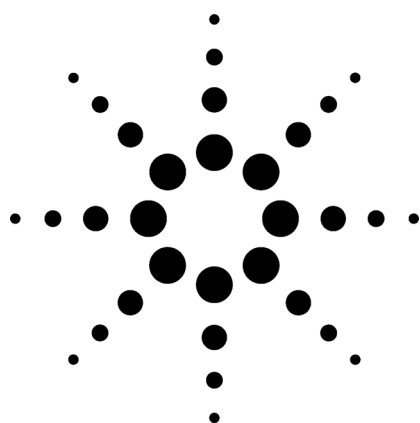


Development of a Screening Analysis by LC Time-Of-Flight MS for Drugs of Abuse

Application



Forensics

Authors

Courtney Milner and Russell Kinghorn
Baseline Separation Technologies Pty Ltd.

Abstract

The screening for drugs of abuse in human samples has relevance in much of today's society; employers, police and prison officials, and forensic pathologists all rely on the accuracy of drug screening. Currently, the most common method of this analysis is a straight-forward immunoassay technique, which although allowing for a rapid turnaround of screening samples, involves a slower confirmatory test of derivatization and detection by gas chromatography/mass spectrometry (GC/MS).

This application note presents the potential for the Agilent Time-of-Flight Mass Spectrometer (LC/MSD TOF) for use as both a screening and a confirmation tool in one analytical run of 30 minutes.

Introduction

History has shown that the human race has a long-standing fascination with the consumption and experimentation of mind and body altering substances. Traditionally, all of these substances occurred naturally. Even today, many drugs of

choice for this use are derived directly from natural substances. One of the most common of these, *cannabis*, is reported to have been tried on at least one occasion by up to 32% of the American population, and amongst the younger generation, 56% of high school senior students. These large numbers and the associated law enforcement issues have resulted in the decriminalization of its private use in many states and territories, but this does not diminish the potential impact on workplace performance or motor skills impairment, which may result in serious accidents. An extensive review of the illicit drug market in 25 major U.S. cities is provided in the Office of National Drug Control Policy Document "Pulse Check" [1].

Over the last 100 years, the physiological effects of many of the current illicit drugs were evaluated and reviewed, resulting in their subsequent banishment from society. During this time, a plethora of new drugs were developed, many finding wide acceptance within the medical community for the treatment of specific ailments. Unfortunately, the undesirable side effects of addiction or long-term abuse were often associated with the use of these drugs. The opiate class of drugs, which provided substantial improvement for the comfort of many patients, is an excellent example of one such class, as they are highly addictive and subject to abuse.



Agilent Technologies

Interest in the analysis of drugs of abuse covers many areas, all with different concerns in the results obtained. Some of the areas of significance include:

- Workplace screening
- Therapeutic monitoring
- Forensic pathology
- Accident investigation
- Crime scene investigation

Today, screening of drugs of abuse is performed through a variety of methods, with the most common lab-based technique being an Enzyme Multiple Immunoassay Test (EMIT), with a confirmatory analysis by GC/MS, if required. This immunoassay technique allows for screening to be performed and reported in as little as 2 hours, yet more commonly a 36–48 hour turnaround time is required. A further disadvantage of the EMIT technique is that it lacks the specificity to identify anything more than the class of drug detected.

The current analytical confirmatory technique of GC/MS was developed in order to achieve the sensitivity and specificity required to accurately determine the exact type and level of the drug compound, within the class indicated by the immunoassay technique. In order to achieve this detection, many of the drugs require derivatization to ensure adequate volatility and/or thermal stability required for GC analysis. See Table 1.

Table 1. National Institute of Drug Analysis Compound Class and Detection Limit Summary

| Compound class | Detection limits (ng/mL) | Confirmation |
|-----------------------|---------------------------------|---------------------|
| Amphetamines | 1000 | EMIT/GC/MS |
| Barbiturates | 300–3000 | EMIT/GC/MS |
| Cocaine | 300 | EMIT/GC/MS |
| Methadone | 300 | EMIT/GC/MS |
| Opiates | 300 | EMIT/GC/MS |
| Phencyclidine | 25 | EMIT/GC/MS |
| Propoxyphene | 300 | EMIT/GC/MS |
| Benzodiazepines | 300 | EMIT/GC/MS |
| Methaqualone | 300 | EMIT/GC/MS |
| Cannabinoids | 50 | EMIT/GC/MS |

Recently published Agilent application notes have shown the potential of LC/MS for the screening analysis and therapeutic monitoring of drugs of abuse using a single quadrupole instrument [2, 3]. Numerous other publications discuss selected drugs of abuse, or drug classes, illustrating the potential for the technique to one day replace GC/MS as either the confirmatory tool or as both the screening and confirmatory tool in one analysis.

Accurate mass measurement, such as that provided by the Agilent LC/MSD TOF, greatly increases the confidence of identification because it inherently limits the possible number of candidate compounds. The better the precision and accuracy of the mass measurement, the fewer the number of compounds theoretically possible for a given accurate mass. This is particularly useful for the analysis of samples from a variety of sources, each with their own potential interferences, such as those encountered with explosives residue analysis.

This application note provides an overview of the power of the Agilent TOF mass spectrometer for the screening and confirmation analysis of drugs of abuse. The TOF mass spectrometer provides accurate mass determinations (<3 ppm) with good linearity, proving its use as an excellent tool for the detection, confirmation, and quantitation of different drug classes. The method used here is not intended to represent one that will determine the lowest possible level of any one particular analyte or class of analytes, but rather is a procedure that could be expanded to cover a wider range of components used in screening analyses.

The compounds studied and their molecular formulas are shown in Table 2.

Table 2. Compounds Included in Study

| Compound | Molecular formula | Drug class |
|---------------------------------------|--|-------------------|
| α -hydroxyalprazolam | C ₁₇ H ₁₃ N ₄ OCl | Benzodiazepine |
| 7-Aminoclonazepam | C ₁₅ H ₁₂ N ₃ OCl | Benzodiazepine |
| Diazepam | C ₁₆ H ₁₃ N ₂ OCl | Benzodiazepine |
| Oxazepam | C ₁₅ H ₁₁ N ₂ O ₂ Cl | Benzodiazepine |
| Temazepam | C ₁₆ H ₁₃ N ₂ O ₂ Cl | Benzodiazepine |
| 7-Aminoflunitrazepam | C ₁₆ H ₁₄ N ₃ OF | Benzodiazepine |
| 7-Aminonitrazepam | C ₁₅ H ₁₃ N ₃ O | Benzodiazepine |
| dl-11-nor-9-carboxy- δ -9-THC | C ₂₁ H ₂₈ O ₄ | Cannabinoid |
| Codeine | C ₁₈ H ₂₁ NO ₃ | Opiate |
| Morphine 3 β -d-glucuronide | C ₂₃ H ₂₇ NO ₉ | Opiate |
| 6-acetylmorphine | C ₁₉ H ₂₁ NO ₄ | Opiate |
| EDDP perchlorate | C ₂₀ H ₂₄ NO ₄ Cl | Opiate |
| (+)-ephedrine | C ₁₀ H ₁₅ NO | Stimulant |
| Fenfluramine | C ₁₂ H ₁₆ NF ₃ | Stimulant |
| dl-MBDB:HCL | C ₁₂ H ₁₈ NO ₂ Cl | Stimulant |
| (\pm) BDB Hydrochloride | C ₁₁ H ₁₆ NO ₂ Cl | Stimulant |
| dl-MDEA | C ₁₂ H ₁₇ NO ₂ | Stimulant |
| dl-MDA | C ₁₀ H ₁₃ NO ₂ | Stimulant |
| dl-MDMA | C ₁₁ H ₁₅ NO ₂ | Stimulant |
| dl-Methamphetamine | C ₁₀ H ₁₅ N | Stimulant |
| dl-Amphetamine | C ₉ H ₁₃ N | Stimulant |
| Phentermine | C ₁₀ H ₁₅ N | Stimulant |
| (+)-Pseudoephedrine | C ₁₀ H ₁₅ NO | Stimulant |
| (-)-Cotinine | C ₁₀ H ₁₂ N ₂ O | Other |
| 4'-Hydroxynordiazepam | C ₁₅ H ₁₁ ClN ₂ O ₂ | Benzodiazepine |
| Nordiazepam | C ₁₅ H ₁₁ N ₂ OCl | Benzodiazepine |
| Flunitrazepam | C ₁₆ H ₁₂ N ₃ O ₃ F | Benzodiazepine |
| Flurazepam | C ₂₁ H ₂₃ N ₃ OClF | Benzodiazepine |
| Desalkylflurazepam | C ₁₅ H ₁₀ N ₂ OClF | Benzodiazepine |
| (-)- δ -9-THC | C ₂₁ H ₃₀ O ₂ | Cannabinoid |
| (\pm)-11-hydroxy- δ -9-THC | C ₂₁ H ₃₀ O ₃ | Cannabinoid |
| Cocaine | C ₁₇ H ₂₁ NO ₄ | Cocaine |
| Benzoylcegonine | C ₁₆ H ₁₉ NO ₄ | Cocaine |
| Buprenorphine | C ₂₉ H ₄₁ NO ₄ | Opiate |
| Morphine | C ₁₇ H ₁₉ NO ₃ | Opiate |
| Normorphine | C ₁₆ H ₁₇ NO ₃ | Opiate |
| Meperidine | C ₁₅ H ₂₁ NO ₂ | Opiate |
| Normeperidine | C ₁₄ H ₁₉ NO ₂ | Opiate |
| dl-Methadone | C ₂₁ H ₂₇ NO | Opiate |
| EMPD | C ₁₉ H ₂₁ N | Opiate |
| Naloxone | C ₁₉ H ₂₁ NO ₄ | Opiate |
| Oxycodone | C ₁₈ H ₂₁ NO ₄ | Opiate |
| LSD | C ₂₀ H ₂₅ N ₃ O | Hallucinogen |
| Iso-LSD | C ₂₀ H ₂₅ N ₃ O | Hallucinogen |
| (\pm)-phenylpropanolamine:HCL | C ₉ H ₁₃ NO:HCl | Stimulant |
| Fluoxetine:HCL | C ₁₇ H ₁₈ F ₃ NO:HCl | Prozac |
| GHB | C ₄ H ₇ O ₃ Na | Other |
| (-)-Nicotine | C ₁₀ H ₁₄ N ₂ | Other |

Methodology

The work undertaken in this study was performed on an Agilent 1100 system consisting of:

Binary pump
Standard auto-sampler
Thermostated column compartment
Diode Array Detector (DAD)
G1969 LC/MSD TOF.

