

Software-assisted, high-throughput identification of main metabolites of pharmaceutical drugs

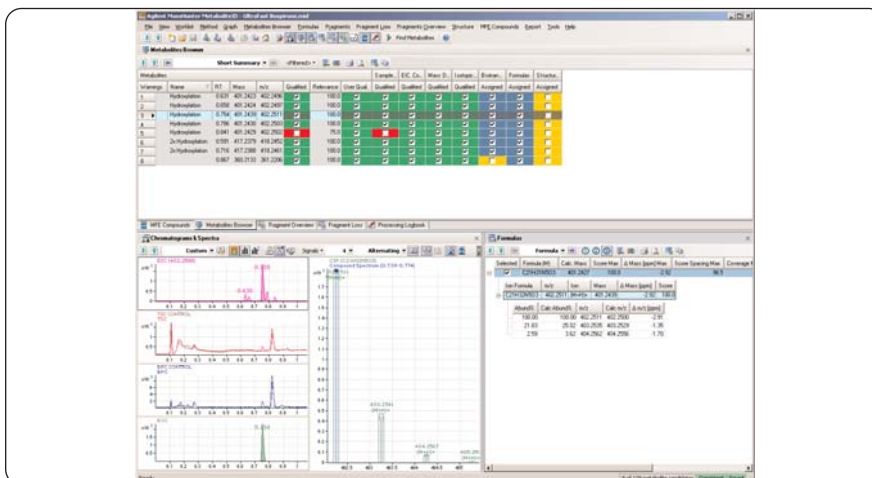
Rapid data acquisition by Agilent 1290 Infinity LC, TOF and Q-TOF instrumentation, and subsequent identification of metabolites by Agilent MassHunter Metabolite Identification software

Application Note

Metabolite identification in drug discovery and drug development

Author

Edgar Naegele
Agilent Technologies
Waldbronn, Germany



Abstract

This Application Note describes:

- Rapid separation of metabolites generated from in-vitro experiments using the Agilent 1290 Infinity LC, system
- Fast acquisition of TOF mass spectra using Agilent 6530 Accurate-Mass Quadrupole Time-of-Flight LC/MS systems
- Fast, software-assisted identification of main metabolites from in-vitro experiments using Agilent MassHunter Metabolite Identification software
- Generation of reports for the identified metabolites using Agilent MassHunter software



Agilent Technologies

Introduction

In modern pharmaceutical drug development it is of crucial importance to analyze the adsorption, distribution, metabolism and excretion (ADME) 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 ADME properties, which requires to work in an high-throughput environment. This Application Note describes the application of the Agilent 1290 Infinity LC system, the Agilent 6530 Q-TOF MS system and the MassHunter Metabolite Identification software for fast, high-throughput identification of main metabolites of new pharmaceutical drug candidate compounds.

Experimental

Equipment

- Agilent 1290 Infinity LC system consisting of 1290 Infinity Binary Pump with integrated degasser, 1290 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 μ m

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 0.33 unit/ μ L

NADPH regeneration system: 1.6 mL NADP solution + 1.6 mL Isocitrate solution + 100 μ L IDH solution.

Incubation mixture: 3.85 μ L substrate + 200 μ L NADPH regeneration system + 746.15 μ L phosphate buffer + 50 μ L S9.

Incubation was carried out at 37 °C for 60 minutes. A 100 μ L aliquot was taken at the beginning (t=0) and at t=60 min. The reaction was stopped by adding 6 μ L perchloric acid and 100 μ L 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.

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: 1.1.0 min
Post time: 1 min.
Injection: Volume 5 μ L, 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

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) and the data file that contained only the parent drug (control sample). 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 reported¹.

Results and discussion

To achieve fast separation of the metabolites on a 50 mm, 1.8 μ m particle size column, a 1 minute gradient was applied by the Agilent 1290 Infinity LC system. The metabolites were generated from the pharmaceutical test compound buspirone in an in-vitro assay. For adequate detection with the time-of-flight mass spectrometer the instrument was operated at a data rate of 5 Hz.

After generation the data was 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-hydroxyl metabolite (figures 3 to 5) and a dihydroxy metabolite (figures 6 to 8) of buspirone are discussed here.

