

# Application Report 286

## Analysis of US EPA Method 632.1 Pesticides Using Ascentis™ C8

US EPA Method 632.1 specifies a procedure for the determination of pesticides in municipal and industrial wastewaters. The suitability of the Ascentis C8 for the separation of the three analytes specified in the method is demonstrated. The structures along with the optimized chromatogram obtained on the Ascentis C8 are presented.

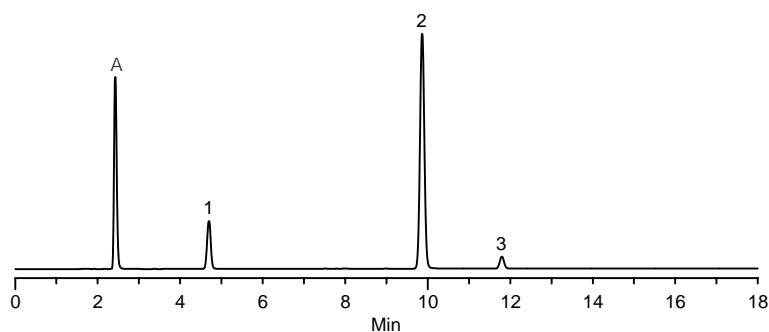
### Key Words

napropamide, 36175, 15299-99-7, propanil, 45639, 709-98-8, vacor, 53558-25-1, pesticide, US EPA Method 632.1, Ascentis C8, 581424-U

Author: Michael D. Buchanan

Acquisition System: "Pikes Peak"  
(Thermo Instrument)

Notebook Reference: 1562-69



G003121

### Conditions

column: Ascentis C8, 15 cm x 4.6 mm I.D., 5 µm particles (581424-U)  
mobile phase A: DI water  
mobile phase B: acetonitrile  
flow rate: 1.0 mL/min.  
temp.: 35 °C  
det.: UV at 254 nm  
injection: 100 µL  
sample: as indicated in 2.2% 90:10 acetonitrile:acetone in DI water

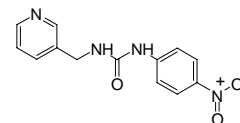
gradient:	Min	%A	%B
	0	65	35
	13	25	75
	18	25	75

### Peak IDs

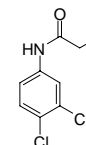
- A. Acetone
1. Vacor (10 µg/mL)
2. Propanil (10 µg/mL)
3. Napropamide (10 µg/mL)

### Structures

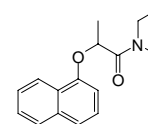
Vacor - G003119



Propanil - G003118



Napropamide - G003117



Ascentis is a trademark of Sigma-Aldrich Co.