

# Fast Analysis of Zidovudine by UHPLC Using a Synchronis C18 1.7 $\mu\text{m}$ Column

M. Dolci, Thermo Fisher Scientific, Runcorn, Cheshire, UK

## Key Words

- Zidovudine
- UHPLC
- USP
- Synchronis C18

## Abstract

Thermo Scientific Synchronis C18 is a highly retentive stationary phase, with a surface area of 320  $\text{m}^2/\text{g}$ .

This application note demonstrates the use of Synchronis™ C18, packed with 1.7  $\mu\text{m}$  particles, for the fast HPLC analysis of zidovudine, allowing an increase in sample throughput. The method of analysis was adapted from the USP monograph, by using an in-house method transfer calculator [1].

## Introduction

Zidovudine, or azidothymidine (tradename: Retrovir) is an antiretroviral drug used for the treatment of HIV/AIDS. It acts by stopping HIV from infecting healthy cells, but it can not help cells already infected with the virus [2]. It was the first HIV treatment to be marketed and it is now a widely used generic drug.

Zidovudine is included in the World Health Organisation's "Essential Drugs List" (list of minimum medical needs for a basic healthcare system) [3].

The United States Pharmacopeia (USP) provides worldwide guidance for the chromatographic analysis of zidovudine [4], which is based on High Performance Liquid Chromatography (HPLC) L1 columns. The implementation of an L1 type column, such as Synchronis C18, packed with 1.7  $\mu\text{m}$  particles, allowed for the zidovudine USP method to be transferred to UHPLC (ultra High Performance Liquid Chromatography).

## Experimental details

The analysis was run on an Thermo Scientific Accela UHPLC system. The data was acquired and processed using Thermo Scientific ChromQuest 5.0 Software.

The separation of zidovudine was achieved on a Synchronis C18 column, a recently developed, highly retentive reversed phase material, specifically designed to target consistent chromatography for a wide range of compounds.



## Sample Preparation

A 1000  $\mu\text{g}/\text{mL}$  of zidovudine standard solution was prepared in methanol; this solution was then diluted to 100  $\mu\text{g}/\text{mL}$  in mobile phase (20:80 (v/v) methanol/water) and used for the analysis.

Thermo Scientific Column	Part Number
Synchronis C18 1.7 $\mu\text{m}$ , 50 x 2.1 mm	97102-052130

## Thermo Scientific Accela HPLC/UHPLC

Column temperature	25 $^{\circ}\text{C}$
Injection volume	0.5 $\mu\text{L}$
Flow rate	0.8 $\text{mL}/\text{min}$
UV detection	265 nm
Mobile phase	20:80 (v/v) methanol/water

Consumables	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

## Results

The original USP analytical conditions, based on a 5 µm, 250 x 4.0 mm column were scaled down using our method transfer calculator to accommodate for the column geometry reduction (as shown in Figure 1). The final run time achieved on the 1.7 µm, 50 x 2.1 mm column was eight times faster than the original run, with zidovudine eluting at 2.42 min (retention time of 26 min on the 5 µm, 250 x 4.0 mm column). Figure 2 shows the chromatograms obtained employing Synchronis C18 1.7 µm, 50 x 2.1 mm column and Synchronis C18 5µm, 250 x 4.0 mm column.. The USP acceptance criteria were met, as demonstrated in Table 1. The statistical assessment is based upon data derived from six replicate injections (see Table 1 for mean values).

**HPLC Method Development Calculator**

**Isocratic Method Transfer**

Enter your dimensions / particle sizes in the green boxes below.

