

Direct quantitation of 77 toxicology and therapeutic drugs in dried blood spots using the fully automated Transcend DSX-1 System

Jingshu Guo, Courtney Patterson, Richard Gibson, Kristine Van Natta, Stephanie Samra
Thermo Fisher Scientific, San Jose, CA, USA,

Abstract

Purpose: Demonstrate a complete and fully automated workflow for dried blood spot analysis across 11 classes of drugs of abuse.

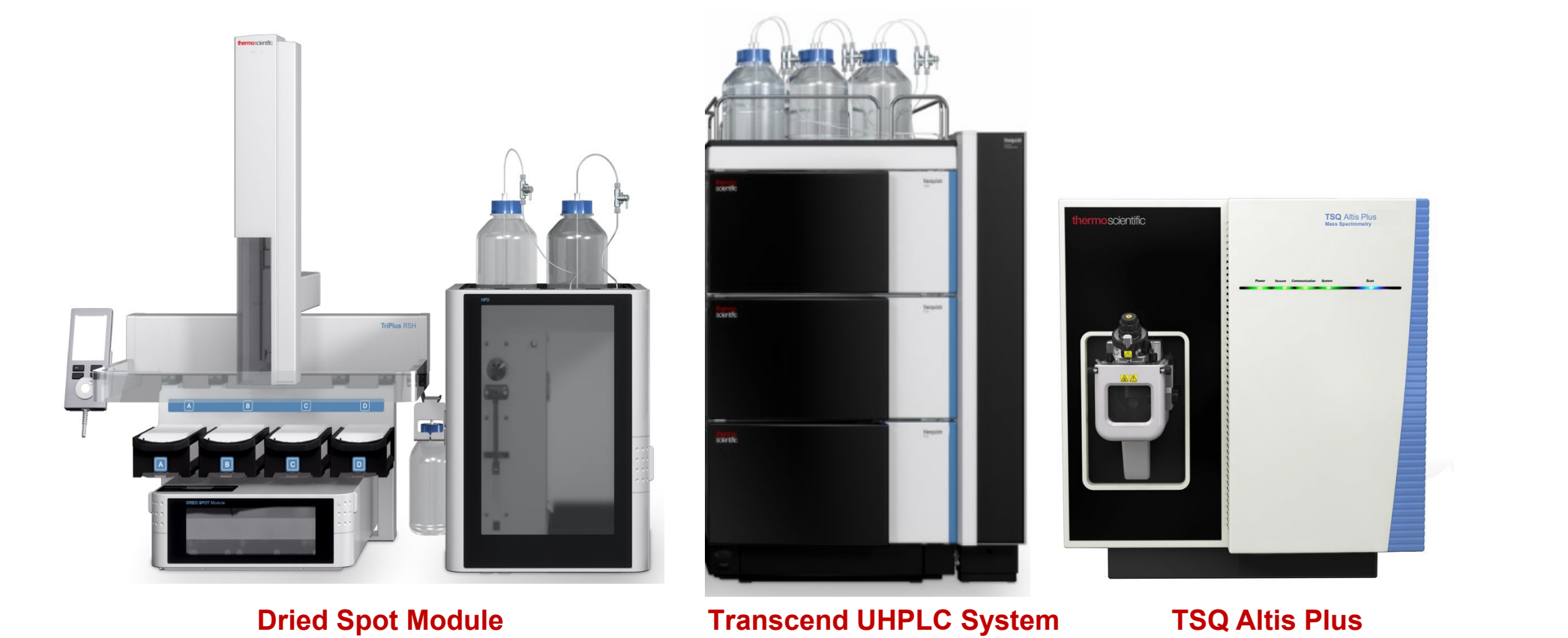
Methods: A 4.3 min analytical method was developed on the Thermo Scientific™ Transcend™ DSX-1 system consisting of a dried matrix spot module coupled with Thermo Scientific™ TurboFlow™ technology for online sample cleanup and a triple quadrupole mass spectrometer to quantify 77 analytes from 6 µL DBS in 364 SRM transitions with retention time scheduling.

Results: All analytes were successfully quantitated at or below ng/mL concentration levels, meeting the screening sensitivity needs of analytical methodologies for routine laboratories, including those monitoring forensic and sport anti-doping samples.

