

Application Report 234

Analysis of the Antiulceratives Using Ascentis™ RP-Amide

This application demonstrates the suitability of the Ascentis RP-Amide for the analysis of the antiulceratives cimetidine, famotidine, nizatidine, and ranitidine.

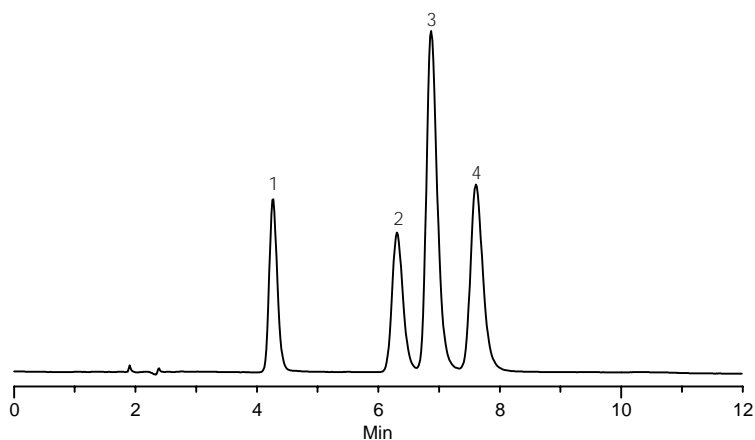
Key Words

cimetidine, 51481-61-9, C4522, famotidine, 76824-35-6, F6889, nizatidine, 76963-41-2, N7035, ranitidine, 66357-35-5, R0633, Ascentis RP-Amide, 565324-U

Author: Hugh M. Cramer

Acquisition System: Cinnabar
Waters 2690

Notebook Reference: 1551



G002920

Conditions

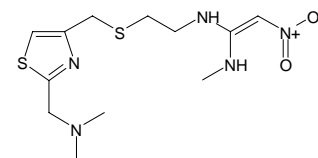
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 95:5, 25 mM potassium phosphate (pH 2.5 with phosphoric acid):acetonitrile
flow rate: 1.0 mL/min.
temp.: 35 °C
det.: UV at 214 nm
injection: 10 µL
sample: as indicated in mobile phase

Peak IDs

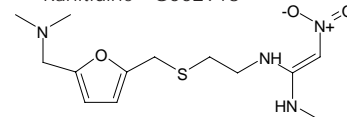
1. Nizatidine (50 µg/mL)
2. Ranitidine (50 µg/mL)
3. Cimetidine (50 µg/mL)
4. Famotidine (50 µg/mL)

Structures

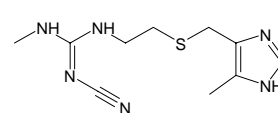
Nizatidine - G002919



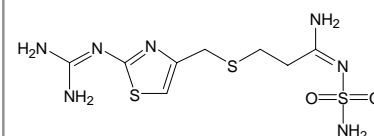
Ranitidine - G002918



Cimetidine - G002917



Famotidine - G002916



Ascentis is a trademark of Sigma-Aldrich Co.