

# Strategies for the Transfer of Liquid Chromatographic Methods Between Different Instruments

Theresa Riley, Maria Grübner, Carsten Paul, Mauro De Pra, Frank Steiner, Thermo Fisher Scientific, Dornierstrasse 4, 82110 Germering, Germany

## ABSTRACT

**Purpose:** To evaluate the effect of different LC instrument parameters on method transfer.

**Methods:** Two pharmaceutical methods and five different LC instruments were utilized.

**Results:** Gradient delay volume, column thermostating, pre-column volume and DAD bandwidth settings should be considered during the transfer of LC methods between different LC instruments.

## INTRODUCTION

A challenging task that frequently occurs in all kinds of analytical industries is the transfer of liquid chromatographic (LC) methods from one instrument to another. This is straightforward in case the transfer is between identical instruments. However, the situation becomes more complicated when instruments of different configurations, generations and/or vendors are used. As all LC hardware components to some extent have influence on the chromatographic results, instrumental differences will also affect the analytical outcome of a transferred LC method. Method robustness as well as the degree of instrumental deviation determine the analytical deviation. Means to counteract these effects depend on the requirements of the operator. For example, if adequate resolution and congruent quantitative results are obtained and sufficing, no effort in adaption is needed. However if in addition retention times need to fit exact specifications, the effort might increase.<sup>1,2</sup> In our study we investigated several strategies to overcome difficulties in method transfer caused by hardware differences between several instrumental platforms. Quaternary and binary systems were considered. Here our focus was on Thermo Scientific™ Ultimate™ 3000 UHPLC Systems, Thermo Scientific™ Vanquish™ UHPLC Systems, Agilent Infinity 1260 and Waters Acuity. We examined strategies to modify system dwell volumes such as different mobile phase mixers and sample loop sizes as well as the adjustable delay volume of the Thermo Scientific™ Vanquish™ Split Sampler. The effect of pre-column volumes on peak shape was demonstrated for strong solvent injections and the impact of detector settings like the bandwidth was shown.

## MATERIALS AND METHODS

### Samples

Sample 1: solution of the active pharmaceutical ingredient (API) acetaminophen (1 mg/mL) and its impurities (according to USP<sup>3</sup>) B, C, D, J and 4-aminophenol (10 µg/mL each) in methanol

Sample 2: EP reference standard mebendazole for system suitability<sup>4</sup> (containing API and impurities A, B, C, D, E, F and G according to EP<sup>5</sup>) in dimethylformamide at 1 mg/mL

### Experimental

Chromatographic methods are given in Table 1 and were derived from the pharmacopeial monographs: USP monograph for acetaminophen<sup>3</sup> and EP monograph for mebendazole<sup>5</sup>. Instruments used in the current study are summarized in Tables 2 and 3. The acetaminophen assay was applied with quaternary systems, the mebendazole method with binary systems. System control and data analysis was performed with Thermo Scientific™ Chromeleon™ 7.2.9 CDS software.

Table 1. Chromatographic conditions

	Acetaminophen	Mebendazole
<b>Column</b>	Thermo Scientific™ Hypersil GOLD™ C8 column, 4.6x100 mm, 3 µm, 175 Å (p/n 25203-104630)	Thermo Scientific™ Hypersil GOLD™ C18 Selectivity LC column, 4.6x100 mm, 3 µm, 175 Å (p/n 25003-104630)
<b>Eluents</b>	A: 1.7 g/L KH <sub>2</sub> PO <sub>4</sub> and 1.8 g/L of Na <sub>2</sub> HPO <sub>4</sub> in water B: Methanol	A: 7.5 g/L Ammonium acetate in water B: Acetonitrile
<b>Gradient</b>	min % A % B 0 99 1 3 99 1 7 19 81 7.1 99 1 12 99 1	min % A % B 0 80 20 15 70 30 20 10 90 25 10 90 25.1 80 20 30 80 20
<b>Flow rate</b>	1 mL/min	1.2 mL/min
<b>Column temp.</b>	35 °C (with eluent preheating)	40 °C (with eluent preheating)
<b>Inj. volume</b>	1 µL	5 µL
<b>Detection</b>	230 nm, 10 Hz data collection rate, 0.5 s response time	250 nm, 10 Hz data collection rate, 0.5 s response time / normal filter time (0.2 s)