Introduction

The dried blood spot (DBS) sampling technique is advantageous over the traditional liquid blood collection due to its minimal invasiveness, smaller sample volume, improved analyte stability, and ease of storage, transportation and tracking, resulting in its increasing usage in fields such as sports anti-doping. Here, we describe a fully automated workflow to rapidly extract and quantify a wide range of drugs of abuse in DBS using the Transcend DSX-1 system (Figure 1) that performs internal standard addition, analyte extraction, 2-dimension LC matrix cleanup, and analyte separation without any manual intervention.

Figure 1. Fully automated Transcend DSX-1 System



Materials and methods

Sample preparation

The certified reference material of each synthetic standard and their isotope-labeled internal standards (IS) were purchased from Cerilliant® (Cerilliant Corporation, Round Rock, TX). K2-EDTA stabilized normal human whole blood was obtained from BioIVT (BioIVT, Westbury, NY) and stored at 4 °C. The analytes were spiked into the whole blood at 0, 0.1, 0.25, 0.5, 1.0, 2.5, 5.0, 10.0, 25.0, 50.0, 100, 200, 300, 400 ng/mL, and 6 µL of the mixture was spotted to the Ahlstrom-Munksjö AutoCollect™ dried blood spot (DBS) cards (Ahlstrom 226 grade paper). The DBS cards were dried at room temperature for at least 3 hours and placed directly onto the cardholder in the dried spot module (DSM).

Automated DBS extraction

The analytes were extracted from DBS cards with a 6 mm clamp via the flow-through desorption (FTD™) technology using the Loading Pump solution A (Figure 2). Internal standards (25 ng/mL in water) were introduced using the built-in IS pump in the DSM module that overfilled a 20 µL IS loop to ensure reproducible IS addition (automated IS addition, AISA™). Every sample spot was photographed with the Intelligent Vision Camera (IVC™) in “Full Spot” sample recognition mode prior to and after each run for sample tracking and traceability. The “Full Spot” mode located the actual positions of the DBS samples and accurately positioned the sample spot in the center of the extraction clamp (Figure 4). The innovative features of Transcend DSX-1 system are summarized in Figure 3.

Online sample cleanup and chromatography.

Automated online cleanup and chromatographic separation were performed on a Transcend TLX-1 system utilizing TurboFlow technology. The Transcend system was controlled by Thermo Scientific™ Aria™ MX software and configured in “Focus mode”. The analysis process and flow path are shown in Figure 4. Analytes and matrix are exacted from the DBS and loaded directly to the TurboFlow column, where the interfering molecules were washed away while analytes are retained. After washing, the analytes were eluted from the TurboFlow column using the high organic eluant stored in the “transfer loop” and refocused on the analytical column. The analyte separation was performed on the analytical column while the TurboFlow column was washed. To prepared for the subsequent analysis, the transfer loop was filled with eluant while the analytical column was washed and equilibrated. The gradient, mobile phases, clamp washes, and columns used are described in Figure 2.

Figure 2. LC conditions for the online sample cleanup and separation controlled by Aria MX software.

