

# Quantitation of drugs of abuse and their metabolites in urine using PaperSpray tandem mass spectrometry for clinical research and forensic toxicology

Authors: Scott A. Borden<sup>1,2</sup>, Armin Saatchi<sup>1</sup>,  
Chris G. Gill<sup>1,2</sup>, Neloni R. Wijeratne<sup>3</sup>

<sup>1</sup>Applied Environmental Research Laboratories  
Vancouver Island University, Nanaimo, BC

<sup>2</sup>University of Victoria, Victoria, BC

<sup>3</sup>Thermo Fisher Scientific, San Jose, CA

Keywords: Illicit drugs, drugs of abuse, opiates, opioids, amphetamines, benzodiazepines, urine, PS-MS/MS, TSQ Fortis MS, VeriSpray PaperSpray ion source, TraceFinder, forensic toxicology, clinical research



TSQ Fortis™ mass spectrometer with the Thermo Scientific™ VeriSpray™ PaperSpray ion source.

## Application benefits

- Development of a robust, reliable quantitative assay for 41 drugs of abuse and their metabolites in urine in a single quantitative method
- No sample preparation, reduced cost per sample, and increased sample throughput

## Goal

To develop a reliable quantitative MS-based urinalysis procedure for drugs of abuse and their metabolites for clinical research or forensic toxicology able to meet established cutoff values using the Thermo Scientific™

## Introduction

Forensic and clinical laboratories routinely quantify drugs of abuse in biological matrices, urine being one of the most common due to ease of procurement. Most often, drugs are quantified using time-consuming chromatographic methods that require significant sample preparation. Therefore, an alternative method that provides comparable high-throughput results is of significant value to forensic and clinical laboratories. PS-MS is a viable alternative to chromatography for the rapid analysis of drugs of abuse in urine, demonstrating sufficient sensitivity, precision, and accuracy to be of clinical and forensic relevance.

Paper spray mass spectrometry (PS-MS) was first described in 2010<sup>1</sup> as a method for the direct analysis of complex samples without sample preparation. PS-MS uses small sampling volumes (<10  $\mu$ L) directly deposited onto a strip of paper.<sup>1</sup> Directly in front of the inlet of the mass spectrometer, solvent is applied to the paper to extract analytes and high voltage is applied to ionize them. A chromatogram of the ion current is collected for a short duration (usually <1 min) and is integrated to quantitate the analyte. Paper spray ionization coupled with triple-quadrupole mass spectrometry allows for rapid and sensitive (low ng/mL) quantitative measurements in a variety of complex samples, such as blood and urine. The VeriSpray PaperSpray system includes a VeriSpray plate loader and magazine that holds up to 10 VeriSpray cartridges, each containing 24 paper strips. This fully automated system allows for high-throughput analysis without sample carryover. Figure 1 shows the VeriSpray system mounted to the TSQ Fortis MS, as well as the removable magazine and VeriSpray cartridges.

In this study, the VeriSpray PaperSpray ion source coupled to a TSQ Fortis triple-quadrupole mass spectrometer was evaluated as a tool for the quantification of drugs of abuse and their metabolites in human urine samples for

applications in clinical research and forensic toxicology. A suite of 41 different drugs and metabolites chosen to represent major drugs classes (e.g., opiates and opioids, amphetamines, fentanyl and its analogs, and benzodiazepines) was quantified in pooled human urine samples, with method validation carried out using matrix-matched quality control samples.

## Experimental

### Sample preparation

A selection of 41 drugs and metabolites from various drug classes (including opioids, amphetamines, benzodiazepines, and novel psychoactive substances) were used to create calibration curve standards in pooled blank human urine samples from five sources. Nine-point calibration curves (n = 6 replicates) were prepared accordingly to meet established cutoff values and encompass the range of concentrations expected in urine samples. Working solutions of drugs and isotopically labeled internal standards were prepared in methanol and spiked into the pooled urine matrix. For analysis, 8  $\mu$ L aliquots of spiked urine were spotted on VeriSpray cartridges and allowed to dry at ambient temperature for 30 minutes.