Instrument Conditions

| Pump | | | |
|-------------------|-----------------------------------|--------------------------------------|---------------------------|
| Time (min) | % Water (0.1% formic acid) | % Methanol (0.1% formic acid) | Flow rate (mL/min) |
| 0 | 90 | 10 | 0.4 |
| 4 | 90 | 10 | 0.4 |
| 22 | 0 | 100 | 0.4 |
| 29 | 0 | 100 | 0.4 |
| 29.6 | 90 | 10 | 0.4 |
| 30 | 90 | 10 | 0.4 |

Post time: 5 minutes
Total run time: 35 minutes
Injection volume: 10 μ L, with needle wash
Column temperature: 30 $^{\circ}$ C
Column: ZORBAX SB-AQ,
150 mm \times 2.1 mm \times 3.5 μ m

MS Detection

| Ionization | ESI Positive |
|----------------------|---------------------|
| Gas temp | 350 $^{\circ}$ C |
| Drying gas | 10 L/min |
| Nebulizer pressure | 40 psig |
| Capillary V (+ve) | 3500 V |
| MS Conditions | |
| Scan m/z range | 100–1000 |
| Fragmentor | 125 V |
| Storage mode | Profile |
| Skimmer | 60 V |
| Oct RF | 200 V |

Reference Mass Introduction with LC-TOF

The Agilent TOF MS uses a reference mass in the generation of reliable high level accurate mass. The electro-spray source for the TOF is a unique dual spray assembly that allows the simultaneous constant introduction of a reference mass component.

The reference mix 1 used in these experiments consists of 2 mL of purine and 0.8 mL of HP-0921. This mixture was prepared in 1 L of 90:10 methanol:water to better represent the mobile phase.

The control software enables the use of the following reference masses:

Positive Ion Detection
121.050873
922.009798

Analysis of Drugs of Abuse by LC-TOF MS

An overwhelming advantage of using TOF MS for the trace level detection of any component is the confirmatory information that is provided through accurate mass. An example of this mass accuracy is shown in Table 3, where the observed masses for each component are detected, and their deviation from the theoretical masses for the adduct are shown.

The ability to closely match the expected mass and the observed mass provides the analyst with a higher level of confidence in the assignment given to a chromatographic peak. In the screening for components such as drugs, which may have a significant impact on the life of a person, this additional confidence is of great importance. This capability also allows the possibility of using this technique as a screening tool for a wide range of components.

Table 3. Theoretical Accurate Mass, Observed Mass and Mass Error

| Compound | Monoisotopic mass | Retention time | Adduct | Observed mass | Adduct accurate mass | Mass error (ppm) |
|---------------------------------------|-------------------|----------------|------------------------------------|---------------|------------------------|------------------|
| α -Hydroxyalprazolam | 324.0778 | 17.76 | [M+H] ⁺ | 325.0852 | 325.0850 | 0.41 |
| 7-Aminoclonazepam | 285.0669 | 13.32 | [M+H] ⁺ | 286.0739 | 286.0741 | -0.93 |
| Diazepam | 284.0716 | 19.15 | [M+H] ⁺ | 285.0796 | 285.0789 | 2.39 |
| Oxazepam | 286.0509 | 17.4 | [M+H] ⁺ | 287.0579 | 287.0581 | -0.98 |
| Temazepam | 300.0666 | 18.2 | [M+H] ⁺ | 301.0741 | 301.0738 | 0.89 |
| 7-Aminoflunitrazepam | 283.1121 | 15.3 | [M+H] ⁺ | 284.1191 | 284.1093 | -0.94 |
| 7-Aminonitrazepam | 251.1059 | 8.82 | [M+H] ⁺ | 252.1134 | 251.1131 | 1.04 |
| dl-11-nor-9-carboxy- δ -9-THC | 344.1988 | 21.38 | [M+H] ⁺ | 345.2061 | 345.2060 | 0.18 |
| Codeine | 299.1521 | 5.5 | [M+H] ⁺ | 300.1592 | 300.1594 | -0.73 |
| Morphine 3 β -d-glucuronide | 461.1686 | 1.7 | [M+H] ⁺ | 462.1764 | 462.1758 | 1.17 |
| 6-Acetylmorphine | 327.1471 | 8.9 | [M+H] ⁺ | 328.1542 | 328.1543 | -0.41 |
| EDDP perchlorate | 377.1394 | 15.42 | [M-O ₄ Cl] ⁺ | 278.1909 | 278.1903 | 2.06 |
| (+)-Ephedrine | 165.1154 | 2.46 | [M+H] ⁺ | 166.1225 | 166.1226 | -0.85 |
| Fenfluramine | 231.1235 | 12.9 | [M+H] ⁺ | 232.1303 | 232.1307 | -1.98 |
| dl-MBDB:HCL | 243.1026 | 10.63 | [M-Cl] ⁺ | 208.1337 | 208.1332 | 2.37 |
| (\pm) BDB hydrochloride | 229.087 | 9.5 | [M-Cl] ⁺ | 194.1181 | 194.1175 | 2.81 |
| dl-MDEA | 207.1259 | 9.6 | [M+H] ⁺ | 208.1332 | 208.1332 | -0.03 |
| dl-MDA | 179.0946 | 4.9 | [M+H] ⁺ | 180.1019 | 180.1191 | -0.03 |
| dl-MDMA | 193.1103 | 6.4 | [M+H] ⁺ | 194.1174 | 194.1175 | -0.08 |
| dl-Methamphetamine | 149.1204 | 3.85 | [M+H] ⁺ | 150.1281 | 150.1277 | 2.49 |
| dl-Amphetamine | 135.1048 | 3.05 | [M+H] ⁺ | 136.1125 | 136.112 | 3.11 |
| Phentermine | 149.1204 | 5.34 | [M+H] ⁺ | 150.1278 | 150.1277 | 0.49 |
| (+)-Pseudoephedrine | 165.1154 | 2.76 | [M+H] ⁺ | 166.1231 | 166.1226 | 2.76 |
| (-)-Cotinine | 176.095 | 2.56 | [M+H] ⁺ | 177.1023 | 177.1022 | 0.34 |
| 4'-Hydroxynordiazepam | 286.0509 | 14.23 | [M+H] ⁺ | 287.0582 | 287.0581 | 0.06 |
| Nordiazepam | 270.056 | 18.1 | [M+H] ⁺ | 271.0634 | 271.0632 | 0.49 |
| Flunitrazepam | 313.0863 | 18.1 | [M+H] ⁺ | 314.0924 | 314.0935 | -3.65 |
| Flurazepam | 387.1514 | 15.23 | [M+H] ⁺ | 388.1591 | 388.1586 | 1.17 |
| Desalkylflurazepam | 288.0466 | 18.2 | [M+H] ⁺ | 289.0535 | 289.0538 | -1.19 |
| (-)- δ -9-THC | 314.2246 | 22.31 | [M+H] ⁺ | 315.2328 | 315.2318 | 2.99 |
| (\pm)-11-hydroxy- δ -9-THC | 330.2195 | 21.07 | [M+H] ⁺ | 331.2267 | 331.2267 | -0.22 |
| Cocaine | 303.1471 | 12.6 | [M+H] ⁺ | 304.1545 | 304.1543 | 0.54 |
| Benzoyllecgonine | 289.1314 | 12.1 | [M+H] ⁺ | 290.1386 | 290.1386 | -0.29 |
| Buprenorphine | 467.3036 | 16.11 | [M+H] ⁺ | 468.3107 | 468.3108 | -0.29 |
| Morphine | 285.1365 | 2.2 | [M+H] ⁺ | 286.1438 | 286.1437 | -0.1 |
| Normorphine | 271.1208 | 2 | [M+H] ⁺ | 272.1286 | 272.1281 | 1.7 |
| Meperidine | 247.1572 | 12.4 | [M+H] ⁺ | 248.1652 | 248.1645 | 2.8 |
| Normeperidine | 233.1416 | 12.5 | [M+H] ⁺ | 234.1493 | 234.1488 | 1.9 |
| dl-Methadone | 309.2093 | 16.72 | [M+H] ⁺ | 310.2166 | 310.2165 | 0.19 |
| EMPD | 263.1674 | 16.4 | [M+H] ⁺ | 264.1754 | 264.1746 | 2.7 |
| Naloxone | 327.1471 | 5.74 | [M+H] ⁺ | 328.1541 | 328.1543 | -0.72 |
| Oxycodone | 315.1471 | 7.2 | [M+H] ⁺ | 316.1547 | 316.1543 | 1.15 |
| LSD | 323.1998 | 14.7 | [M+H] ⁺ | 324.2073 | 324.207 | 0.8 |
| Iso-LSD | 323.1998 | 14.55 | [M+H] ⁺ | 324.2078 | 324.207 | 2.34 |
| (\pm)-Phenylpropanolamine:HCL | 187.0764 | 1.95 | [M-Cl] ⁺ | 152.1069 | 152.1069 | -0.59 |
| Fluoxetine:HCL | 345.1107 | 16.3 | [M-Cl] ⁺ | 310.1412 | 310.1413 | 1.85 |
| GHB | 126.0293 | 1.72 | [M+H] ⁺ | 127.0369 | 127.0365 | 2.63 |
| (-)-Nicotine | 162.1157 | 1.6 | [M+H] ⁺ | 163.1233 | 163.1229 | 1.99 |
| 2-oxo-3-hydroxy-LSD | 355.1896 | n.d. | | | Not detected in +veESI | |

A greater than two-fold increase in sensitivity for many components is seen with the narrowing of the mass-extraction window. Figure 1 shows the reduction in noise that is observed with the extraction of a smaller mass range for flunitrazepam, commonly known as Rohypnol, a date-rape drug. The ability of TOF-MS to accurately determine the presence of components such as Rohypnol at low levels may assist with investigations into reported abuse of the illicit substance, and prove to be a critical factor in confirmation when dealing with complex matrices.

