

# Application Report 145

## Analysis of Non-steroidal Anti-inflammatory Drugs (NSAIDs) Using Ascentis™ RP-Amide

This application demonstrates the suitability of Ascentis RP-Amide for the efficient separation of NSAIDs ketoprofen, naproxen and ibuprofen using Ascentis RP-Amide.

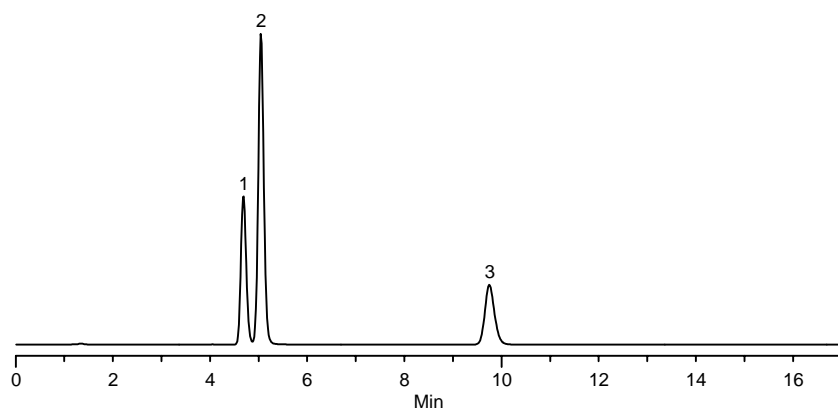
### Key Words

Ascentis RP-Amide, 565324-U, anti-inflammatory, NSAIDs, ketoprofen, 2207-15-4, K1751, naproxen, 26159-34-2, N5160, ibuprofen, 15687-27-1, I4883

Author: Jacinth A.M. McKenzie

Acquisition System: Waters

Notebook Reference: 1550-41



G002621

### Conditions

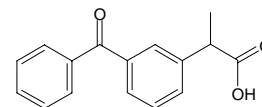
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)  
mobile phase: 45:55, 10 mM dibasic potassium phosphate (pH 3.0):acetonitrile  
flow rate: 1.0 mL/min.  
temp.: 35 °C  
det.: UV, 230 nm  
injection: 5 µL  
sample: as indicated in 10 mM dibasic potassium phosphate (pH 3.0)

### Peak IDs

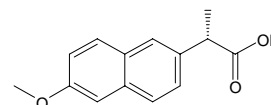
1. Ketoprofen (250 µg/mL)
2. Naproxen (50 µg/mL)
3. Ibuprofen (250 µg/mL)

### Structures

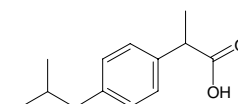
Ketoprofen - G002618



Naproxen - G002619



Ibuprofen - G002620



Ascentis is a trademark of Sigma-Aldrich Co.