

# Application Report 265

## Separation of Analgesics Using Ascentis™ RP-Amide

This application demonstrates the separation of analgesics, oxyphenbutazone, ketoprofen, naproxen, phenylbutazone, flunixin and meclofenamic acid using an Ascentis RP-Amide column.

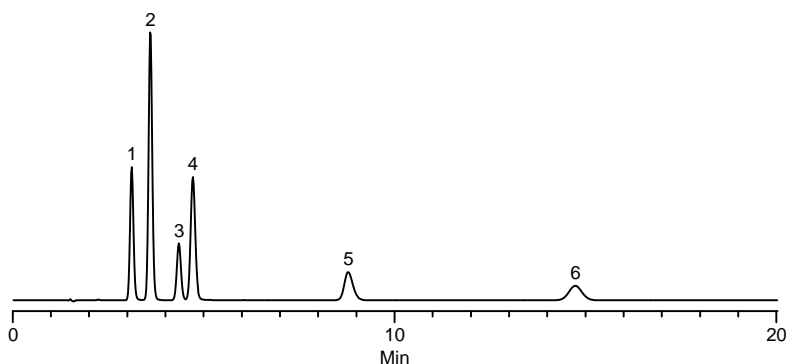
### Key Words

Ascentis RP-Amide, 565324-U, analgesics, oxyphenbutazone hydrate 156025 7081-38-1, ketoprofen 47,190-9 22161-81-5, naproxen N5160 26159-34-2, phenylbutazone P8386 50-33-9, flunixin meglumine F0429 38677-85-9, meclofenamic acid M4531 6385-02-0.

Author: Jacinth A.M. McKenzie

Acquisition System: Mt Hoyt  
Jay\_Truesdail, export id # 1188

Notebook Reference: 1550



G002997

### Conditions

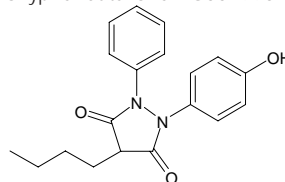
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)  
mobile phase: 30:70, 50 mM potassium phosphate (pH 3.0):methanol  
flow rate: 1.0 mL/min.  
temp: 35 °C  
det: UV at 254 nm  
injection: 10 µL  
sample: as indicated in 50 mM potassium phosphate (pH 3.0)

### Peak IDs

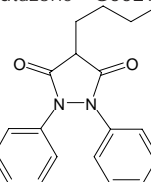
1. Oxyphenbutazone (50 µg/mL)
2. Ketoprofen (50 µg/mL)
3. Naproxen (50 µg/mL)
4. Phenylbutazone (50 µg/mL)
5. Flunixin (50 µg/mL)
6. Meclofenamic acid (50 µg/mL)

### Structures

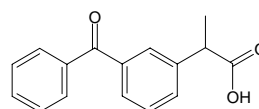
Oxyphenbutazone - G002995



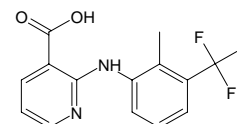
Phenylbutazone - G002994



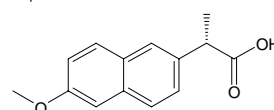
Ketoprofen - G002619



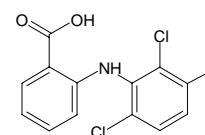
Flunixin - G002993



Naproxen - G002620



Meclofenamic acid - G002996



Ascentis is a trademark of Sigma-Aldrich Co.