Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes



Orbitrap quantitation: lab of the future

Close your uncertainty gap with the selectivity and confidence of HRAM





Introduction

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Thermo Scientific[™] Orbitrap[™] mass spectrometer technology is disrupting the use of mass spectrometry (MS) for qualitative, and more importantly for quantitative analyses, in applied markets. Why is this occurring?

For many years, liquid chromatography-triple quadrupole mass spectrometry (LC-QQQ MS) technology has been the standard for quantitation of many types of compounds, from pesticide residues in food and feed, to testosterone in plasma and urine for research applications. Due to their selectivity and sensitivity, quadrupole-based mass analyzers provide reliable, cost-efficient quantitation and identification of target analytes.

For targeted analyses, nominal-mass-resolution quadrupole-based instruments are operated in either the selected ion monitoring (SIM) or selected reaction monitoring (SRM) data acquisition modes. During SIM and SRM data acquisition, the scope of analysis is limited to the target ions selected and therefore suffers from non-detection of analytes that are present. Undetected compounds and unknown metabolites can be a source of uncertainty and concern.

Set-up of QQQ instruments require careful tuning and optimization of data acquisition parameters for each target analyte and ongoing monitoring of data acquisition time windows to ensure analyte detection. These tasks require time and effort, and limit the number of possible measurements that can be made during an analytical run.

Time-of-flight (TOF) technology has also been used for quantitation, but its dynamic range is limited due to detector saturation effects. Moreover as shown in Figure 1, TOF instruments do not have adequate resolution, especially at the lower masses measured in small molecule analyses, to provide adequate selectively. Poor selectivity often results in false identification (false negatives and false positives) of an analyte when it is not present and the inability to quantitate the analytes of interest.¹

1. Rajski, Ł.; Gómez-Ramos Mdel, M.; Fernández-Alba, A. R. Large pesticide multiresidue screening method by liquid chromatography-Orbitrap mass spectrometry in full scan mode applied to fruit and vegetables. J. Chromatogr. A. 2014, Sep 19;1360:119-27.



Quantitation: historical perspective



Introduction

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

High resolving power: quantitative capability for the lab of the future

Included in approximately 5000 peer-reviewed publications, Orbitrap mass spectrometer technology is well established as a powerful LC-MS platform. Its high-resolution, accurate-mass (HRAM) measurements, together with its high dynamic range, have enabled rigorous characterization of complex mixtures.²

Why is high resolution important? High resolving power is particularly important for experiments involving complex mixtures, such as biological, environmental, and food samples generated from a matrix, since these will contain a significant number of background ions in addition to the analytes of interest. In such cases, high resolving power will make the difference between detecting and not detecting analytes at low concentrations due to the masking effect of isobaric interferences. In other words, accurate quantitation relies on high selectivity, which is the ability to resolve compounds of interest from background interferences.

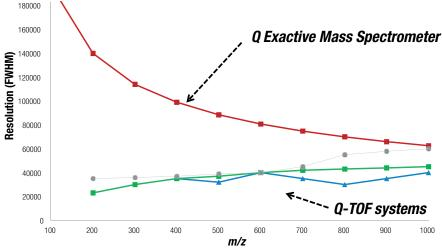


Figure 1. Comparison of Q Exactive mass spectrometer and Q-TOF resolution.

2. The Thermo Scientific Exactive Mass Spectrometer. The ultimate screening machine and overview of scientific literature. Thermo Scientific Literature Review: XX63548_E 06/12M



Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

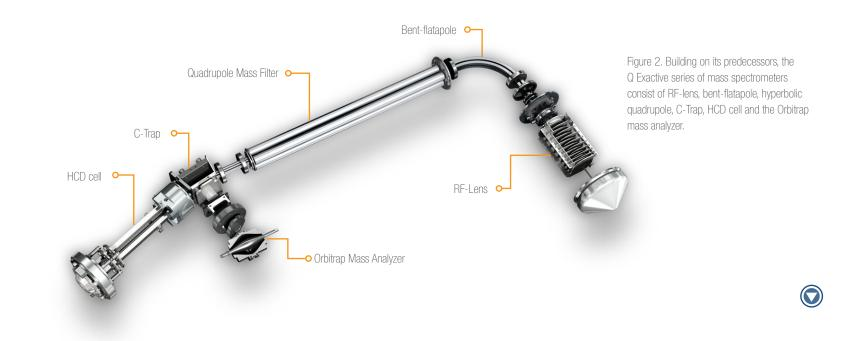
Conclusion

Application Notes

Orbitrap mass spectrometer technology has revolutionized the field of mass spectrometry. The <u>Thermo Scientific</u>[™]Q Exactive[™] series of quadrupole-Orbitrap mass spectrometers are the next step in that revolution, combining high-performance quadrupole precursor selection with HRAM detection. The Thermo Scientific Q Exactive mass spectrometer and <u>Thermo Scientific Q Exactive Focus mass spectrometer</u> are benchtop LC-MS systems composed of six unique sections: RF-lens, bent-flatapole, hyperbolic quadrupole mass filter, C-Trap, HCD cell, and the Orbitrap mass analyzer (Figure 2).

