

Application Report 232

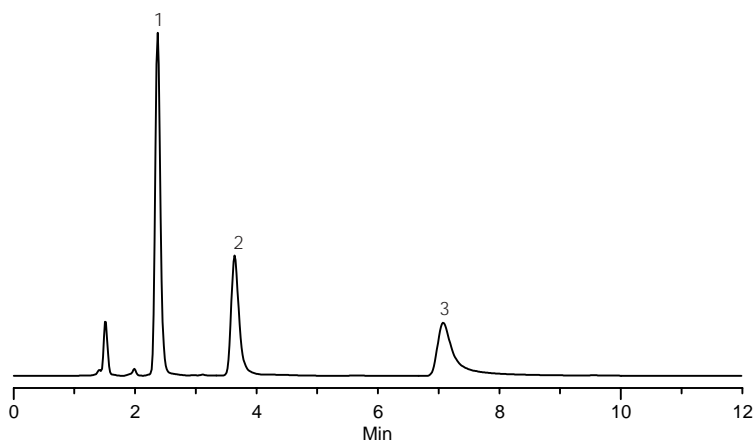
Analysis of the Angiotensins Using Ascentis™ RP-Amide

This application demonstrates the suitability of the Ascentis RP-Amide for the analysis of angiotensins I, II, and III.

Key Words

Angiotensin I, 70937-97-2, A9650, Angiotensin II, 68521-88-0, A9525, Angiotensin III, 100900-06-9, WA13840, Ascentis RP-Amide, 565324-U

Author: Hugh M. Cramer
Acquisition System: Mt Hoyt
Waters 2690
Notebook Reference: 1551



G002914

Conditions

column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 77.5:22.5, 5 mM ammonium phosphate (pH 7.0 with ammonium hydroxide):acetonitrile
flow rate: 1.0 mL/min.
temp.: 35 °C
det.: UV at 215 nm
injection: 10 µL
sample: as indicated in mobile phase

Peak IDs

1. Angiotensin II (100 µg/mL)
2. Angiotensin III (100 µg/mL)
3. Angiotensin I (100 µg/mL)

Structures

Angiotensin II
DRVYIHPF

Angiotensin III
RVYIHPF

Angiotensin I
DRVYIHPFHL

Ascentis is a trademark of Sigma-Aldrich Co.