

Application Report 231

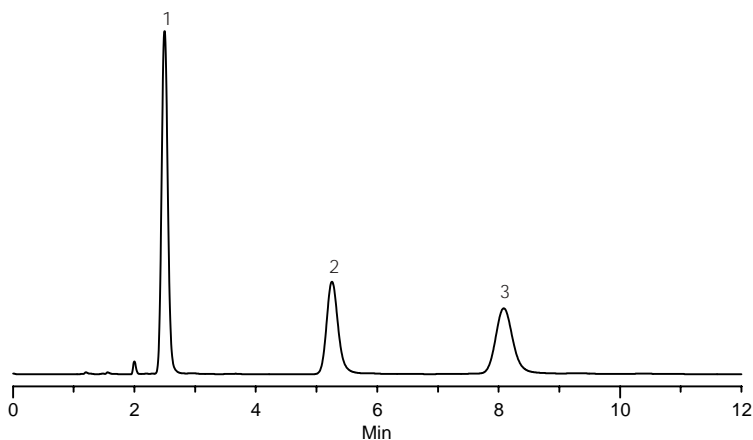
Analysis of the Angiotensins Using Ascentis™ C18

This application demonstrates the suitability of the Ascentis C18 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 C18, 581324-U

Author: Hugh M. Cramer
Acquisition System: Mt McKinley Thermo
Notebook Reference: 1551



G002913

Conditions

column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-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.