lons generated in the heated electrospray (H-ESI II) ion source enter the RF-lens through a heated capillary. The RF-lens acts as an ion funnel that focuses and transmits ions to the "bent-flatapole" where they are redirected by 90 degrees. This reduces instrument footprint and allows neutrals to be removed before the ions stream into the mass analyzer. Next, ions enter the hyperbolic quadrupole, which allows for the selection of ions in a broad or narrow mass window down to 0.4 amu. Ions then are captured in the C-Trap, an ion storage device where the user can choose to (1) send ions to the Orbitrap mass analyzer for full-scan detection (MS-level) or (2) fragment the ions inside the nitrogen-filled higher collisional dissociation (HCD) cell, send them back to the C-Trap, and then to the Orbitrap mass analyzer for fragment (MS/MS) detection. Mass spectra are acquired via image current detection. The vacuum inside the Orbitrap mass analyzer is maintained within the 10-10 mbar range.

Q Exactive series of mass spectrometers







Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Comprised of two outer electrodes and a central electrode, the Orbitrap mass analyzer acts both as an analyzer and a detector (Figure 3). When ions enter the Orbitrap analyzer they are captured via "electrodynamic squeezing." Once captured, the ions oscillate around the central electrode and in between the two outer electrodes. Different ions oscillate at different frequencies resulting in mass separation. Mass spectra are acquired via image current detection by measuring the oscillation frequencies induced by ions on the outer electrodes. In this way, the Orbitrap mass analyzer

is a Fourier Transform (FT) mass analyzer analog of well-known FT- ion cyclotron resonance (ICR) technology, with the advantage of allowing smaller instrument size and easier instrument operation.

Orbitrap mass analyzer

The Orbitrap mass analyzer's uniquely high resolution is due to three factors: 1) nanometer-range accuracy Orbitrap electrodes, 2) high voltage supplies, and 3) mass-to-charge measurements delivered as a function of oscillation frequency.

Figure 3. Orbitrap mass analyzer technology provides uniquely high resolution.







Q Exactive series of mass spectrometers

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Close your uncertainty gap: identify, quantitate, and confirm with selectivity and confidence

Orbitrap mass spectrometer technology is rapidly growing in popularity because it uniquely provides quantitative and qualitative (Quan/Qual) HRAM analytical capability in a <u>single platform</u> and often in a single analytical run (Table 1). Accurate mass assignment and high resolution in both MS and MS/MS modes increase selectivity and reduce uncertainty,

benefiting most applications imaginable. Full-scan HRAM data acquisition captures all sample data, enabling identification of "unexpected" compounds and retrospective data analysis without need to re-run samples.

Application/Need	HRAM	QQQ
High Sensitivity	•	•
Confirm Analyte Structure	•	•
Analysis of "Unknowns"	•	•
Retrospective Search for New Compounds	•	•
Low Cost	•	•
Established Methods	•	•

Table 1. Comparison of HRAM and QQQ mass spectrometer capabilities.



Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Q Exactive mass spectrometers can be operated in several scan modes, including: full-scan MS (full MS), all-ion fragmentation (AIF) in the HCD collision cell, SIM for single ion detection, and parallel reaction monitoring (PRM) for detecting all fragment ions in parallel. Fragments generated in-source or in the multipole collision cell by AIF provide additional information that complements full-scan data.

AlF fragments all ions in a single fragmentation event without precursor ion isolation and records all fragment ions in a single mixed spectrum. MS/MS (MS²) data can be collected using data-dependent acquisition (DDA) where precursor ions selected using predetermined rules are subjected to MS/MS, or by using data-independent acquisition (DIA)

where a full scan is followed by a number of wide isolation MS² scans, which together cover the same isolation range as the preceding full scan. The benefit of DIA is the ability to capture both all data and MS² HR data scans without need for additional sample injections, enabling the acquisition of full information for all compounds and their fragments.

The ability to perform fast polarity switching allows acquisition of both positive and negative scans within a single acquisition method, eliminating the need for extra sample injections, which is of particular importance when only limited amounts of sample are available. Positive and negative polarity switching can be performed on the chromatographic timescale and does not require instrument recalibration.

Powerful quantitative capability



Powerful quantitative capability

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Sunitinib

100

0.0125 to 5,000 ng/mL

Concentration range in plasma extract:

Area: 15,121 0.0125ng/mL m/z 399.2191

MEW= 10ppm

ESI +- HR full scan

m/z 200 to 1,000 R: 70,000 at mz 200 Area

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

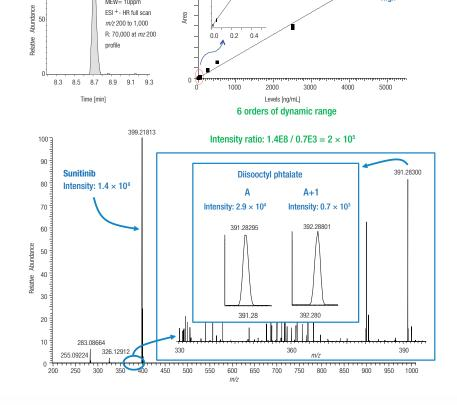
Pharmaceutical Discovery

Conclusion

Application Notes

The Q Exactive series of mass spectrometers provide up to six orders of magnitude of linear quantitative dynamic range (Figure 4) and up to five orders of intra-scan dynamic range (Figure 5) for accurate quantitation in low-level components in complex matrices.* In addition, these mass spectrometers are easily calibrated in just a few minutes and maintain that calibration for days at a time.

