

Application Report 168

Analysis of Anticholinergics Using Ascentis™ RP-Amide

This application demonstrates the suitability of Ascentis RP-Amide for the efficient separation of anticholinergics scopolamine and methyl scopolamine by HPLC.

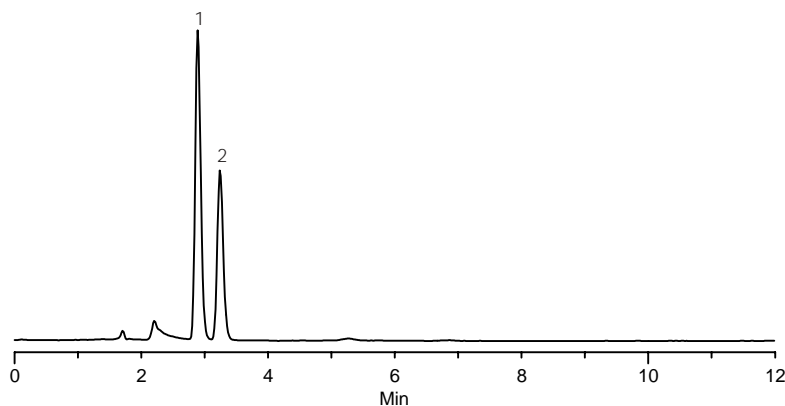
Key Words

Ascentis RP-Amide, methyl scopolamine, methscopolamine, 155-41-9, S8502, scopolamine, 55-16-3, S1013, 565324-U, dramamine

Author: Hugh Cramer

Acquisition System: Waters Alliance
2690 Mt Hoyt

Notebook Reference: 1551



G002663

Conditions

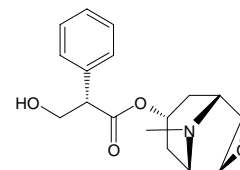
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 85:15, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):acetonitrile
flow rate: 1 mL/min.
temp.: 35 °C
det.: UV, 220 nm
injection: 10 µL
sample: 50 µg/mL each in mobile phase

Peak IDs

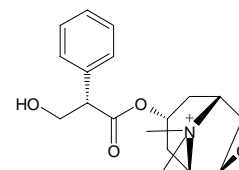
1. Scopolamine
2. Methyl scopolamine

Structures

Scopolamine - G002353



Methyl scopolamine - G002352



Ascentis is a trademark of Sigma-Aldrich Co.