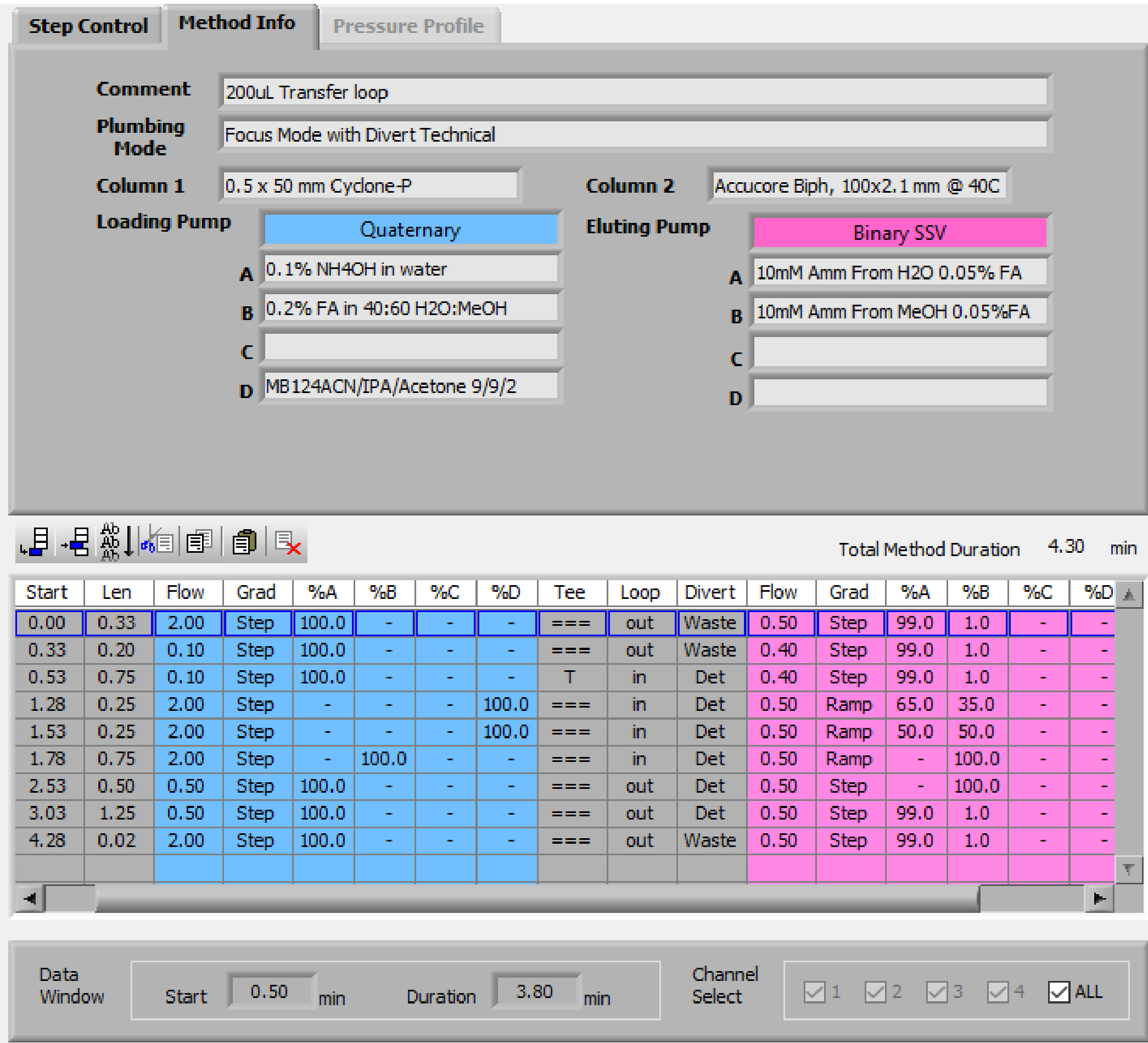