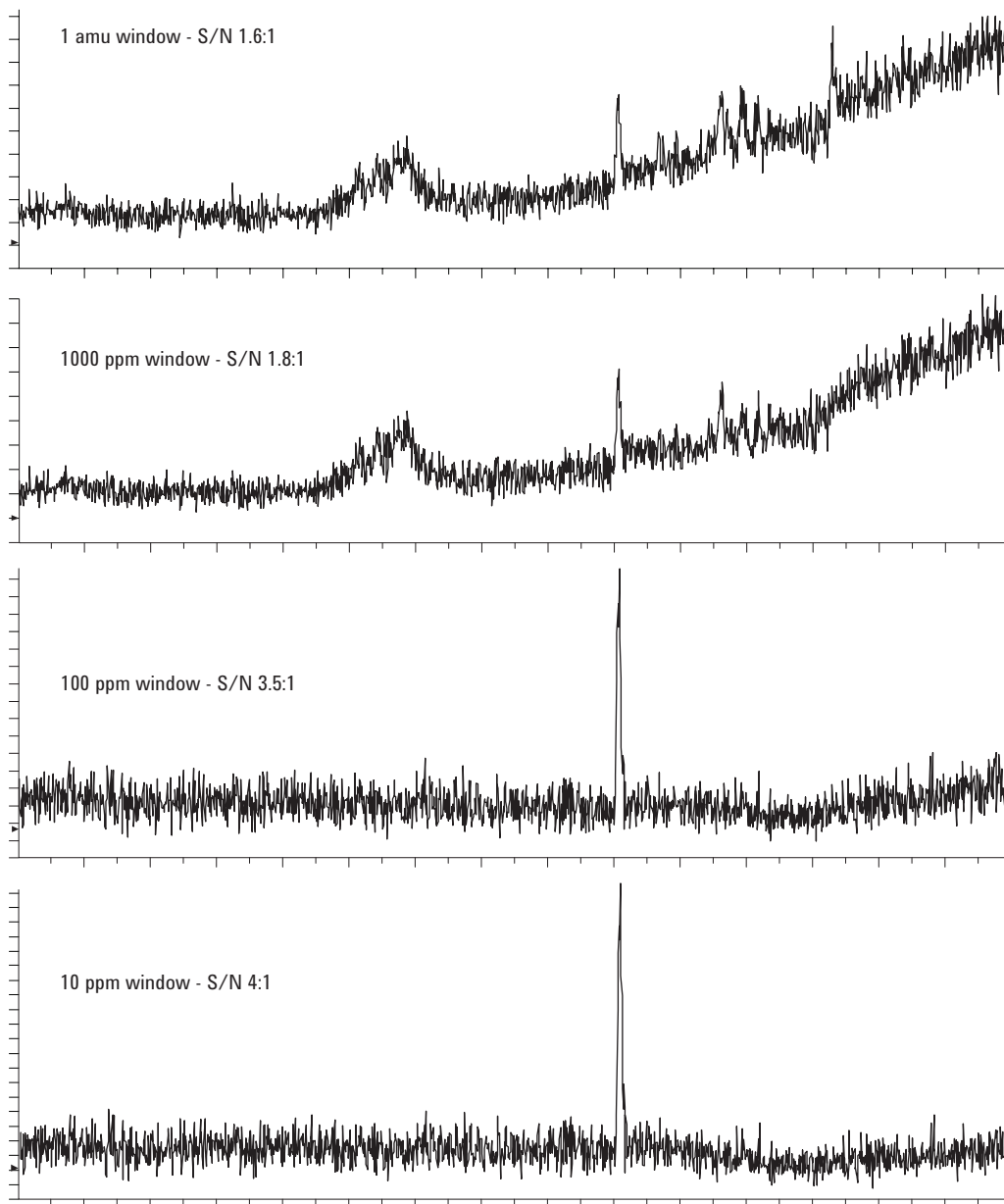


Figure 1. Effect of extracted ion range on noise – 1 ng/mL flunitrazepam.

TOF Linearity

TOF-MS has traditionally been considered as unsuitable for quantitation due to the use of time-to-digital conversion of data. The Agilent TOF MS uses analog-to-digital conversion, allowing for far better quantitative data than the alternative technology of time-to-digital conversion. Several of the components analyzed by TOF were tested for

linearity as part of this study. Figures 2–5 show the linearity of four selected components, most displaying linearity over three orders of magnitude from 1 ng/mL to 1000 ng/mL. However, some components, such as δ -9-THC and fluoxetine, only exhibit a narrower linear range, a result of their ionization behavior (Figures 6 and 7). Nevertheless, regression values of over 0.999 were seen for each of these components.

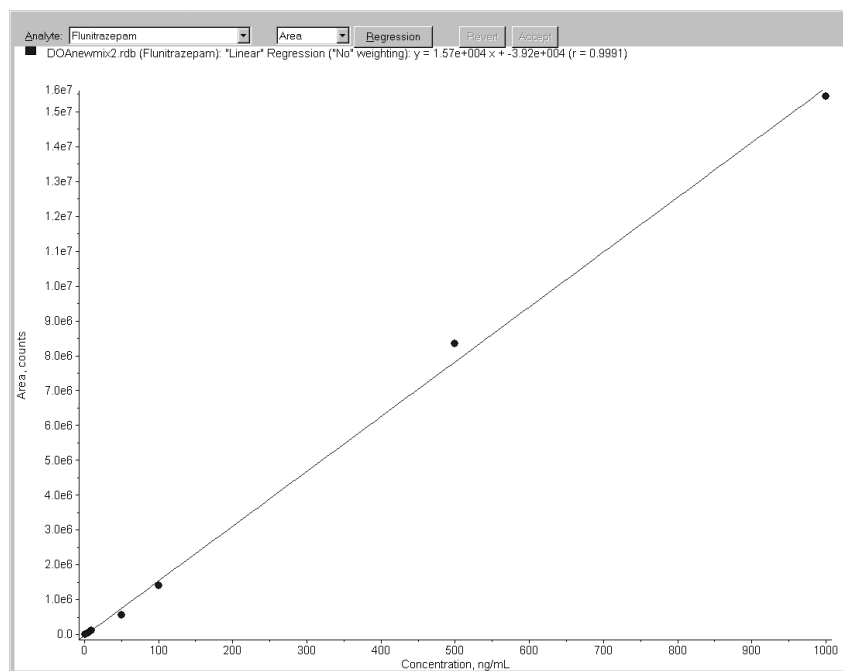


Figure 2. Calibration curve for flunitrazepam from 1 ng/mL to 1000 ng/mL with TOF-MS.

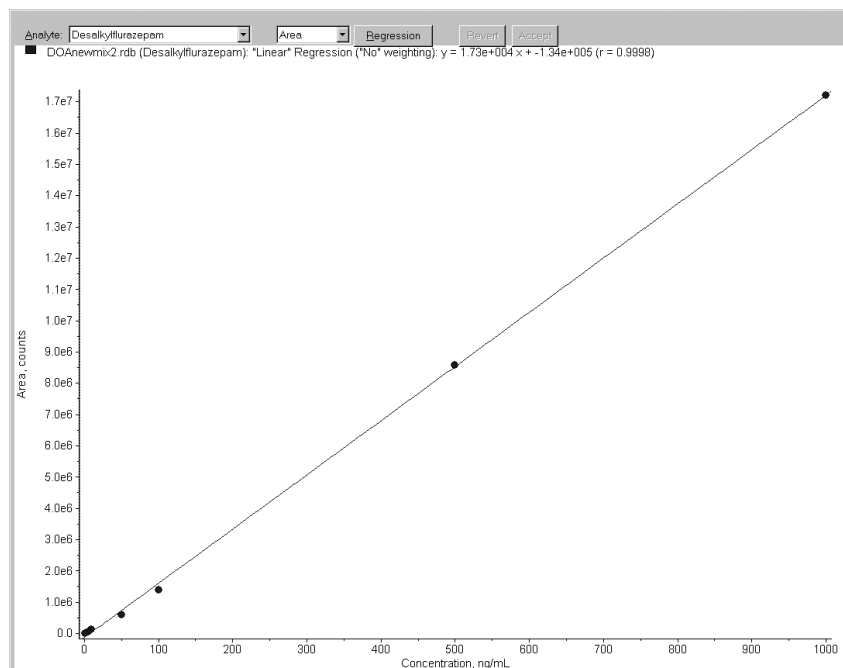


Figure 3. Calibration curve for desalkylflurazepam from 1 ng/mL to 1000 ng/mL with TOF-MS.

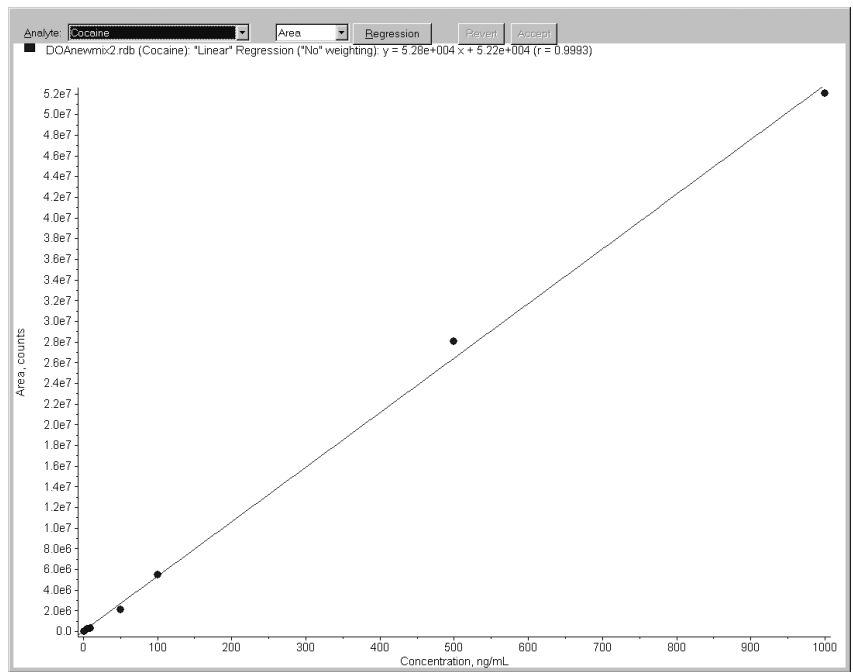


Figure 4. Calibration curve for cocaine from 1 ng/mL to 1000 ng/mL with TOF-MS.

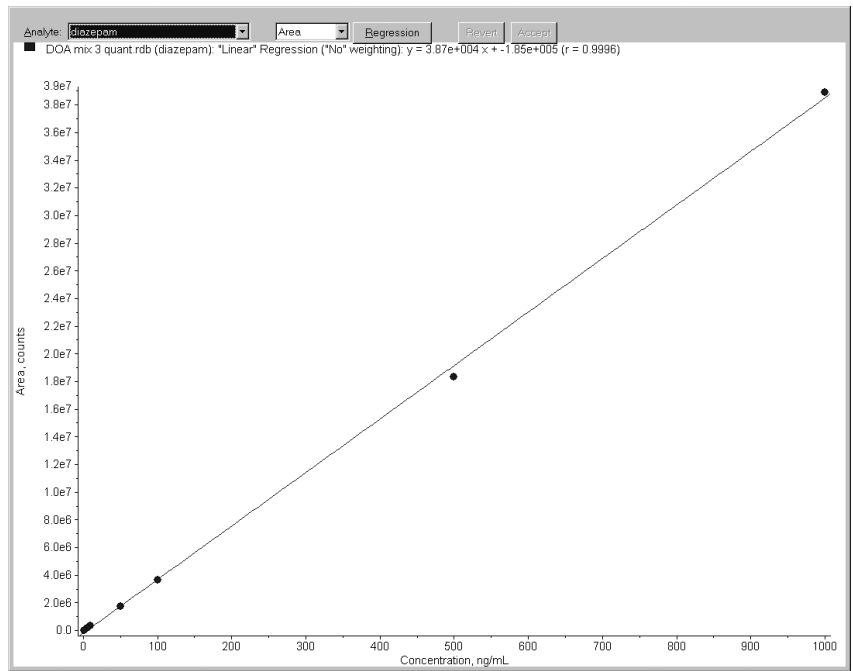


Figure 5. Calibration curve for diazepam from 1 ng/mL to 1000 ng/mL with TOF-MS.