Figure 1. (A) VeriSpray PaperSpray system mounted to TSQ Fortis triple quadrupole MS, (B) magazine, and (C) VeriSpray cartridge

## PaperSpray and MS conditions

The TSQ Fortis triple quadrupole mass spectrometer was used for all analyses and was coupled with the VeriSpray PaperSpray system for ionization. PaperSpray solvents used (sample rewet and spray solvents) were acetonitrile/water 90/10 with 0.1% acetic acid and applied according to the settings in Table 1. Mass spectrometry parameters were optimized using a spray voltage of +3.8 kV, a cycle time of 1.2 s, and Q1 and Q3 resolution at 0.7 Da FWHM. Source parameters are outlined in Table 2. The optimum RF lens settings and collision energies for the product ions were determined by direct infusion of standards into the mass spectrometer using the heated electrospray ionization source. Optimized SRM transitions for all target analytes and internal standards are listed in Table 3. The paper tip to MS inlet distance was set to 5.0 mm.

**Table 1. VeriSpray solvent application parameters.**

Each rewetting and solvent dispense is 10 µL.

Rewetting dispense	Delay (s)
1	1
Solvent dispense	Delay (s)
1	1
2	1
3	1
4	1
5	3
6	3
7	5
8	5
9	5
10	5
11	7
12	7
13	7
14	7
15	7

**Table 2A. TSQ Fortis parameters**

Ion source parameter	Value
Spray voltage	Time dependent
Positive ion	3800 V
Sweep gas	0 Arb
Ion transfer tube temperature	300 °C
Q1 resolution (FWHM)	0.7
Q3 resolution (FWHM)	0.7
CID gas	2 mTorr

**Table 2B. Time-dependent spray voltage settings**

Time (min)	Voltage (V)
0	0
0.1	3800
1.1	0

## Method validation

Method validation was done according to the Scientific Working Group for Forensic Toxicology's (SWGTOX) standard practices for method validation in forensic toxicology.<sup>2</sup> Low, medium, and high QC samples were prepared such that the low QC sample was near the cutoff value, the high QC sample was at 80% of the highest calibrator, and the medium QC was at the midpoint of the high and low QC samples. The QC samples were analyzed daily (n = 5 replicates), over a period of five days to determine within-run and between-run precision (%CV) and accuracy (%Bias) values:

$$\text{Within-run \%CV} = \frac{\text{Standard deviation of a single day of samples}}{\text{Mean calculated value of a single day of samples}} \times 100 \quad (1)$$

$$\text{Between-run \%CV} = \frac{\text{Standard deviation of grand mean}}{\text{Grand mean}} \times 100 \quad (2)$$

$$\text{\%Bias} = \frac{\text{Grand mean of calculated concentration} - \text{nominal concentration}}{\text{Nominal concentration}} \times 100 \quad (3)$$

## Data acquisition and analysis

Thermo Scientific™ TraceFinder™ software version 4.1 SP5 optimized for clinical research was used for automated data acquisition and processing. All linear regressions used a 1/x weighting. Lower limits of quantitation (LLOQs) were determined from the lowest calibration standard that met the following criteria:  $R^2 \geq 0.98$ ,  $\%CV \leq 15$ , and  $\%Bias \leq \pm 20$ , and a S/N value > 4. Noise was defined as the average of 10 replicate measurements of pooled human urine.

