

Application Report 233

Analysis of the Antiulceratives Using Ascentis™ C18

This application demonstrates the suitability of the Ascentis C18 for the analysis of the antiulceratives cimetidine, famotidine, nizatidine, and ranitidine.

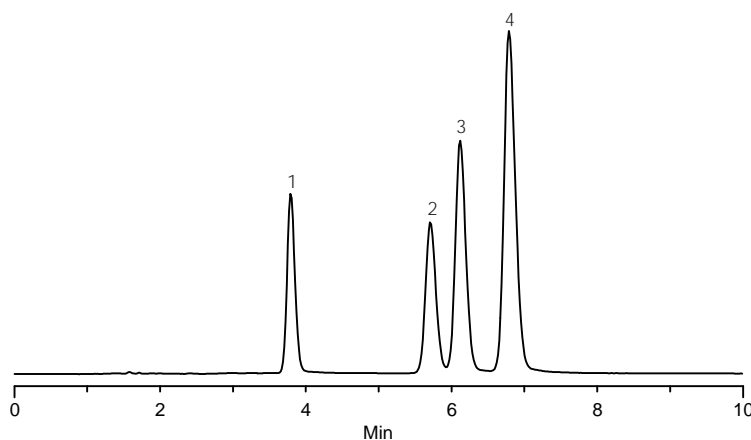
Key Words

cimetidine, 51481-61-9, C4522, famotidine, 76824-35-6, F6889, nizatidine, 76963-41-2, N7035, ranitidine, 66357-35-5, R0633, Ascentis C18, 581324-U

Author: Hugh M. Cramer

Acquisition System: Cinnabar
Waters 2690

Notebook Reference: 1551



G002915

Conditions

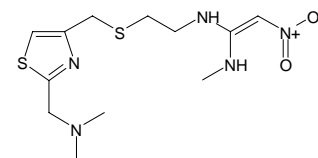
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)
mobile phase: 87.5:12.5, 25 mM potassium phosphate (pH 2.5 with phosphoric acid):acetonitrile
flow rate: 1.0 mL/min.
temp.: 35 °C
det.: UV at 214 nm
injection: 10 µL
sample: as indicated in mobile phase

Peak IDs

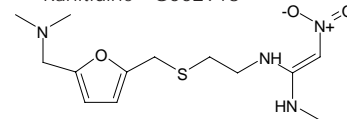
1. Nizatidine (50 µg/mL)
2. Ranitidine (50 µg/mL)
3. Famotidine (50 µg/mL)
4. Cimetidine (50 µg/mL)

Structures

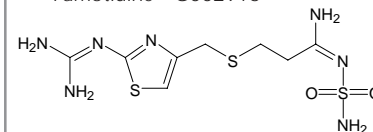
Nizatidine - G002919



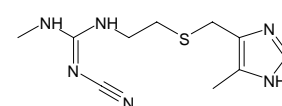
Ranitidine - G002918



Famotidine - G002916



Cimetidine - G002917



Ascentis is a trademark of Sigma-Aldrich Co.