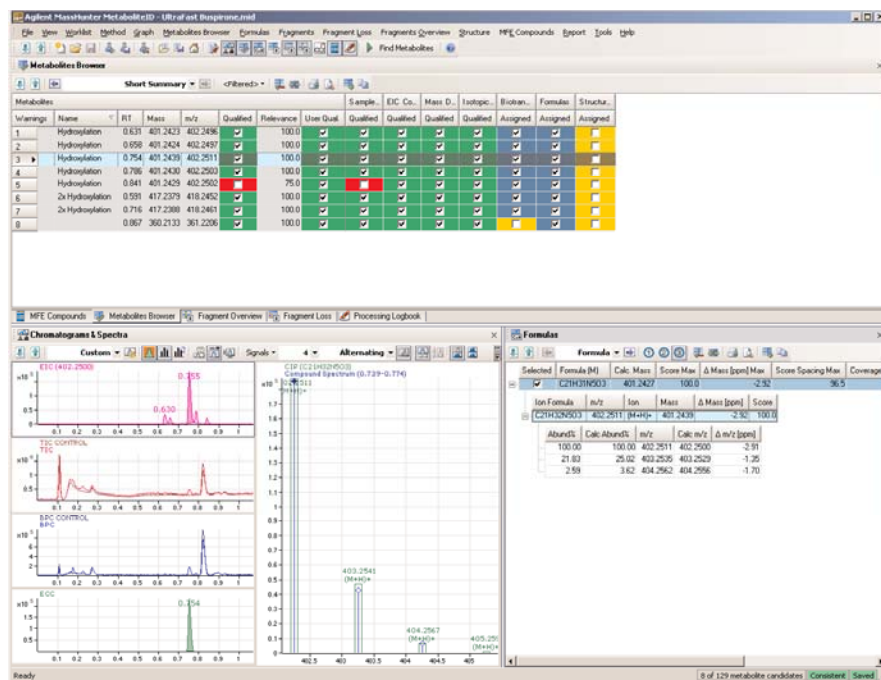


Figure 1
Result table showing an at-a-glance summary of buspirone metabolite analysis with overall identified metabolites, extracted ion chromatograms (EIC), extracted compound chromatograms (ECC), isotopic pattern analysis and calculated formulas.

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 by Score, User = Qualified by User.

The calculation of the formula was outlined in the detailed formula report (figure 7).

The EIC of m/z 418.24 showed about five significant peaks for possible dihydroxylated metabolites of buspirone with the selected peak at 0.71 minutes (figure 7A). The ECC showed the extracted MFE compound for the molecular mass of 417.2388 at retention time 0.71 identical to the EIC (figure 7B). The measured and calculated isotopic pattern of this compound is shown in figure 7C.

Metabolite Information						
Name	2x Hydroxylation		BioXF Name	2x Hydroxylation		
Formula	C21H31N5O4		Mass	417.2388		
m/z	418.2461		RT	0.716		
<u>Formula Summary</u>						
Selected	Score	Formula	Ion Formula	Mass	Calc. Mass	Δ Mass [ppm]
TRUE	100.0	C21H31N5O4	C21H32N5O4	417.2388	417.2376	-2.87
<u>Formula Details</u>						
Formula (M)	Selected					
C21H31N5O4	TRUE					
Species	m/z					
(M+H)+	418.2461					
Formula Results						
Ion Formula	Score	Mass	Δ Mass [mDa]	Δ Mass [ppm]	DBE	
C21H32N5O4	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.

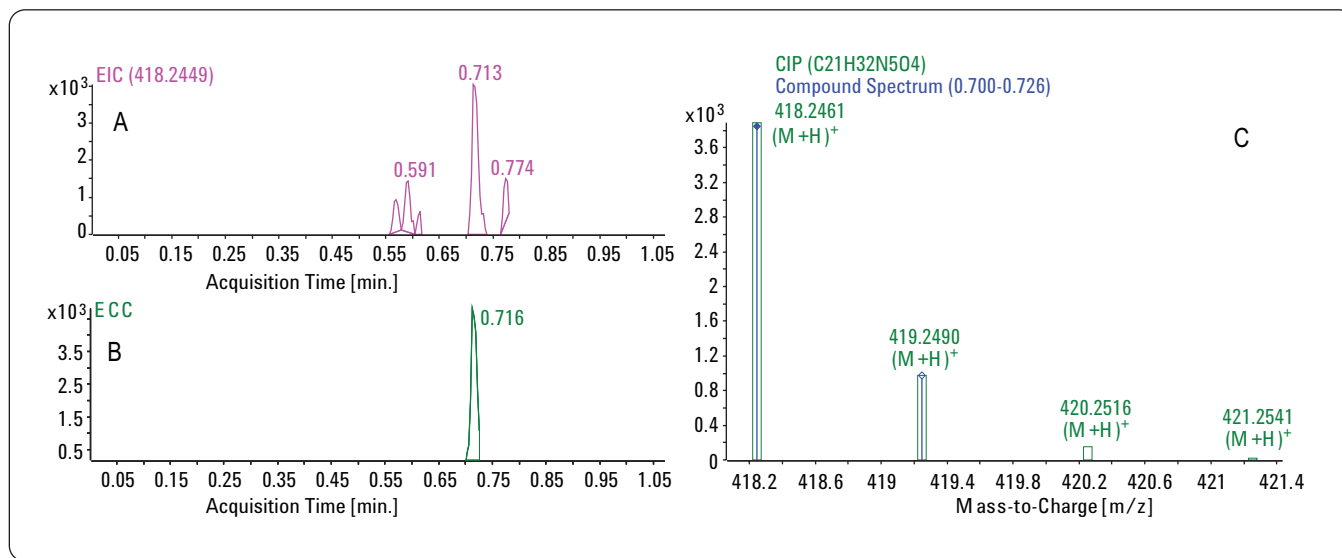


Figure 8
Detailed metabolite report for dihydroxy metabolite buspirone at retention time 0.71 min:
A) Extracted Ion Chromatograms (EIC) of compounds with mass 418.24
B) Extracted Compound Chromatogram (ECC) of dihydroxy metabolite of buspirone at retention time 0.71 min
C) Measured and calculated isotopic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

Conclusion

This Application Note demonstrated the use of the Agilent 1290 Infinity LC system with an Agilent 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.

References

1. E. Naegele, F. Wolf, U. Nassal, R. Jäger, H. Lehmann, F. Kuhlmann, K. Subramanian, "An interwoven, multi-algorithm approach for computerassisted identification of drug metabolites", *Agilent Technologies Application Note, publication number 5989-7375EN*, **2007**.

www.agilent.com/chem/metid

For Research Use Only. Not for use in
diagnostic procedures.

This information is subject to change without
notice.

© Agilent Technologies, Inc., 2009
Published May 1, 2009
Publication Number 5989-9924EN



Agilent Technologies