Hiah



0.2 04

0.0

Figure 4. The Q Exactive mass spectrometer provides up to six orders of linear dynamic range in high-resolution full-scan acquisition mode for Sunitinib. The concentration range in plasma extract was 0.0125 to 5,000 ng/mL. MEW refers to mass extraction window.

Figure 5. The Q Exactive series of mass spectrometers provide up to five orders of intra scan dynamic range for Sunitinib. MA refers to mass accuracy.

* Courtesy of Dr. Bertrand Rochat (ASMS 2015), CHUV, Lausanne, Switzerland.





Powerful quantitative capability

Choice of quantitation experiments

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Using the Q Exactive mass spectrometer's powerful modes of operation, there are generally three types of targeted quantitation experiments, and one type of non-targeted quantitation experiment, that are performed.

Targeted quantitation

1) Full-scan MS with data-dependent tandem mass spectrometry (full MS/ddMS² with inclusion list). In this workflow, the user defines a list of target parent ions and a retention time window for each. During the analysis, data acquisition is performed in full-scan mode. If any of the parent ions in the target list are detected within the defined retention time window, the instrument automatically performs an MS/MS scan, and then resumes full-scan operation. Data obtained in the full-scan mode are used for detection and quantitation, whereas data obtained from the MS/MS scans are used for identification with confirmation. Characteristics of the full MS/ddMS² quantitation workflow include:

- Peak areas of the extracted ion chromatograms of the parent ions of interest are used for quantitation
- Selectivity is based on high resolution
- Enables multi-residue methods in complex matrices
- Minimal method development required for food safety, clinical research, forensic toxicology, or environmental applications

- 2) Targeted SIM with data-dependent tandem mass spectrometry (t-SIM/ddMS²). Characteristics include:
 - Peak areas of the extracted ion chromatograms of the parent ions of interest are used for quantitation
 - Scheduled target list (retention time, m/z)
 - Most useful where ultimate sensitivity is needed
 - Minimum MS method development needed (the user must define parent ions and mass isolation width)
- 3) Parallel reaction monitoring (PRM). Characteristics include:
 - Peak areas of the extracted ion chromatograms of the precursor ion selected fragment ions are used for quantitation
 - Post-acquisition processing of the fragment extracted ion chromatograms because all fragments are recorded
 - Scheduled target list (retention time, *m/z*, collision energy)
 - Provides the most sensitivity and selectivity for quantitation of compounds in very complex matrices

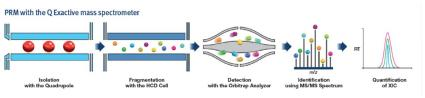


Figure 6. The Q Exactive analyzer provides parallel detection of all target product ions in one coordinated HRAM analysis.







Powerful quantitative capability

Choice of quantitation experiments

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Non-targeted quantitation

Full-scan MS with DIA delivers reproducible quantitation with complete full-scan MS/MS qualitative information for unknown analysis. In the DIA approach, a number of fragmentation events with wide isolation windows are used to cover the entire range of precursor ions of interest. Compared with AIF, DIA can improve selectivity because the product ions are derived from a smaller range of precursor ions. Furthermore, it improves the sensitivity of detection of fragment ions because a higher number of analyte precursor ions are delivered to the C-Trap. The DIA approach allows for absolute quantitation through the use of reference standards, as well as relative quantitation when using <u>Thermo Scientific</u>[™] <u>TraceFinder</u>[™] <u>software</u> to compare different samples from one data set in one workflow. If needed, this workflow allows general unknown screening as an extension to analyses done previously with several options for suspect identification.



Growing body of publications

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Q Exactive mass spectrometers are now used extensively for the quantitation of small molecule analytes in a broad range of clinical research, forensic toxicology, pharmaceutical discovery, environmental analysis, and food testing applications. The result is a steadily-increasing collection of peer-reviewed publications highlighting the benefits of Orbitrap mass spectrometer-based quantitation (Figure 6). In the overview of scientific peer-review publications that follows, three important themes emerge:

 Easier environmental method development: Compared to QQQ MS multi-residue methods that involve optimization of hundreds of SIM or SRM transitions in numerous time windows, Orbitrap mass spectrometer-based quantitation methods do not require special tuning and optimization nor comprehensive MS knowledge to set up.

- Greater confidence due to less possibility of false detection of the analytes of interest: The higher resolving power provided by the Orbitrap mass analyzer of 70,000 (FWHM) at *m/z* 200 or greater, provides better analytical selectivity for low molecular weight compounds in complex matrices.
- Unique ability to detect non-target compounds and to analyze data retrospectively, without re-running samples: Full-scan HRAM data can be acquired across the full mass range of interest.

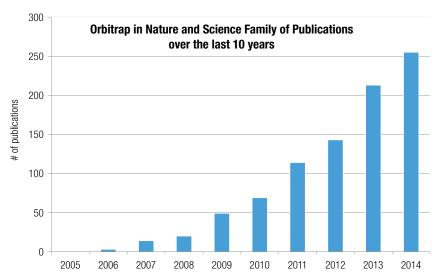


Figure 6. Growing collection of scientific peer-reviewed publications highlighting Orbitrap mass spectrometerbased quantitation.



Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes



Food safety testing

- Targeted screening and quantitation of food contaminants: workflow
- Pesticides in fruits and vegetables
- Synthetic hormones and other veterinary drugs in animal products
- Mycotoxins in dairy products
- Multiclass contaminants in food
- Marine biotoxins
- Dyes in wine
- Herbal medicines and dietary supplements





Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

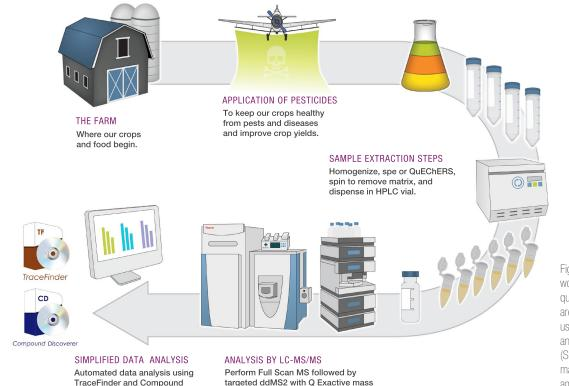
Application Notes

Targeted screening and quantitation of food contaminants: workflow

Ensuring the safety of food supplies is of utmost importance for public health and safety. Screening and quantitating foodstuffs for known and unknown contaminants, such as pesticides, herbicides, antibiotics, and adulterants in complex matrices must be fast and simple. Orbitrap mass spectrometer-based HRAM analysis is ideally suited for screening and quantitating hundreds or thousands of contaminants in a single analysis. These analyses are performed using variations of the workflow shown in Figure 7.

Didyouknow?

Screening, identification, and quantiation of food contaminants is faster and more certain with Thermo Scientific[™] High-Resolution Accurate-Mass MS/MS Spectral Libraries. To learn more: <u>click here</u>



spectrometer or perform Full Scan High

Speed Screening with Exactive Plus,

Discoverer software.

Figure 7. Q Exactive mass spectrometer workflow for targeted screening and quantitation of food contaminants. Samples are processed and contaminants are extracted using various methods such as QuEChERS and online and offline solid phase extraction (SPE), and then are analyzed on the Q Exactive mass spectrometer. Data is scrutinized and the amount of contaminate is determined using automated software.



Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Del Mar Gómez-Ramos et al. applied the Q Exactive mass spectrometer to LC-MS analysis of 139 pesticide residues in QuEChERS extracts of tomato, pepper, orange and green tea.³ An analysis of 100 real samples was performed to evaluate the mass spectrometer's identification and quantitation capability and the results were compared to a QQQ MS/MS analysis of the same samples. When the authors analyzed the samples using QQQ MS/MS, the results were consistent with those produced by the Q Exactive mass spectrometer—the same pesticides were found and false positives and false negatives were not reported. The authors noted that compared to the QQQ instrument, the Q Exactive mass spectrometer increased selectivity and, in full-scan mode, permitted retrospective analysis of data.

Pesticides in fruits and vegetables

Wang et al. applied the Q Exactive mass spectrometer to LC-MS analysis of 166 pesticide residues in fruits and vegetables.⁴ Pesticides were extracted using the QuEChERS method. Full-scan data (full MS) were used for quantitation. Data-dependent MS/MS (ddMS²) product-ion spectra were used for confirmation. Matrix-matched standard calibration curves with isotopically labeled standards, or chemical analogues as internal standards, were used for quantitation. The authors evaluated overall recovery, intermediate precision, and measurement uncertainty. Approximately 90.3 to 91.5% of the pesticides had good recoveries between 81 and 110%, 92.1 to 97.6% had intermediate precision of \leq 20%, and 89.7 to 95.2% had measurement uncertainty of \leq 40%. Confirmation of targets was based on a mass accuracy \leq 5 ppm and LC retention time tolerance within ±2.5%. The authors concluded that for quantitation and confirmation of pesticide residues in fresh fruits and vegetables, the Q Exactive mass spectrometer demonstrated good performance.

• Attain selectivity surpassing QQQ instruments

• Analyze full-scan data retrospectively

3. Del Mar Gómez-Ramos, M.; Rajski, Ł.; Heinzen, H.; Fernández-Alba, A. R. Liquid chromatography Orbitrap mass spectrometry with simultaneous full scan and tandem MS/MS for highly selective pesticide residue analysis. Anal. Bioanal. Chem. 2015 Aug; 407(21):6317-26.



4. Wang, J.; Chow, W.; Leung, D.; Chang, J. Application of ultrahigh-performance liquid chromatography and electrospray ionization quadrupole orbitrap high-resolution mass spectrometry for determination of 166 pesticides in fruits and vegetables. J. Agric. Food Chem. 2012, Dec 12;60(49):12088-104.



Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Wang et al. went on to apply the Q Exactive mass spectrometer to the determination of 451 pesticide residues in ten fruit and vegetable matrices.⁵ As before, the QuEChERS method was used to extract target pesticides. Full-scan data were used for quantitation and full MS/dd-MS² generated product ion spectra were used to identify compounds. To achieve optimal method accuracy, quantitation was performed using matrix-matched standard calibration curves along with isotopically labeled standards, or a chemical analogue, as internal standards. The method was validated based on overall recovery, intermediate precision, and measurement uncertainty. In the ten matrices studied, 94.5% of the pesticides in fruits and 90.7% in vegetables had recoveries between 81 and 110%; 99.3% of the pesticides in fruits and 99.1% of the pesticides

Pesticides in fruits and vegetables

in vegetables had an intermediate precision of \leq 20%; and 97.8% of the pesticides in fruits and 96.4% of the pesticides in vegetables showed measurement uncertainty of \leq 50%. The measurement uncertainty met the recommended default value of the European Commission SANCO/12495/2011 method validation and quality control procedures for pesticide residues analysis in food and feed. The authors concluded that the method demonstrated acceptable performance for the quantitation of pesticide residues in fruits and vegetables. Further, full MS/dd-MS² with library matching demonstrated the potential to improve pesticide identification in routine practice.

5. Wang, J.; Chow, W.; Chang, J.; Wong, J. W. Ultrahigh-performance liquid chromatography electrospray ionization Q-Orbitrap mass spectrometry for the analysis of 451 pesticide residues in fruits and vegetables: method development and validation. J. Agric. Food Chem. 2014, Oct 22;62(42):10375-91.



Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Zomer et al. developed and validated a Q Exactive mass spectrometerbased method using DIA for the simultaneous quantitation, identification, and qualitative screening of pesticides in fruits and vegetables.⁶ Thermo Scientific TraceFinder software was used to process data. Both quantitative and qualitative performance was evaluated. The goal of the quantitative assessment was to determine whether non-targeted fullscan MS can replace QQQ MS. Quantitative performance was assessed by spiking 184 pesticide compounds in lettuce and orange matrices at 10 and 50 ng/g and guantitative validation was performed according SANCO/12571/2013. The goal of the qualitative assessment was to determine the ability of the method to detect the presence or absence of pesticides in an automated fashion. Qualitative performance was tested by analyzing nine additional matrices (apple, French bean, broccoli, carrot, celery, grape, leek, nectarine, and tomato) spiked with the same 184 compounds at 10, 50, and 200 ng/g. Reproducibility of qualitative performance was tested by repeating the analysis on new extracts, about 4 weeks later.

Data-independent acquisition provided a fully non-targeted approach for data acquisition. As described by the authors, a full-scan acquisition event without fragmentation at resolving power 70,000 was followed by five consecutive DIA fragmentation events at resolving power 35,000. According to the authors, the advantages of using a full-scan technique are that it is not necessary to decide or know beforehand which compounds should be targeted and the ease with which the number of compounds in a method can be increased beyond the practical maximum number of compounds typical of a QQQ MS method.

Pesticides in fruits and vegetables

The quantitative validation demonstrated that the majority of the compounds met the criteria for trueness and precision set forth in SANCO/12571/2013. For the gualitative validation of the untargeted screening capabilities of the method, an overall detection rate of 92% was achieved at 10 ng/g, which increased to 98% at 200 ng/g. A screening detection limit (SDL) of 10 ng/g was achieved for 134 of the pesticides. For 39 of the compounds, the SDL was 50 ng/g. For two pesticides, the SDL was 200 ng/g. For the other nine compounds no SDL could be established. The recommended ion ratio identification criteria were met for 93% of the detected pesticide/matrix/concentration combinations. Based on these results, the authors concluded that the method can be used to combine, in one measurement, the quantification and identification of pesticides typically detected using QQQ MS/MS with qualitative screening to find a range of less frequently detected compounds. Qualitative screening uses the same data with another data-processing approach involving automatic detection by the TraceFinder software.

- Detect both target and untargeted compounds in full scan mode
- Analyze more compounds per run than practical using QQQ instruments
- Quantify, identify, and screen in a single run

6. Zomer, P. and Mol H. G. J. Simultaneous quantitative determination, identification and qualitative screening of pesticides in fruits and vegetables using LC-Q-Orbitrap[™]_-MS. Food Addit. Contam: Part A. 2015 Sep 15:1-9. [Epub ahead of print].



Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Because complex matrices present a challenge to fast and accurate screening and quantitation of pesticides in food, **Yang et al**. studied the relationship between matrix effects and LC separation and elution of pesticides and matrix components using an Orbitrap mass spectrometer.⁷ The study used two LC columns containing different adsorbents. The 108 samples were prepared in solvent and five different sample matrices (avocado, spinach, orange, hazelnut, and honey) using calibration standards of 381 pesticides at three dilution levels of 1×, 1/10×, and 1/100×.

Principal component analysis and slope ratios of calibration curves demonstrated that the $1/100 \times$ sample dilution could minimize ion suppression (matrix effects) for most of the pesticides analyzed. If a pesticide coeluting with matrix components had a peak intensity of 25 times or higher, the suppression for the pesticide persisted to $1/100 \times$ dilution.

