

# Application Report 215

## Analysis of the Pesticides Aldicarb, Carbofuran, and Atrazine Using Ascentis™ RP-Amide

This application demonstrates the suitability of the Ascentis RP-Amide for the analysis of the pesticides aldicarb, carbofuran, and atrazine.

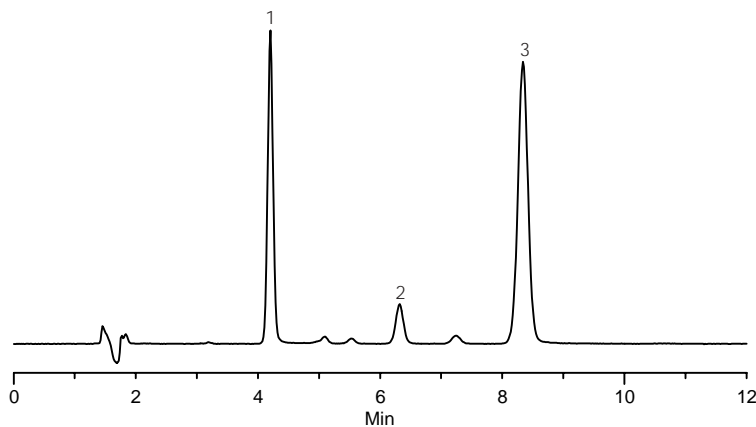
### Key Words

aldicarb, 116-06-3, 33386, carbofuran, 1563-66-2, 45370, atrazine, 1912-24-9, 45330, Ascentis RP-Amide, 565324-U

Author: Hugh M. Cramer

Acquisition System: Cinnabar  
Waters 2690

Notebook Reference: 1551



G002869

### Conditions

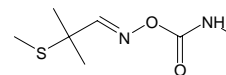
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)  
mobile phase: 60:40, water:acetonitrile  
flow rate: 1.0 mL/min.  
temp.: 35 °C  
det.: UV at 254 nm  
injection: 10 µL  
sample: as indicated in mobile phase

### Peak IDs

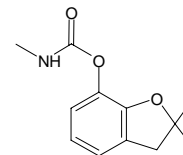
1. Aldicarb (50 µg/mL)
2. Carbofuran (50 µg/mL)
3. Atrazine (50 µg/mL)

### Structures

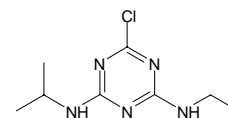
Aldicarb - G002865



Carbofuran - G002866



Atrazine - G002867



Ascentis is a trademark of Sigma-Aldrich Co.