

# Quantitation of Bath Salts/Cathinones in Urine by LC-MS/MS

Kristine Van Natta, Marta Kozak, Thermo Fisher Scientific, San Jose, CA

## Key Words

Cathinone, bath salts, designer drugs, MDPV, methylone, mephedrone, ethylone, butylone, naphyrone, methedrone, TSQ Quantum Ultra, liquid/liquid extraction, forensic toxicology

## Goal

To develop an LC-MS/MS method for the analysis of the three Schedule I cathinones (MDPV, methylone and mephedrone), as well as other substituted cathinones (methedrone, ethylone, butylone and naphyrone) in urine with LOQs of 1 ng/mL for forensic toxicology.

## Introduction

Substituted cathinones, sometimes known as “bath salts,” have become the latest abused designer drugs. Based on cathinone, a substance found in the African *Catha edulis* (khat) plant, substituted cathinones are stimulants with amphetamine- and cocaine-like effects. As with many designer drug classes, variations on base structure abound (Figure 1). On October 21, 2011 the United States Drug Enforcement Agency (US DEA) listed three of the most common substituted cathinones: methylenedioxy-pyrovalerone (MDPV), methylone, and mephedrone, as Schedule I drugs, thereby making them illegal. As these drugs are not detected by current ELISA drug screening tests, new methods are needed to detect and quantify them.

## Experimental

### Sample Preparation

Deuterated internal standards were available for all compounds except methedrone and naphyrone. Butylone- $d_3$  was used as internal standard for methedrone and MDPV- $d_8$  was used for naphyrone.

Sample preparation was a liquid-liquid extraction (LLE). First, 200  $\mu$ L of urine and 10  $\mu$ L of internal standard mix solution (2  $\mu$ g/mL of each deuterated IS) were basified with 100  $\mu$ L of 1 N NaOH. Extraction was performed by adding 1 mL of ethylacetate/hexane (1:1), mixing, and centrifuging. Then, an 800  $\mu$ L aliquot of the resulting supernatant was transferred to a clean test tube containing 20  $\mu$ L of DMSO to prevent complete evaporation of solvent. Analytes have low molecular weight, are slightly volatile, and will evaporate if left too long in the evaporator. The supernatant was evaporated at 37  $^{\circ}$ C under nitrogen for 15 minutes. Samples were diluted with 200  $\mu$ L of 5% methanol and transferred to an HPLC vial equipped with a limited-volume insert. Finally, 20  $\mu$ L was injected into the LC-MS system.

### Liquid Chromatography

Chromatographic separations were performed under gradient conditions using a Thermo Scientific™ Accela™ 1250 pump and Accela Open autosampler. The analytical column was a Thermo Scientific™ Hypersil GOLD™ column (50  $\times$  2.1 mm, 1.9  $\mu$ m particle size). The column was maintained at room temperature. The injection volume was 20  $\mu$ L. Mobile phases A and B consisted of 10 mM ammonium formate with 0.1% formic acid in water and methanol, respectively. Mobile phase C was acetonitrile/1-propanol/acetone (45:45:10). All mobile phases were Fisher Chemical™ brand solvents. A shallow gradient at a flow rate of 500  $\mu$ L/min was used to separate isomeric ethylone and butylone. The total run time was 5 minutes.

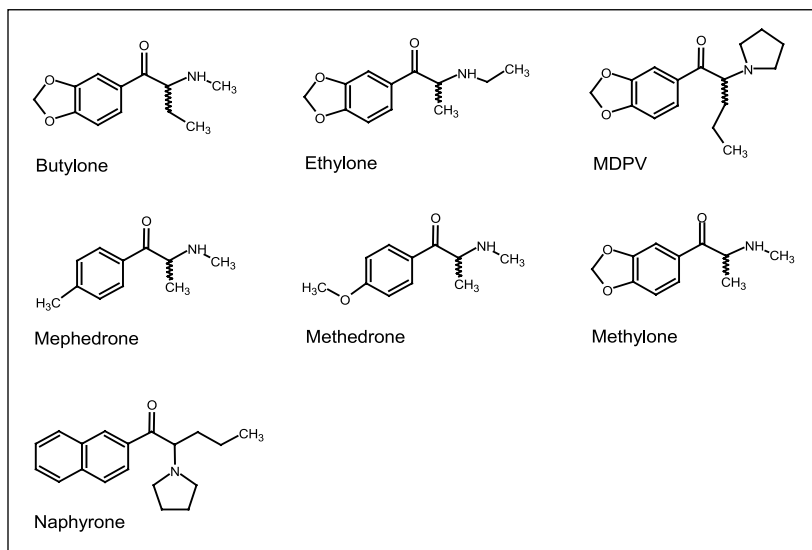


Figure 1. Structures of substituted cathinones (bath salts)

## Mass Spectrometry

MS analysis was carried out using a Thermo Scientific™ TSQ Quantum Ultra™ triple-stage quadrupole mass spectrometer equipped with a heated electrospray ionization (HESI-II) probe. Two selected-reaction monitoring (SRM) transitions were monitored for each analyte and each deuterated internal standard to provide ion ratio confirmations (IRC). Data acquisition and processing were performed using Thermo Scientific™ TraceFinder™ software.

