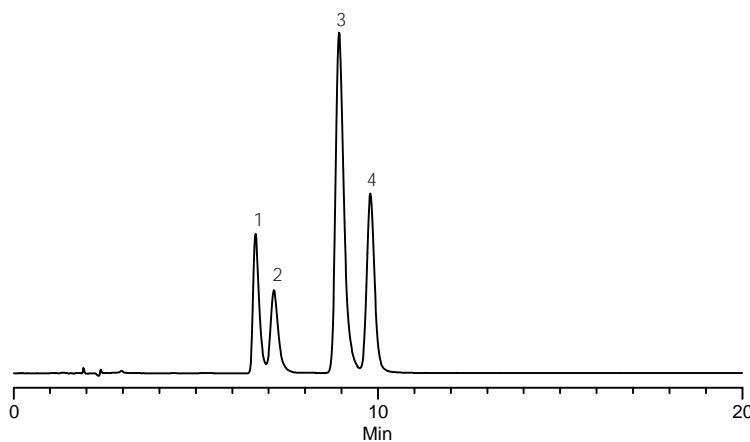
Key Words

acetaminophen, 103-90-2, A7085, pseudoephedrine, 345-78-8, 294616, codeine, 76-57-3, C1653, doxylamine, 562-10-7, D3775, Ascentis RP-Amide, 565324-U

Author: Hugh M. Cramer

Acquisition System: Cinnabar
Waters 2690

Notebook Reference: 1551



G002924

Conditions

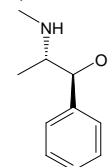
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 95:5, 25 mM potassium phosphate (pH 2.5 with phosphoric acid):
acetonitrile
flow rate: 1.0 mL/min.
temp.: 35 °C
det.: UV at 215 nm
injection: 10 µL
sample: as indicated in buffer

Peak IDs

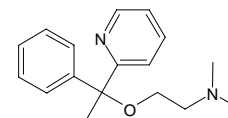
1. Pseudoephedrine (50 µg/mL)
2. Doxylamine (50 µg/mL)
3. Codeine (50 µg/mL)
4. Acetaminophen (50 µg/mL)

Structures

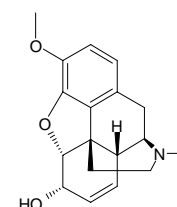
Pseudoephedrine - G002922



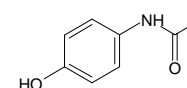
Doxylamine - G002923



Codeine - G002350



Acetaminophen - G002257



Ascentis is a trademark of Sigma-Aldrich Co.