**Table 3. Optimized SRM transitions using positive polarity**

Compound	Precursor (m/z)	Product (m/z)	Collision energy (V)	RF lens (V)
N-phenyl-1-(2-phenylethyl)-4-piperidinamine (4-ANPP)	281.2	188.1	17	146
	281.2	105.1	31	146
4-ANPP-D <sub>5</sub>	286.2	188.3	17	96
	286.2	105.0	32	96
Acetylfentanyl	323.2	188.2	23	97
	323.2	105.1	36	97
Acetylfentanyl- <sup>13</sup> C <sub>6</sub>	329.3	188.1	23	100
	329.3	105.1	36	100
Acetylmorphine	328.2	211.1	26	114
	328.2	165.1	38	114
Acetylmorphine-D <sub>6</sub>	334.2	211.1	27	128
	334.2	165.1	38	128
Acrylfentanyl	335.2	188.2	22	104
	335.2	105.1	36	104
Acrylfentanyl-D <sub>5</sub>	340.3	188.1	22	102
	340.3	105.1	37	102
Alprazolam	309.1	281.1	25	111
	309.1	205.1	41	111
Alprazolam-D <sub>5</sub>	314.2	286.1	26	111
	314.2	210.1	42	111
Aminoclonazepam	286.1	222.1	25	113
	286.1	250.1	20	113
Aminoflunitrazepam	284.2	135.2	27	102
	284.2	227.1	25	102
Amphetamine	136.1	91.2	17	73
	136.1	65.1	36	73
Amphetamine-D <sub>11</sub>	147.2	98.1	19	79
	147.2	70.1	38	79
Benzoylecgonine	290.1	168.2	19	97
	290.1	77.1	47	97
Benzoylecgonine-D <sub>8</sub>	298.2	171.2	20	93
	298.2	82.2	48	93
Butyrylfentanyl	351.2	188.2	23	104
	351.2	105.1	38	104
Carfentanil	395.2	335.2	18	100
	395.2	363.2	13	100
Carfentanil-D <sub>5</sub>	400.3	340.2	18	110
	400.3	368.3	14	110
Codeine	300.2	215.1	25	108
	300.2	165.1	41	108
Codeine-D <sub>3</sub>	303.2	215.1	26	123
	303.2	165.1	41	123
Cyclopropylfentanyl	349.2	188.2	23	98
	349.2	105.1	37	98
Desalkylflurazepam	289.1	140.1	30	119
	289.1	226.1	28	119
Desalkylflurazepam-D <sub>4</sub>	293.1	140.0	30	123
	293.1	230.2	29	123
Diazepam	285.1	193.0	32	117
	285.1	154.1	28	117
Ecgonine methyl ester	200.1	82.1	26	80
	200.1	119.0	25	80
Ecgonine methyl ester-D <sub>3</sub>	203.1	85.2	26	80
	203.1	119.0	25	80
2-Ethylidene-1,5-dimethyl-3,3-diphenylpyrrolidine (EDDP)	278.2	234.2	31	95
	278.2	249.2	24	95
EDDP-D <sub>3</sub>	281.2	234.1	31	98
	281.2	249.2	25	98
Ethylpentylone	250.2	202.1	18	81
	250.2	189.1	24	81
Ethylpentylone-D <sub>5</sub>	255.2	207.1	19	82
	255.2	194.1	24	82
Fentanyl	337.2	188.2	23	101
	337.2	105.1	37	101
Fentanyl-D <sub>5</sub>	342.3	188.2	23	100
	342.3	105.1	37	100
4-Fluoroisobutyrylfentanyl (FIBF)	369.2	188.1	24	108
	369.2	105.1	39	108
Fluorofentanyl	355.2	188.1	23	107
	355.2	105.1	37	107
Fluorofentanyl-D <sub>3</sub>	358.2	188.1	24	108
	358.2	105.1	38	108
Furanylfentanyl	375.2	188.2	22	106
	375.2	105.1	38	106
Furanylfentanyl-D <sub>5</sub>	380.2	188.1	23	102
	380.2	105.1	38	102
HO-Alprazolam	325.1	216.1	39	117
	325.1	205.1	45	117
HO-Alprazolam-D <sub>5</sub>	330.2	210.1	45	127
	330.2	221.1	40	127
Hydrocodone	300.2	171.1	39	114
	300.2	241.1	27	114
Hydrocodone-D <sub>3</sub>	303.2	171.1	39	117
	303.2	241.1	27	117