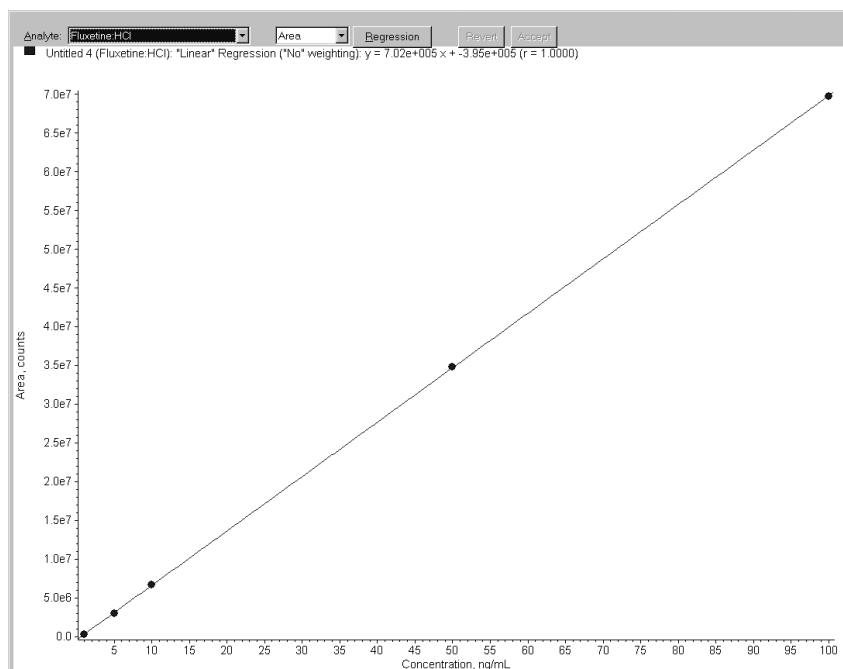


Figure 6. Calibration curve for (-)- δ -9-THC from 1 ng/mL to 100 ng/mL with TOF-MS.

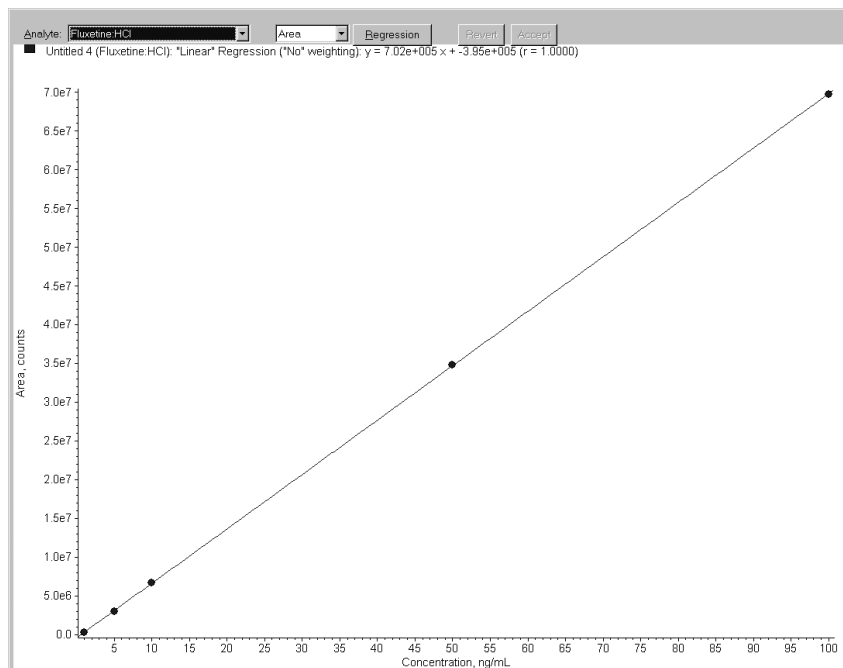
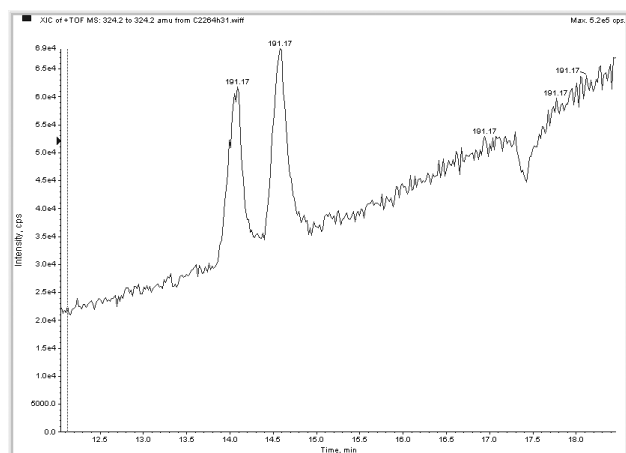
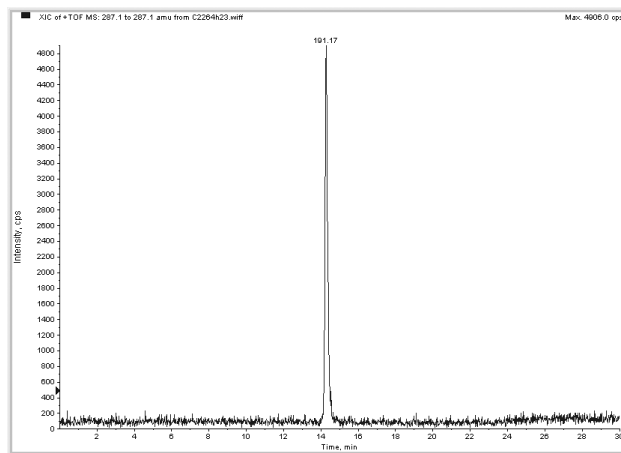


Figure 7. Calibration curve for fluoxetine from 1 ng/mL to 100 ng/mL with TOF-MS.

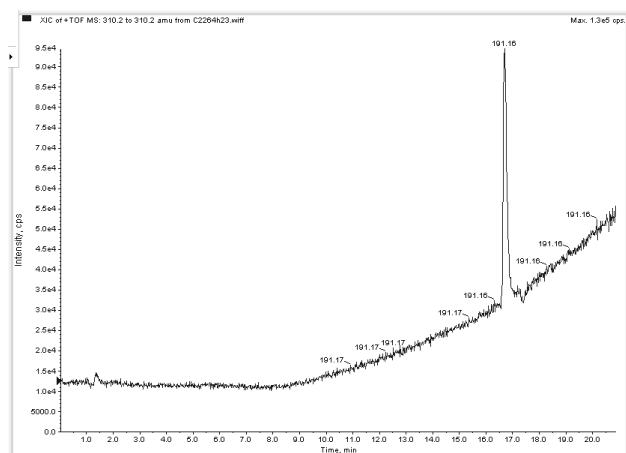
Chromatogram examples for four components at 1 ng/mL are shown below in Figure 8 with 10 ppm extraction windows.



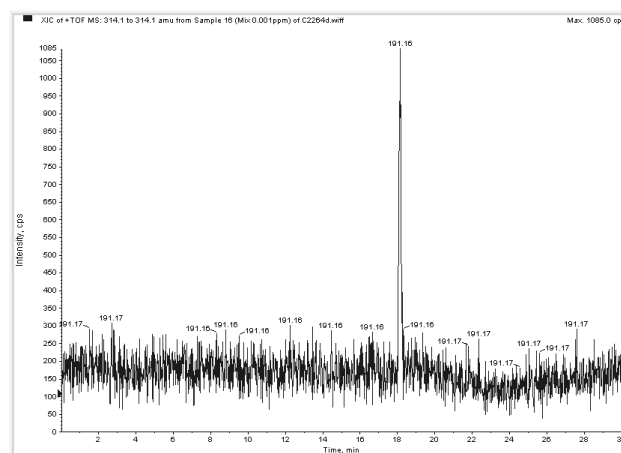
1 ng/mL iso-LSD and LSD (respectively)



1 ng/mL 4'-hydroxynordiazepam



1 ng/mL dl methadone



1 ng/mL flunitrazepam (Rohypnol)

Figure 8. Ten ppm extraction window of 1 ng/mL solutions of four drugs of abuse.

Table 4 summarizes the limits of detection (LOD) for each of the components analyzed under this method. Note that while the method has not been optimized for any one component, it is designed to provide a broad screening tool for the analysis of drugs of abuse.

