

# HPLC Analysis of Zileuton

Monica Dolci, Thermo Fisher Scientific, Runcorn, Cheshire, UK

## Key Words

- Zileuton
- HPLC

## Abstract

This application note demonstrates the use of the Thermo Scientific Synchronis C18 column for the analysis of zileuton.

## Introduction

Zileuton is a 5-lipoxygenase inhibitor [1] which inhibits the immunoregulators leukotrienes. Leukotrienes are involved in defence mechanisms such as inflammations. Controlling their biosynthesis can lead to the treatment of asthma and other allergic conditions. Zileuton is used for the maintenance treatment of asthma.

The original immediate-release formulation of zileuton, known as ZYFLO, was introduced in 1996 by Abbott Laboratories. The structure of this compound is shown in Figure 1.

Zileuton is a moderately polar compound, with a log P of 1.4, and therefore its retention under reversed phase LC requires a retentive stationary phase.

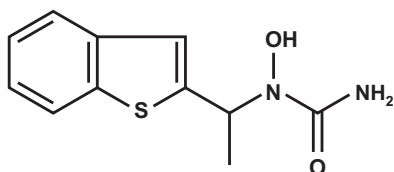


Figure 1: Structure of zileuton

Synchronis™ C18 columns are based on highly pure 100 Å silica, with a surface area of 320 m<sup>2</sup>/g and a carbon load of 16%. This ensures good retention of analytes with a range of hydrophobicities.

This application note shows an efficient and reproducible method for the analysis of zileuton, with excellent peak shape.



## Experimental Detail

Chemicals and Reagents	Part Number
Fisher Scientific HPLC grade water	W/0106/17
Fisher Scientific HPLC grade methanol	M/4056/17
Thermo Scientific 2 mL clear vial and Si/PTFE seal	60180-600

## Sample Preparation

A 1000 µg/mL standard solution of zileuton was prepared in mobile phase; this solution was then diluted to 20 µg/mL in mobile phase and used for analysis.

Separation Conditions	Part Number
Instrumentation:	Thermo Scientific Accela HPLC/UHPLC system
Column:	Synchronis C18 5 µm, 50 x 2.1mm 97105-052130
Injection volume:	1 µL
Flow rate:	0.2 mL/min
UV detection:	227 nm
Mobile phase:	30:70 (v/v) acetonitrile/water

## Results

Zileuton was retained on the Synchronis C18 column using an isocratic method. Figure 2 shows the chromatogram obtained employing Synchronis C18 5  $\mu\text{m}$ , 50 x 2.1 mm column, with zileuton eluting at 3.49 min (retention factor of 3.36).

The following chromatographic parameters were monitored: retention time ( $t_R$ ), peak area, tailing factor and efficiency. Mean and % RSD values for the above parameters were based on data derived from six replicate injections, and are reported in Table 1.

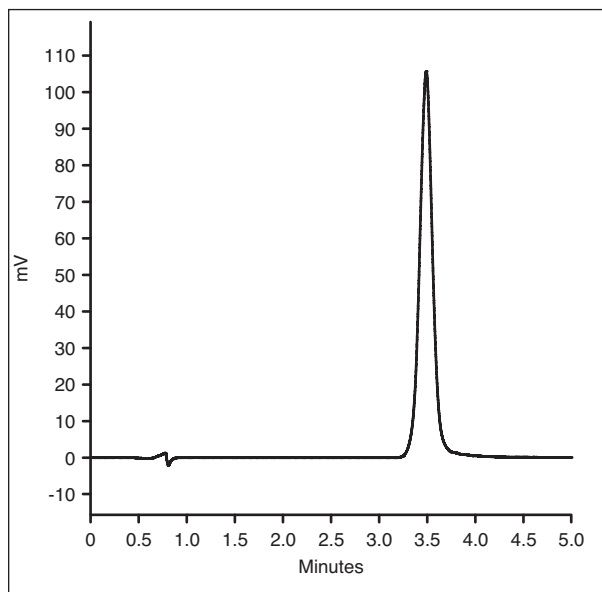


Figure 2: Chromatogram of zileuton, separated on a Synchronis C18 5  $\mu\text{m}$ , 50 x 2.1 mm column

Zileuton	$t_R$ (min)	Tailing Factor	Efficiency (Plates/m)	Peak Area
Mean	3.49	1.01	63646	10121320
% RSD	0.02	0.39	0.24	0.06

Table 1: Average and Method Precision (%RSD) of chromatographic parameters, for the analysis of zileuton on a Synchronis C18 5  $\mu\text{m}$ , 50 x 2.1 mm column (data calculated from six replicate injections)

## Conclusions

The use of a Synchronis C18 column allowed the successful retention of zileuton. Synchronis C18 columns are an excellent choice for the reversed-phase separation of zileuton, affording highly reproducible analysis.

## References

[1] F.J. Alvarez and R.T. Slade, *Pharmaceutical Research*, Vol 9, N 11, 1992.

*In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.*

**North America  
USA and Canada**  
+1 800 332 3331

**Europe  
France**  
+33 (0)1 60 92 48 34

**Germany**  
+49 (0) 2423 9431 -20  
or -21

**United Kingdom**  
+44 1928 534110

**Asia  
Japan**  
+81 3 5826 1615

**China**  
+86-21-68654588  
or +86-10-84193588  
800-810-5118

**India**  
+91-22-6742 9494

**Thermo Fisher  
Scientific Australia  
Pty Ltd**  
1300 735 292 (free call  
domestic)

**Thermo Fisher  
Scientific New  
Zealand Ltd**  
0800 933 966 (free call  
domestic)

**All Other Enquiries**  
+44 (0) 1928 534 050

## Technical Support

**North America**  
800 332 3331

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

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

© 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.

ANCCSZILEUTSYNC 0112