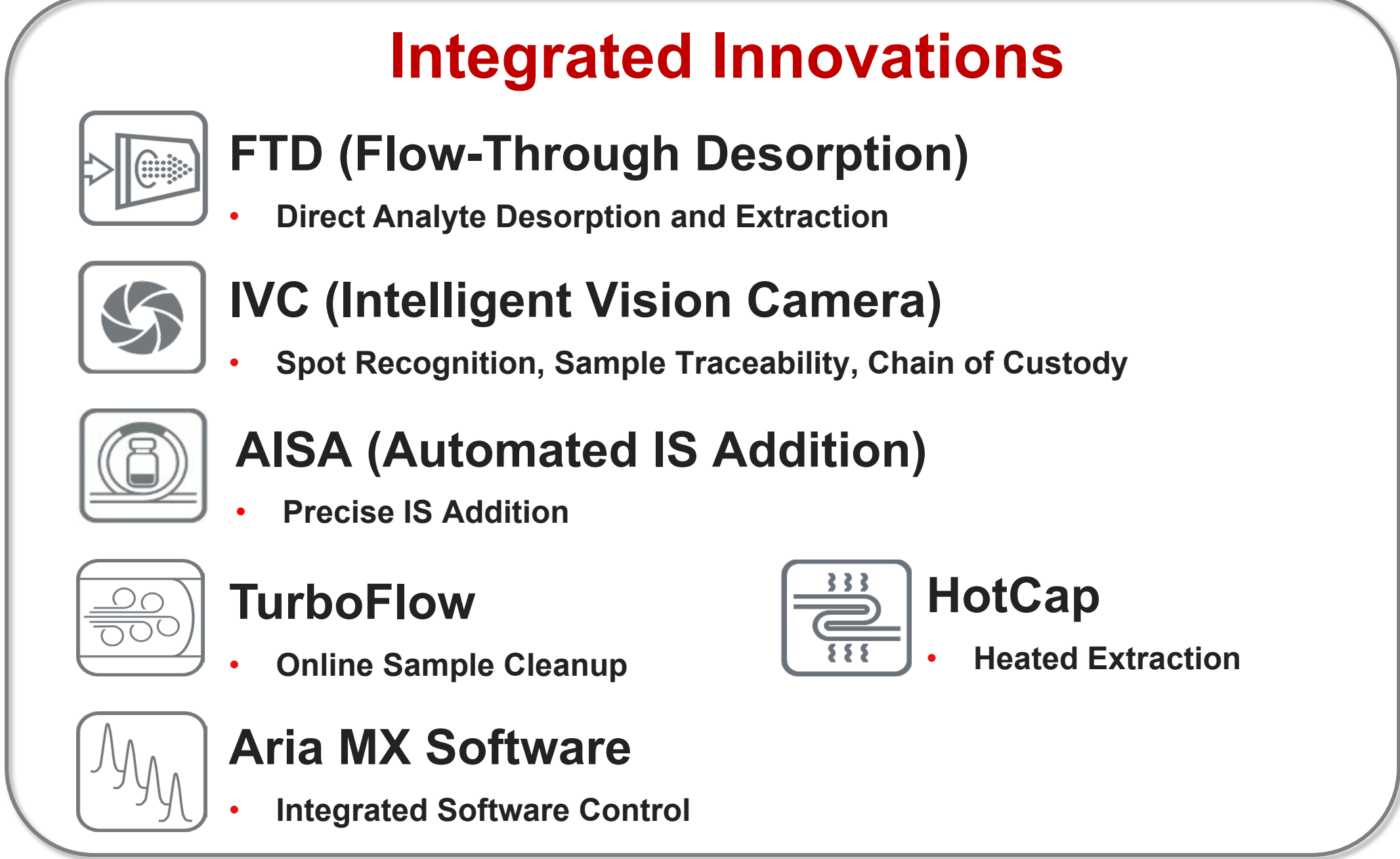