Table 2. Utilized quaternary systems

	Agilent 1260 Quaternary	UltiMate 3000 SD Quaternary	Vanquish Flex Quaternary
<b>Pump</b>	Quaternary pump (G1311B)	Standard quaternary pump LPG-3400SD (p/n 5040.0031)	Quaternary pump F (p/n VF-P20-A)
<b>Sampler</b>	High Performance Autosampler (G1367E) with thermostat module (G1330B)	Wellplate Autosampler WPS-3000TRS (p/n 5840.0020) with 7 µL eluent preheater (p/n 6722.0540)	Split Sampler FT (p/n VF-A10-A)
<b>Column Compartment</b>	TCC with 6 µL heat exchanger (G1316A)	TCC-3000SD (p/n 5730.0010)	Column Compartment H (p/n VH-C10-A)
<b>Detector</b>	Diode array detector DAD VL (G1315D)	Diode array detector DAD-3000 (p/n 5082.0010)	Diode array detector DAD FG (p/n VF-D11-A)
<b>Flow Cell</b>	standard: 10 mm, 13 µL (G1315-60022)	analytical: 10 mm, 13 µL (p/n 6082.0100)	standard bio: 10 mm, 13 µL (p/n 6083.0540)

Table 3. Utilized binary systems

	Acuity	Vanquish Horizon
<b>Pump</b>	Binary Solvent Manager	Binary Pump H (p/n VH-P10-A)
<b>Sampler</b>	Sample Manager	Split Sampler FT (p/n VF-A10-A)
<b>Sample loop</b>	10 µL	default 25 µL (V=50 µL, p/n 6850.1911)
<b>Column Compartment</b>	High Temperature Column Heater	Column Compartment H (p/n VH-C10-A)
<b>Detector</b>	Tunable Ultraviolet Detector	Variable Wavelength Detector F (p/n VF-D40-A)
<b>Flow Cell</b>	analytical (10 mm, 500 nL)	semi-micro (7 mm, 2.5 µL, p/n 6077.0360)

## RESULTS

### Gradient delay volume adaption

The gradient delay volume (GDV) of a LC system is defined as the volume between the point of gradient mixing and the column entry. Contributors are pump, sampler and capillary volumes. As the GDV delays the arrival of a particular solvent composition at the column it has a strong impact on elution times. Thus during method transfer GDV adaptations are frequently applied to compensate retention time differences between the sending and receiving LC system.<sup>2</sup> For the transfer of the acetaminophen assay from a Thermo Scientific™ UltiMate™ 3000 SD system to a Thermo Scientific™ Vanquish™ Flex system only a fine-tuning of the GDV was required, which can be accomplished by the adaption of the idle volume of the metering device in the Vanquish autosampler. Figure 1 shows the working principle of that device and Figure 2 the overlaid chromatograms before and after adaption.

If the range of the idle volume (up to 100 µL) is not sufficient for the GDV adjustment the next level is the replacement of the sample loop by a higher volume one. Figure 3 depicts an example for the transfer from an Agilent 1260 system to a Vanquish Flex system by exchanging the default loop (25 µL, V=50 µL) by the 100 µL loop (V=130 µL) and fine tune by the metering device. However, for GDV differences of major amount a change of the mobile phase mixer should be considered. In Figure 4 that approach is shown for the transfer from an UltiMate 3000 SD instrument to an Agilent 1260 system. Substitution of the 350 µL static mixer by the 750 µL mixer distinctly overcompensated the actual GDV difference. Thus a gradient prestart was applied to match retention times, meaning that the gradient program started at a negative time (-0.27 min) and the injection was conducted at 0 min. All three strategies achieved very good results for all compounds, which eluted during the gradient, without any detrimental effects on the chromatographic performance. However, the GDV does not affect isocratically eluted peaks like the 4-aminophenol peak in Figures 2-4. Mismatches here might be caused by slight deviations in mobile phase proportioning or column thermostating.

