

Software-assisted, high-throughput identification of main metabolites of pharmaceutical drugs from time-of-flight mass spectrometry data

Edgar Naegele, Agilent Technologies, Waldbronn, Germany



IMSC 2009
Poster PMM 093

Introduction

In modern pharmaceutical drug development it is of crucial importance to analyze the metabolism properties of possible new drug candidates as quickly as possible in order to make decisions about further investments in the development of a special compound. To find compounds with the correct properties it is essential to screen a large number of compounds for their metabolism properties, which requires to work in an high-throughput environment. This work describes the application of an ultra high performance liquid chromatography (UHPLC) device connected to a time-of-flight mass spectrometer and a special metabolite identification software for ultra fast, high throughput identification of main metabolites of new pharmaceutical drug candidates.

Experimental

Equipment:

- Agilent 1290 Infinity LC system consisting of 1290 Infinity Binary Pump with integrated degasser, 1290 Infinity High Performance Autosampler with thermostat, and 1290 Infinity Thermostatted Column compartment
 - Agilent 6530 Accurate-Mass Q-TOF LC/MS system
 - Agilent MassHunter Metabolite Identification (MetID) software
 - Column: ZORBAX SB-C18, 2.1 x 50 mm, 1.8 µm
- LC method:**
Solvent A: Water + 0.1 % formic acid
Solvent B: ACN + 0.1 % formic acid
Flow: 0.8 mL/min
Gradient 0 min, 5 %B; 0.10 min, 5 %B; 1.10 min, 75 %B. Stop time: 1.10 min. Post time: 1 min.
Injection: Volume 5 µL, sample cooler at 4°C, needle wash in 50 % methanol for 5 s, injection loop to bypass at 0.1 min with flush out factor 16
Column: Temperature 60 °C

TOF MS method:

Source: ESI positive. Capillary: 3500 V. Dry gas: 12 L/min. Nebulizer: 55 psi. Gas temp.: 350 °C. Skimmer: 65 V. Fragmentor: 200 V. Mass range: 100-1000 m/z. Acquisition rate: 5 spectra/s. Reference masses: 121.0508 and 922.0080.

Experimental

Sample preparation:

The following stock solutions were used:

- 20 mg/mL microsomal S9 preparation
- 0.1 mg/mL buspirone in water
- 1.6 mg NADP in 1.6 mL 0.1 M phosphate buffer, pH 7.4
- 50 mM isocitrate/MgCl₂ (203 mg MgCl₂·6H₂O + 258.1 mg isocitrate in 20 mL H₂O)
- Isocitrate dehydrogenase (IDH) 0.33 unit/µL
- NADPH regeneration system: 1.6 mL NADP solution + 1.6 mL Isocitrate solution + 100 µL IDH solution.
- Incubation mixture: 3.85 µL substrate + 200 µL NADPH regeneration system + 746.15 µL phosphate buffer + 50 µL S9.

Incubation was carried out at 37 °C for 60 minutes. A 100 µL aliquot was taken at the beginning (t=0) and at t=60 min. The reaction was stopped by adding 6 µL perchloric acid and 100 µL acetonitrile followed by centrifugation for 15 min at 14,000 rpm. The supernatant was evaporated to dryness using a SpeedVac concentrator and reconstituted with water containing 0.1 % formic acid for LC/MS analysis. The incubation sample stopped at 0 min was used as control.

Data analysis method in the MetID software:

The first step in the analysis comprised a comparison between the data file that contained the metabolite compounds (metabolite sample, t=60) and the data file that contained only the parent drug (control sample, t=0). All detectable mass signals were extracted from the MS level data using the Molecular Feature Extraction (MFE) algorithm. Related compound isotope masses and adduct masses were grouped together into discrete molecular features, and chemical noise was removed. The compounds lists of the metabolized sample and the control were then compared. All new compounds or those that increased twofold in the metabolized sample were considered potential metabolites and were subjected to further analysis by different algorithms. The algorithms can identify and qualify new metabolites, or just qualify metabolites found by another algorithm. In this high throughput experiment all algorithms' results were weighted equally and combined into a final identification relevance score. Metabolites were qualified when their final score was above the stringently defined relevance threshold. The results from all algorithms were collated in a results table, which could be inspected at-a-glance and reported.