Figure 3. Summary of innovative features of the Transcend DSX-1 system.



Materials and methods (cont’)

Mass spectrometry

Analyte detection was performed using a Thermo Scientific™ TSQ Altis™ Plus mass spectrometer equipped with a heated electrospray ionization probe (HESI) and operated in the Selected Reaction Monitoring (SRM) mode. The MS parameters are shown in Table 1. The SRM transitions were imported from Thermo Scientific™ Tox Explorer™ Collection (Thermo Scientific™ TSQ Quantis Plus™ platform)² and mzCloud Mass Spectral Library (<https://www.mzcloud.org/>). Analytes and their isotope-labeled internal standards were monitored in a total of 364 SRM transitions with retention time scheduling (tR ± 0.1 min). The dwell time per transition and the number of transitions per cycle are shown in Figure 4.

Data analysis

Post-acquisition data analysis was carried out using Thermo Scientific™ TraceFinder™ software (v. 5.1).

Table 1. MS source parameters and SRM properties.

Capillary voltage	4000 (+) / 2500 (-)	Cycle Time (s)	0.5
Sheath Gas (Arb)	50	Q1 Resolution (FWHM)	0.7
Aux Gas (Arb)	15	Q3 Resolution (FWHM)	1.2
Sweep Gas (Arb)	1	Source Fragmentation	5
Ion Transfer Tube Temp. (°C)	320	Chromatographic Peak Width (sec.)	6
Vaporizer Temp. (°C)	350	CID Gas (mTorr)	1.5

Results

A total of 77 therapeutic and clinical toxicology drugs from 11 classes, including anticonvulsants, antidepressants, antihistamines, antipsychotics, benzodiazepines, cocaine, dissociative, opioids, and stimulants, are quantified in a single injection from DBS cards using a rapid automated method on a Transcend DSX-1 system. DSX-1 combines a dried spot autosampler for direct analyte extraction with Transcend TLX-1 system for online sample separation using the TurboFlow technology. The method only takes 4.3 minutes from analyte extraction to MS detection. The overlay of the extracted chromatograms of the analytes is shown in Figure 4. Analyte carryover was estimated to be below 0.5% by measuring analyte concentrations in a blank sample after the highest calibration sample.