Table 3 (continued). Optimized SRM transitions using positive polarity

Compound	Precursor (m/z)	Product (m/z)	Collision energy (V)	RF lens (V)
Hydromorphone	286.1	185.1	30	116
	286.1	157.1	41	116
Hydromorphone-D <sub>3</sub>	289.2	185.1	30	134
	289.2	157.1	42	134
Isobutyrylfentanyl	351.2	188.2	24	101
	351.2	105.1	38	101
Ketamine	238.1	125.0	27	80
	238.1	207.0	14	80
3,4-Methylenedioxy-amphetamine (MDA)	180.1	105.1	22	83
	180.1	135.1	19	83
MDA-D <sub>5</sub>	180.1	133.1	18	83
	185.1	110.1	23	102
3,4-Methylenedioxy-methamphetamine (MDMA)	185.1	138.1	19	102
	185.1	137.1	20	102
MDMA-D <sub>5</sub>	194.1	163.1	13	75
	194.1	135.1	21	75
Meperidine	248.2	220.1	21	100
	248.2	174.2	21	100
Meperidine-D <sub>4</sub>	252.2	224.2	22	101
	252.2	178.2	21	101
Methadone	310.2	265.2	15	72
	310.2	105.0	27	72
Methadone-D <sub>9</sub>	319.3	268.2	16	77
	319.3	105.1	28	77
Methamphetamine	150.1	91.1	19	78
	150.1	119.1	11	78
Methamphetamine-D <sub>11</sub>	161.2	97.1	19	80
	161.2	127.2	12	80
Morphine	286.1	152.1	55	109
	286.1	201.1	26	109
Morphine-D <sub>6</sub>	292.2	153.1	43	134
	292.2	201.1	26	134

Compound	Precursor (m/z)	Product (m/z)	Collision energy (V)	RF lens (V)
Methamphetamine-D <sub>11</sub>	161.2	97.1	19	80
	161.2	127.2	12	80
Morphine	286.1	152.1	55	109
	286.1	201.1	26	109
Morphine-D <sub>6</sub>	292.2	153.1	43	134
	292.2	201.1	26	134
Normeperidine	234.2	160.1	16	99
	234.2	56.1	23	99
Normeperidine-D <sub>4</sub>	238.2	164.2	16	93
	238.2	58.1	23	93
Ocfentanil	371.2	188.2	22	101
	371.2	105.1	38	101
Phencyclidine (PCP)	244.3	86.2	13	68
	244.3	91.1	31	68
Phencyclidine-D <sub>5</sub>	249.2	86.2	13	70
	249.2	96.1	32	70
Sufentanil	387.2	238.2	19	92
	387.2	111.0	36	92
Sufentanil-D <sub>5</sub>	392.2	238.2	20	97
	392.2	111.0	36	97
Temazepam	301.1	255.1	22	103
	301.1	177.0	38	103
Temazepam-D <sub>5</sub>	306.1	260.2	24	112
	306.1	177.0	38	112
U-47700	329.2	284.1	17	87
	329.2	173.0	32	87
Valeryl fentanyl	365.3	188.2	24	102
	365.3	105.1	39	102
Zolpidem	308.2	235.2	35	108
	308.2	263.2	26	108
Zolpidem-D <sub>7</sub>	315.2	242.2	36	104
	315.2	270.2	27	104

## Results and discussion

Calibrations with good linearity ( $R^2 > 0.98$ ) were achieved for all 41 drugs and metabolites and used for the quantitation of quality control samples prepared in pooled blank urine. The calibration models, LLOQs, and cutoff values are outlined in Table 4. Method validation was carried out according to SWGTOX guidelines; acceptable results are indicated when precision (%CV) values do not exceed 20% and accuracy (%Bias) values do not exceed  $\pm 20\%$ .

Typical results from representative compounds of the major drug classes are outlined in Table 5. Determined LLOQ values are below or near established cutoff values for forensic or clinical toxicology. Total analysis time for dried urine spots, including extraction (ca. 1.1 min) followed by mass spectrometric analysis (1.2 min), was approximately 2.3 minutes.

