

HPLC-UV Method for the Determination of Telmisartan Using a Core Enhanced Technology Accucore Column

Eilidh MacRitchie, Thermo Fisher Scientific, Runcorn, Cheshire, UK

Key Words

- Telmisartan
- Accucore RP-MS
- Hypertension
- Solid Core
- Core Enhanced Technology
- High speed

Abstract

This application note demonstrates the use of the Thermo Scientific Accucore RP-MS column for the determination of telmisartan by HPLC-UV.

Introduction

Accucore™ HPLC columns use Core Enhanced Technology™ 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 optimised 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.

Telmisartan is an angiotension II receptor antagonist used for the treatment of hypertension.

This application note demonstrates the successful analysis of telmisartan using an Accucore RP-MS column.



Experimental Details

Chemicals and Reagents	Part Number
Fisher Scientific HPLC grade water	W/0106/17
Fisher Scientific HPLC grade ammonium acetate	A/3446/50
Fisher Scientific HPLC grade acetonitrile	A/0626/17
Telmisartan purchased from Sigma Aldrich	

Sample Handling Equipment

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

Separation Conditions	Part Number
Instrumentation:	Thermo Scientific Accela UHPLC system
Column:	Accucore RP-MS 2.6 µm, 50 x 2.1 mm 17626-052130
Mobile phase:	70:30 (v/v) 20 mM ammonium acetate/acetonitrile
Flow rate:	0.4 mL/min
Column temperature:	25 °C
Injection details:	1 µL partial loop
Injection wash solvent:	water
UV detector wavelength:	230 nm
Backpressure:	150 bar

Solutions

Working standard contained 20 µg/mL of telmisartan in water.

Results

The analysis was performed on an Accucore RP-MS 2.6 μm , 50 x 2.1 mm column. As shown in Figure 1, telmisartan was analyzed in less than 2 minutes. Table 1 shows the results from six replicate injections.

	Telmisartan
Retention time (minutes)	1.15
%RSD on retention time	0.2
Asymmetry	1.31
%RSD on asymmetry	0.7

Table 1: Retention time and asymmetry results for telmisartan

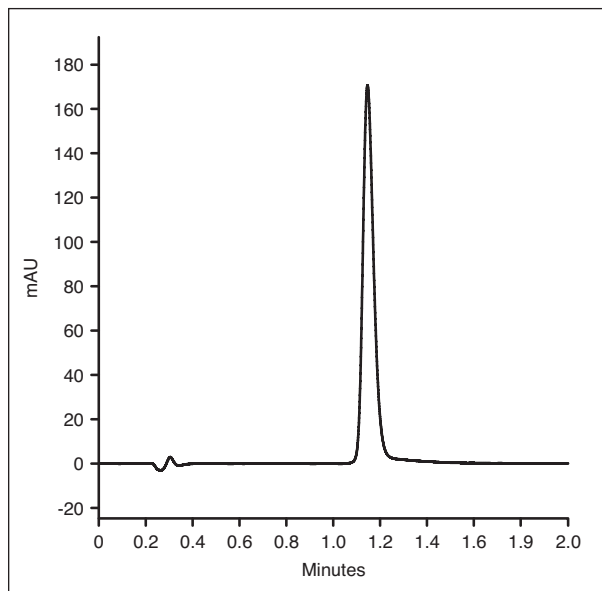


Figure 1: Chromatogram of telmisartan analyzed using an Accucore RP-MS 2.6 μm , 50 x 2.1 mm column

Conclusions

Replicate injections of telmisartan showed that Accucore RP-MS produced stable and reproducible results. This demonstrates that Accucore RP-MS is an excellent choice of column for the rapid analysis of telmisartan.

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

© 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.

ANCCSCETELMIS 1211