Results and Discussion

After generation the data were loaded into the MetID software and analyzed using a common method. The result was displayed by the MetID software in an at-a-glance table, in which the result for each metabolite could be examined in more detail (figure 1). From the results table a summary report was generated, which showed the available information for each metabolite (figure 2). The more extensive report contained the detailed results for each metabolite. As example the result for a mono hydroxy metabolite (figures 3 to 5) and a dihydroxy metabolite (figures 6 to 8) of buspirone are discussed here.

| Name | Mass | RT | Rel. | Qual. | User | SC | IPM | EIC | MDF | Form. | BioXF |
|------------------|----------|------|--------|-------|------|----|-----|-----|-----|-------|-------|
| 2x Hydroxylation | 417.2379 | 0.59 | 100.00 | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ |
| Hydroxylation | 401.2423 | 0.63 | 100.00 | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ |
| Hydroxylation | 401.2424 | 0.66 | 100.00 | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ |
| 2x Hydroxylation | 417.2388 | 0.72 | 100.00 | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ |
| Hydroxylation | 401.2439 | 0.75 | 100.00 | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ |
| Hydroxylation | 401.2430 | 0.79 | 100.00 | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ |
| Buspirone | 385.2478 | 0.82 | - | - | - | - | ✓ | ✓ | ✓ | ✓ | - |
| Hydroxylation | 401.2429 | 0.84 | 75.00 | × | ✓ | × | ✓ | ✓ | ✓ | ✓ | ✓ |

Figure 2) Summary result report, including qualified metabolites sorted by their retention times (RT), with their metabolite names and relative score, molecular mass and the passed flag for individual algorithm results. SC=Sample-control comparison, IPM = Isotopic Pattern Matching, EIC = Extracted Ion Chromatogram, MDF = Mass Defect Filter, Form. = Calculated Formula, BioXF = Assigned Biotransformation, Qual. = Qualified, User = Qualified by User.

| Metabolite Information | | | | | | | | | | | |
|-----------------------------------|---|--------------------|--------------------|-------------------|----------------------|------------|--------|--|--|--|--|
| Name | Hydroxylation | BioXF Name | Hydroxylation | | | | | | | | |
| Formula | C ₂₁ H ₃₁ N ₅ O ₃ | Mass | 401.2439 | | | | | | | | |
| m/z | 402.2511 | Species | (M+H) ⁺ | | | | | | | | |
| RT | 0.754 | Sample Type | MetaboliteSample | | | | | | | | |
| MFE Compound Search | | | | | | | | | | | |
| Mass | m/z | Species | RT | Start Time | End Time | Volume | Height | | | | |
| 401.2439 | 402.2511 | (M+H) ⁺ | 0.754 | 0.739 | 0.774 | 192448 | 187344 | | | | |
| EIC Compound Search | | | | | | | | | | | |
| Mass | m/z | Species | RT | Start Time | End Time | Area | Area % | | | | |
| 401.2427 | 402.2500 | (M+H) ⁺ | 0.755 | 0.739 | 0.774 | 149323 | 100.00 | | | | |
| Sample Comparison Results | | | | | | | | | | | |
| Qualified | Changed | Resp. Ratio | Corr. RT | Normalized Height | | | | | | | |
| IS | Now | | | | | | | | | | |
| Isotopic Pattern Matching Results | | | | | | | | | | | |
| Qualified | Score | Delta m/z | | | | | | | | | |
| IS | 95.91 | 0.00 | | | | | | | | | |
| Mass Defect Filter Results | | | | | | | | | | | |
| Qualified | Delta Mass [mDa] | | | | | | | | | | |
| IS | -3.91 | | | | | | | | | | |
| Formula Results | | | | | | | | | | | |
| Assigned | Neutral Formula | Calc. Mass | Delta Mass [mDa] | Delta Mass [ppm] | Calculation Base | | | | | | |
| IS | C ₂₁ H ₃₁ N ₅ O ₃ | 401.2427 | -1.17 | -2.92 | MetCompoundMSpectrum | | | | | | |
| Biotransformation Results | | | | | | | | | | | |
| Assigned | Name | Phase | Offset Formula | Delta Mass [mDa] | Delta Mass [ppm] | Calc. Mass | | | | | |
| IS | Hydroxylation | I | +O | 1.17 | 2.92 | 401.2427 | | | | | |