**Table 4. Compound calibration models (9 levels, 6 replicates),  $R^2$  values, lower limits of quantitation (LLOQ) and cutoff values**

Compound	Internal standard	[Internal standard] (ng/mL)	Calibration range (ng/mL)	$R^2$	LLOQ (ng/mL)	Cutoff (ng/mL)
4-ANPP	4-ANPP-D <sub>5</sub>	25	0.5 – 600	0.9956	1	2
Acetylfentanyl	Acetylfentanyl- <sup>13</sup> C <sub>6</sub>	25	0.5 – 600	0.9966	1	2
Acetylmorphine	Acetylmorphine-D <sub>6</sub>	100	7.5 – 900	0.9880	7.5	10
Acrylfentanyl	Acrylfentanyl-D <sub>5</sub>	25	0.5 – 600	0.9946	1	2
Alprazolam	Alprazolam-D <sub>5</sub>	100	7.5 – 1800	0.9968	3	10
Aminoclonazepam	Alprazolam-D <sub>5</sub>	100	7.5 – 1800	0.9803	30	10
Aminoflunitrazepam	Alprazolam-D <sub>5</sub>	100	7.5 – 1800	0.9809	75	5
Amphetamine	Amphetamine-D <sub>11</sub>	500	5 – 6000	0.9963	25	50
Benzoylcegonine	Benzoylcegonine-D <sub>8</sub>	500	5 – 6000	0.9964	10	50
Butyrylfentanyl	Acrylfentanyl-D <sub>5</sub>	25	0.5 – 600	0.9967	1	2
Carfentanil	Carfentanil-D <sub>5</sub>	25	0.5 – 600	0.9970	1	2
Codeine	Codeine-D <sub>3</sub>	500	5 – 6000	0.9937	250	20
Cyclopropylfentanyl	Furanylfentanyl-D <sub>5</sub>	25	0.5 – 600	0.9913	1	2
Desalkylflurazepam	Desalkylflurazepam-D <sub>4</sub>	100	7.5 – 1800	0.9906	7.5	25
Diazepam	Alprazolam-D <sub>5</sub>	100	1.5 – 1800	0.9904	7.5	20
Ecgonine methyl ester	Ecgonine methyl ester-D <sub>3</sub>	500	5 – 6000	0.9953	25	50
EDDP	EDDP-D <sub>3</sub>	100	1.5 – 1800	0.9976	1.5	10
Ethylpentylone	Ethylpentylone-D <sub>5</sub>	100	1.5 – 1800	0.9974	3	30
Fentanyl	Fentanyl-D <sub>5</sub>	25	0.5 – 600	0.9956	1	2
FIBF	FIBF-D <sub>3</sub>	25	0.5 – 600	0.9936	0.5	2
Fluorofentanyl	Fluorofentanyl-D <sub>3</sub>	25	0.5 – 600	0.9983	0.5	2
Furanylfentanyl	Furanylfentanyl-D <sub>5</sub>	25	0.5 – 600	0.9966	1	2
HO-Alprazolam	HO-Alprazolam-D <sub>5</sub>	100	1.5 – 1800	0.9921	75	10
Hydrocodone	Hydrocodone-D <sub>3</sub>	500	5 – 6000	0.9975	10	20
Hydromorphone	Hydromorphone-D <sub>3</sub>	500	5 – 6000	0.9955	10	20
Isobutyrylfentanyl	Acrylfentanyl-D <sub>5</sub>	25	0.5 – 600	0.9967	1	2
Ketamine	Ethylpentylone-D <sub>5</sub>	100	1.5 – 1800	0.9967	1.5	30
MDA	MDA-D <sub>5</sub>	500	5 – 6000	0.9971	25	100
MDMA	MDMA-D <sub>5</sub>	500	5 – 6000	0.9971	5	100
Meperidine	Meperidine-D <sub>4</sub>	100	1.5 – 1800	0.9973	1.5	10
Methadone	Methadone-D <sub>9</sub>	100	1.5 – 1800	0.9970	1.5	10
Methamphetamine	Methamphetamine-D <sub>11</sub>	500	5 – 6000	0.9979	10	50
Morphine	Morphine-D <sub>6</sub>	500	25 – 3000	0.9800	250	20
Normeperidine	Normeperidine-D <sub>4</sub>	100	1.5 – 1800	0.9930	1.5	10
Ocfentanil	Fluorofentanyl-D <sub>3</sub>	25	0.5 – 600	0.9963	1	2
Phenylcyclidine	Phenylcyclidine-D <sub>5</sub>	100	1.5 – 1800	0.9947	1.5	10
Sufentanil	Sufentanil-D <sub>5</sub>	25	0.5 – 600	0.9952	1	2
Temazepam	Temazepam-D <sub>5</sub>	100	1.5 – 1800	0.9907	7.5	20
U-47700	Fluorofentanyl-D <sub>3</sub>	25	0.5 – 600	0.9932	2.5	2
Valerylfentanyl	Fluorofentanyl-D <sub>3</sub>	25	0.5 – 600	0.9847	10	2
Zolpidem	Zolpidem-D <sub>7</sub>	100	1.5 – 1800	0.9972	1.5	10

