

# Application Report 115

## Analysis of Tranquilizer Fentanyl Citrate Using Ascentis™ C18

This application demonstrates the suitability of Ascentis C18 for the efficient elution of the tranquilizer fentanyl by HPLC.

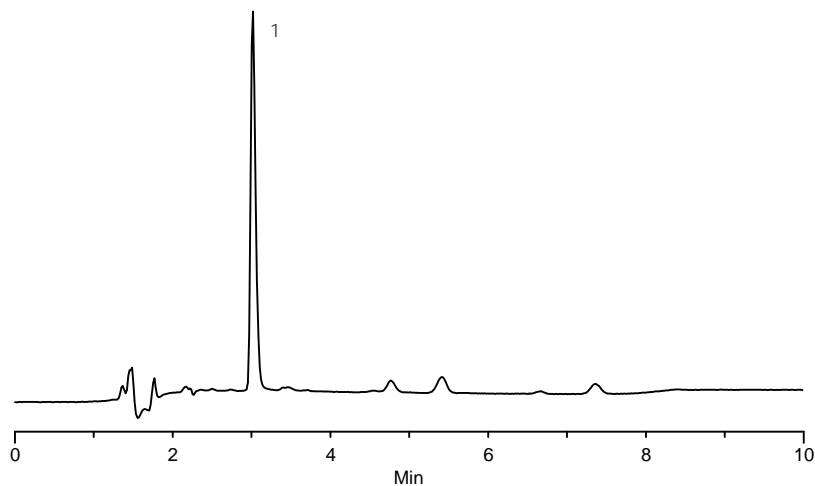
### Key Words

Ascentis C18, fentanyl, 990-73-8, F5886, 581324-U, tranquilizer

Author: Hugh Cramer

Acquisition System: Waters Alliance  
2690 ID 9371

Notebook Reference: 1495-09



G002435

### Conditions

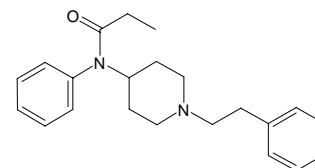
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)  
mobile phase: 67:33, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):acetonitrile  
flow rate: 1 mL/min.  
temp.: 35 °C  
det.: UV at 215 nm  
injection: 10 µL  
sample: 20 µg/mL in mobile phase

### Peak IDs

1. Fentanyl

### Structures

Fentanyl - G002436



Ascentis is a trademark of Sigma-Aldrich Co.