

# Analysis of Non-Steroidal Anti-Inflammatory Drugs Using a Highly Pure, High Surface Area C18 HPLC Column

Jamil Ali, Thermo Fisher Scientific, Runcorn, Cheshire, UK

## Key Words

Synchronis C18, diclofenac, naproxen, aspirin, ibuprofen, sulindac, piroxicam, fenoprofen, NSAIDs, non-steroidal anti-inflammatory drugs

## Abstract

This application note demonstrates the use of the Thermo Scientific™ Synchronis™ C18 HPLC column for the analysis of non-steroidal anti-inflammatory drugs. Synchronis C18 columns provide a fast simple method with good hydrophobic retention, excellent peak shape and high resolution.

## Introduction

One of the key goals for the chromatographer is to achieve a consistent, reproducible separation. The selection of a highly reproducible HPLC column is essential if this goal is to be attained. The Synchronis column range has been engineered to provide exceptional reproducibility due to its highly pure, high surface area silica, dense bonding and double endcapping, all controlled and characterized through the use of rigorous testing.

Non-steroidal Anti-Inflammatory Drugs (NSAIDs) are medications used in reducing inflammation, relieve pain (analgesic) and to lower temperature (fever). NSAIDs are the most prescribed medications for treating conditions such as pain, arthritis, fever and migraine.<sup>1,2</sup> Separating non-steroidal anti-inflammatory drugs with good resolution can be problematic in liquid chromatography. In this application the Synchronis C18 phase was employed to achieve the separation of seven most commonly used NSAIDs.



## Experimental Details

### Consumables

Fisher Scientific™ HPLC grade water	W/0106/17
Fisher Scientific HPLC grade acetonitrile	A/0626/17
Fisher Scientific HPLC grade methanol	M/4056/17
Thermo Scientific Autosampler vial kit	A4954-010

Separation Conditions	Part Number
Instrumentation:	Synchronis C18, 3 $\mu$ m, 100 x 3.0 mm Thermo Scientific HPLC system
Column Temperature:	30 °C
Injection volume:	10 $\mu$ L
Flow rate:	0.60 mL/min
UV detection:	260 nm
Mobile phase A:	0.1% formic acid in water
Mobile phase B:	0.1% formic acid in acetonitrile
Gradient:	50-75% B in 5 minutes
Run time:	7 minutes
Pressure:	207 bar
Wash solvent:	90:10 (v/v) water: acetonitrile

### Sample Preparation

Stock solutions of diclofenac, naproxen, aspirin, ibuprofen, sulindac and fenoprofen were prepared in methanol at a concentration of 1.0mg/mL. Piroxicam was prepared in a methanol:water (1:1 v/v) solution at 1.0mg/mL. The stock solutions were mixed and diluted with water to yield a final working solution containing each compound at the following concentration:

Compound	Concentration ( $\mu$ g/mL)
<b>Naproxen</b>	15
<b>Ibuprofen</b>	100
<b>Diclofenac</b>	30
<b>Aspirin</b>	30
<b>Sulindac</b>	30
<b>Piroxicam</b>	30
<b>Fenoprofen</b>	30

### Results

Analysis was carried out on a 3  $\mu$ m Synchronis C18 column as illustrated in Figure 1. The Synchronis column showed excellent resolving power and provided a very simple method for the separation of all seven compounds. The high surface area silica, dense bonding and double endcapping assist in minimizing secondary silanol interactions and as a result Synchronis columns provide good peak shapes and high resolving power. The reproducibility of the method was assessed by doing six consecutive runs. The method was shown to be very robust with consistent retention times. The chromatograms are shown overlaid in Figure 2. The relative standard deviation of the retention from the six consecutive runs was calculated to be less than 0.12% for all seven compounds. (Table 1)

