# Application Note: ANCCSCETARTDIHY

# Analysis of Artesunate and Dihydroartemisinin Using a Core Enhanced Technology Accucore HPLC Column

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

## **Key Words**

- Accucore RP-MS
- Fused core
- Superficially porous
- Core Enhanced Technology
- Artesunate
- Dihydroartemisinin
- Malaria

# Abstract

This application note demonstrates the use of the Thermo Scientific Accucore RP-MS HPLC column for the fast analysis of artesunate and its active metabolite dihydroartemisinin.

## Introduction

Accucore<sup>™</sup> HPLC columns use Core Enhanced Technology<sup>™</sup> to facilitate fast and high efficiency separations. The 2.6 µm diameter particles are not totally porous, but rather have a solid core and a porous outer layer. The optimized phase bonding creates a series of high coverage, robust phases. Accucore RP-MS uses an optimized alkyl chain length for more effective coverage of the silica surface. This coverage results in a significant reduction in secondary interactions and thus highly efficient peaks with very low tailing. The tightly controlled 2.6 µm diameter of Accucore particles results in much lower backpressures than typically seen with sub-2 µm materials.

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<sub>MAX</sub> of artesunate is 0.09+/-0.04 µg/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, in addition to this the speed of analysis that can be obtained using Accucore RP-MS will also be demonstrated



## **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

NSC Mass Spec Certified 2 mL clear vial with blue	MSCERT4000-34W
bonded PTFE silicone cap	

#### Sample Preparation

Working standard contained 1 mg/mL of artesunate and 2 mg/mL of  $\alpha$  and  $\beta$  isomers of dihydroartemisinin in acetonitrile.

#### **Separation Conditions**

Accela 600 HPLC System	
Accucore RP-MS 2.6 μm,     17626-052130       50 x 2.1 mm     17626-052130	
185 bar	
30 °C	
1 µL	
0.6 mL/min	
210 nm	
60:40 (v/v) water + 0.1% formic acid/ acetonitrile + 0.1% formic acid	



# Results

The analysis was carried out on an Accucore RP-MS 2.6  $\mu$ m, 50 x 2.1 mm column. Artesunate and its active metabolites,  $\alpha$  and  $\beta$  dihydroartemisinin are eluted in under 2 minutes. The low peak intensity (Figure 1) is due to the lack of UV chromophore indicating that the method would be suited to MS detection.

Replicate injections of the test mix showed that Accucore RP-MS produced reproducible retention and peak shape (Table 1).

	α- Dihydroartemisinin	$\beta$ -Dihydroartemisinin	Artesunate
Peak position	1а	1b	2
Average As	1.26	1.16	1.25
Average Rs	0.00	7.38	3.15
%RSD t <sub>R</sub>	0.43	0.46	0.44

Table 1: Method precision (%RSD) for artesunate and  $\alpha$  and  $\beta$  dihydroartemisinin (data calculated from six replicate injections)

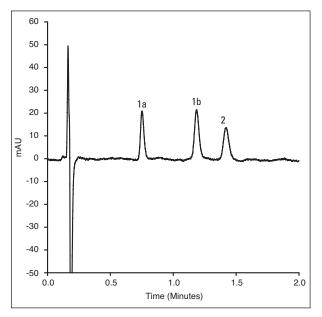


Figure 1: Chromatogram for  $\alpha$  and  $\beta$  dihydroartemisinin (1a), (1b) and artesunate (2) separated on an Accucore RP-MS 2.6  $\mu m$ , 50 x 2.1 mm column

# Conclusions

Accucore RP-MS successfully separated artesunate and dihydroartemisinin in less than 2 minutes. Accucore RP-MS columns are an excellent choice for fast analysis with a MS compatible mobile phase.

## References

 Artesunate: Final text for revision of The International Pharmacopoeia (December 2009) World Health Organisation (2009). Working document QAS/09.340/Final December 2009.
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

+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

**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.