**Table 5. Method validation results for representative compounds of each drug class tested at three different QC levels**

Compound	Low QC				Medium QC				High QC			
	[QC] (ng/mL)	Within run (%CV)	Between run (%CV)	% Bias	[QC] (ng/mL)	Within run (%CV)	Between run (%CV)	% Bias	[QC] (ng/mL)	Within run (%CV)	Between run (%CV)	% Bias
Amphetamine	60	8.3	7.8	-0.1	2430	6.0	4.0	-4.8	4800	6.6	3.7	-0.7
Alprazolam	30	8.0	4.7	2.8	729	7.1	4.8	-2.2	1440	6.2	5.0	-0.4
Benzoylcegonine	100	5.0	4.1	9.4	2430	3.1	2.6	-4.9	4800	3.5	3.3	-0.4
Carfentanil	10	6.5	4.9	13.4	243	4.8	3.5	-5.2	480	6.0	4.2	0.8
EDDP	18	5.1	3.8	9.1	729	2.7	2.0	-4.5	1440	5.9	4.0	0.6
Ethylpentylone	30	4.3	3.4	10.3	729	4.1	2.8	-3.0	1440	4.4	3.5	1.7
Fluorofentanyl	6	3.5	3.3	16.5	243	4.4	3.2	-4.4	480	3.2	2.4	-1.4
Meperidine	18	11.6	6.9	15.0	729	6.3	3.8	-4.2	1440	3.5	2.4	0.9
Methadone	18	7.0	4.9	10.6	729	6.7	4.7	-3.8	1440	5.6	3.5	-0.1
Methamphetamine	60	4.2	2.9	4.2	2430	4.8	2.8	-4.9	4800	4.7	2.8	-0.4
Phencyclidine	18	10.8	6.0	12.7	729	5.0	3.8	-5.7	1440	6.2	3.6	0.0
Zolpidem	18	7.6	5.0	14.5	729	4.2	3.2	-4.9	1440	3.1	2.5	1.8

## Conclusion

The VeriSpray system was successfully applied for the accurate and precise quantitation of various drugs of abuse and their metabolites in human urine samples for use in clinical research or forensic toxicology. A total analysis time of ~2 minutes along with removing the requirement of sample pre-treatment/preparation and separation enables every clinical research and/or forensic toxicology laboratory to address more samples with increased confidence, and in turn, address their business and scientific goals.

## References

1. Liu, J.; Wang, H.; Manicke, N.E.; Lin, J-M.; Cooks, R.G.; Ouyang, Z. Development, characterization, and application of paper spray ionization. *Analytical Chemistry*. **2010**, *82*(6), 2463–71.
2. SWGTOX. Scientific Working Group for Forensic Toxicology (SWGTOX) standard practices for method validation in forensic toxicology. *Journal of Analytical Toxicology*. **2013**, *37*, 452–74.

Find out more at [thermofisher.com/verispray](http://thermofisher.com/verispray)