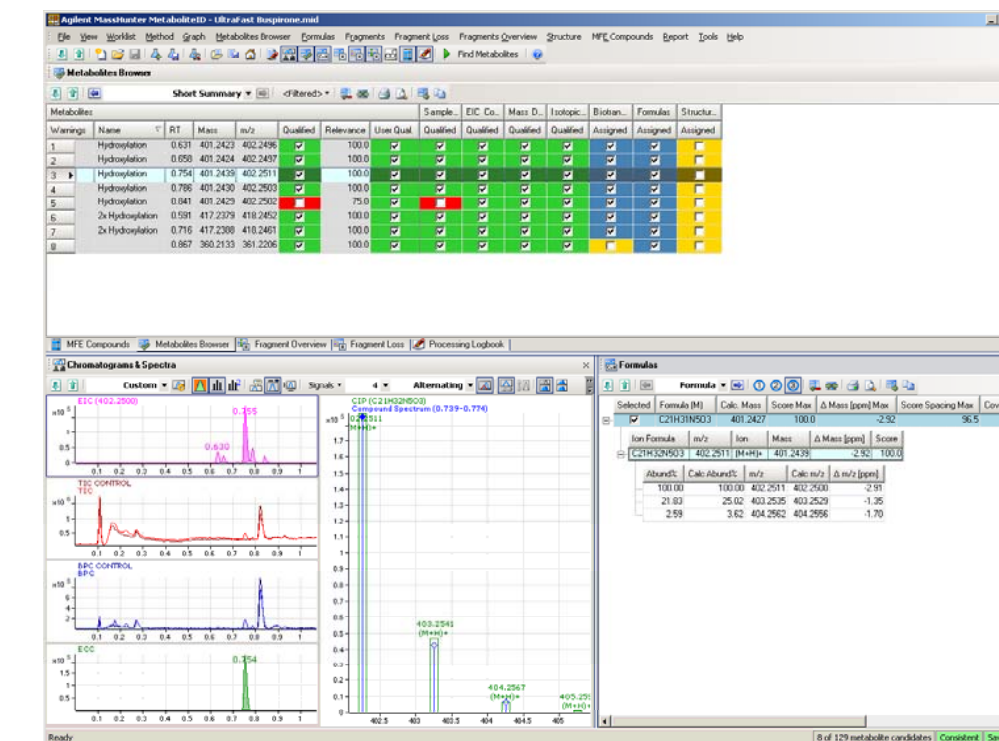


Figure 1) Result table showing an at-a-glance summary of buspirone metabolite analysis with overall identified metabolites, extracted ion chromatograms (EIC), extracted compound chromatograms (ECC), isotopic pattern analysis and calculated formulae.