Table 4. LOD for Components by LC/MSD TOF

| Component | Accurate mass | LOD (ng/mL) |
|-----------------------------|------------------------|--------------------|
| α-Hydroxyalprazolam | 325.085 | 5 |
| 7-Aminoclonazepam | 286.0741 | 1 |
| Diazepam | 285.0789 | 1 |
| Oxazepam | 287.0581 | 5 |
| Temazepam | 301.0738 | 5 |
| 7-Aminoflunitrazepam | 284.1093 | 5 |
| 7-Aminonitrazepam | 251.1131 | 1 |
| dl-11-nor-9-carboxy-δ-9-THC | 345.206 | 5 |
| Codeine | 300.1594 | 50 |
| Morphine 3β-d-glucuronide | 462.1758 | 5 |
| 6-Acetylmorphine | 328.1543 | 10 |
| EDDP perchlorate | 278.1903 | 1 |
| (+)-Ephedrine | 166.1226 | 20 |
| Fenfluramine | 232.1307 | 1 |
| dl-MBDB:HCL | 208.1332 | 5 |
| (±) BDB Hydrochloride | 194.1175 | 10 |
| dl-MDEA | 208.1332 | 10 |
| dl-MDA | 180.1191 | 50 |
| dl-MDMA | 194.1175 | 50 |
| dl-Methamphetamine | 150.1277 | 20 |
| dl-Amphetamine | 136.112 | 50 |
| Phentermine | 150.1277 | 20 |
| (+)-Pseudoephedrine | 166.1226 | 20 |
| (-)-Cotinine | 177.1022 | 20 |
| 4'-Hydroxynordiazepam | 287.0581 | 0.5 |
| Nordiazepam | 271.0632 | 0.5 |
| Flunitrazepam | 314.0935 | 1 |
| Flurazepam | 388.1586 | 1 |
| Desalkylflurazepam | 289.0538 | 1 |
| (-)-δ-9-THC | 315.2318 | 1 |
| (±)-11-Hydroxy-δ-9-THC | 331.2267 | 1 |
| Cocaine | 304.1543 | 1 |
| Benzoyllecgonine | 290.1386 | 5 |
| Buprenorphine | 468.3108 | 1 |
| Morphine | 286.1437 | 10 |
| Normorphine | 272.1281 | 10 |
| Meperidine | 248.1645 | 1 |
| Normeperidine | 234.1488 | 1 |
| dl-Methadone | 310.2165 | 1 |
| EMPD | 264.1746 | 1 |
| Naloxone | 328.1543 | 5 |
| Oxycodone | 316.1543 | 20 |
| LSD | 324.207 | 1 |
| Iso-LSD | 324.207 | 1 |
| (±)-Phenylpropanolamine:HCL | 152.1069 | 1 |
| Fluoxetine:HCL | 310.1413 | 1 |
| GHB | 127.0365 | 20 |
| (-)-Nicotine | 163.1229 | 5 |
| 2-oxo-3-hydroxy-LSD | Not detected in +veESI | ND |

Detection of Drugs of Abuse in Bodily Fluids

Urine is the matrix of choice for the detection of drugs of abuse in areas such as workplace screening, and was therefore chosen to evaluate the LC/TOF MS method developed. To present a “worst case scenario”, neat urine was spiked for this analysis and run directly with no sample clean-up. This would not normally be done; however, it was used as an illustration of the method’s ability to provide a quick screening result without cleanup.

A further discussion of a solid phase extraction (SPE) sample preparation method that may be considered can be found in Agilent Technologies application note 5989-2260EN [4].

In the first instance, neat and spiked urine was scanned for both cocaine and benzoylecgonine, a metabolite of cocaine (Figures 9 and 10). Recoveries of the spiked samples (100 ng/mL – representing a level lower than the traditional EMIT screen) were both approximately 100%.

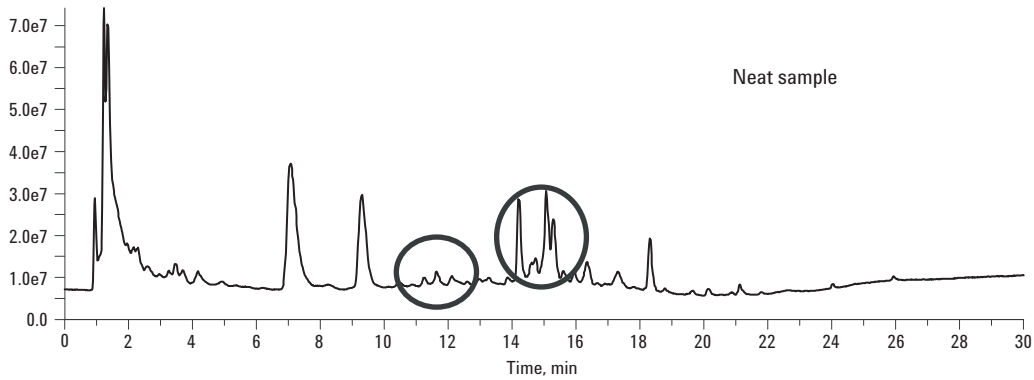


Figure 9. TIC of Blank Urine – Injected neat – circle shows expected retention time of analytes of interest.

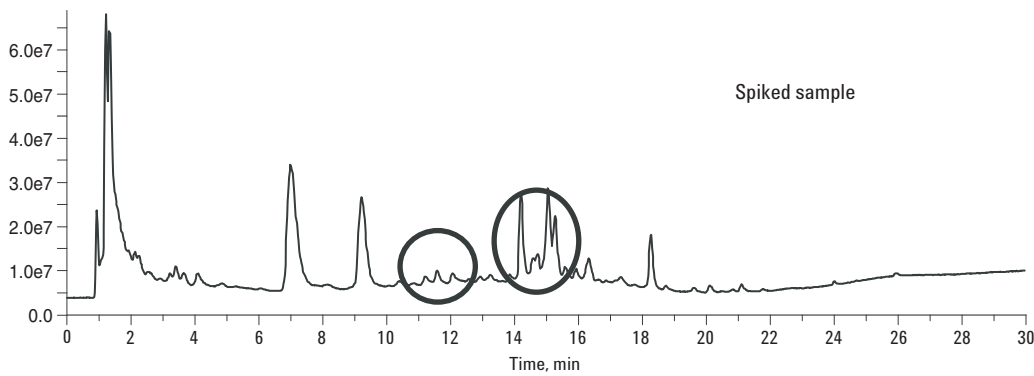


Figure 10. TIC of Neat Urine with 100 ng/mL spike of cocaine, benzoylecgonine, flunitrazepam, and 7-aminoflunitrazepam.

It can be seen from Figures 11 and 12 that the ability to narrow the mass extraction window greatly reduces the noise for a given mass, and with retention time information can provide a high level of confidence in the assignment of a component.

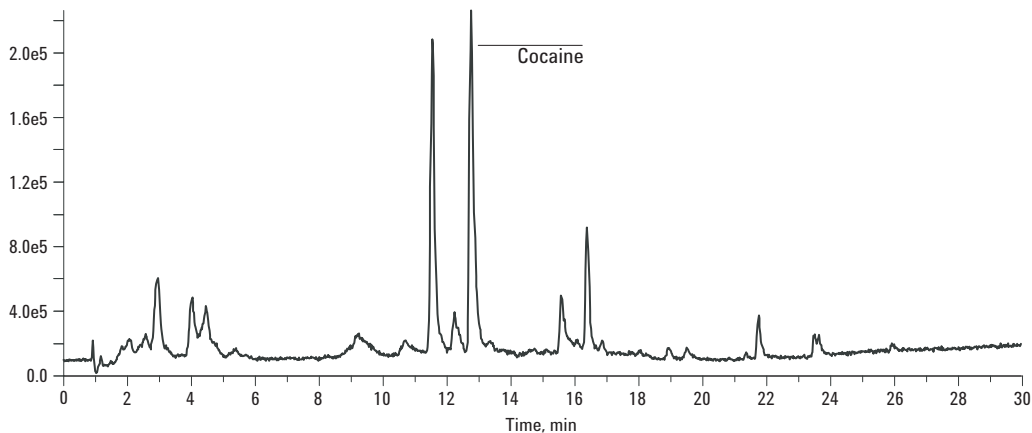


Figure 11. One amu extraction window of scanned target mass 304.1543 – cocaine.

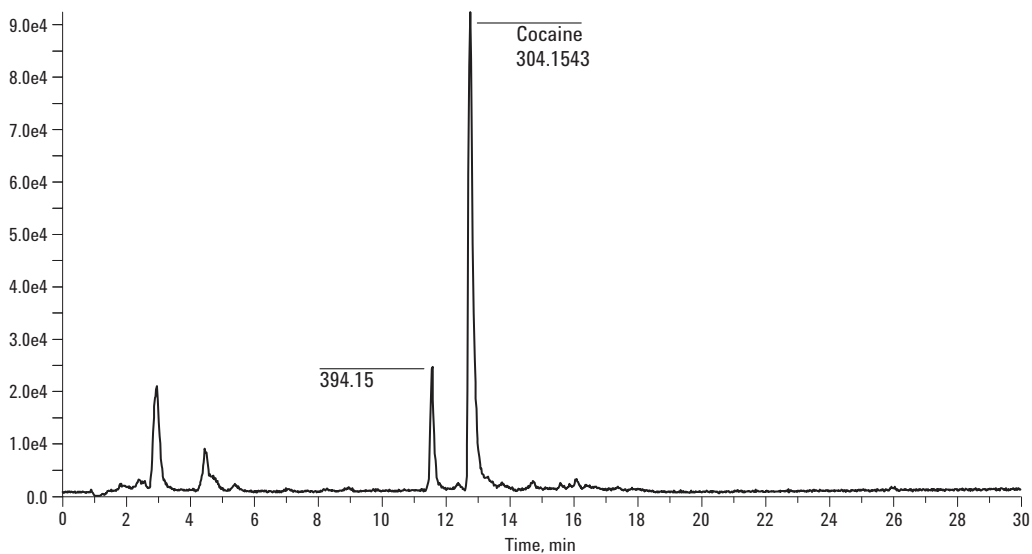


Figure 12. Ten ppm extraction window of scanned target mass 304.1543 – cocaine.

In the instance with Figure 12, the larger peak at approximately 12.9 minutes shows an excellent match with cocaine (mass error of ~ 0.2 ppm), while the earlier peak at 11.5 minutes has a mass of 394.15 and a fragmentation product in the extraction window (Figures 13 and 14).

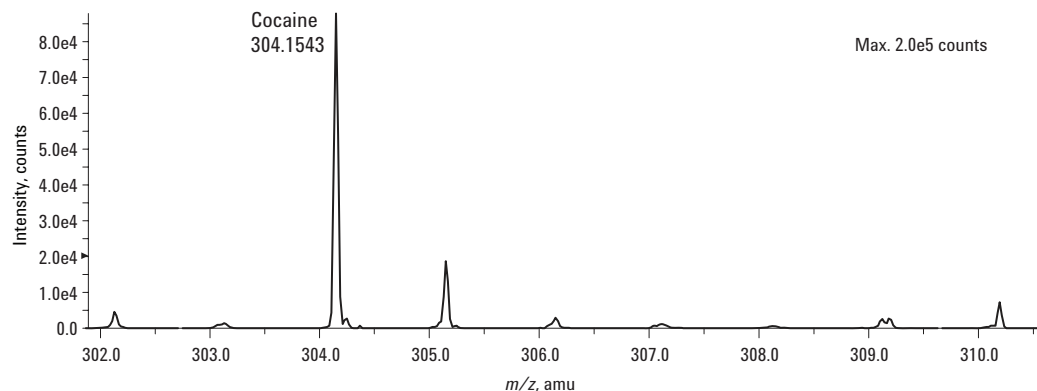


Figure 13. Confirmation of mass of cocaine.

