

# Application Report 223

## Analysis of the Herbicide Triazines Using Ascentis™ C18

This application demonstrates the suitability of the Ascentis C18 for the analysis of the herbicide triazines simazine, cyanazine, and atrazine.

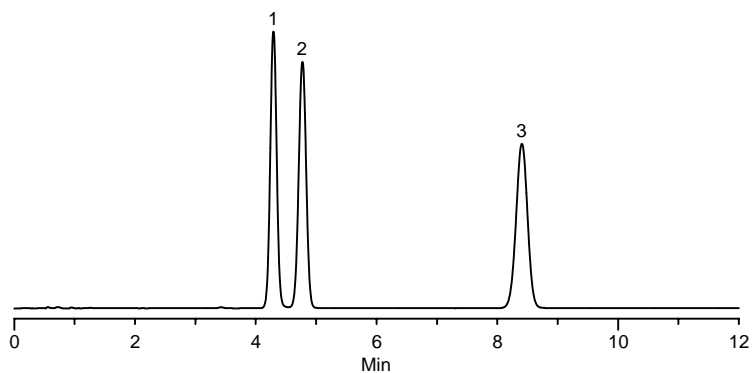
### Key Words

simazine, 122-34-9, cyanazine, 21725-46-2, 48592, atrazine, 1912-24-9, 45330, Ascentis C18, 581324-U

Author: Hugh M. Cramer

Acquisition System: Mt McKinley  
Thermo

Notebook Reference: 1551



G002892

### Conditions

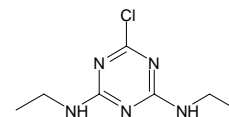
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)  
mobile phase: 70:30, water:acetonitrile  
flow rate: 2.0 mL/min.  
temp.: 35 °C  
det.: UV at 214 nm  
injection: 10 µL  
sample: as indicated in mobile phase

### Peak IDs

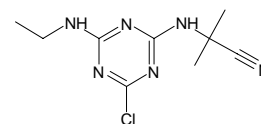
1. Simazine (50 µg/mL)
2. Cyanazine (50 µg/mL)
3. Atrazine (50 µg/mL)

### Structures

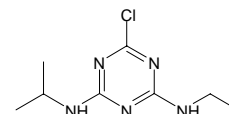
Simazine - G001005



Cyanazine - G002891



Atrazine - G002867



Ascentis is a trademark of Sigma-Aldrich Co.