| Metabolite Information | | | | | | | | | | | |
|---|---|---|---|--------------|------------|--------------|--|--|--|--|--|
| Name | Hydroxylation | BioXF Name | Hydroxylation | | | | | | | | |
| Formula | C ₂₁ H ₃₁ N ₅ O ₃ | Mass | 401.2439 | | | | | | | | |
| m/z | 402.2511 | RT | 0.754 | | | | | | | | |
| Formula Summary | | | | | | | | | | | |
| Selected | Score | Formula | Ion Formula | Mass | Calc. Mass | Δ Mass [ppm] | | | | | |
| TRUE | 100.0 | C ₂₁ H ₃₁ N ₅ O ₃ | C ₂₁ H ₃₂ N ₅ O ₃ | 401.2439 | 401.2427 | -2.92 | | | | | |
| Formula Details | | | | | | | | | | | |
| Formula (M) | Selected | | | | | | | | | | |
| C ₂₁ H ₃₁ N ₅ O ₃ | TRUE | | | | | | | | | | |
| Species | m/z | | | | | | | | | | |
| (M+H) ⁺ | 402.2511 | | | | | | | | | | |
| Formula Results | | | | | | | | | | | |
| Ion Formula | Score | Mass | Δ Mass [mDa] | Δ Mass [ppm] | DBE | | | | | | |
| C ₂₁ H ₃₂ N ₅ O ₃ | 100.0 | 401.2439 | -1.17 | -2.92 | 9 | | | | | | |
| Isotopic Peak Information | | | | | | | | | | | |
| Abund % | Calc Abund% | m/z | Calc m/z | Δ m/z [ppm] | | | | | | | |
| 100.00 | 100.00 | 402.2511 | 402.2500 | -2.91 | | | | | | | |
| 21.83 | 25.02 | 403.2535 | 403.2529 | -1.35 | | | | | | | |
| 2.59 | 3.62 | 404.2562 | 404.2556 | -1.70 | | | | | | | |

Figure 4) Detailed metabolite report about the formula including isotopic pattern, calculated for the buspirone hydroxy metabolite at retention time 0.75 min..

Figure 3) Detailed metabolite report for the buspirone hydroxy metabolite at retention time 0.75 min. This part of the report gives detailed information about the identified metabolite and the identifying algorithms. Other detailed information about formula (figure 4), chromatograms and isotopic pattern (figure 5) are also available.

Results and Discussion

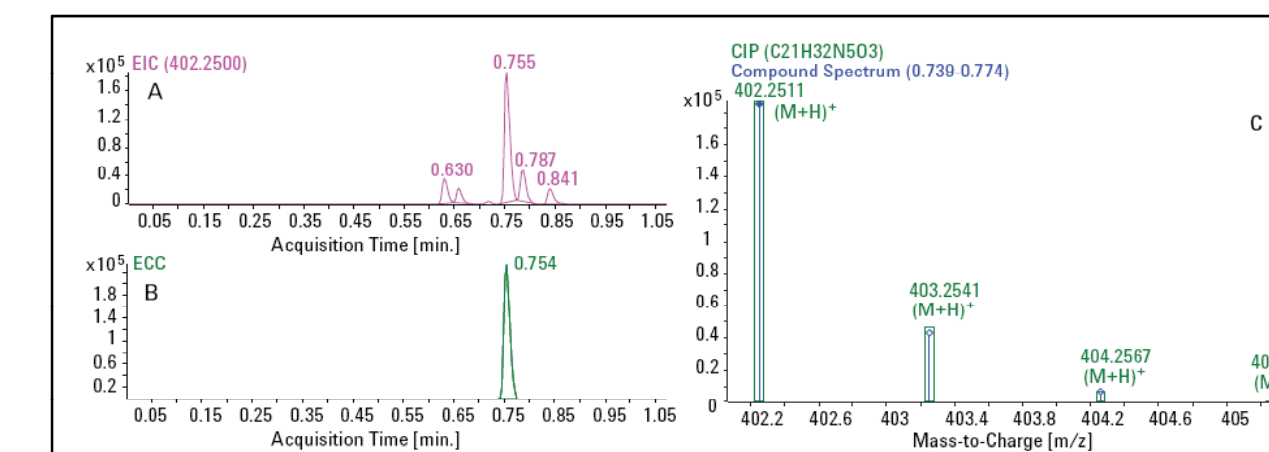


Figure 5) Detailed metabolite report for buspirone hydroxy metabolite at retention time 0.75 min: A) Extracted Ion Chromatograms (EIC) of compounds with mass 402.25 B) Extracted Compound Chromatogram (ECC) of buspirone hydroxy metabolite at retention time 0.75 min C) Measured isotopic pattern of buspirone hydroxy metabolite at retention time 0.75 min (blue lines) and calculated isotopic pattern (CIP, green box).