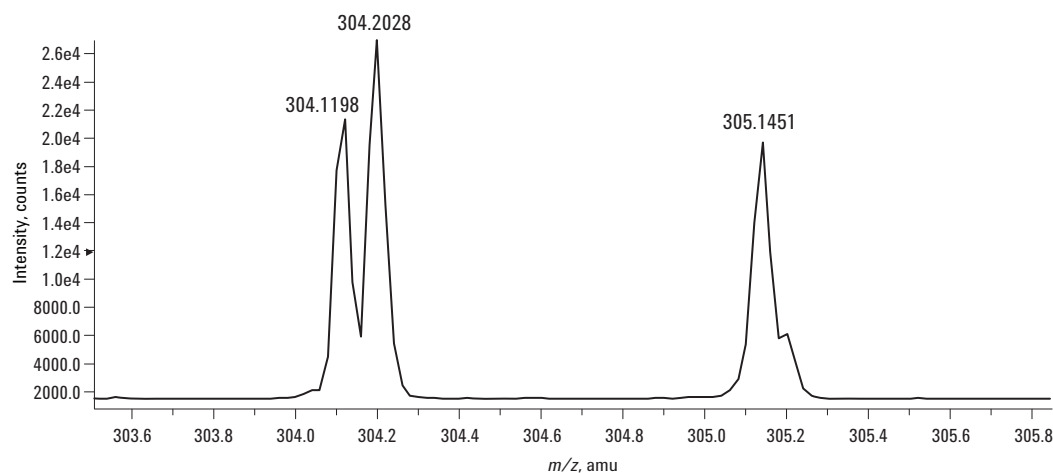


Figure 14. Mass spectrum 304 – 305 amu for peak at 11.9 minutes – no match of mass for cocaine.

A second drug that was of particular interest was flunitrazepam, another substance used as a date rape drug (Rohypnol), and one of its metabolites, 7-aminoflunitrazepam. Again, these components were spiked at 100 ng/mL into neat urine and injected directly. The recoveries achieved for these two components were approximately 40%, which is likely a result of ion suppression in the source. However, due to the excellent detection limits possible, even with up to 60% suppression of the signal, a clear peak can be seen for both compounds at 100 ng/mL (a level well below the EMIT screening reporting limits) when extracting a narrow mass window (Figures 15 and 16).

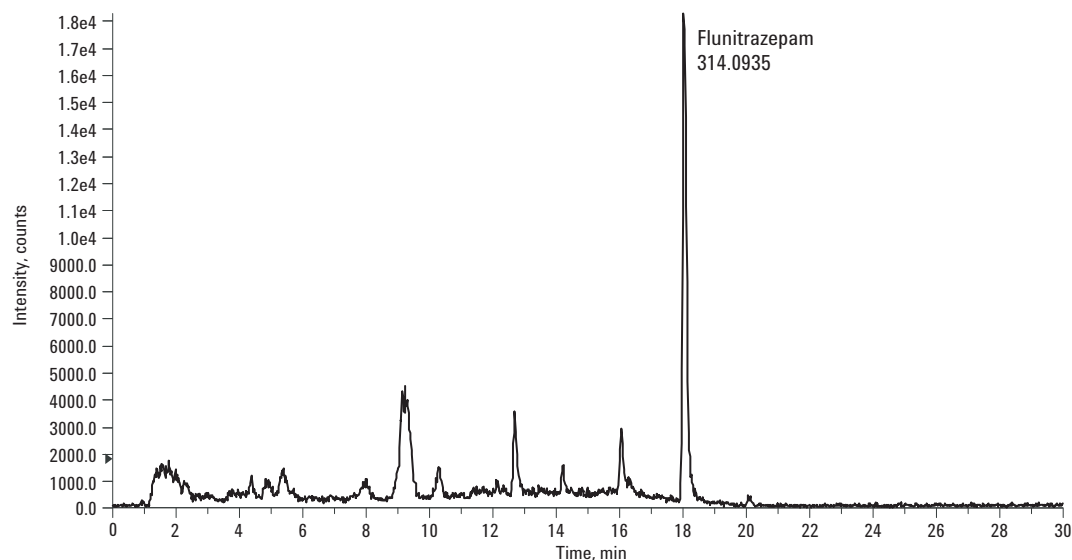


Figure 15. Ten ppm extraction window of scanned target mass 314.0935, flunitrazepam in neat urine. Retention time helps confirm presence at 18.1 minutes.

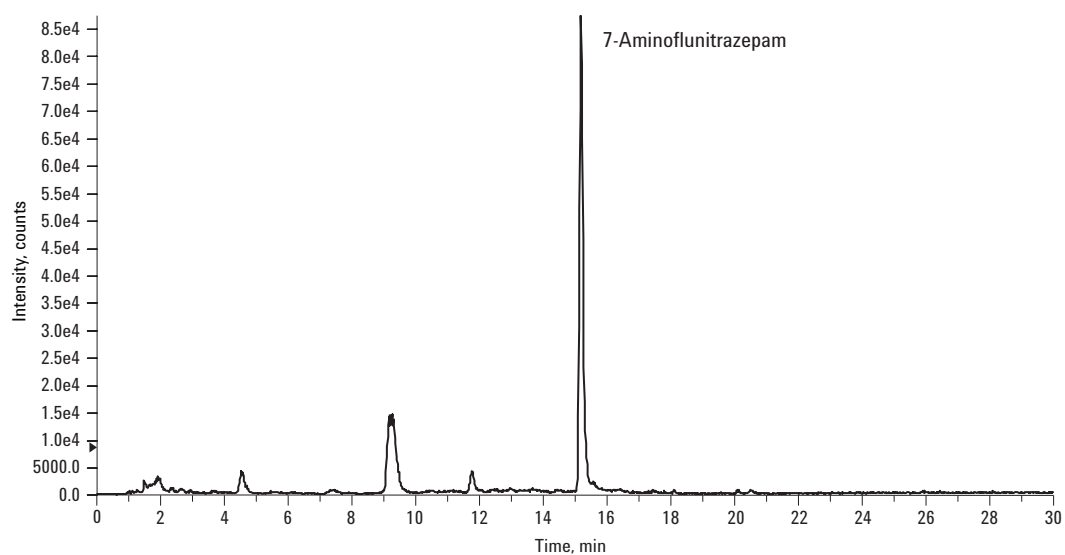


Figure 16. Ten ppm extraction window of scanned target mass 284.1191, 7-aminoflunitrazepam in neat urine. Retention time helps confirm presence at 15.3 minutes.

The second common matrix encountered in the screening of drugs of abuse is blood and plasma. To test the method when analyzing plasma, a sample was spiked with desalkylflurazepam. Sample preparation was again kept to a minimum, with a simple acetonitrile precipitation performed on the sample prior to injection. The effect of the mass extraction window on the detection of peaks, seen in Figures 17 to 20, shows the removal of interferences from the spiked plasma sample.

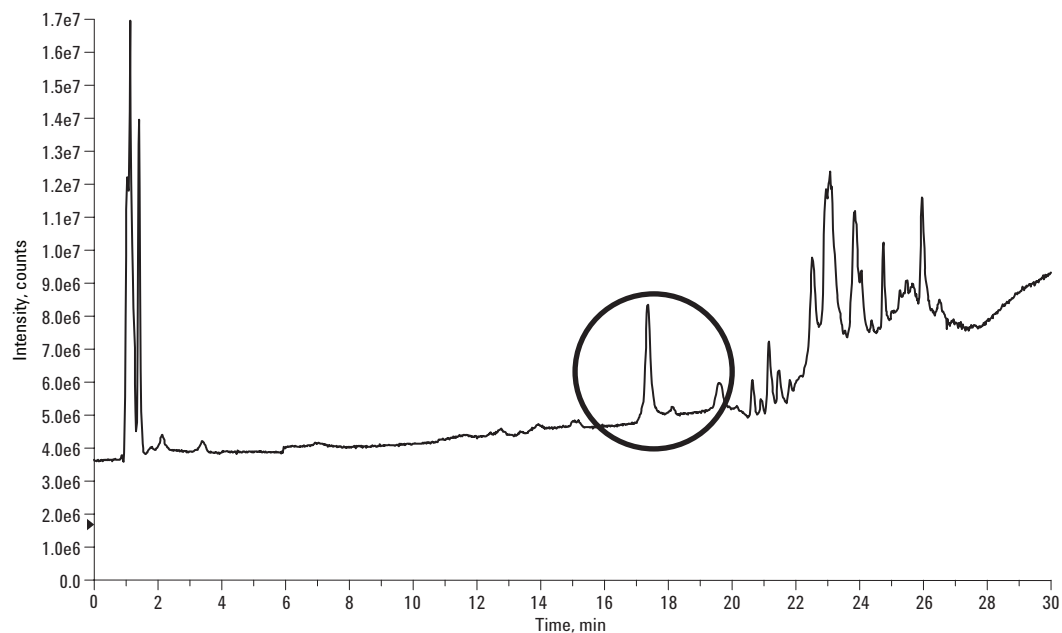


Figure 17. TIC of unspiked plasma.

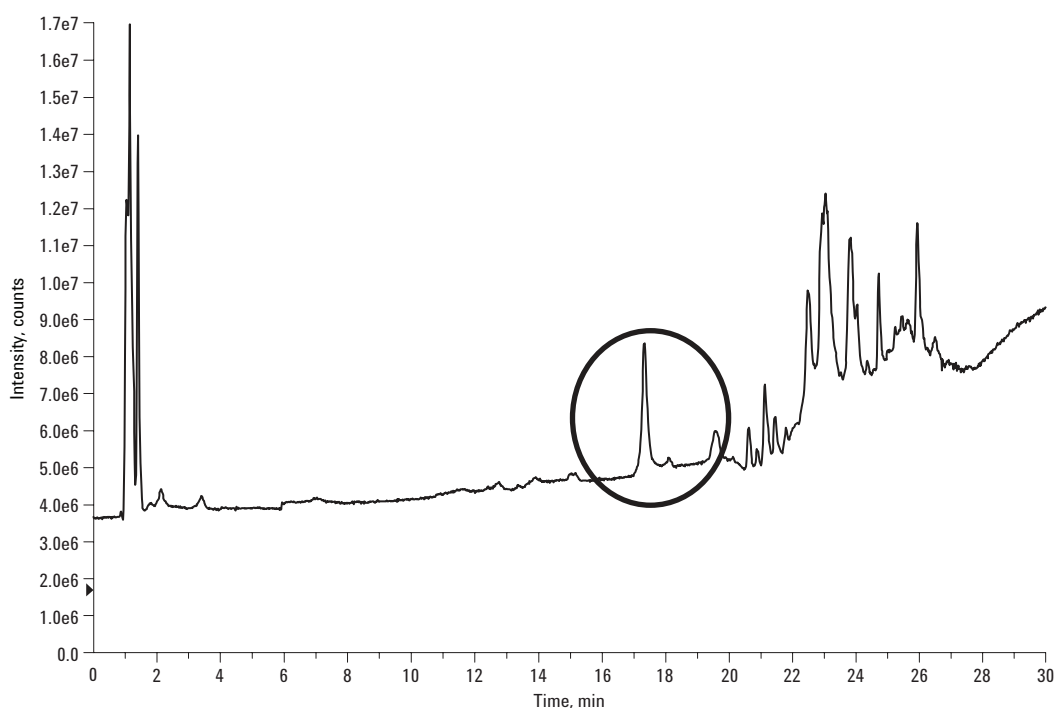


Figure 18. TIC of plasma spiked with 200 ng/mL of desalkylflurazepam.

