

Application Report 214

Analysis of the Pesticides Aldicarb, Carbofuran, and Atrazine Using Ascentis™ C18

This application demonstrates the suitability of the Ascentis C18 for the analysis of the pesticides aldicarb, carbofuran, and atrazine.

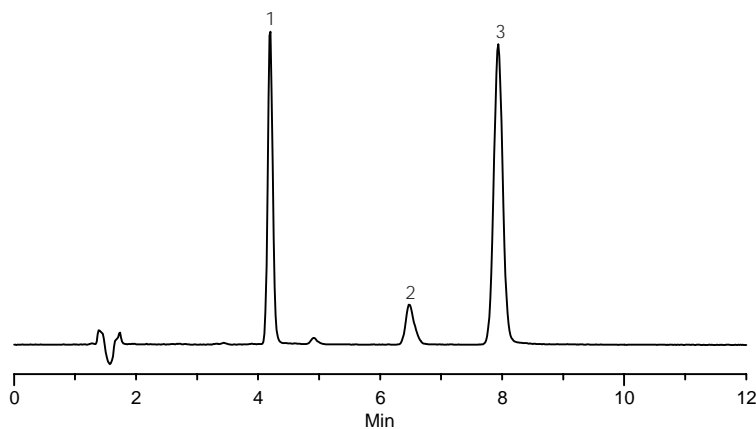
Key Words

aldicarb, 116-06-3, 33386, carbofuran, 1563-66-2, 45370, atrazine, 1912-24-9, 45330, Ascentis C18, 581324-U

Author: Hugh M. Cramer

Acquisition System: Cinnabar
Waters 2690

Notebook Reference: 1551



G002868

Conditions

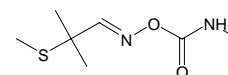
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)
mobile phase: 60:40, water:acetonitrile
flow rate: 1.0 mL/min.
temp.: 35 °C
det.: UV at 254 nm
injection: 10 µL
sample: as indicated in mobile phase

Peak IDs

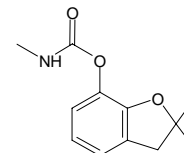
1. Aldicarb (50 µg/mL)
2. Carbofuran (50 µg/mL)
3. Atrazine (50 µg/mL)

Structures

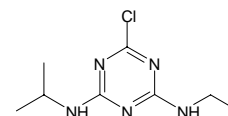
Aldicarb - G002865



Carbofuran - G002866



Atrazine - G002867



Ascentis is a trademark of Sigma-Aldrich Co.