

Application Report 193

Analysis of Ascorbic Acid and Dehydroascorbic Acid Using Ascentis™ RP-Amide

This application demonstrates the suitability of Ascentis RP-Amide for the efficient separation of ascorbic acid and dehydroascorbic acid by HPLC.

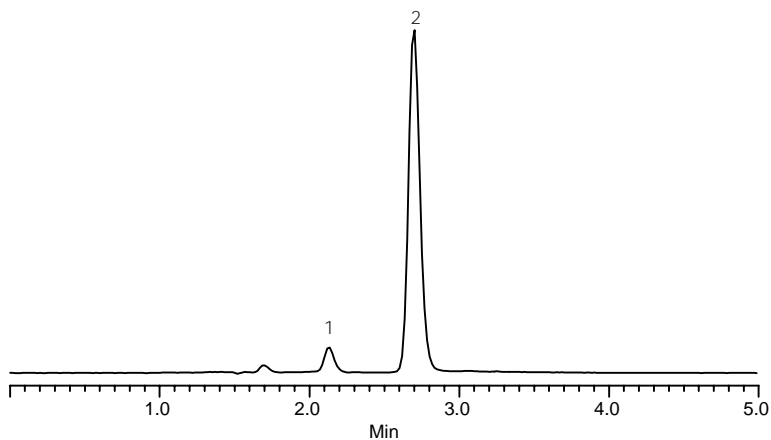
Key Words

Ascentis RP-Amide, 565324-U, ascorbic acid, A4403, 50-81-7, dehydroascorbic acid, 261556, 490-83-5

Author: Hugh Cramer

Acquisition System: Waters Alliance
2690 Cinnabar

Notebook Reference: 1551



G002722

Conditions

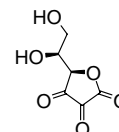
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 25 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid)
flow rate: 1 mL/min.
temp.: 35 °C
det.: UV, 230 nm
injection: 10 µL
sample: as listed each in mobile phase

Peak IDs

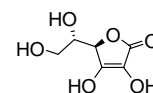
1. Dehydroascorbic acid (100 µg/mL)
2. Ascorbic acid (20 µg/mL)

Structures

Dehydroascorbic acid - G002719



Ascorbic acid - G002720



Ascentis is a trademark of Sigma-Aldrich Co.