## **ThermoFisher** SCIENTIFIC

# Quantitation of an Oral Fluid Drug Panel Including THC Using High Resolution Accurate-Mass (HRAM) Orbitrap Mass Spectrometry

# Courtney Patterson (1) and Kerry Hassell (1)

(1)Thermo Fisher Scientific, San Jose, California, United States

### Abstract

**Purpose**: The objective of this work was to accurately confirm and quantitate 31 drugs in oral fluid required by SAMHSA and the National Safety Council (NSC) by liquid chromatography and high-resolution, accurate mass (HRAM) Orbitrap mass spectrometry.

Methods: Human oral fluid samples were spiked with the 31 drugs of abuse at nine different concentration levels, extracted using DPX INTip<sup>™</sup> SPE, separated chromatographically, and detected on an Orbitrap mass spectrometer.

**Results:** All drugs achieved lower LOQ's than the cutoffs suggested by the new SAMHSA guidelines. Each of the drugs were also confirmed below those guideline levels using library search, isotopic pattern, and fragment matching.

#### Data Analysis

Data was acquired and processed with Thermo Scientific<sup>™</sup> TraceFinder<sup>™</sup> software, version 5.2 which utilizes its Compound Database that stores information including molecular formula, exact mass, retention time and fragment ions for all compounds of interest. A mass window of 5ppm was used as the precursor ion criteria of HRAM data.

### Results

The combined extracted ion chromatogram of each drug is depicted in Figure 2. This 7-minute method is able to separate each of the isomers and elute each compound in under 6-minutes.

#### Table 1. Criteria assigned in TraceFinder software for limits.

Parameter	Criteria
Limit of Quantitation (LOQ)	Back-calculated concentration on calibration curve within 20%.
Upper Limit of Linearity (ULOL)	Highest calibrator that achieves linearity
Limit of Identification (LOI)	<u>Orbitrap</u> : IP = passing isotopic pattern score (70) FI = presence of diagnostic fragment ions LS = passing library score (70)

 Table 2. Criteria assigned in TraceFinder software for Orbitrap

Table 3. Calibration and confirmation results of the 31 analytes in oral fluid. LOQ, ULOL, and LOI are in ng/mL.

	LOQ	ULOL	LOI	SAMHSA 2023	NSC Tier 1 Cutoff
6-MAM	0.5	1000	0.5	2	1
7-aminoclonazepam	0.5	1000	0.5	-	1
Alprazolam	0.5	1000	0.5	-	1
Amphetamine	1	1000	1	25	20
Benzoylecgonine	2	1000	2	8	8
Buprenorphine	0.5	1000	0.5	-	2
Carisoprodol	10	5000	10	-	500
Clonazepam	1	1000	1	-	1
Cocaethylene	1	1000	1	-	8
Cocaine	2	1000	2	8	8

2

0.5

0.5

5

5

25

2

0.5

1

0.5

1

15

15

15

25

25

25

15

0.5

20

20

500

10

20

10

10

1000

1000

1000

1000

1000

1000

1000

1000

5000

1000

1000

1000

1000

0.5

0.5

25

0.5

0.5

### Introduction

As the clinical and forensic communities move towards oral fluid matrix for ease of collection and roadside testing, it is important to be able to test for a wide range of analytes and achieve the required sensitivity. With the new SAMHSA guidelines providing LOQ levels for a list of drugs, the extraction protocol and instrumentation need to be sensitive enough to accomplish these cut-offs. Including tetrahydrocannabinol (THC) into the assay provides challenges in the extraction procedure as most drugs of abuse are basic and THC is neutral. This extraction workflow, which can extract THC alongside other drugs of abuse, coupled with the Orbitrap<sup>™</sup> mass spectrometer generates high-resolution accurate mass data that offers improved sensitivity, selectivity, and accuracy for the detection and quantitation of drugs of abuse in oral fluid.

### **Materials and methods**

#### **Sample Preparation**

Nine calibration levels ranging from 0.5 to 1,000 ng/mL (1.25 to 5,000 ng/mL for Meprobamate and Carisoprodol) were prepared by spiking stock solution of the 31 target analytes into human oral fluid. Samples were diluted with Quantisal<sup>™</sup> Buffer 1:3 (oral fluid: buffer) and 500 µL were aliquoted for extraction. Each sample was spiked with 125 µL of internal standard stock. The samples were then extracted using DPX INTip SCX/WAX SPE. The tips were conditioned with 800 µL x2 of 50% MeOH and then the samples were aspirated with the tips 4-5x. The tips were then washed with 500 µL x3 of 30% MeOH. Finally, the analytes were eluted with 500 µL x2 of 48% ACN, 48% MeOH, 4% Ammonium Hydroxide (v/v/v). Samples were dried down at 50°C for 25 minutes. The samples were reconstituted in 20 µL of MeOH + 0.1% Formic Acid and 80 µL of

Figure 2. Combined extracted ion chromatogram of 31 drugs in oral fluid



#### **Recovery Study**

A brief recovery study was performed to test the amount of analyte recovered from a pre-extraction spike compared to post-extraction spike. Figure 3 highlights the percent recovered per each compound.