The effect and optimization of mass spectrometer parameters on the analysis of pesticide residues in complex food matrices is of great interest. **Rajski et al**. evaluated of the impact of Orbitrap mass spectrometer operating parameters such as resolution on analytical performance.⁸ Analyses were performed on QuEChERS extracts of tomato, pepper, orange, and green tea. The extracts were spiked with 170 pesticides at concentrations of 10 μ g/kg, 50 μ g/kg, 100 μ g/kg, and 500 μ g/kg and were diluted 5-fold prior to UHPLC-MS analysis in the full-scan mode. Three resolution settings—17,500, 35,000, and 70,000—were tested at each concentration level.

Pesticides in fruits and vegetables

Using a resolution setting of 17,500 with 5 ppm of mass tolerance, the percent pesticides detected at 10 µg/kg ranged from 91% in tomato to 83% in green tea. The percent pesticides detected increased when higher resolution settings were used. Analysis at 35,000 resolution produced better results—peak areas were more reproducible and more pesticides were detected—and was sufficient for analysis of the tomato, pepper and orange matrices. Green tea, the most problematic matrix, required analysis at 70,000 resolution. The rates of compounds detected at 70,000 resolution ranged from 98% in tomato to 88% in green tea, thus the authors determined the resolution of 70,000 to be best with the smallest percentage of false negatives at low concentrations. False negative detects were mainly due to a lack of sensitivity for a particular compound, combined with ion suppression effects of the matrix.

Reproducibility improved at resolution levels of 35,000 or higher. Linearity was evaluated from 2–100 ng/mL (10–500 µg/kg in the sample). Unlike other high-resolution MS technologies such as TOF, the Orbitrap mass spectrometer did not experience poor linearity due to detector saturation. The results obtained in this study were comparable those obtained using QQQ MS technology.

• Detect more pesticides using higher-resolution settings (70,000)

• Avoid detector saturation and poor linearity of Q-TOF technology

7. Yang, P.; Chang, J. S.; Wong, J. W.; Zhang, K.; Krynitsky, A. J.; Bromirski, M.; Wang, J. Effect of Sample Dilution on Matrix Effects in Pesticide Analysis of Several Matrices by Liquid Chromatography-High-Resolution Mass Spectrometry. J. J. Agric. Food Chem. 2015, Jun 3;63(21):5169-77.





8. Rajski, Ł.; Gómez-Ramos Mdel, M.; Fernández-Alba, A. R. Large pesticide multiresidue screening method by liquid chromatography-Orbitrap mass spectrometry in full scan mode applied to fruit and vegetables. J. Chromatogr. A. 2014, Sep 19;1360:119-27.



Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Synthetic hormones and other veterinary drugs in animal products

Because the synthetic steroid hormones used to treat animals have been found to affect cancer risk, consumers are concerned about exposure via the animal products they eat. However, the complexity of food matrices makes it difficult to confidently detect and quantify these contaminants at low concentrations. **Kumar et al**. studied the application of the Q Exactive mass spectrometer in four acquisition modes: full MS/AIF, full MS/tMS², full MS/ddMS², and t-SIM/ddMS², to the analysis of synthetic hormones.⁹

Finding it most suitable, the authors developed a t-SIM/ddMS² confirmation method for the analysis of eight synthetic hormones (trenbolone, 17α ethinylestradiol, zeranol, stanozolol, dienestrol, diethylstilbestrol, hexestrol, taleranol) and one naturally occurring hormone (zearalenone) in animal urine. Quadrupole precursor ion selection (SIM) acted as a powerful filter to reduce ion suppression. The method was validated according to the European Commission Decision 2002/657/EC for analysis of residues of veterinary medicinal products. The decision limit (CC α) and the detection capability (CC β) ranged from 0.11 µg/L to 0.69 µg/L, and 0.29 µg/L and 0.90 µg/L, respectively. Overall, the results suggested that the Q Exactive mass spectrometer provides sensitivity similar to that of QQQ instruments, with enhanced selectivity.

Kaufmann et al. applied the Q Exactive mass spectrometer to the quantitation of over 100 compounds belonging to a variety of veterinary drug classes in milk.¹⁰ Instead of traditional extraction and clean up approaches such as solid phase extraction (SPE), QuEChERS and ultra-filtration, the authors introduced a new technique—salting out supported liquid extraction (SOSLE)—to enhance extraction efficiency and sample clean-up of polar analytes. The method was validated based on European Commission Decision 2002/957/EC as it applies to quantitative veterinary drug analysis methods. The authors concluded that the Q Exactive mass spectrometer provided good quantitative data and superior compound confirmation. The SOSLE technique produced extracts of equal or superior cleanliness and with higher average recoveries than those obtained using QuEChERS or SPE. The largest improvement was obtained for polar analytes as such penicillines, quinolones, and tetracyclines.

• Reduce ion suppression with quadruple precursor ion selection

• Attain sensitivity equal to QQQ instruments, with improved selectivity

9. Kumar, P.; Rúbies, A.; Centrich, F.; Granados, M.; Cortés-Francisco, N.; Caixach, J.; Companyó, R. Targeted analysis with benchtop quadrupole-orbitrap hybrid mass spectrometer: application to determination of synthetic hormones in animal urine. Anal. Chim. Acta. 2013, May 30;780:65-73.