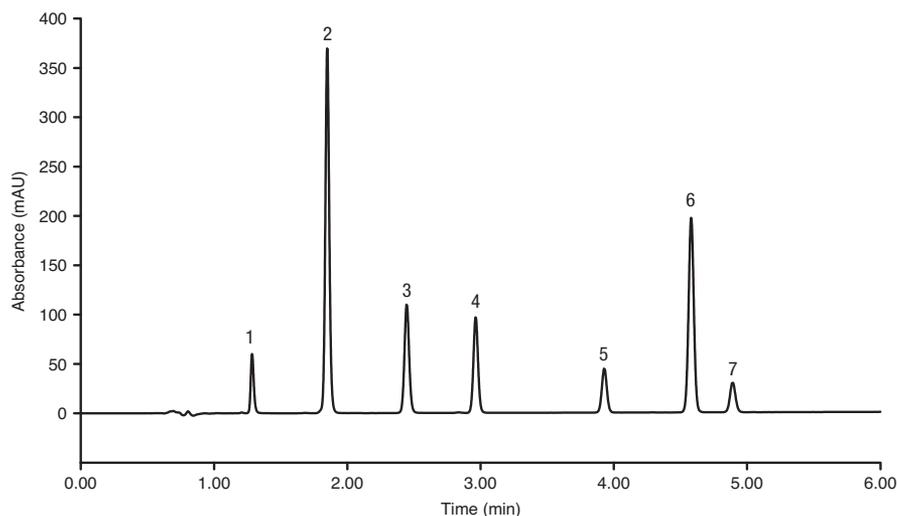


Figure 1: Separation of non-steroidal anti-inflammatory drugs

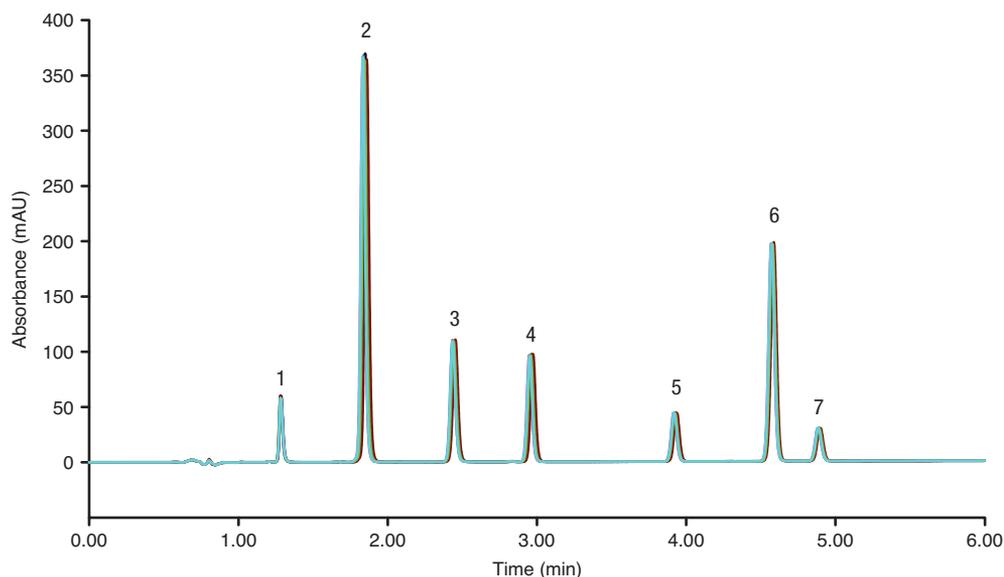


Figure 2: Overlaid chromatograms showing method robustness

Peak ID	Retention Time (min)	Asymmetry (EP)	Resolution(USP)	T <sub>R</sub> RSD (%) n=6
1	1.28	1.21	12.05	0.03
2	1.86	1.08	9.97	0.11
3	2.45	1.08	8.13	0.06
4	2.97	1.04	14.52	0.07
5	3.94	1.02	9.19	0.04
6	4.59	1.02	4.18	0.03
7	4.90	1.05	n.a.	0.02

Table 1: Results summary

## Conclusion

The Synchronis C18 column successfully separated all seven non-steroidal anti-inflammatory drugs with a relatively short retention time and a very simple method. This application demonstrates that Synchronis C18 columns

- retain and resolve non-steroidal anti-inflammatory drugs
- give excellent peak shape
- provide excellent resolution
- give consistent and highly reproducible methods

The method overall proved to be robust, simple and highly reproducible.

For Research Purposes Only

[thermoscientific.com/chromatography](http://thermoscientific.com/chromatography)

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific Inc. products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. 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

**Thermo**  
 SCIENTIFIC

Part of Thermo Fisher Scientific