Figure 3. Recovery study of the 31 drugs of abuse in oral fluid showing percent recovered



#### confirmation.

0.10 0.09 0.08 0.07 0.06 0.05 0.04

Parameter	Criteria
Isotopic Pattern	< 10 ppm mass deviation, < 20% intensity deviation, fit > 70%
Fragment Ion	At least 2 fragments with < 10 ppm mass deviation in $MS^2$ spectra
mzVault HRAM Library	Reverse search with > 70% match of ddMS2 spectra

Figure 4. Extracted ion chromatograms of A) THC and B) fentanyl at LOQ along with their respective calibration curves



	Oxazepam	0.5	1000	0.5	-			
14 3	Oxycodone	0.5	1000	0.5	15			
	Oxymorphone	0.5	1000	0.5	15			
	РСР	0.5	1000	0.5	10			
	Temazepam	0.5	1000	0.5	-			
	ТНС	1	1000	1	2			
	Tramadol	2	1000	2	-			
	Zolpidem	1	1000	1	-			
	THC Troubleshooting							
5.1 5.2	Throughout this study, some initial challenge themselves in regards to detecting THC in or							

Codeine

Diazepam

Fentanyl

Hydrocodone

Hydromorphone

Lorazepam

MDA

**MDMA** 

Meprobamate-Na

Methadone

Methamphetamine

Morphine

Nordiazepam

THC Troubleshooting Throughout this study, some initial challenges presented themselves in regards to detecting THC in oral fluids. One of the most important factors to consider in extracting THC is that it is a neutral compound and sticky to glass vials and other consumables. To combat this, silanized glass tubes and vial inserts were used to help prevent THC from sticking. It is also imperative to add MeOH to every step of the extraction process. This means adding enough volume of MeOH diluted internal standard to the sample (roughly 30%) and a wash solvent that contains 30% MeOH. This will improve the binding of THC to the SPE resin prior to its elution. When reconstituting, add the MeOH portion of the reconstitution solvent and vortex before adding the remaining aqueous portion.

### Conclusions

This fast and quantitative method was developed around 31 drugs

#### H2O + 0.1% Formic Acid.

#### Liquid Chromatography

Analytes were separated with the Thermo Scientific<sup>™</sup> Vanquish<sup>™</sup> Horizon ultra-high performance liquid chromatography (UHPLC) system by a 7-min gradient (Figure 1) using a Thermo Scientific<sup>™</sup> Accucore<sup>™</sup> Vanquish<sup>™</sup> Biphenyl column (2.6 µm, 50 x 2.1 mm). Mobile phases consisted of 0.1% formic acid in both water (mobile phase A) and methanol (mobile phase B). 5-µL of each standard were injected in triplicate.

# Figure 1. UHPLC Chromatographic gradient used for data acquisition.



#### Mass Spectrometry

Targeted analysis and quantitation were performed a Thermo Scientific<sup>™</sup> Orbitrap Exploris<sup>™</sup> 120 HRAM mass spectrometer. Fullscan, targeted, data-dependent MS2 (ddMS2) mode was used with an inclusion list for the targeted compounds. Resolutions of 60,000 (FWHM at m/z 200) for full scan and 15,000 for MS2 were employed. An isolation window of m/z 1.5 and compound specific collision energies were applied to generate rich HRAM MS2 spectra.

#### Quantitation

The limit of quantitation (LOQ), upper limit of linearity (ULOL), and limit of identification (LOI) were evaluated for each of the 31 analytes. Table 1 shows the criteria for each of these limits. For confirming each drug, isotopic pattern, fragment ion matching, and library search were employed. Table 2 defines these specific identification parameters.

Each of the 31 drugs of abuse achieved lower LOQ's and LOI's than the recommended confirmation cutoffs in the new 2023 SAMHSA guidelines. Additionally, each of the compounds had lower LOQ's and LOI's than the National Safety Council's Tier 1 drug cutoffs except for Diazepam which achieved an LOI of 2 ng/mL. More work will need to go into fine tuning to achieve the 1 ng/mL cutoff.

Figure 4 depicts the extracted ion chromatograms of two of the analytes, THC and fentanyl, at their respective LOQ concentrations with their corresponding calibration curves.



of abuse specified by SAMHSA and the National Safety Council. A complete workflow was presented that involved sample preparation using DPX INTip SPE. Linearity was achieved from LOQ's as low as 0.5 ng/mL up to 1,000 ng/mL which exemplifies the sensitivity of these instruments and extraction procedure. This method is fast and highly flexible for oral fluid testing in toxicology and clinical research labs as additional compounds may be added barring validation. This extraction and LCMS protocol sufficiently passes the SAMHSA guidelines and could be used by labs that want to be certified in oral fluid testing.

### **Acknowledgements**

Many thanks to Nick Chestara, Bill Kemnitzer, and Madison Kilpatrick from DPX Technologies for their continued support and expertise in developing the extraction protocol for this project.

### **Trademarks/licensing**

© 2024 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 the use of these products in any manner that might infringe on the intellectual property rights of others.