Application Note: ANCCSMETFORMIN

Analysis of Metformin on a Perfluorophenyl Stationary Phase by HPLC/UV

W. Faulkner, Thermo Fisher Scientific, Runcorn, Cheshire, UK

Introduction

Metformin

Key Words

- Melamine
- Polar compounds
- Perfluorophenyl phase,
- Hypersil GOLD PFP
- Metformin (500 mg) tablets

Metformin [3-(diaminomethylidene)-1,1dimethylguanidine] is an orally administered, antihyperglycaemic drug used for the treatment of Type 2 (non-insulin dependent) diabetes mellitus. Metformin improves glycaemic control by both enhancing insulin sensitivity, and, by decreasing intestinal absorption of glucose.

As a direct consequence of its high polarity $(\log P = -2.64)$, metformin is difficult to retain on conventional stationary phases (e.g., C18 and C8) employed in reversed-phase chromatography [1]. Moreover, the strong interaction with polar media (e.g., silica and cyano) may mean that the biguanide drug cannot be eluted easily during its examination by normal phase chromatography [2, 3].

In view of the potential shortcomings associated with both of the above, alternative chromatographic approaches may be adopted for the separation of metformin and its related compounds (e.g., cyanoguanidine, melamine and 1-methylbiguanide), including ion-pair chromatography [4,5], hydrophilic interaction liquid chromatography (HILIC) [6,7], aqueous normal phase (ANP) chromatography [8,9] and ionexchange (IEX) chromatography [10,11]. In fact, the US pharmacopeial procedure for the inspection of metformin specifies that separations must be performed in an ionexchange manner using USP L9 column packings. However, ion-exchange techniques can suffer from some practical constraints. In particular, when used with MS, both the precipitation of salt and the suppression of ionisation of the analyte(s) under investigation can be observed.

Despite the existence of a variety of chemically bonded phases for the retention of hydrophilic molecules, perfluorinated phases (e.g., pentafluorophenyl) have attracted particular attention [12, 13]. Perfluorinated and fluorinated stationary phases have been found to exhibit novel selectivity for several classes of compounds and, in many instances, provide appealing alternatives to traditional C18 and C8 phases.

The purpose of this chromatographic study is to demonstrate the effectiveness of a perfluorophenyl stationary phase for the retention of metformin and its related compound melamine under typical reversed-phase conditions.



Experimental details

Chemicals and reagents

Fisher Scientific Ammonium acetate (AR grade)	A/3440/50	
Fisher Scientific Ammonia solution (HPLC grade, SG – 0.88, 35 %)	A/3295/PB05	
Fisher Scientific Acetonitrile (HPLC grade)	A/0626/17	
Fisher Scientific Water (HPLC gradient grade)	W/0106/17	
Metformin (500 mg) tablets		
Metformin Hydrochloride (≥ 98 %)		
Melamine (≥ 99 %)		

Sample handling equipment

Fisher Scientific Finnipipette F2 pipettor kit 10 μL – 100 μL, 100 μL – 1000 μL, 1 mL – 10 mL	PMP-020-220F
Fisher Scientific Finnitip pipette tips, 10 µL	PMP-107-110W
Fisher Scientific Finnitip pipette tips, 200 µL	PMP-107-600F
Fisher Scientific Finnitip pipette tips, 1000 µL	PMP-103-206K
Fisher Scientific Finnitip pipette tips, 10 mL	PMP-107-040R
Fisher Scientific Syringe filters (PVDF filter membrane, 30 mm 0.D., 0.45 µm pores)	HCA-442-013Q
Thermo Scientific Borosilicate glass vials (2 mL, 12 mm x 32 mm) with 8 mm black screw cap fitted with a silicone/PTFE seal	60180-600

Sample Preparation

Analytical standards

Primary analytical standards of both metformin hydrochloride and melamine were prepared separately by the dissolution of approximately 0.01 g (weighed accurately) of reference material in water (10.0 mL). The concentration of each analyte was approximately 1000 µg/mL.

Thereafter, a mixed working standard was prepared by combining 1 part of each primary standard with 98 parts of mobile phase (90:10 (v/v) NH₄OAc -MeCN). The concentration of each analyte in the working standard was approximately 10 µg/mL



Extraction of Metformin (500 mg) tablets

A number (x 4) of metformin tablets was ground to a powder using a mortar and pestle. Thereafter, a portion (approximately 120 mg) of the powder, containing a nominal quantity of about 100 mg of metformin hydrochloride, was transferred into a conical flask and subsequently brought into contact with water (100 mL). Following agitation of the mixture for 30 minutes at room temperature, a portion of the slurry (10.0 mL) was removed and allowed to pass through a 0.45 μ m PVDF filter membrane. After discarding the first few millilitres of solution, the resultant, clear liquid was collected. Finally, the aqueous extract was diluted 100-fold by mixing 1 part of aqueous extract with 99 parts of mobile phase (90:10 (v/v) NH₄OAc – MeCN).