## Validation

Standard calibration curves were prepared by fortifying pooled blank human urine with analytes. Quality control (QC) samples were prepared in a similar manner at low (LQC), middle (MQC), and high (HQC) concentrations. Intrarun variability and robustness were determined by processing six replicates of each QC level along with a calibration curve, as outlined in the Sample Preparation section, on three different days. Matrix effects were investigated by comparing peak areas of analyte at 10 ng/mL and internal standard prepared in twelve different lots of urine to those of a sample prepared in water.

## Results and Discussion

MDPV, methyone, mephedrone, methedrone, ethylone, and butylone were all linear from 1–1000 ng/mL. Figure 2 shows representative calibration curves for all compounds tested. Figure 3 shows representative chromatograms at 1 ng/mL for all compounds. Interassay quality control statistics shown in Table 1 demonstrate the method to be reproducible across the calibration range for the above compounds. Limited matrix effects were seen for the above compounds. These effects were largely mediated by deuterated internal standards. The absolute recoveries of all cathinones tested in various lots of urine, compared to a sample prepared in water, ranged from 85% to 132%. Relative recoveries ranged from 107% to 124%. Precision across all lots also improved when deuterated internal standards were used. Table 2 shows average statistics for all lots showing improvement in both precision and accuracy when internal standards were used.

