

# Application Report 208

## Analysis of the Pesticides Carbaryl, Diuron, and Propachlor Using Ascentis™ C18

This application demonstrates the suitability of the Ascentis C18 for the analysis of the pesticides carbaryl, diuron, and propachlor.

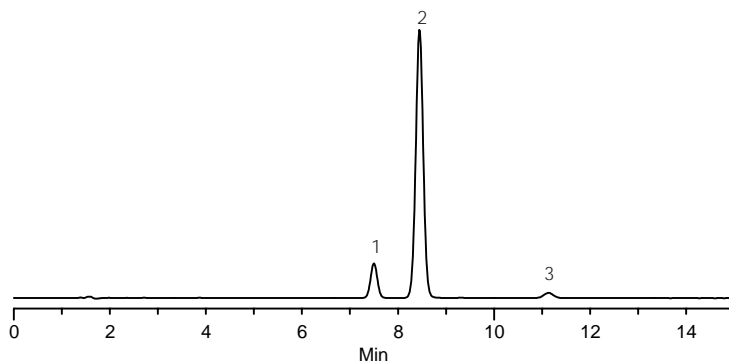
### Key Words

carbaryl, 63-25-2, 45367, diuron, 330-54-1, 45463, propachlor, 1918-16-7, 442765, Ascentis C18, 581324-U

Author: Hugh M. Cramer

Acquisition System: Cinnabar  
Waters 2690

Notebook Reference: 1551



G002854

### Conditions

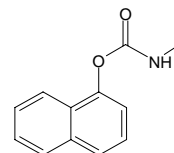
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)  
mobile phase: 60:40, water:acetonitrile  
flow rate: 1.0 mL/min.  
temp.: 35 °C  
det.: UV at 254 nm  
injection: 10 µL  
sample: as indicated in mobile phase

### Peak IDs

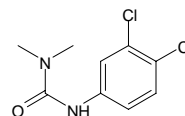
1. Carbaryl (50 µg/mL)
2. Diuron (50 µg/mL)
3. Propachlor (50 µg/mL)

### Structures

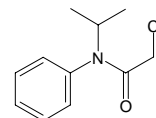
Carbaryl - G002851



Diuron - G002852



Propachlor - G002853



Ascentis is a trademark of Sigma-Aldrich Co.