

# High-sensitivity, high-throughput pesticide analysis by micro-flow LC-MS

Runsheng Zheng<sup>1</sup>, Christopher Pynn<sup>1</sup>, Alec Valenta<sup>2</sup>, Katherine Lovejoy<sup>1</sup>, Ece Aydin<sup>1</sup>, Kenneth Matuszak<sup>2</sup>, Wim Decrop<sup>1</sup>, Martin Samonig<sup>1</sup>

<sup>1</sup>Thermo Fisher Scientific, Germany; <sup>2</sup>Thermo Fisher Scientific, USA

## Introduction

**Purpose:** Demonstrate the benefits of using micro-flow (50  $\mu\text{L}/\text{min}$ , 1.0 mm i.d.) for pesticide analysis vs analytical flow (300  $\mu\text{L}/\text{min}$ , 2.1 mm) without impacting sample throughput.

**Methods:** A novel LC-MS configuration comprised of a Thermo Scientific™ Vanquish™ Neo UHPLC system coupled to an Thermo Scientific™ Orbitrap Exploris™ 240 mass spectrometer affording a 15-min micro-flow separation of 50 pesticide standards at 50  $\mu\text{L}/\text{min}$  using a 1.0 mm i.d. x 100 mm Thermo Scientific™ Hypersil GOLD™ HPLC column. Results were compared to an analytical flow configuration using a Thermo Scientific™ Vanquish™ Flex Binary UHPLC system at 300  $\mu\text{L}/\text{min}$  using a 2.1 mm i.d. column with identical chemistry.

**Results:** This study highlights the advantages of micro-flow LC-MS which affords a 2 to 4-fold increase in peak area across fifty pesticides compared with analytical flow LC-MS (2.1 mm i.d., 300  $\mu\text{L}/\text{min}$ ). The sensitivity gain using micro-flow results from the combined benefits of reduced column i.d. and increased ionization efficiency, yielding higher peak heights and larger peak areas without impacting in peak widths (FWHM). The micro-flow method also provided low retention time variation without sacrificing sample throughput. More, the reduced flow rate resulted in a 6-fold reduction in solvent usage. Taken together, these data suggest that microflow LC-MS provides a more sensitive, cost efficient, and environmentally friendly alternative to analytical flow for pesticide analysis.

## Materials and methods

### Sample Preparation

A pesticide standard containing 50 compounds (Cat. No. 31975, RESTEK GmbH, Germany) was diluted to 100 pg/ $\mu\text{L}$  using a solvent mixture of 12.5% acetonitrile, 0.1% formic acid (v/v). The dilution was performed in a Thermo Scientific™ SureSTART™ high recovery vial (P/N: 6PSV9-V1) which was then sealed with a talcum-free screw cap (P/N: 6PSC9STB1).

### LC-MS Method

**Table 1** presents the LC parameters for both methods. The required hardware components of the Vanquish Neo UHPLC system are listed in **Table 2**. The Vanquish Neo UHPLC system was configured in the micro-flow Direct Injection workflow using 50  $\mu\text{m}$  I.D. capillaries. The hardware components of the Vanquish Flex Binary UHPLC system are listed in **Table 3**.

The Vanquish Flex binary pump was equipped with a 35  $\mu\text{L}$  mixer set for low gradient delay volume (GDV) applications. Data were acquired on an Orbitrap Exploris 240 MS in full scan mode (MS1). More detail can be found in **Table 4**.

### Data Analysis

Data analysis was conducted using Skyline<sup>1</sup> software and R script<sup>2</sup>.