| Metabolite Information | | | | | | | | | | | |
|-----------------------------------|---|--------------------|--------------------|-------------------|----------------------|------------|--------|--|--|--|--|
| Name | 2x Hydroxylation | BioXF Name | 2x Hydroxylation | | | | | | | | |
| Formula | C ₂₁ H ₃₁ N ₅ O ₄ | Mass | 417.2388 | | | | | | | | |
| m/z | 418.2461 | Species | (M+H) ⁺ | | | | | | | | |
| RT | 0.716 | Sample Type | MetaboliteSample | | | | | | | | |
| MFE Compound Search | | | | | | | | | | | |
| Mass | m/z | Species | RT | Start Time | End Time | Volume | Height | | | | |
| 417.2388 | 418.2461 | (M+H) ⁺ | 0.716 | 0.700 | 0.726 | 3865 | 3869 | | | | |
| EIC Compound Search | | | | | | | | | | | |
| Mass | m/z | Species | RT | Start Time | End Time | Area | Area % | | | | |
| 417.2376 | 418.2449 | (M+H) ⁺ | 0.713 | 0.703 | 0.739 | 3483 | 100.00 | | | | |
| Sample Comparison Results | | | | | | | | | | | |
| Qualified | Changed | Resp. Ratio | Corr. RT | Normalized Height | | | | | | | |
| IS | Now | | | | | | | | | | |
| Isotopic Pattern Matching Results | | | | | | | | | | | |
| Qualified | Score | Delta m/z | | | | | | | | | |
| IS | 91.50 | 0.00 | | | | | | | | | |
| Mass Defect Filter Results | | | | | | | | | | | |
| Qualified | Delta Mass [mDa] | | | | | | | | | | |
| IS | -6.37 | | | | | | | | | | |
| Formula Results | | | | | | | | | | | |
| Assigned | Neutral Formula | Calc. Mass | Delta Mass [mDa] | Delta Mass [ppm] | Calculation Base | | | | | | |
| IS | C ₂₁ H ₃₁ N ₅ O ₄ | 417.2376 | -1.20 | -2.87 | MetCompoundMSpectrum | | | | | | |
| Biotransformation Results | | | | | | | | | | | |
| Assigned | Name | Phase | Offset Formula | Delta Mass [mDa] | Delta Mass [ppm] | Calc. Mass | | | | | |
| IS | 2x Hydroxylation | I | +O2 | 1.20 | 2.87 | 417.2376 | | | | | |

Figure 6) Detailed metabolite report for dihydroxy metabolite of buspirone at retention time 0.71 min. This part of the report gives detailed information about the identified metabolite and the identifying algorithms. Other detailed information about formula (see figure 7), chromatograms and isotopic pattern (see figure 8) are also available.

Conclusions

This work demonstrates the use of the Agilent 1290 Infinity LC system with the Agilent 6530 accurate mass Q-TOF LC/MS system for fast separation and accurate mass measurement of compounds in an in-vitro metabolite sample under high-throughput conditions. The metabolite compounds were separated in a run time below one minute and the width of the peaks extracted by the Metabolite ID software were below one second (FWHH). The major metabolites were identified quickly by means of the Agilent Metabolite identification software. A summary report as well as detailed reports for each metabolite were generated.