Existing Method		New Method	
Parameter	Value	Parameter	Value
Column geometry	5 µm	Column geometry	1.7 µm
Particle Size	250 mm	Particle size	50 mm
Column length	4.0 mm	Column length	2.1 mm
Column ID		Column ID	
Method conditions		Method conditions	
Flow rate	1 ml/min	Flow Rate	0.81 ml/min
Injection volume	10 µl	Injection volume	0.95 µl
Analysis Time	30 min	Analysis Time	2 min
Pressure	54 bar	Pressure	275 bar
		Adjusted Flow Rate	1.7 ml/min
			1.7 min
			339 bar

**Comparisons**

Existing Method vs. New Column: Analysis Time (13x faster), Pressure (5x higher)

Existing Method vs. Variable Flow: Analysis Time (13x faster), Pressure (6x higher)

Time Saved: 28 min, 28.3 min

Solvent Saved: 28.3 ml

Figure 1: Isocratic method transfer calculator used in this application note to scale down the original USP chromatographic conditions [1]

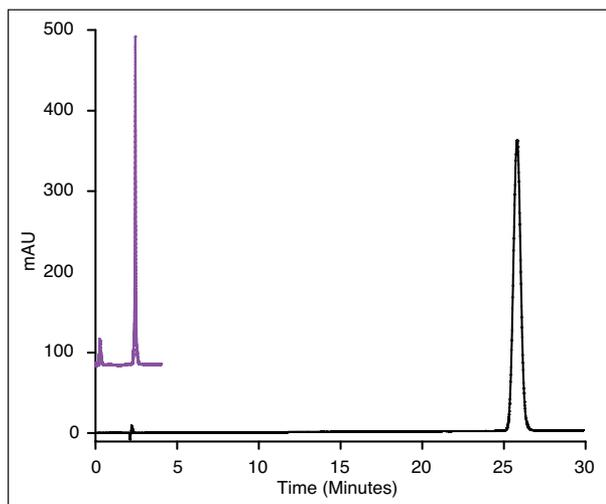


Figure 2: Chromatograms of 100 µg/mL of zidovudine separated on a Synchronis C18 1.7 µm, 50 x 2.1 mm column (top chromatogram), and on a Synchronis C18 5 µm, 250 x 4.0 mm column (bottom chromatogram)

## Conclusions

The use of a Synchronis C18, packed with 1.7 µm particles column, allowed to successfully scale down the USP method for the analysis of zidovudine, in order to significantly increase sample throughput. The analytical results passed the specifications stated in the USP monograph. Synchronis C18 1.7 µm columns are therefore an excellent choice for the high speed, high efficiency analysis of zidovudine, allowing a high sample throughput.

## References

- <http://www.separatedbyexperience.com/products/IsocraticMethod.aspx>
- W. V. Caufield; J. T. Stewart; *Chromatographia* (2001); Vol 54; pp 561-568.
- [http://whqlibdoc.who.int/hq/2005/a87017\\_eng.pdf](http://whqlibdoc.who.int/hq/2005/a87017_eng.pdf). World Health Organization. March 2005 Retrieved 2006-03-12
- [http://www.pharmacopeia.cn/v29240/usp29nf24s0\\_m89510.html](http://www.pharmacopeia.cn/v29240/usp29nf24s0_m89510.html)

	T <sub>r</sub> (min)	Tailing Factor	Efficiency	Peak Area
Average	2.416	0.970	118567	242096
% RSD	0.05	0.30	0.78	0.79
USP acceptance criteria	≤ 2.0 (% RSD)	≤ 2.0	≤ 2.0 (% RSD)	≤ 2.0 (% RSD)

Table 1: Average and Method Precision (%RSD) of chromatographic parameters, derived from the analysis of zidovudine on a Synchronis C18 1.7 µm, 50 x 2.1 mm column (data calculated from six replicate injections) USP acceptance criteria are also reported.

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

**Switzerland**  
+41 56 618 41 11

**United Kingdom**  
+44 1928 534110

**Asia  
Japan**  
+81 3 5826 1615

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

**India**  
1800 22 8374 (toll-free)  
+91 22 6716 2200

**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)

© 2011 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.

ANCCSSYNZIDOV 0811