**Table 1. LC Methods and consumables**

	Analytical flow	Micro-flow
LC	Vanquish Flex UHPLC system	Vanquish Neo UHPLC system
Column	Hypersil GOLD 100 x 2.1 mm, 1.9 $\mu\text{m}$ (25002-102130)	Hypersil GOLD 100 x 1.0 mm, 1.9 $\mu\text{m}$ (25302-101030)
Mobile Phases	A) Water/MeOH (95/5, v/v, %), 5 mM ammonium formate, 0.1% FA B) Water/MeOH (5/95, v/v, %), 5 mM ammonium formate, 0.1% FA	
Washes	Rear seal wash: water/methanol (90/10, v/v, %) Needle wash: water/methanol (80/20, v/v, %)	Rear seal wash: water/isopropanol (25/75, v/v, %) with 0.1% FA Strong needle wash: 100% acetonitrile with 0.1% FA Weak needle wash: 100% water with 0.1% FA
Temperature	50 °C, Still Air Mode	
Injection Amount	1 $\mu\text{L}$ (100 pg/pesticide)	
Gradient	0.0 - 1.0 min: 2%B; 1.0 - 2.0 min: 2-50% B; 2.0 - 9.0 min: 50-99% B; 9.0 - 11.0 min: 99% B; 11.0 - 11.1 min: 2% B; 11.1 - 14.0 min: 2%B	
Flow Rate	300 $\mu\text{L}/\text{min}$	50 $\mu\text{L}/\text{min}$