Calibration curves were built using a weighting factor of 1/x from a lower limit of quantification of 0.1 ng/mL to an upper limit of quantification of 400 ng/mL. All calibration curves achieved R² values greater than 0.98. The LOQ values were defined with % RSD and % CV < 15, | % Diff | < 20, and relative ion ratio < %20, and are reported in Table 2. The extracted chromatograms of norfentanyl and its internal standards at the LOQ level, and its calibrations curved with 0.1 to 2.5 ng/mL zoomed-in, are shown as examples in Figure 5. The LOQ values are all in the low ng/mL levels, which largely meet the screening sensitivity needs of analytical methodologies in routine toxicology laboratories.

Figure 4. Representative combined chromatogram for all analytes monitored in this method. Dwell time of each SRM transition and the number of transitions per cycle (0.5 s) during the chromatographic separation (left inserts). For each cycle, the minimum dwell time for any transition was >2.5 ms. Images of a 6 µL dried blood spot before and after extraction (right inserts). The Intelligent Vision Camera (IVC) recognizes the actual location of the blood spot and positions the clamp to the center of the sample spot.

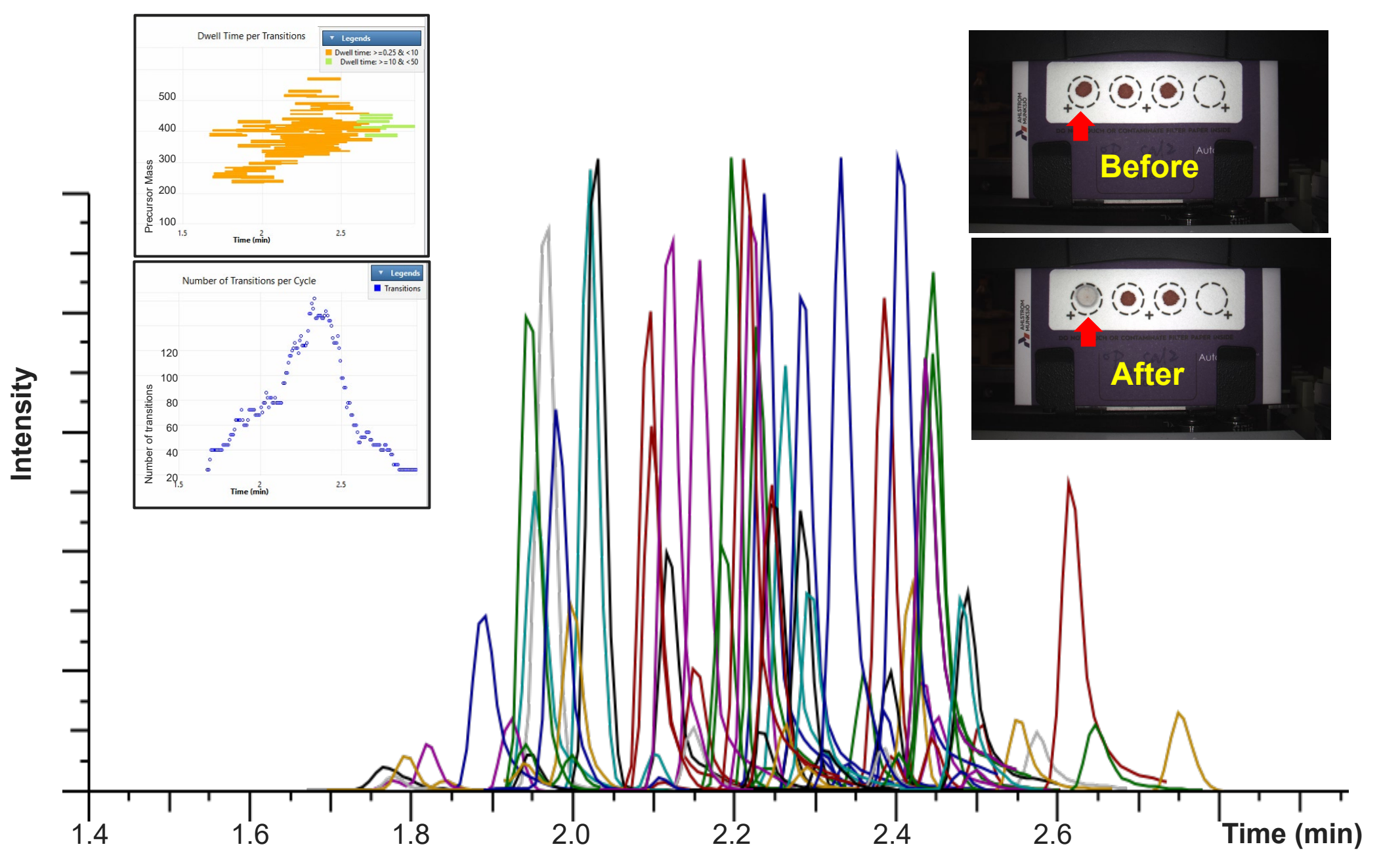


Figure 5. Representative quantification results of norfentanyl and its internal standard in DBS, demonstrating peak area, peak height, number of scans across the peak, and the calibration curve of 0.1 to 400 ng/mL with 0.1 to 2.5 ng/mL zoomed-in shown in the insert figure.

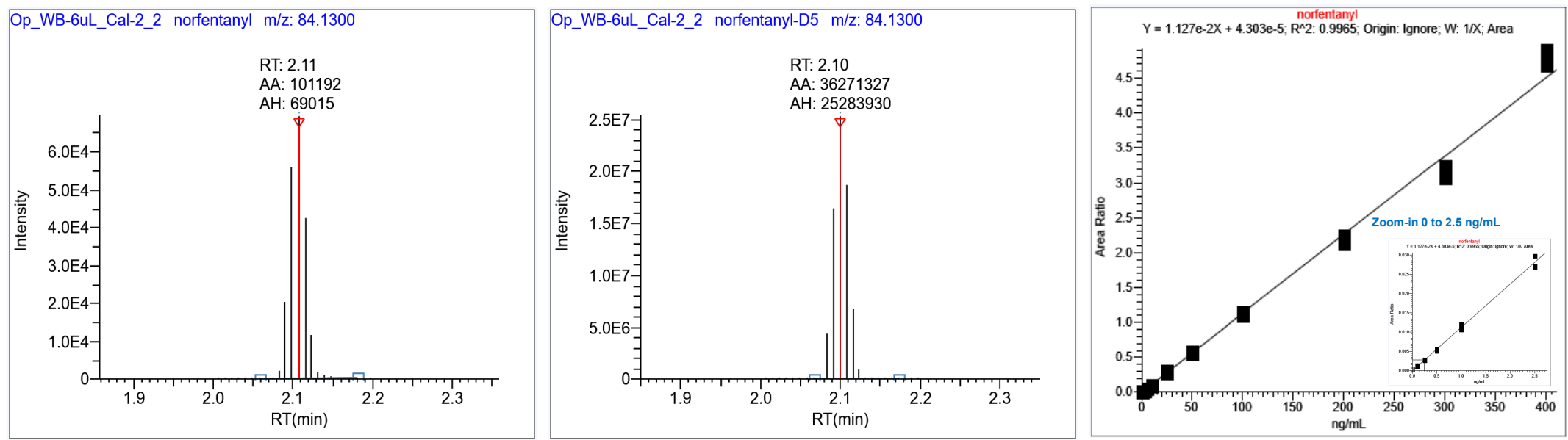


Table 2. Calibration results of the 77 analytes in 6 µL DBS. LOQ is defined as the lowest concentration with % RSD and % CV < 15, | % Diff | < 20, and relative ion ratio < %20 (N = 3).

