

Application Report 224

Analysis of the Herbicide Triazine Using Ascentis™ RP-Amide

This application demonstrates the suitability of the Ascentis RP-Amide for the analysis of the herbicides simazine, cyanazine, and atrazine.

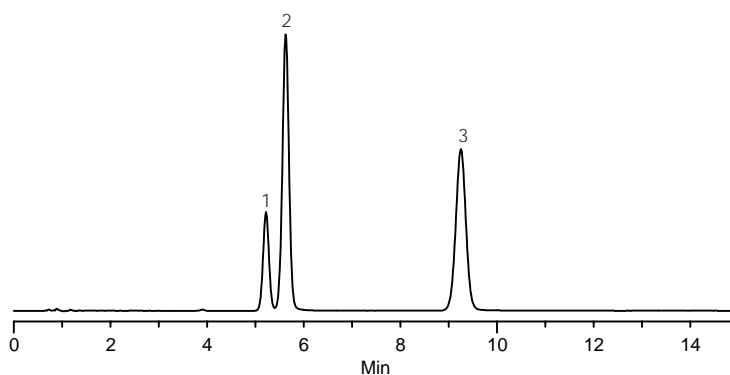
Key Words

simazine, 122-34-9, cyanazine, 21725-46-2, 48592, atrazine, 1912-24-9, 45330, Ascentis RP-Amide, 565324-U

Author: Hugh M. Cramer

Acquisition System: : Mt Hoyt
Waters 2690

Notebook Reference: 1551



G002893

Conditions

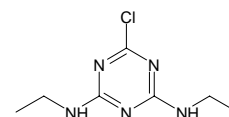
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 70:30, water:acetonitrile
flow rate: 2.0 mL/min.
temp.: 35 °C
det.: UV at 214 nm
injection: 10 µL
sample: as indicated in mobile phase

Peak IDs

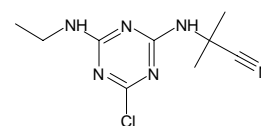
1. Simazine (50 µg/mL)
2. Cyanazine (50 µg/mL)
3. Atrazine (50 µg/mL)

Structures

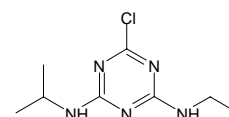
Simazine - G001005



Cyanazine - G002891



Atrazine - G002867



Ascentis is a trademark of Sigma-Aldrich Co.