Spray needle	100 $\mu\text{m}$ I.D.	50 $\mu\text{m}$ I.D.

**Table 2. Micro-flow UHPLC system**

Module	Part number
Vanquish Neo UHPLC system	VN-S10-A-01
Vanquish Display, Column Compartment N	6036.1180, VN-C10-A-01

**Table 3. Analytical Flow UHPLC system**

Module	Part number
Binary Pump F with 35 $\mu\text{L}$ mixer set	VF-P10-A and 6044.3870
Split Sampler F, Column Compartment, and System Base	VF-A10-A, VH-C10-A, VH-S01-A

**Table 4. Orbitrap Exploris 240 parameters**

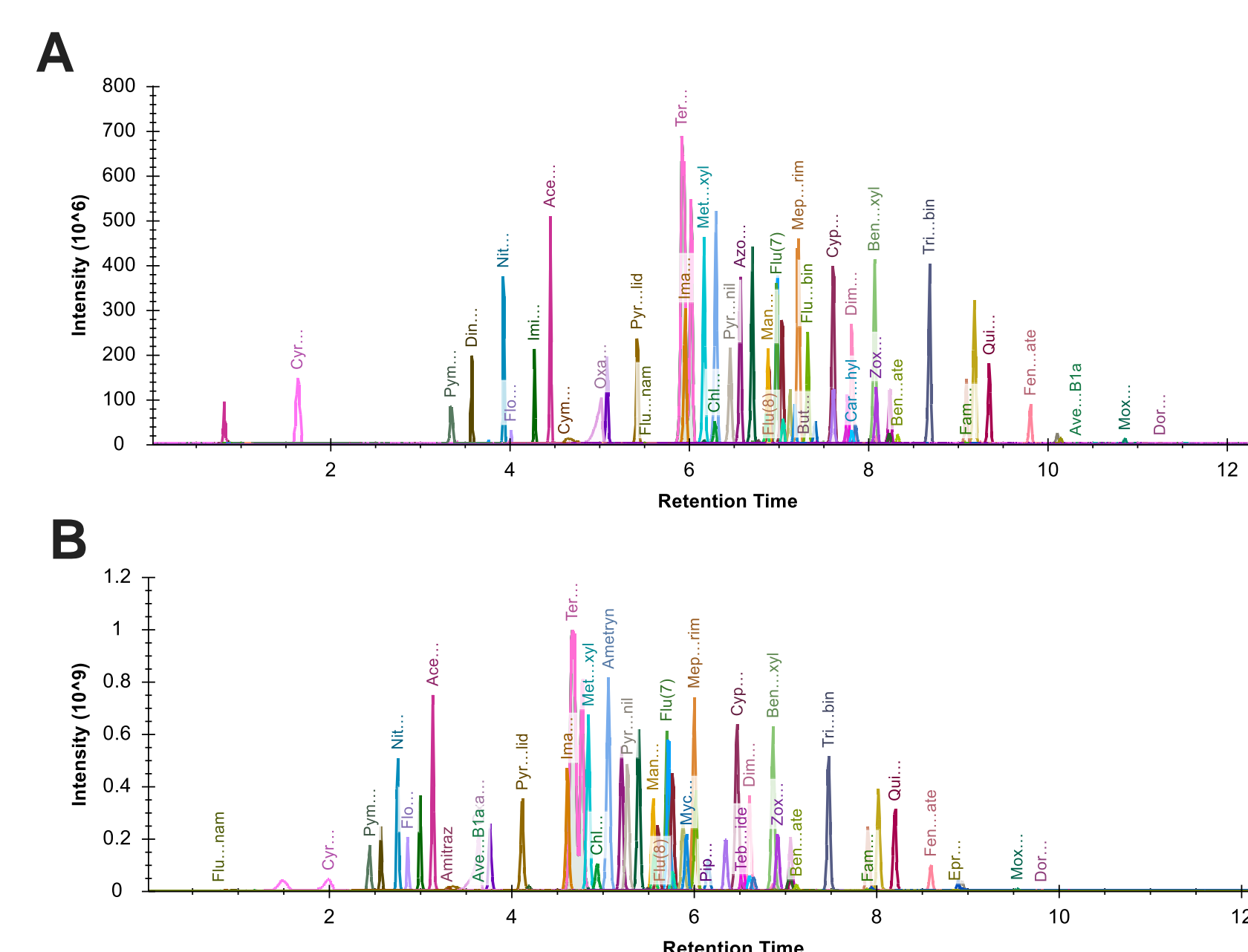
Parameter	Analytical Flow	Micro Flow
Spray Voltage (static, positive)		3300V
Sheath Gas		30
Auxiliary Gas		10
Sweep Gas		0
Ion Transfer Tube Temperature		325
Vaporizer Temperature		350
Acquisition Mode		Full Scan (MS1)
Polarity		Positive
RF Lens (%)		60
Orbitrap Resolution		60,000
Scan Range (m/z)		110-1100
AGC target		Custom (300%)
Maximum Injection Time		Auto
Data Type		Profile
Sweep cone	On	Off
Sweep gas	1	0