Drug Classes	Compound	t _R (min)	IS	R ²	LOQ (ng/mL)	
Opioids	6-Acetyl mor phine	1.95	6-Acetyl mor phine- ² H ₆	0.9973	0.25	
	Buprenorphine	2.39	Buprenorphine- ² H ₄	0.9918	1.00	
	Codeine	1.94	Codeine- ² H ₆	0.9972	0.10	
	Fentanyl	2.34	Fentanyl- ² H ₅	0.9980	0.25	
	Hydrocodone	2.00	Hydrocodone- ² H ₃	0.9979	0.25	
	Hydromorphone	1.80	Hydromorphone- ² H ₃	0.9963	0.25	
	Methadone	2.49	Methadone- ² H ₃	0.9957	0.25	
	Morphine	1.77	Morphine- ² H ₆	0.9971	0.25	
	Norbuprenorphine	2.24	Norbuprenorphine- ² H ₃	0.9938	2.50	
	Norfentanyl	2.10	Norfentanyl- ² H ₄	0.9965	0.10	
	O-Desmethyltramadol	1.95	O-Desmethyltramadol- ² H ₆	0.9977	0.25	
	Oxycodone	1.98	Oxycodone- ² H ₃	0.9959	0.25	
	Oxymorphone	1.79	Oxymorphone- ² H ₃	0.9982	0.25	
	Tramadol	2.13	Tramadol- ¹³ C ₇ H ₉	0.9973	0.10	
Anticonvulsants	10-OH-carbazepine	2.25	10-OH-carbazepine- ¹³ C ₆	0.9960	1.00	
	Carbamazepine	2.45	Carbamazepine-epo- ¹³ C ₆	0.9942	0.50	
	Gabapentin	1.87	Gabapentin- ² H ₁₀	0.9913	10.00	
	Lamotrigine	2.10	Methamphetamine- ² H ₅	0.9971	0.10	
Antidepressants	Levetiracetam	1.92	Levetiracetam- ² H ₆	0.9972	0.25	
	Primidone	2.12	Primidone- ² H ₅	0.9981	0.50	
	Topiramate	2.28	Topiramate- ² H ₁₂	0.9951	5.00	
	Amitriptyline	2.45	Amitriptyline- ² H ₃	0.9950	1.00	
	Bupropion	2.18	Bupropion- ² H ₉	0.9966	0.25	
	Citalopram	2.29	Citalopram- ² H ₆	0.9960	0.50	
	Clomipramine	2.50	Clomipramine- ² H ₃	0.9935	0.10	
	Desipramine	2.41	Desipramine- ² H ₃	0.9961	1.00	
	Doxepin	2.33	Doxepin- ² H ₃	0.9963	0.50	
	Fluoxetine	2.30	Fluoxetine- ² H ₃	0.9956	0.25	
	Imipramine	2.43	Imipramine- ² H ₃	0.9962	0.10	
	Mirtazapine	2.27	Mirtazapine- ² H ₃	0.9983	0.25	
	Nortriptyline	2.43	Nortriptyline- ² H ₃	0.9957	1.00	
	O-Desmethylvenlafaxine	2.03	O-Desmethylvenlafaxine- ² H ₆	0.9971	0.25	
	Paroxetine	2.40	Paroxetine- ² H ₆	0.9960	2.50	
	Sertraline	2.48	Sertraline- ² H ₃	0.9939	10.00	
	Trazodone	2.47	Trazodone- ² H ₆	0.9980	0.10	
	Venlafaxine	2.22	Venlafaxine- ² H ₆	0.9959	0.25	
Antihistamines	Brompheniramine	2.28	Risperidone- ² H ₄	0.9957	0.10	
	Chlorophenylpiperazine	2.12	Chlorophenylpiperazine- ² H ₆	0.9983	0.50	
	Chlorpheniramine	2.24	Chlorpheniramine- ² H ₆	0.9985	0.25	
	Diphenhydramine	2.30	Diphenhydramine- ² H ₃	0.9975	0.10	
	Doxylamine	2.23	Doxylamine- ² H ₃	0.9976	0.25	
	Hydroxyzine	2.41	Hydroxyzine- ² H ₆	0.9952	0.10	
	Methorphan	2.39	Methorphan- ² H ₃	0.9938	1.00	
	Norchlorcyclizine	2.40	Paroxetine- ² H ₆	0.9912	25.00	
	Promethazine	2.40	Promethazine- ² H ₃	0.9964	1.00	
	Cocaine	Benzoylcocaine	2.16	Benzoylcocaine- ² H ₆	0.9983	2.50
	Cocacethylene	2.25	Cocacethylene- ² H ₃	0.9972	0.25	
