

# Application Report 97

## Analysis of Antiinflammatory Drugs Using Ascentis™ C18

This application demonstrates the suitability of Ascentis C18 for the efficient separation of indomethacin and piroxicam by HPLC.

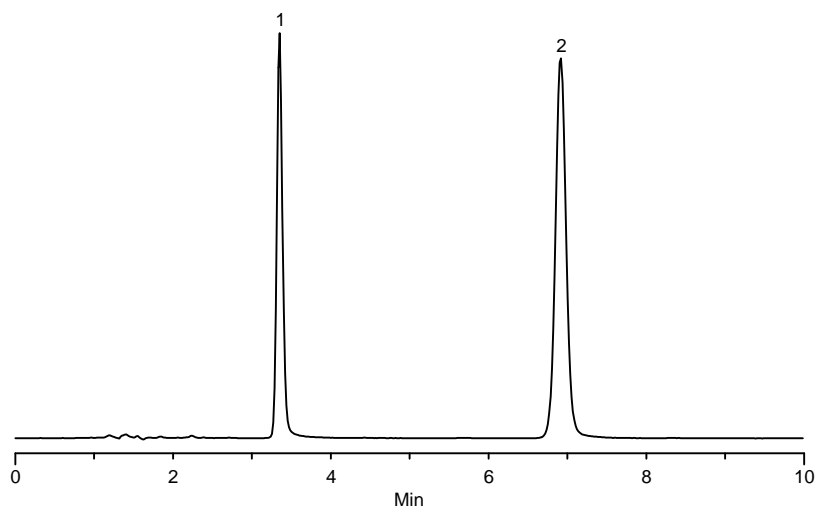
### Key Words

Ascentis C18, indomethacin, 53-86-1, I7378, piroxicam, 36322-90-4, P5654, 581324-U

Author: Hugh Cramer

Acquisition System: Waters Alliance  
2690 ID 9371

Notebook Reference: 1486-69



G002391

### Conditions

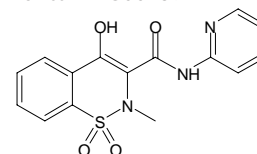
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)  
mobile phase: 45:55, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):acetonitrile  
flow rate: 1 mL/min.  
temp.: 35 °C  
det.: UV at 220 nm  
injection: 10 µL  
sample: 25 µg/mL each in mobile phase

### Peak IDs

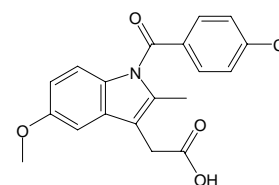
1. Piroxicam
2. Indomethacin

### Structures

Piroxicam - G002392



Indomethacin - G002393



Ascentis is a trademark of Sigma-Aldrich Co.