## Results

### Micro-flow chromatographic performance

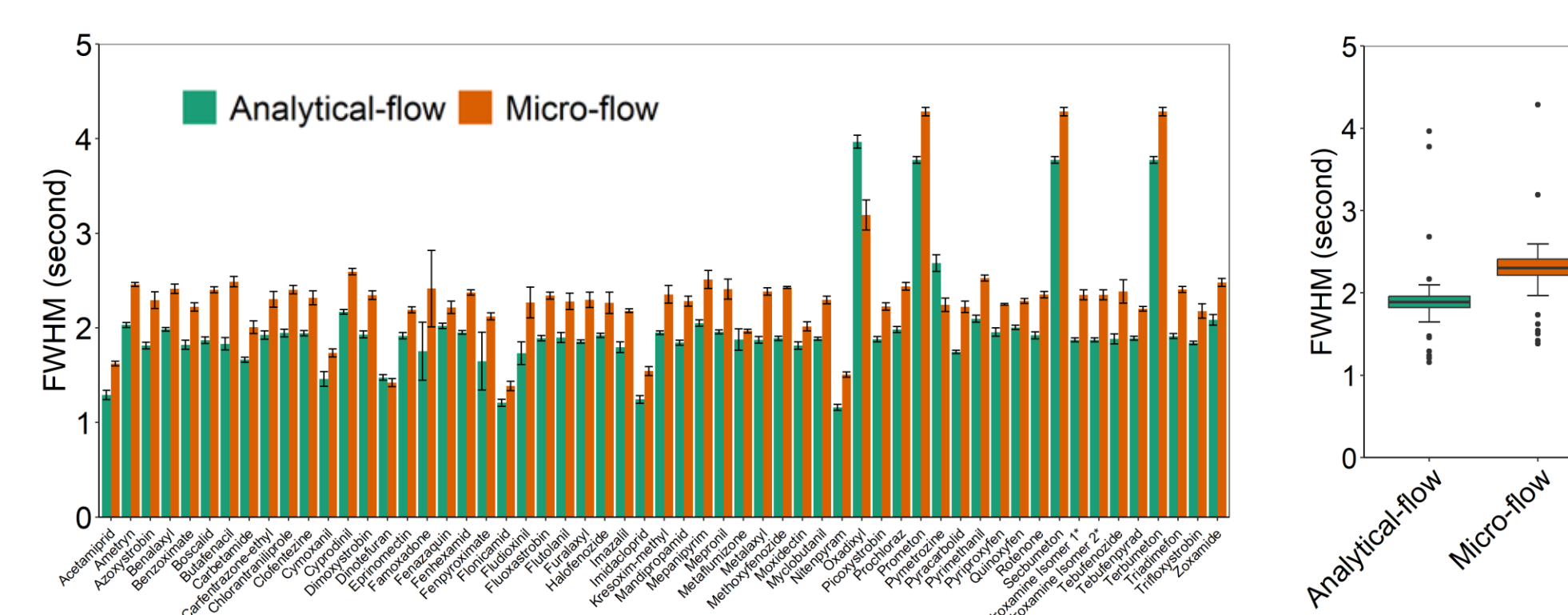
One concern of running low-flow applications is a reduction in sample throughput. Due to a low GDV (<2  $\mu\text{L}$ ) in the micro-flow configuration, the Vanquish Neo UHPLC system using a 1.0 mm x 100 mm column at 50  $\mu\text{L}/\text{min}$  delivers a comparable chromatogram