	Cocaine	2.19	Cocaine- ² H ₃	0.9967	0.10	
Dissociatives	Ketamine	2.18	Ketamine- ² H ₄	0.9991	0.10	
	Norketamine	2.13	Norketamine- ² H ₄	0.9992	0.25	
	Phencyclidine	2.38	Phencyclidine- ² H ₅	0.9971	1.00	
Benzodiazepine	7-Aminoclonazepam	2.24	7-Aminoclonazepam- ² H ₄	0.9976	0.25	
	Alprazolam	2.68	Alprazolam- ² H ₅	0.9982	0.25	
	Clonazepam	2.52	Clozapine- ² H ₄	0.9950	5.00	
	Diazepam	2.75	Diazepam- ² H ₅	0.9979	0.25	
	Lorazepam	2.47	Lorazepam- ² H ₄	0.9987	0.50	
	Nordiazepam	2.59	Nordiazepam- ² H ₅	0.9987	0.50	
	Oxazepam	2.50	Oxazepam- ² H ₅	0.9991	0.50	
	Temazepam	2.64	Temazepam- ² H ₅	0.9990	0.25	
	Zolpidem	2.43	Zolpidem- ² H ₆	0.9994	0.10	
	α-Hydroxyalprazolam	2.57	α-Hydroxyalprazolam- ² H ₅	0.9988	0.50	
	Antipsychotics	9-Hydroxyrisperidone	2.27	9-Hydroxyrisperidone- ² H ₄	0.9981	0.25
	Chlorpromazine	2.50	Chlorpromazine- ² H ₃	0.9982	0.50	
	Clozapine	2.32	Clozapine- ² H ₄	0.9965	0.50	
	Olanzapine	2.12	Benzoylcocaine- ² H ₆	0.9857	2.50	
	Quetiapine	2.45	Quetiapine- ² H ₆	0.9974	0.25	
	Risperidone	2.38	Risperidone- ² H ₄	0.9962	0.10	
Stimulants	Amphetamine	1.91	Amphetamine- ² H ₅	0.9978	1.00	
	Methamphetamine	1.97	Methamphetamine- ² H ₅	0.9974	0.10	
	MDA	1.98	MDA- ² H ₅	0.9972	0.50	
	MDMA	2.02	MDMA- ² H ₅	0.9968	0.25	
Miscellaneous	Carisoprodol	2.34	Carisoprodol- ² H ₇	0.9954	1.00	
	Cyclobenzaprine	2.43	Cyclobenzaprine- ² H ₃	0.9970	2.50	
	Meprobamate	2.16	Meprobamate- ² H ₃	0.9985	0.50	
	Zolpidem carboxylic acid	2.23	Zolpidem carboxylic acid- ² H ₄	0.9962	0.50	
	Zopiclone	2.31	Zopiclone- ² H ₄	0.9974	0.50	
OTC	Acetaminophen	1.79	Acetaminophen- ² H ₄	0.9990	5.00	

Conclusions

A comprehensive, 4.3 min LC-MS-based method was set up to extract and reliably quantify 77 drugs of abuse across 11 classes from DBS using a fully automated and integrated Transcend DSX-1 system complete with sample tracking and IS addition.

References

- Wu, I., et. al. Thermo Fisher Scientific, 2021. TN000625
- Van Natta, K. et. al. Thermo Fisher Scientific, 2022. TN001256

Acknowledgements

We thank Edwin Gortz and Peter Ringeling from Spark Holland™ (The Netherlands) for the technical consultant of the dried matrix spot module.

Trademarks/licensing

© 2023 Thermo Fisher Scientific Inc. All rights reserved. Cerilliant is a registered trademark of Cerilliant Corporation, BioIVT of BioIVT, LLC, and AutoCollect for Ahlstrom. FTD, IVC, and AISA are trademarks of Spark Holland. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage the use of these products in any manner that might infringe on the intellectual property rights of others.