

Analysis of Artesunate and Dihydroartemisinin Using a Synchronis C18 HPLC Column

Joanne Gartland, Thermo Fisher Scientific, Runcorn, Cheshire, UK

Key Words

- Synchronis C18
- Artesunate
- Dihydroartemisinin
- Malaria

Abstract

This application note demonstrates the use of the Thermo Scientific Synchronis C18 HPLC column for the analysis of artesunate and its active metabolite dihydroartemisinin.

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.

Artesunate is obtained as an extract from the plant *Artemisia annua* and is commonly used for the treatment of malaria. Artesunate (active metabolite dihydroartemisinin) and artesunate-based combination therapy (ACT) is recommended by the World Health Organization (WHO) for the treatment of severe and multidrug resistant malaria. The analytical method for artesunate described in the WHO monograph, uses a mobile phase which includes potassium phosphate, which is incompatible with mass spectrometry [1]. Considering the C_{MAX} of artesunate is $0.09 \pm 0.04 \mu\text{g/mL}$ [2] and that artesunate lacks an intensive chromophore for UV absorption, MS detection is required to give the required sensitivity. Within this application note we will demonstrate an alternative method using an MS-compatible mobile phase.



Experimental Details

Chemicals and Reagents	Part Number
Fisher Scientific HPLC grade water	W/0106/17
Fisher Scientific HPLC grade acetonitrile	A/0626/17

Sample Handling Equipment	Part Number
NSC Mass Spec Certified 2 mL clear vial with blue bonded PTFE silicone cap	MSCERT4000-34W

Sample Preparation
Working standard contained 1 mg/mL of artesunate and 2 mg/mL of α and β isomers of dihydroartemisinin in acetonitrile.

Separation Conditions	Part Number	
Instrumentation:	Accela 600 HPLC System	
Column:	Synchronis™ C18 5 μm , 100 x 2.1 mm	97105-102130
Measured pressure:	50 bar	
Column temperature:	30 °C	
Injection volume:	1 μL	
Flow rate:	0.2 mL/min	
UV detection:	210 nm	
Mobile phase:	40:60 (v/v) water + 0.1% formic acid/ acetonitrile + 0.1% formic acid	

Results

The analysis was carried out on a Synchronis C18 5 μ m, 100 x 2.1 mm column. As shown on Figure 1, artesunate and its active metabolites, α and β dihydroartemisinin are eluted in less than 5 minutes with outstanding peak symmetry. The low peak intensity (Figure 1) is due to the lack of a UV chromophore indicating that the method would be suited to MS detection.

Replicate injections of the test mix showed that Synchronis C18 produced highly reproducible retention and peak shape (Table 1).

	α -Dihydroartemisinin	β -Dihydroartemisinin	Artesunate
Peak position	1a	1b	2
Average As	1.05	1.05	1.05
Average Rs	0.00	4.49	2.52
%RSD t_R	0.06	0.11	0.10

Table 1: Method precision (%RSD) for artesunate and α and β dihydroartemisinin (data calculated from six replicate injections)

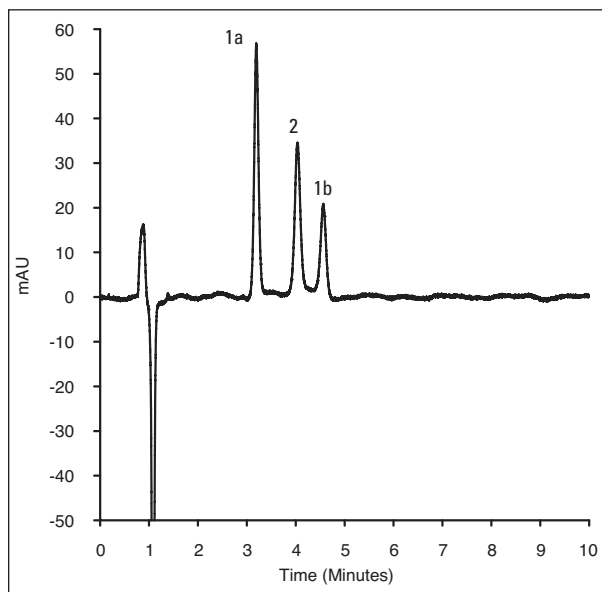


Figure 1: Chromatogram for α and β dihydroartemisinin (1a), (1b) and artesunate (2) separated on a Synchronis C18 5 μ m, 100 x 2.1 mm column

Conclusions

Synchronis C18 successfully separated artesunate and dihydroartemisinin in less than 5 minutes. Synchronis C18 columns are an excellent choice for fast highly reproducible analysis with a MS compatible mobile phase.

References

- [1] Artesunate: Final text for revision of The International Pharmacopoeia (December 2009) World Health Organisation (2009). Working document QAS/09.340/Final December 2009.
- [2] Am. J. Trop. Med. Hyg., 58(3), 1998, pp.365-368

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

© 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.

ANCCSSYNCARTDIH 1011