to analytical flow with a 2.1 mm x 10 cm column at 300  $\mu\text{L}/\text{min}$  (**Figure 1**).

**Figure 1. Extracted ion chromatograms of analytical (A) and micro-flow (B) methods.**



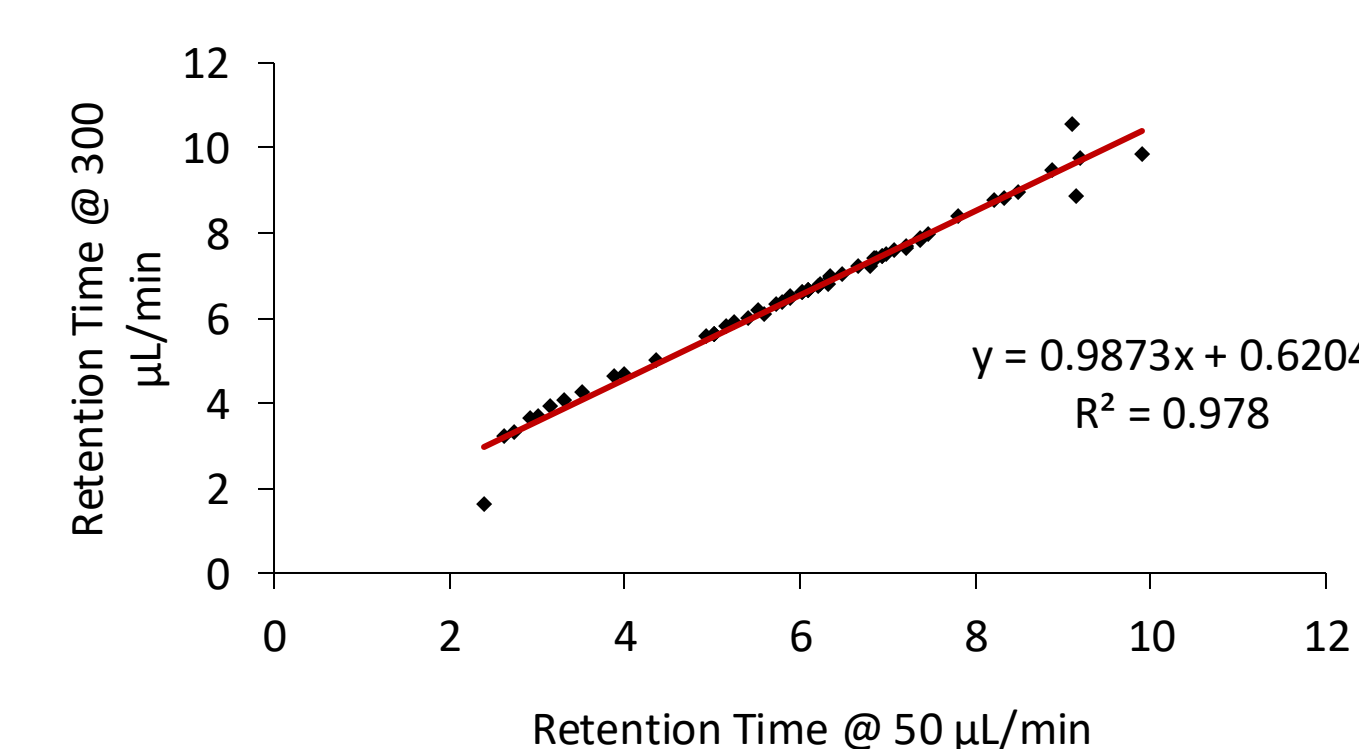
In addition, micro-flow yields similar peak FWHM, illustrating no significant loss in separation efficiency downscaling (**Figure 2**).