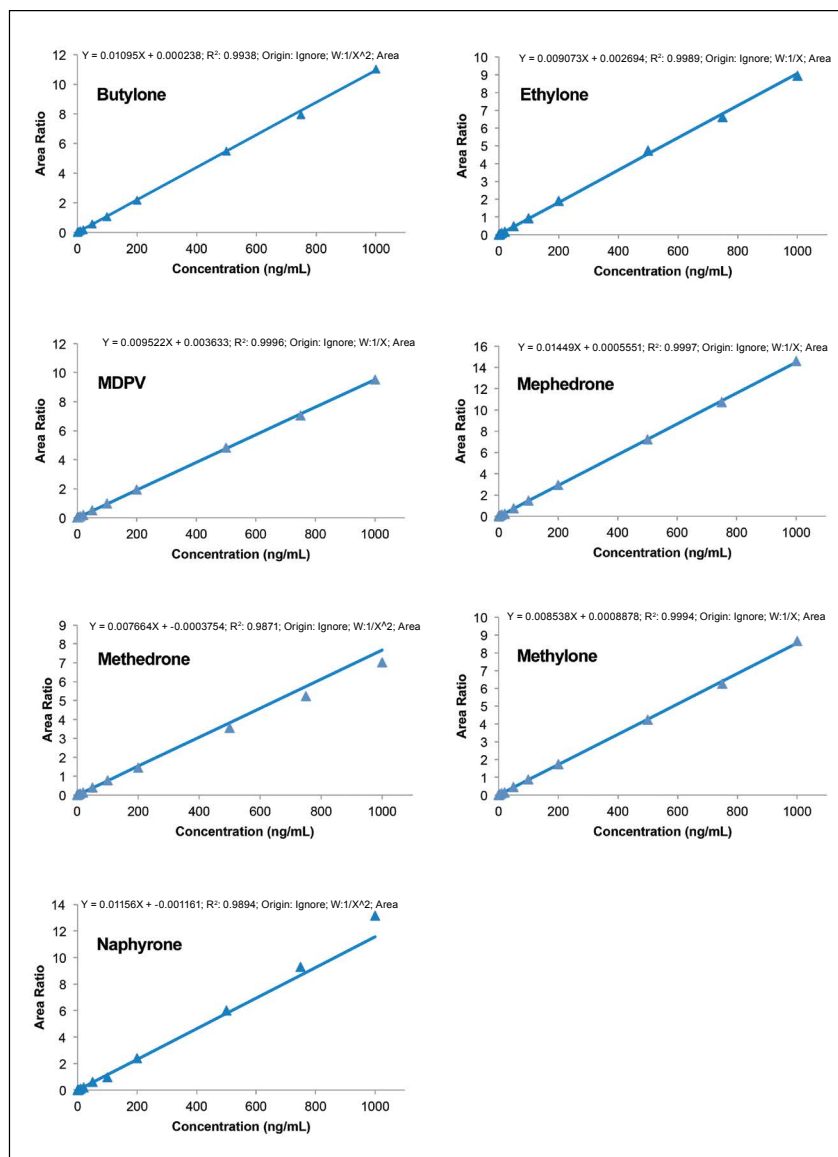


Figure 2. Representative calibrations curves for cathinones in urine

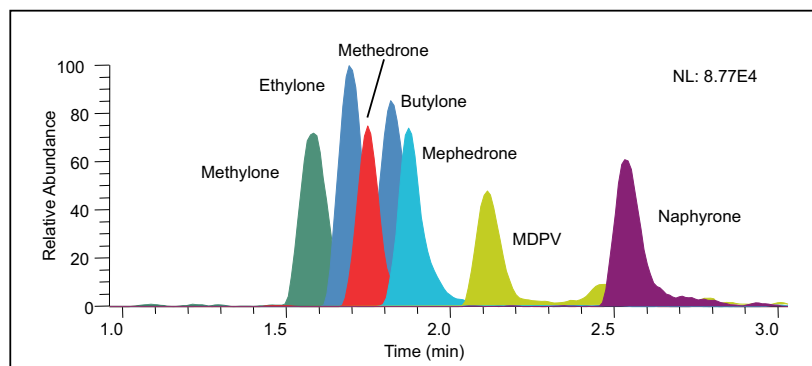


Figure 3. Representative chromatogram of cathinones at 1 ng/mL

Table 1. Interassay QC results

n = 18	LQC		MQC		HQC	
	%Bias	%CV	%Bias	%CV	%Bias	%CV
butylone	-5.80%	5.13%	1.40%	4.81%	-2.86%	3.76%
ethylone	-7.36%	5.93%	6.21%	1.72%	0.777%	1.71%
MDPV	-8.32%	5.48%	5.48%	2.89%	-0.907%	2.92%
mephedrone	-3.23%	2.79%	7.89%	2.64%	0.978%	2.00%
methedrone	-0.0565%	6.21%	8.09%	2.85%	-2.44%	1.79%
methylone	-3.95%	4.86%	6.06%	2.55%	0.394%	2.07%
naphyrone	-40.2%	59.7%	-18.2%	10.9%	-10.3%	8.68%

Table 2. Average imprecision and bias across all lots of urine

	Absolute Recovery		Relative Recovery	
	Imprecision	Bias	Imprecision	Bias
butylone	4.2%	12%	2.5%	13%
butylone- $d_3$	4.0%	-1.2%	na	na
ethylone	4.8%	16%	1.9%	18%
ethylone- $d_5$	4.9%	-1.5%	na	na
MDPV	3.0%	25%	1.6%	17%
MDPV- $d_8$	2.9%	7.1%	na	na
mephedrone	4.7%	21%	2.2%	16%
mephedrone- $d_3$	5.4%	4.3%	na	na
methedrone <sup>1</sup>	5.4%	18%	2.7%	19%
methylone	6.3%	15%	1.4%	17%
methylone- $d_3$	6.0%	-1.9%	na	na
naphyrone <sup>2</sup>	17%	49%	16%	39%

<sup>1</sup> Butylone- $d_3$  used as IS

<sup>2</sup> MDPV- $d_8$  used as IS

Although naphyrone was detected at 1 ng/mL, it showed more variability than the other compounds and a greater matrix effect from lot to lot. Absolute recoveries for naphyrone ranged from 113% to 207% while relative recoveries using MDPV- $d_8$  as internal standard ranged from 111% to 191%. All available internal standards were tried, and MDPV- $d_8$  showed the best results. A lack of a deuterated analog for naphyrone does not allow for matrix effect corrections and negatively affects method precision. In this assay, naphyrone should be considered qualitative.

## Conclusion

We achieved our goal of a 1 ng/mL LOQ for the three DEA-regulated cathinones, MDPV, mephedrone, and methylone, as well as methylone, ethylone, and butylone in urine for forensic toxicology. Naphyrone, which shows greater variability, can be detected down to 1 ng/mL in a qualitative manner. Deuterated internal standards are essential for rigorous quantitation of these compounds.

For Forensic Toxicology use only.

## [www.thermofisher.com](http://www.thermofisher.com)

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

<b>Africa</b> +43 1 333 50 34 0	<b>Denmark</b> +45 70 23 62 60	<b>Japan</b> +81 45 453 9100	<b>Singapore</b> +65 6289 1190
<b>Australia</b> +61 3 9757 4300	<b>Europe-Other</b> +43 1 333 50 34 0	<b>Latin America</b> +1 561 688 8700	<b>Spain</b> +34 914 845 965
<b>Austria</b> +43 810 282 206	<b>Finland</b> +358 9 3291 0200	<b>Middle East</b> +43 1 333 50 34 0	<b>Sweden</b> +46 8 556 468 00
<b>Belgium</b> +32 53 73 42 41	<b>France</b> +33 1 60 92 48 00	<b>Netherlands</b> +31 76 579 55 55	<b>Switzerland</b> +41 61 716 77 00
<b>Canada</b> +1 800 530 8447	<b>Germany</b> +49 6103 408 1014	<b>New Zealand</b> +64 9 980 6700	<b>UK</b> +44 1442 233555
<b>China</b> 800 810 5118 (free call domestic) 400 650 5118	<b>India</b> +91 22 6742 9494	<b>Norway</b> +46 8 556 468 00	<b>USA</b> +1 800 532 4752
	<b>Italy</b> +39 02 950 591	<b>Russia/CIS</b> +43 1 333 50 34 0	

**Thermo**  
SCIENTIFIC

Part of Thermo Fisher Scientific