Figure 1. Vanquish autosampler metering device

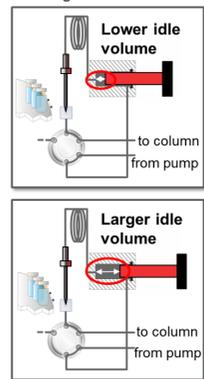


Figure 2. Acetaminophen assay before (A) and after (B) GDV adaption by the idle volume of the metering device in the Vanquish autosampler

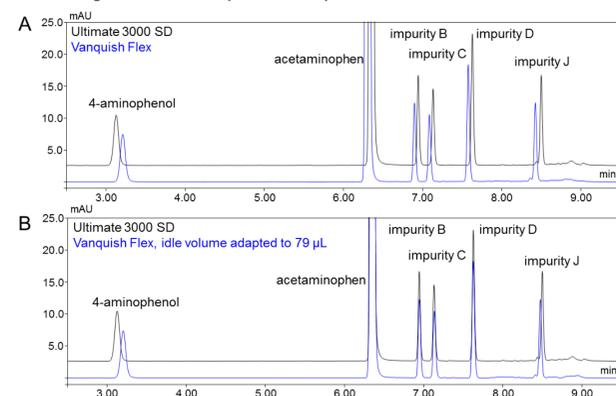


Figure 3. Acetaminophen assay before (A) and after (B) GDV adaption by sample loop and idle volume of the metering device in the Vanquish autosampler

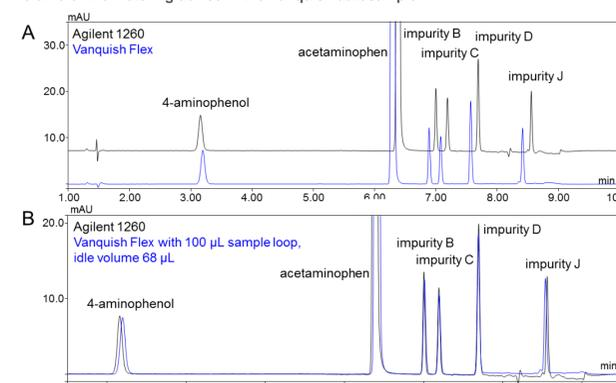


Figure 4. Acetaminophen assay before and after GDV adaption by static mixer exchange (A) and after applying a gradient prestart

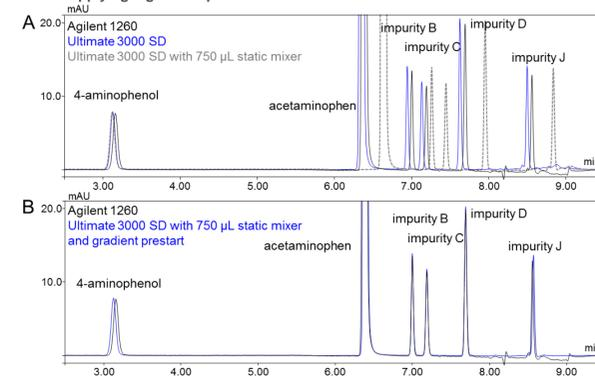
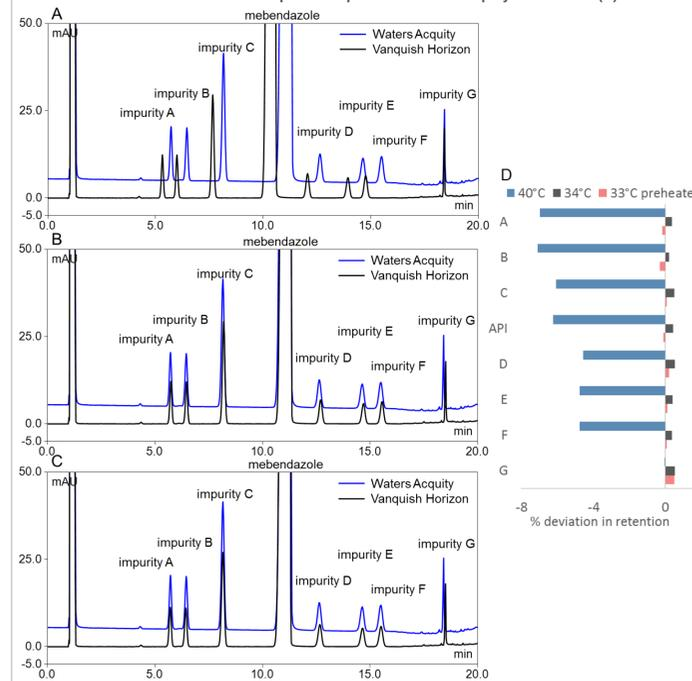


