

Application Report 238

Analysis of Antiviral Nevirapine Using Ascentis™ RP-Amide

This application demonstrates the suitability of the Ascentis RP-Amide for the analysis of nevirapine. The structure along with the optimized chromatogram obtained on the Ascentis RP-Amide are presented.

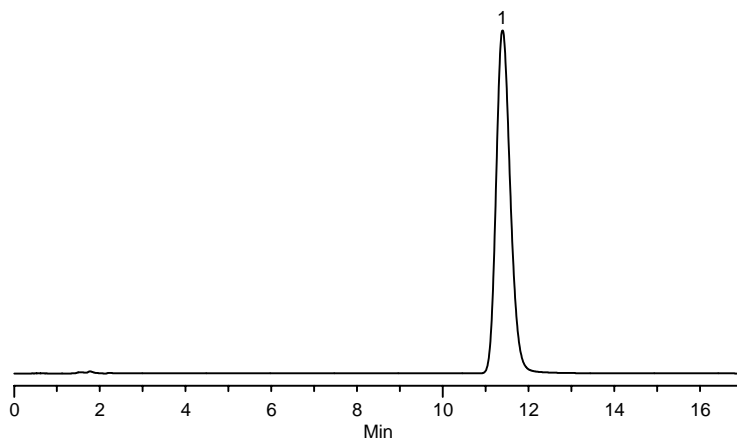
Key Words

Nevirapine, 129618-40-2, Ascentis RP-Amide, 565305-U

Author: Michael D. Buchanan

Acquisition System: "Barkaboom Mountain" (Shimadzu Class-VP)

Notebook Reference: 1562-05



G002928

Conditions

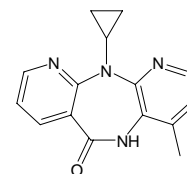
column: Ascentis RP-Amide, 15 cm x 2.1 mm I.D., 5 µm particles (565305-U)
mobile phase: 80:20, 25 mM ammonium phosphate, pH 5.0 with sodium hydroxide:acetonitrile
flow rate: 0.2 mL/min.
temp.: 35 °C
det.: UV at 220 nm
injection: 10 µL
sample: as indicated in mobile phase

Peak IDs

1. Nevirapine (32.5 µg/mL)

Structures

Nevirapine - G002929



Ascentis is a trademark of Sigma-Aldrich Co.