| Metabolite Information | | | | | | | | | | | |
|---|---|---|---|--------------|------------|--------------|--|--|--|--|--|
| Name | 2x Hydroxylation | BioXF Name | 2x Hydroxylation | | | | | | | | |
| Formula | C ₂₁ H ₃₁ N ₅ O ₄ | Mass | 417.2388 | | | | | | | | |
| m/z | 418.2461 | RT | 0.716 | | | | | | | | |
| Formula Summary | | | | | | | | | | | |
| Selected | Score | Formula | Ion Formula | Mass | Calc. Mass | Δ Mass [ppm] | | | | | |
| TRUE | 100.0 | C ₂₁ H ₃₁ N ₅ O ₄ | C ₂₁ H ₃₂ N ₅ O ₄ | 417.2388 | 417.2376 | -2.87 | | | | | |
| Formula Details | | | | | | | | | | | |
| Formula (M) | Selected | | | | | | | | | | |
| C ₂₁ H ₃₁ N ₅ O ₄ | TRUE | | | | | | | | | | |
| Species | m/z | | | | | | | | | | |
| (M+H) ⁺ | 418.2461 | | | | | | | | | | |
| Formula Results | | | | | | | | | | | |
| Ion Formula | Score | Mass | Δ Mass [mDa] | Δ Mass [ppm] | DBE | | | | | | |
| C ₂₁ H ₃₂ N ₅ O ₄ | 100.0 | 417.2388 | -1.20 | -2.87 | 9 | | | | | | |
| Isotopic Peak Information | | | | | | | | | | | |
| Abund % | Calc Abund% | m/z | Calc m/z | Δ m/z [ppm] | | | | | | | |
| 100.00 | 100.00 | 418.2461 | 418.2449 | -2.86 | | | | | | | |
| 23.90 | 25.06 | 419.2488 | 419.2478 | -2.35 | | | | | | | |

Figure 7) Detailed metabolite report about the formula, including isotopic pattern, calculated for dihydroxy metabolite of buspirone at retention time 0.71 min.

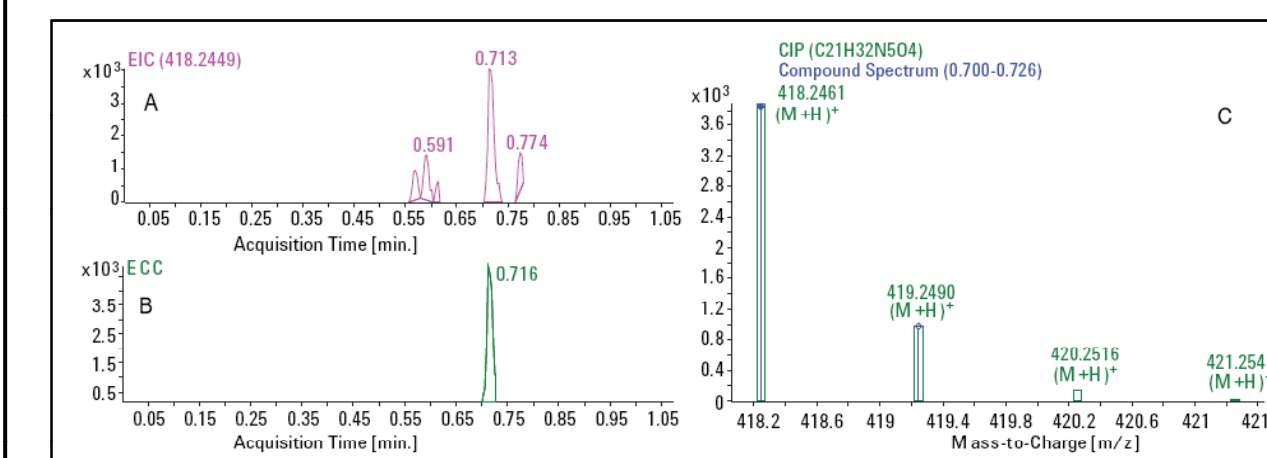


Figure 8) Detailed metabolite report for dihydroxy metabolite of buspirone at retention time 0.71 min: A) Extracted Ion Chromatograms (EIC) of compounds with mass 418.24 B) Extracted Compound Chromatogram (ECC) of dihydroxy metabolite of buspirone at retention time 0.71 min C) Measured and calculated isotopic pattern of dihydroxy buspirone metabolite at retention time 0.71 min