

Application Report 307

Analysis of Analgesics Using Ascentis™ C8

This application demonstrates the suitability of the Ascentis C8 for the analysis of the analgesics buprenorphine and ketorolac. Structures along with the optimized chromatogram obtained on the Ascentis C8 are presented.

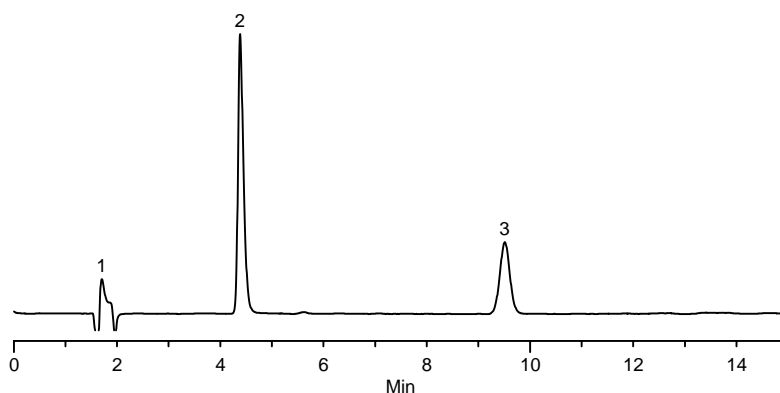
Key Words

buprenorphine, 53152-21-9, B7536, ketorolac, 74103-07-4, K1136, Ascentis C8, 581424-U

Author: Hugh M. Cramer

Acquisition System: Hitachi Mt Fuji

Notebook Reference: 1558



G003180

Conditions

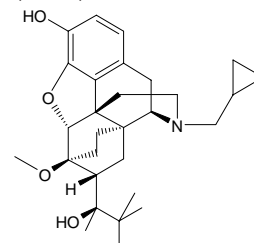
column: Ascentis C8, 15 cm x 4.6 mm I.D., 5 µm particles (581424-U)
mobile phase: 65:35, 10 mM ammonium phosphate monobasic (pH 2.50 with phosphoric acid):acetonitrile
flow rate: 1.0 mL/min.
temp.: 35 °C
det.: UV at 220 nm
injection: 10 µL
sample: as indicated in mobile phase

Peak IDs

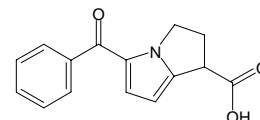
1. Buprenorphine (50 µg/mL)
2. Ketorolac (50 µg/mL)

Structures

Buprenorphine - G002237



Ketorolac - G003179



Ascentis is a trademark of Sigma-Aldrich Co.