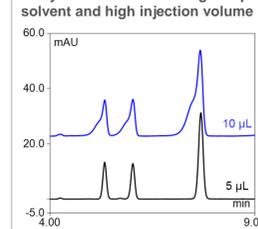
Figure 5. Mebendazole analysis with Waters Acuity system set to 40 °C and Vanquish Horizon system set to 40 °C (A), 34 °C (B) and 40 °C with active eluent preheating set to 33 °C; % deviation of retention times of Vanquish compared to Waters Acuity is shown in (D)



### Column thermostating

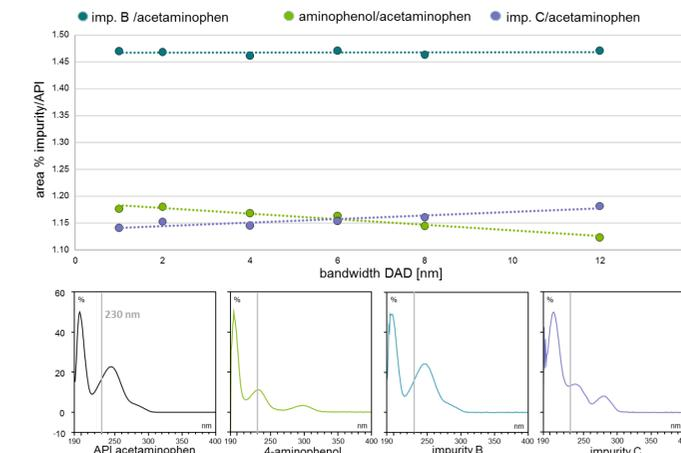
Very distinct effects of unequal column thermostating were observed during the transfer of a mebendazole analysis from a Waters Acuity system to a Vanquish Horizon system. Large differences in retention time, which were observed when both column thermostats were nominally set to 40 °C, were eliminated by reducing the temperature to 34 °C at the Vanquish system or by reducing only the temperature of the active solvent preheater to 33 °C (Figure 5), indicating a more efficient column thermostating of the Vanquish system.

Figure 6. Peak fronting in mebendazole analysis as result of strong sample solvent and high injection volume



As the peak area ratio is frequently used for relative quantification, we compared the outcome of several bandwidth settings of a Vanquish DAD. Figure 7 shows very similar UV spectra for the API and impurity B. Thus their area ratio is constant. In contrast, the spectra of 4-aminophenol and impurity C deviate from the API. As a consequence the relative quantification is affected by the bandwidth setting (in different directions for both compounds). Thus, detector settings should be carefully evaluated during method transfer, particularly when different vendor instruments are used.

Figure 7. Peak area ratios of impurities related to the API in dependence of detector bandwidth



## CONCLUSIONS

- Several strategies for GDV difference compensation during method transfer were evaluated and resulted in straightforward retention time matches of sending and receiving LC system.
- Column thermostating is a critical parameter in method transfer and solvent mismatches of sample and mobile phase should be avoided.
- The bandwidth settings of diode array detectors impact the relative peak ratios of compounds with different UV spectra and should be carefully evaluated during method transfer.

## REFERENCES (if necessary)

- Paul, C.; Grübner, M. et al.: Thermo Scientific White Paper 72711: An instrument parameter guide for successful (U)HPLC method transfer, 2018.
- Grübner, M., Paul, C., Steiner F.: Thermo Scientific Application Note 72717: Method transfer of a USP derived acetaminophen assay from an Agilent 1260 Infinity system to an UltiMate 3000 SD system and a Vanquish Flex UHPLC system, 2018.
- The United States Pharmacopeial Convention, United States Pharmacopeia USP41-NF36 1S, Acetaminophen monograph, 2018.
- European Directorate for the Quality of Medicines & HealthCare; European Pharmacopoeia (Ph. Eur.); 7, Allée Kastner CS 30026, F-67081 Strasbourg (France)
- European Directorate for the Quality of Medicines & HealthCare; European Pharmacopoeia (Ph. Eur.) Online, 9th edition 2018 (9.5), monograph 0845: Mebendazole

## TRADEMARKS/LICENSING

© 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

PO72936-EN 02192

**ThermoFisher**  
SCIENTIFIC