10. Kaufmann, A.; Butcher, P.; Maden, K.; Walker, S.; Widmer, M. Multi-residue quantification of veterinary drugs in milk with a novel extraction and clean up technique: salting out supported liquid extraction (SOSLE). Anal. Chim. Acta. 2014, Apr 11;820:56-68.

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Mycotoxin contamination can occur when molds grow on dairy products or when dairy cattle eat contaminated feeds. **Jia et al**. developed a Q Exactive mass spectrometer-based UHPLC-MS method for simultaneous analysis of 58 mycotoxins in dairy products.¹¹ The QuEChERS sample preparation method was used. The authors validated the method using the guidelines specified in European Commission Decision 2002/657/EC and 401/2006/EC. Extraction recoveries ranged from 86.6 to 113.7%, with a coefficient of variation of < 6.2%. All target compounds were detected within the range of 0.001 to 100 µg/kg, with a correlation coefficient of > 0.99. The limits of detection (LOD) ranged from 0.001 to 0.92 µg/kg. The decision limit (CC α) values were comparable to those of previously reported QQQ methods. Repeatability was less than 6.4%. Three consecutive on-column injections of 1 µg/kg were made to detect any decrease in mass accuracy. No significant decrease was observed and the maximum mass deviation ranged from 0.2 to 2.5 ppm, demonstrating the wide dynamic range of Q-Exactive mass spectrometer at resolution 70,000 (FWHM). Compared with other methods, the method increased sensitivity and mass accuracy by more than five times. In summary, the authors found the method useful for fast screening of mycotoxins in dairy products.

- Realize a wider dynamic range of detection
- Increase sensitivity and mass accuracy by a factor of five

11. Jia, W.; Chu, X.; Ling, Y.; Huang, J.; Chang, J. Multi-mycotoxin analysis in dairy products by liquid chromatography coupled to quadrupole orbitrap mass spectrometry. J. Chromatogr. A. 2014, Jun 6;1345:107-14.



Mycotoxins in dairy products

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Analytical methods for screening and quantitating the many contaminants found in foods must be reliable and meet regulatory requirements for recoveries and limits of quantification (LOQ). **Dzuman et al**. presented a reliable and sensitive method that used HPLC separation in combination with HRAM Orbitrap mass spectrometer-based detection for the determination of 323 pesticide residues, 55 mycotoxins, and 11 plant toxins.¹² The efficiency of a special core-shell HPLC column with relatively high particle size (2.6 mm) with lower operational pressures was also examined. The authors validated the method in three sample matrices, leek, wheat, and tea, which differed in the type and amount of components causing matrix effects. QuEChERs was used for sample preparation and extraction of target analytes. A HRAM MS/MS spectral library containing the spectrum of fragment ions for each analyte was created to facilitate identification and confirmation of the target compounds.

Multiclass contaminants in food

The core-shell analytical column demonstrated good separation efficiency and robustness, yielding retention-time RSDs of less than 0.3% after 2000 injections. Method LOQs for the target analytes were less than 10 µg/kg for 82%, 81%, and 61% for the leek, wheat, and tea matrices, respectively. For the majority of the target analytes, recoveries were 70 to 120%, a range acceptable according to SANCO/12571/2013 analytical quality control and validation procedures for pesticide residues analysis in food and feed. The exception was the highly polar mycotoxin deoxynivalenol-3-glucoside with recoveries of 35%, 47%, and 42% for leek, wheat, and tea matrices, respectively. Calibration curve linearity expressed as coefficients of determination ranged from of 0.9661 to 1.000. Repeatability at the LOQ ranged from 0.25 to 13.51% RSD.

- Screen and quantify multiclass contaminants, reliably and in accordance with regulatory requirements
- Identify and confirm target compounds

12. Dzuman, Z.; Zachariasova, M.; Veprikova, Z.; Godula, M.; Hajslova J. Multi-analyte high performance liquid chromatography coupled to high resolution tandem mass spectrometry method for control of pesticide residues, mycotoxins, and pyrrolizidine alkaloids. Anal. Chim. Acta. 2015, Mar 10;863:29-40.



Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

The method was validated according to the European Commission Decision. Ensuring the safety of baby food is of utmost importance, but screening and quantitating hundreds of known and unknown contaminants such as pesticides and antibiotics can be challenging. **Jia et al**. developed an UHPLC-ESI Q Exactive mass spectrometer method for simultaneous analysis of 333 pesticides and veterinary drug residues in baby food.¹³ QuEChERs was used for sample preparation and extraction.

The method was validated according to the European Commission Decision 2002/657/EC and SANCO/12571/2013. QuEChERs extraction recoveries ranged of 79.8 to 110.7%, with coefficients of variation < 8.3%. The 333 compounds ranged from 0.1 to 1000 µg/kg in concentration, with a correlation coefficient > 0.99. The LODs ranged from 0.01 to 5.35 µg/kg. The LOQs were in the range of 0.01 to 9.27 µg/kg. After successfully screening 93 commercial baby food samples for pesticide and veterinary drug residues, the authors concluded the method is appropriate for rapid screening of foods. In particular, tilmicosin, fenbendazole, tylosin tartrate, and thiabendazole were detected in some samples.