Instrumentation

Thermo Scientific HPLC system equipped with a PDA detector. Data were acquired and processed using ChromQuest 4.2 software.

Chromatographic conditions

Column - Thermo Scienti 150 mm x 4.6 mm	fic Hypersil GOLD PFP 5µm	PN: 25405-154630
Mobile phase	NH ₄ OAc 20 mM, pH - 7.98	/ MeCN (90:10 v/v)
Flow rate	1.0 mL/min	
Column temperature	30 °C	
UV detection	233 nm	
Injection volume	5 µL	
Run time	5 minutes	
Syringe flush	mobile phase	

Results

Under the isocratic conditions adopted for this analysis, adequate retention and separation of these polar molecules can be accomplished in less than 5 minutes. The chromatographic performance of the pentafluorophenyl (PFP) phase during the analysis of a mixture of metformin and melamine is summarised in Table 1. Chromatographic data are matched with excellent precision. Comparative chromatograms of both a standard solution of the polar analytes and a solution derived from an aqueous extract of an amount of powdered 'Metformin 500 mg' tablets are shown in Figure 1.

This qualitative approach precludes any accurate determination of the content of metformin (and melamine) in the tablets. Despite the fact that a detailed mechanism accurately describing the retentive behaviour of the perfluorophenyl phase is still to be elucidated, it has been proposed that the interaction between the carbon-fluorine dipole (on the aromatic PFP ring) and the amino groups in the metformin molecule is primarily responsible for the observation of enhanced retention.

	Retention time (min)		Efficiency (USP Plates/m)		Tailing factor (USP)		Resolution (USP)	
	Mean	%RSD	Mean	%RSD	Mean	%RSD	Mean	%RSD
Melamine	2.101	0.09	71916	0.64	1.32	1.18	-	-
Metformin	3.322	0.23	51755	0.50	1.66	1.02	10.54	0.24

Table 1. Chromatographic performance of Hypersil GOLD® PFP (data derived from 10 replicate injections)



Figure 1. Comparative chromatographic profiles: Upper trace – aqueous extract of metformin (500 mg) tablets; lower trace – standard solution of metformin and melamine; 1. melamine, 2. metformin

Conclusions

Perfluorophenyl bonded phases may be used for the successful retention of the hydrophilic molecules metformin and melamine. Moreover, good retention and separation can be achieved under typical reversed-phase conditions in less than five minutes.

References

- Tache, F., David, V., Farca, A., Medvedovici, A., J. of Microchem., 2001, 68, 13.
- AbuRuz, S., Millership, J., McElnay, J., J. of Chromatography B, 2003, 798 (2), 203
- Chen, X., Gu, Q., Qiu, F., Zhong, D., J. of Chromatography B, 2004, 802, 377
- 4. Ranetti, M. et al, Farmacia, 2009, 57 (6), 728
- David, V., A., Medvedovici, Albu, F., Journal of Liquid Chromatography and related technologies, 2005, 28 (1), 81
- 6. Liu, A., Coleman, S.P., J. of Chromatography B, 2009, 877 (29), 3695
- 7. Huttunen, K.M. et al, J. of Pharmaceutical and Biomedical analysis, 2009, 50 (3), 469
- Pesek, J., Matyska, M.T., LC/GC North America, May 2007
- 9. Pesek, J., Matyska, M.T., 'Conference on Small Molecule Science', Chapel Hill, North Carolina, July 2007
- 10. Bonfigli, A.R. et al, *Therapeutic Drug Monitoring*, 1999, 21 (3), 330
- 11. Rogers, L.A. et al, Journal of Liquid Chromatography and related technologies, 2009, 32 (15), 2246
- 12. Przybyciel, M., 'Recent developments in LC column technology', June 2003, p.2
- 13. Przybyciel, M., LC/GC, LC column technology supplement, April 2006, p. 49

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

Switzerland +41 56 618 41 11

United Kingdom +44 1928 534110

Asia Japan +81 3 5826 1615

China +86-21-68654588 or +86-10-8419358 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 demostic)

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.

