

# Application Report 320

## Analysis of US EPA Method 632 Carbamate and Urea Pesticides Using Ascentis™ C8

US EPA Method 632 specifies a procedure for the determination of carbamate and urea pesticides in municipal and industrial wastewaters. The suitability of the Ascentis C8 for the separation of the nineteen analytes specified in the method is demonstrated. The structures along with the optimized chromatogram obtained on the Ascentis C8 are presented.

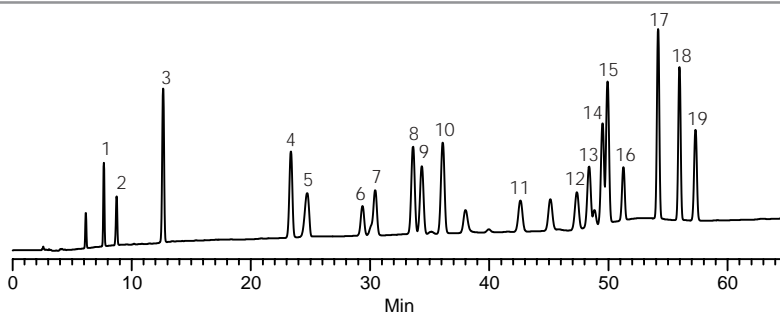
### Key Words

aminocarb, 45322, 2032-59-9, barban, 33390, 101-27-9, carbaryl, 45367, 63-25-2, carbofuran, 45370, 1563-66-2, chlorpropham, 45393, 101-21-3, diuron, 45463, 330-54-1, fenuron, 45494, 101-42-8, fluomethuron, 45502, 2164-17-2, linuron, 36141, 330-55-2, methiocarb, 36152, 2032-65-7, methomyl, 36159, 16752-77-5, mexacarbate, 315-18-4, monuron, 36174, 150-68-5, neburon, 36176, 555-37-3, oxamyl, 36184, 23135-22-0, propham, 45641, 122-42-9, propoxur, 45644, 114-26-1, siduron, 34373, 1982-49-6, swep, 45666, 1918-18-9, carbamate and urea pesticide, US EPA Method 632, Ascentis C8, 581425-U

Author: Michael D. Buchanan

Acquisition System: "Pikes Peak"  
(Thermo Instrument)

Notebook Reference: 1562-72



G003239

### Conditions

column: Ascentis C8, 25 cm x 4.6 mm I.D., 5 µm particles (581425-U)  
mobile phase A: DI water  
mobile phase B: acetonitrile  
flow rate: 1.0 mL/min.  
temp: 30 °C  
det: UV at 210 nm  
injection: 10 µL  
sample: as indicated in 16% acetonitrile in DI water  
gradient:

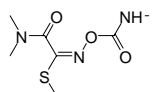
Min	%A	%B
0	85	15
4	75	25
44	55	45
50	40	60
65	40	60

### Peak IDs

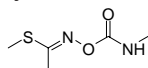
1. Oxamyl (10 µg/mL)
2. Methomyl (10 µg/mL)
3. Fenuron (10 µg/mL)
4. Monuron (10 µg/mL)
5. Aminocarb (10 µg/mL)
6. Propoxur (10 µg/mL)
7. Carbofuran (10 µg/mL)
8. Fluomethuron (10 µg/mL)
9. Carbaryl (10 µg/mL)
10. Diuron (10 µg/mL)
11. Propham (10 µg/mL)
12. Siduron (10 µg/mL)
13. Methiocarb (10 µg/mL)
14. Linuron (10 µg/mL)
15. Swep (10 µg/mL)
16. Mexacarbate (10 µg/mL)
17. Chlorpropham (10 µg/mL)
18. Barban (10 µg/mL)
19. Neburon (10 µg/mL)

### Structures

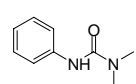
Oxamyl - G003189



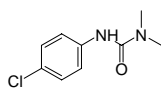
Methomyl - G003186



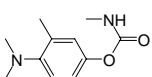
Fenuron - G002858



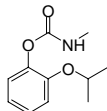
Monuron - G002857



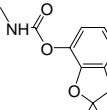
Aminocarb - G003181



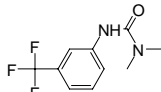
Propoxur - G003191



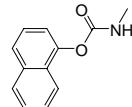
Carbofuran - G002866



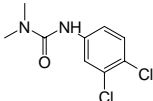
Fluomethuron - G003184



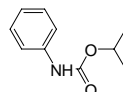
Carbaryl - G002851



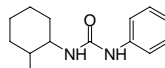
Diuron - G002852



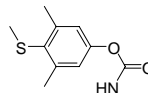
Propham - G003190



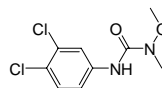
Siduron - G003192



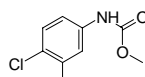
Methiocarb - G003185



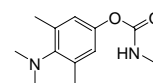
Linuron - G002856



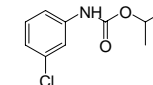
Swep - G003193



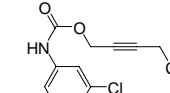
Mexacarbate - G003187



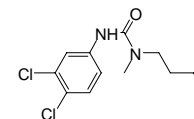
Chlorpropham - G003183



Barban - G003182



Neburon - G003188



Ascentis is a trademark of Sigma-Aldrich Co.