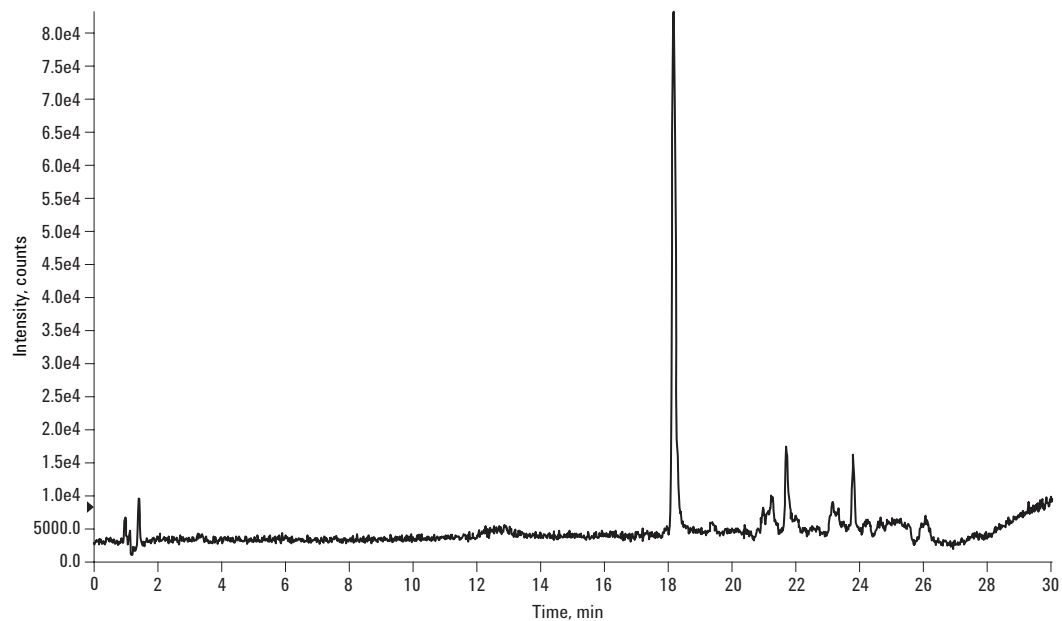


Figure 19. Mass extraction window (0.1 amu) of spiked plasma sample (346 ppm mass window).

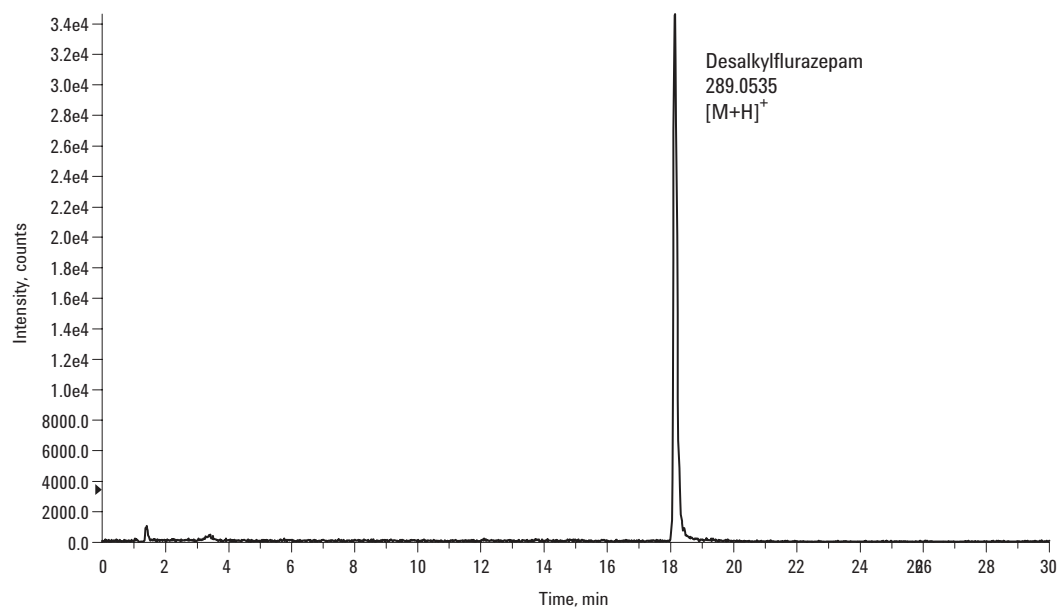


Figure 20. Ten ppm (0.003 amu) mass extraction window of spike plasma – note minimal noise.

Analysis of Coronial Samples – Using a Screener Database

The instrumentation and software provides the user the ability to create a screener database for all components they wish to automatically screen for. The minimum requirement for this database is the empirical formula and name for the component of interest, although the inclusion of a retention time will assist with confidence in the confirmation and reduce analysis time.

Several samples were acquired from the local coronial office to test the procedure that was developed. These samples were provided as butyl chloride extracts of blood samples obtained from deceased persons, for screening using the developed method.

The coronial samples supplied were screened using a database created from the 48 components analyzed under this method.

Sample Preparation

Samples were obtained from 1-mL blood volumes, liquid-liquid extracted with 6–8 mL of butyl chloride following centrifugation. Organic layer evaporated to dryness and then reconstituted in 100- μ L mobile phase for a final 10-fold concentration.

Sample 1

Sample 1 was known to contain amphetamine, codeine, diazepam, and nordiazepam from the previous analysis performed at the coronial office. In addition to the four previously reported components, the screen also indicated the presence of

- Nicotine
- Cotinine
- Acetylmorphine
- Ephedrine
- Methamphetamine
- Pseudoephedrine

The total ion chromatogram (TIC) for this sample is shown in Figure 21.

An excerpt of the screen report is shown in Figure 22 with the details for cotinine. For each component included in the compound database, the screening report displays the extracted ion chromatogram, spectra of detected peak, and enlarged spectra of the target mass. This is accompanied by a summary table with the mass and retention time error. In this instance, an excellent match is seen for both retention time and mass.

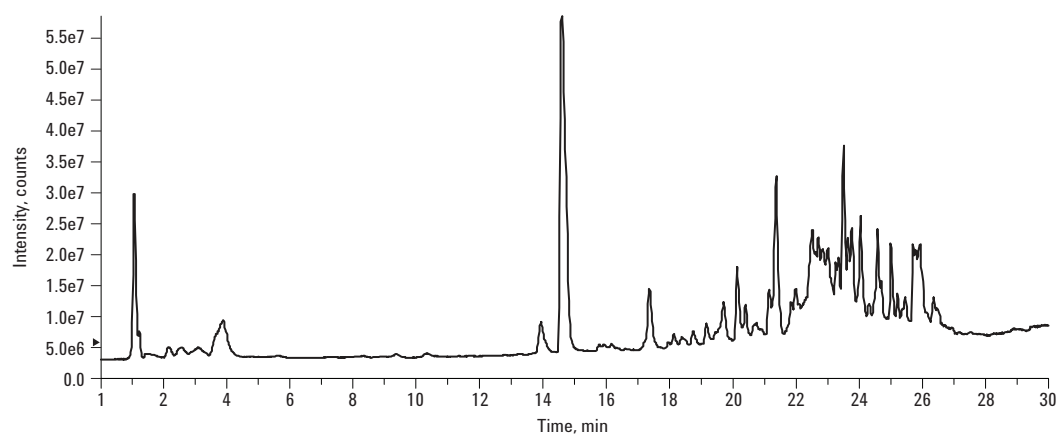
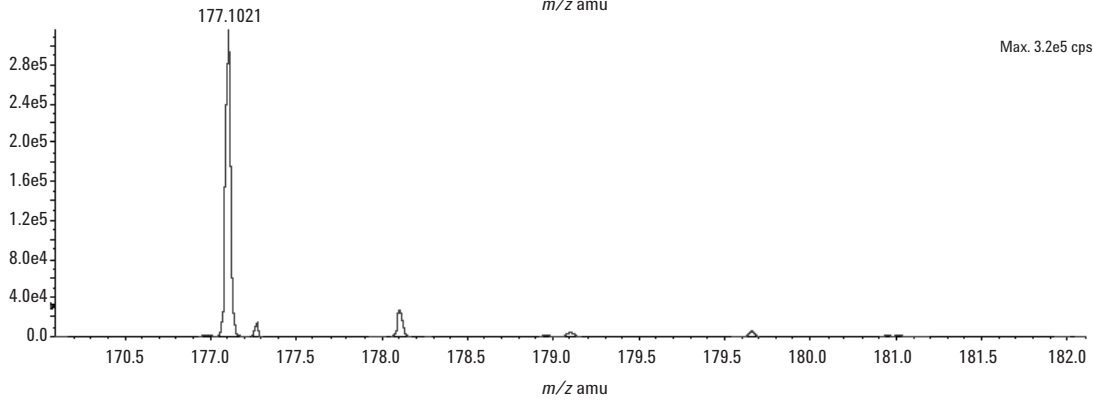
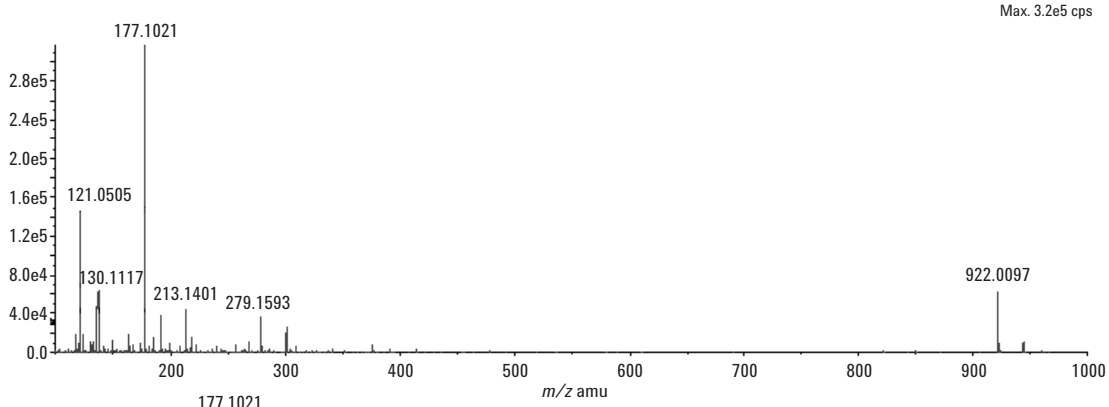
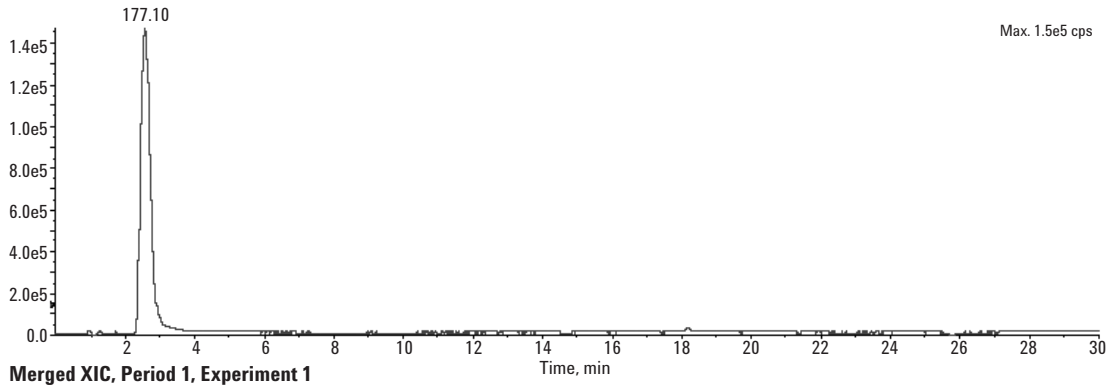