**Figure 2. Pesticide peak width evaluation**



Benefitting from the proportional downscaling of the chromatography conditions (flow rate and column i.d.), the micro-flow separation maintains the elution order of most of the analytes, simplifying method transfer (**Figure 3**).

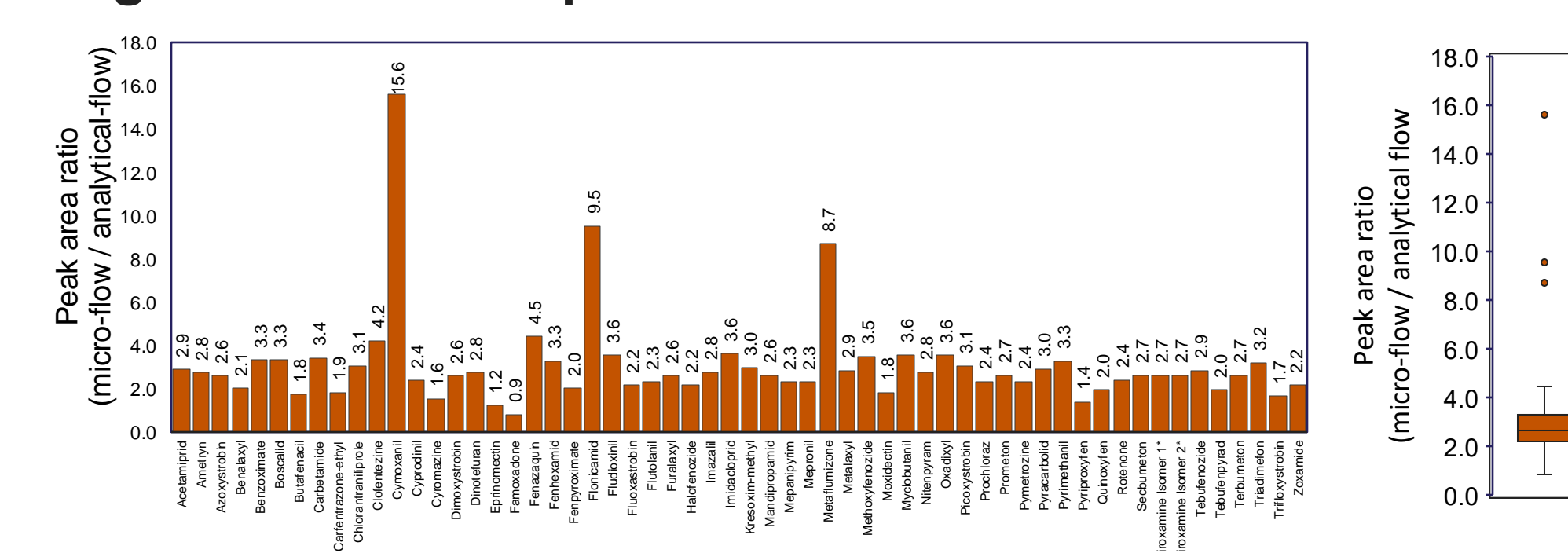
**Figure 3. Retention time comparison**



### Boosted LC-MS micro-flow sensitivity

Using optimized LC-MS conditions micro-flow increases peak intensities and areas up to 16-fold with a median increase of 2.5-fold (**Figure 4**).

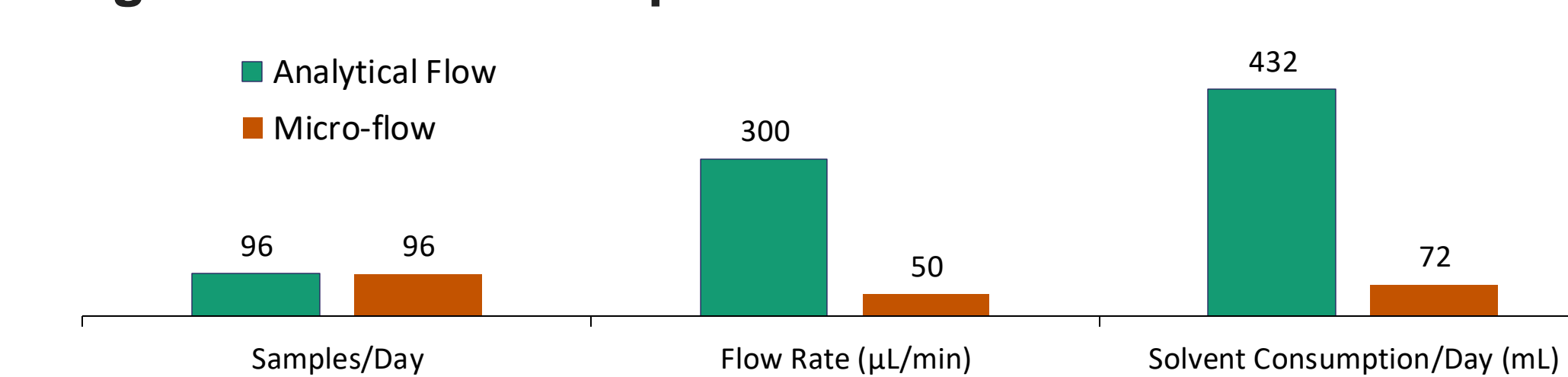
**Figure 4. Pesticide peak area evaluation**



### Economic and environmental impact of downscaling

For a 15-min gradient, micro-flow reduced mobile phase consumption by 6-fold, from 4.5 mL to 0.75 mL per run. Solvent volume savings were ~360 mL/day or 10.8 L/month per instrument, yielding a decrease in cost and environmental footprint (**Figure 5**).

**Figure 5. Method comparison**



## Conclusions

This work demonstrates the feasibility and benefits of analyzing pesticides using micro-flow LC-MS with a Vanquish Neo UHPLC system including:

- Comparable RTs and FWHMs to analytical flow
- Up to 16-fold (median at 2.5-fold) increased peak areas
- A 6-fold reduction in solvent consumption

## References

1. Birgit Schilling, etc. Mol Cell Proteomics. 2012. 11 (5), page 202-214.
2. R Core Team (2021). R: A language and environment for statistical computing. R Foundation for Statistical Computing, Vienna, Austria. URL <https://www.R-project.org/>.

## Acknowledgements

We would like to thank Maryline Carvalho for the technical support and fruitful discussion.

## Trademarks/licensing

© 2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.