# Analysis of Ibuprofen and Valerophenone Using a Thermo Scientific Accucore XL C18 4 µm HPLC Column

Eilidh MacRitchie, Thermo Fisher Scientific, Runcorn, Cheshire, UK

# **Key Words**

Accucore XL C18, solid core, core enhanced technology, ibuprofen, valerophenone

#### **Abstract**

This application note compares the performance of the solid core Thermo Scientific Accucore XL C18 4  $\mu$ m HPLC column with that of a fully porous 5  $\mu$ m traditional HPLC column for the analysis of ibuprofen and valerophenone using a isocratic method based on the USP¹.

#### Introduction

Based on Core Enhanced Technology<sup>TM</sup> using 4 μm solid core particles, Accucore<sup>TM</sup> XL HPLC columns allow users of conventional HPLC methods to enjoy performance far beyond that of columns packed with 5 μm or even 3 μm fully porous particles. Very high separation efficiencies using standard HPLC instruments and conditions provide increased peak resolution and lower limits of detection. An ultra-stable packed bed results in exceptionally robust columns that demonstrate excellent retention and response reproducibility.

Ibuprofen is a non-steroidal anti-inflammatory drug (NSAID) and is a commonly used painkiller. It is particularly useful for relieving the symptoms of inflammation. This application note demonstrates improved chromatographic performance in the analysis of ibuprofen and valerophenone on an Accucore XL HPLC column in relation to a fully porous traditional HPLC column under the same experimental conditions.



## **Experimental Details**

| Consumables   | Part Number |
|---|-------------|
| Fisher Scientific HPLC grade water  | W/0106/17   |
| Fisher Scientific HPLC grade acetonitrile   | A/0626/17   |
| Fisher Scientific HPLC grade orthophosphoric acid   | 0/0515/PB08 |
| Thermo Scientific Borosilicate glass vials (2 mL, 12 mm x 32 mm) with 8 mm black screw cap fitted with a silicone/PTFE seal | 60180-600   |



#### **Sample Preparation**

A mixed working standard containing 50 µg/mL each of ibuprofen and valerophenone was prepared in water.

| Separation Conditions | Part Num  | nber |
|-----------------------|---|------|
| Instrumentation:      | Thermo Scientific Dionex UltiMate 3000 RSLC system                                  |      |
| Columns:              | Accucore XL C18 4 μm, 150 x 4.6 mm 74104-154<br>Fully porous C18 5 μm, 150 x 4.6 mm | 1630 |
| Mobile phase:         | water with phosphoric acid, pH 2.5/ acetonitrile (66.3:33.7 v/v)                    |      |
| Flow rate:            | 2 mL/min  |      |
| Column temperature:   | 30 °C   |      |
| UV detection:         | 214 nm  |      |
| Injection volume:     | 5 μL  |      |

| Data Processing |   |
|-----------------|---|
| Software:       | Thermo Scientific Dionex Chromeleon 7.0<br>Chromatography Data System |

## **Results**

Analysis of ibuprofen and valerophenone was performed on a Accucore XL C18 4  $\mu$ m HPLC column and a 5  $\mu$ m fully porous C18 column using the same experimental conditions. As shown in Figure 1, ibuprofen and valerophenone were well resolved with the USP criteria of resolution of not less than 2 being met on both columns.

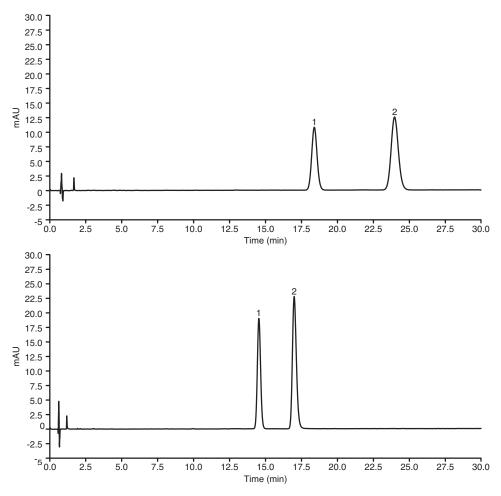


Figure 1: Chromatogram of valerophenone (1) and ibuprofen (2) analyzed using an Accucore XL C18 4  $\mu$ m, 150 x 4.6 mm column (bottom trace) compared to a fully porous 5  $\mu$ m C18, 150 x 4.6 mm column (top trace)

Efficiency for both compounds improve by greater than 70% when using the Accucore XL HPLC column compared to the fully porous column (Table 1). Signal to noise ratio also increased by 112% on average, improving sensitivity.

| Compound      | Plates (USP) |              | Signal to Noise Ratio |              |
|---------------|--------------|--------------|-----------------------|--------------|
|               | Accucore XL  | Fully Porous | Accucore XL           | Fully Porous |
| Valerophenone | 19532        | 11218        | 908                   | 462          |
| Ibuprofen     | 18274        | 10538        | 1202                  | 534          |

Table 1: Efficiency and signal to noise ratio data for valerophenone and ibuprofen

The backpressure for the Accucore XL C18 4  $\mu m$  HPLC column was measured at 312 bar and the 5  $\mu m$  fully porous column backpressure was measured at 239 bar. The increase in performance is gained with a small increase in backpressure and is still within the operating limits of a conventional HPLC system. In addition six replicate injections illustrated excellent reproducibility for both ibuprofen and valerophenone (Table 2).

| Compound      | Accucore XL         |          | Fully Porous        |          |
|---------------|---------------------|----------|---------------------|----------|
|               | t <sub>R</sub> /min | %RSD n=6 | t <sub>R</sub> /min | %RSD n=6 |
| Valerophenone | 14.55               | 0.14     | 18.40               | 0.08     |
| Ibuprofen     | 17.01               | 0.09     | 23.97               | 0.09     |

Table 2: Retention time and precision data calculated from six replicate injections

#### Conclusion

The use of an Accucore XL C18 4 µm HPLC column gave significant performance improvement over a conventional 5 µm fully porous column under the same chromatographic conditions with no changes in system configuration.

Efficiency improved by greater than 73% and signal to noise ratio by 112% on average. Excellent reproducibility was also demonstrated, illustrating that the Accucore XL C18 4  $\mu$ m HPLC column is an ideal choice for the analysis of ibuprofen and valerophenone.

# Reference

1. USP-32, Ibuprofen, Chromatographic purity

#### thermoscientific.com/accucore

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

USA and Canada +1 800 332 3331 France +33 (0)1 60 92 48 34 Germany +49 (0) 2423 9431 20 or 21 United Kingdom +44 (0)1928 534110 Japan +81 3 5826 1615 China +86 21 68654588 +86 10 84193588 +86 20 83145199 800 810 5118 India +91 22 6742 9494 +91 27 1766 2352 Australia 1 300 735 292 (free call domestic) New Zealand 0800 933 966 (free call domestic) All Other Enquiries +44 (0) 1928 534 050 Technical Support North America +1 800 332 3331

**Outside North America** +44 (0) 1928 534 440