Figure 21. TIC of Coronial Sample 1.



| Formula | Compound name | Mass | Peak RT (min) | Peak area | Description | |
|--------------------|--------------------|-----------|---------------|-------------|-------------|-----------------------|
| $C_{10}H_{12}N_2O$ | (-)-cotinine | 176.09496 | 2.57 | 2.78259 E7 | - | |
| Species | Abundance (counts) | Ion mass | Measured mass | Error (mDa) | Error (ppm) | Ret. time error (min) |
| $[M + H]^+$ | 316885.14 | 177.10224 | 177.10215 | -0.00009 | -0.51 | 0.01 |

Figure 22. Excerpt from screener report of coronial Sample 1 showing confirmation of the presence of cotinine.

The results obtained from the analysis of Sample 1 suggest the deceased was a smoker, with the presence of both the nicotine and cotinine in the sample.

Sample 2

Sample 2 was known to contain citralopram, codeine, doxylamine, and tramadol from the previous analysis at the coronial office. In addition to the four previously reported compounds, screening with the database further showed the presence of

- Diazepam
- 6-acetylmorphine
- MDMA
- Methamphetamine
- Cotinine
- Meperidine
- nicotine

A TIC of the sample is shown in Figure 23. It could again be inferred that this patient was a smoker.

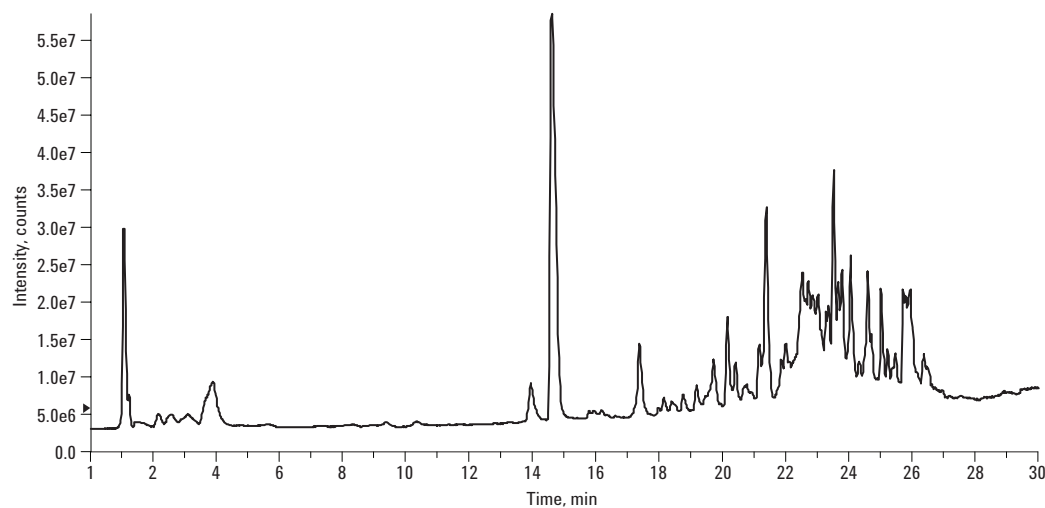


Figure 23. TIC of coronial Sample 2.

An excerpt of the screener report, in this instance for the confirmation of meperidine, is shown in Figure 24. Again, an excellent match to both retention time and mass can be seen.

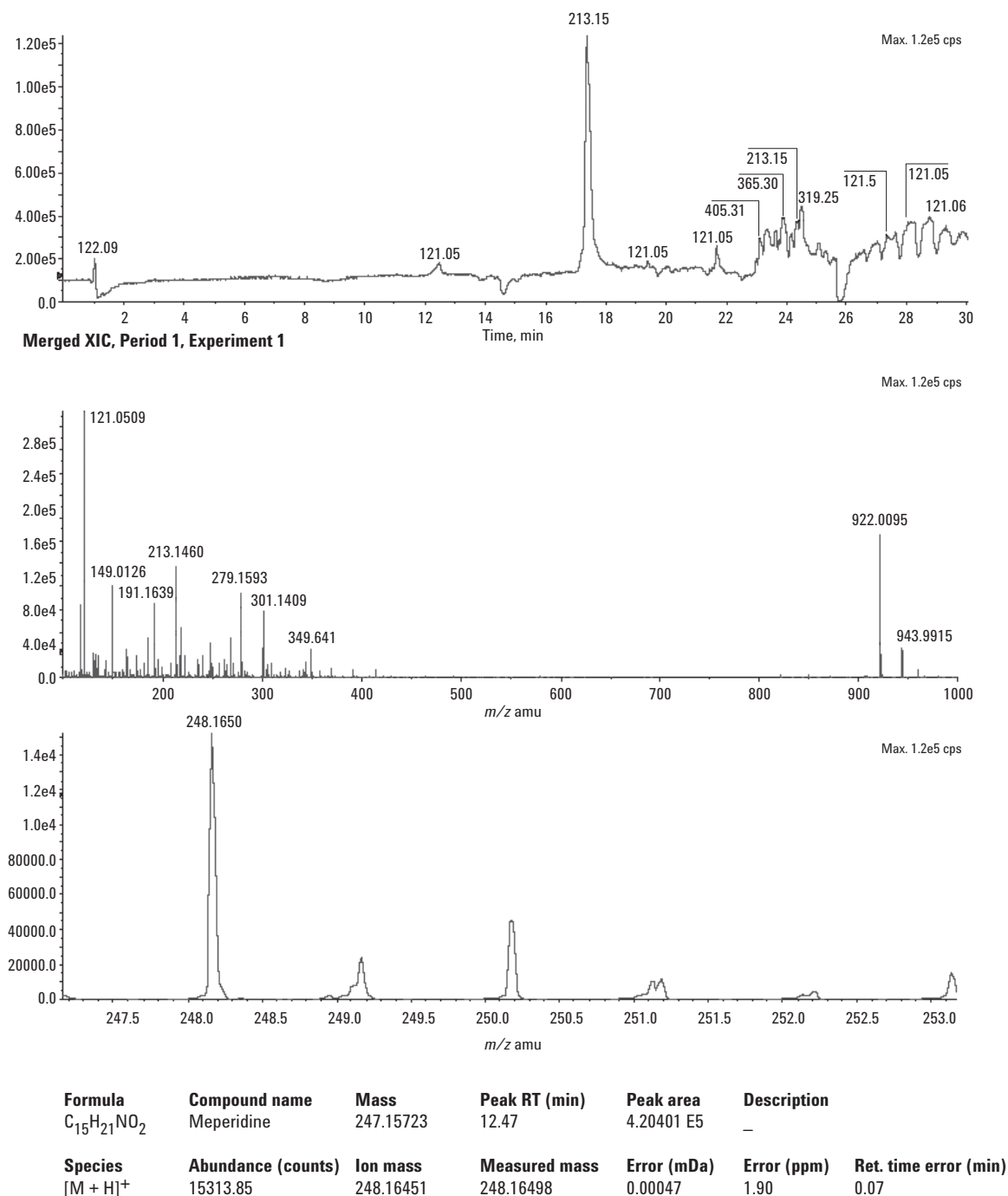


Figure 24. Excerpt from screener report indicating presence of meperidine.

Conclusions

The analysis of drugs of abuse is important in many different areas of our society, from law enforcement to medical monitoring. Current analytical techniques use a two-step screening and confirmation procedure to achieve the required specificity and sensitivity required. This application note has investigated 48 of the more common drugs of abuse and their applicability for determination through LC-TOF MS. It is not intended to be a comprehensive study of all possible components, but provides an excellent launching pad for the inclusion of the full gamut of possibilities

This application note shows the potential of the Agilent LC-TOF-MS as a single tool for both screening and confirmatory analysis, with quantitative information, often at levels below those currently analyzed for today. As a final example of the power of this technique, real-life coronial samples were evaluated under a screening protocol, with an additional seven components other than those previously reported by the coronial office detected.

References

1. Executive Office of the President – Office of National Drug Control Policy “Pulse Check – Drug Markets and Chronic Users in 25 of Americas Largest Cities” January 2004.
2. Kolbjorn Zahlse, Trond Aamo, and Jerry Zweigenbaum, “Screening Drugs of Abuse by LC/MS”, Agilent Technologies, publication 5989-1541EN www.agilent.com/chem
3. Kolbjorn Zahlse, Trond Aamo, and Jerry Zweigenbaum, “Therapeutic Drug Monitoring by LC/MSD – Clozapine, an example”, Agilent Technologies, publication 5989-1267EN www.agilent.com/chem
4. Jerry Zweigenbaum, “Quantitative Analysis of Opiates in Urine Using Accurate Mass LC/MSD TOF”, Agilent Technologies, publication 5989-2206EN www.agilent.com/chem

For More Information

For more information on our products and services, visit our Web site at www.agilent.com/chem.

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc. 2005

Printed in the USA
August 24, 2006
5989-3157EN