More than 300 compounds belonging to several classes of veterinary drugs and pesticides have been found in animal feed. Because medium and high-resolution mass spectrometers provide advantages in multi-residue analysis, **Gómez-Pérez et al**. compared the performance

of medium-resolution (MRMS) TOF and high-resolution Orbitrap mass spectrometers for the analysis of toxic compounds in 18 different chicken, hen, rabbit, and horse feed samples.¹⁴ Sample cleanup procedure was evaluated and several validation parameters were established including matrix effect, linearity, recovery and sensitivity.

The authors obtained better results using the Orbitrap mass spectrometer with sensitivity of 1 to 12.5 µg/kg (below MRL), recovery values of 60–125%, and fewer compounds experienced signal suppression or enhancement. The TOF LOQ values ranged from 5 to 100 µg/kg. Sulfadiazine, trimethoprim, robenidine and monensin sodium veterinary drugs, and the pesticide chlorpyrifos, were identified when the method was applied to the feed samples, demonstrating its applicability as a quantitative method regardless of the type of feed. Other advantages provided by the Orbitrap mass spectrometer important for routine analysis were short analysis time (14 minutes), ability to perform fast screening using Thermo Scientific[™] ToxID[™] automated screening software (which processed more than 450 compounds in less than 5 minutes), and lack of need for a lock mass.

- Obtain better sensitivity and recovery for multi-residue analysis
- Reduce signal suppression and enhancement compared to Q-TOF technology
- Rapidly screen and save time using ToxID software

13. Jia, W.; Chu, X.; Ling, Y.; Huang, J.; Chang, J. High-throughput screening of pesticide and veterinary drug residues in baby food by liquid chromatography coupled to quadrupole Orbitrap mass spectrometry. J. Chromatogr. A. 2014, Jun 20;1347:122-8.



14. Gómez-Pérez, M. L.; Romero-González, R.; Martínez Vidal, J. L.; Garrido Frenich, A. Analysis of veterinary drug and pesticide residues in animal feed by high-resolution mass spectrometry: comparison between time-of-flight and Orbitrap. Food Addit. Contam. Part A Chem. Anal. Control Expo. Risk. Assess. 2015, Mar 18:1-10._

Multiclass contaminants in food



Marine biotoxins

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

During algal blooms, biotoxins produced by certain types of algae can become concentrated in filter feeders such as shellfish. Though the biotoxins don't harm the shellfish, but they can accumulate to levels that cause serious illness or even death in humans and other mammals when eaten. Rúbies et al. developed and validated a high-throughput method to analyze lipophilic marine biotoxins (okadaic acid, dinophysistoxins, azaspiracids, pectenotoxins, vessotoxins, spirolids) in fresh and canned bivalves.¹⁵ The method coupled QuEChERS sample cleanup and extraction with LC-MS analysis using the Q Exactive mass spectrometer operating in tandem MS mode, with resolution set to 70,000 (FWHM) at m/z 200. Separation of analytes was performed in about ten minutes in gradient elution mode with a BEH C18 column and mobile phases based on 6.7 mM ammonia aqueous and acetonitrile mixtures. For each analyte, the molecular ion and one or two product ions were acquired, with a mass accuracy better than 5 ppm. Quantitation was performed using surrogate matrix matched standards, with eprinomectin as the internal standard.

The QuEChERS sample clean up and extraction procedure yielded high absolute recoveries. High-resolution MS/MS data acquisition was powerful in avoiding matrix interferences. Accurate mass data for both the molecular ion and the selected fragments could be obtained for each target analyte, enabling confirmation of compound identity and avoidance of false positives. The authors concluded that the method is straightforward, reliable, and suitable for routine confirmatory quantitative analysis of lipophilic marine biotoxins in fresh and canned bivalves at regulated levels. The method meets the requirements of the EU food safety regulations and is in routine use in a public health laboratory.

• Avoid matrix interferences with high-resolution

• Confirm compound identity and avoid false positives using accurate mass

15. Rúbies, A.; Muñoz, E.; Gibert, D.; Cortés-Francisco, N.; Granados, M.; Caixach, J.; Centrich, F. New method for the analysis of lipophilic marine biotoxins in fresh and canned bivalves by liquid chromatography coupled to high resolution mass spectrometry: a quick, easy, cheap, efficient, rugged, safe approach. J. Chromatogr. A. 2015, Mar 20;1386:62-73.



Marine biotoxins

 $(\bigtriangleup$

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Likewise, **Domènech et al**. developed and validated a Q Exactive mass spectrometer-based LC-MS method for the quantitation and confirmation of lipophilic marine biotoxins in mussels.¹⁶ Compounds representative of each lipophilic toxin group were analyzed: Okadaic acid (OA), yessotoxin, azaspiracid-1, gymnodimine, 13-desmethyl spirolide C, pectenotoxin-2 and Brevetoxin B. Identification and confirmation criteria were established. Fragment and isotope ions and ion ratios were evaluated for use in confirmation. The authors found that both fragment ion and isotope ion ratios can be used to confirm a positive result, but for each compound one or the other can be more suitable.

Accuracy (trueness and precision), linearity, calibration curve check, LOQ and specificity were used as method validation parameters and the validation was performed at 0.5 times the European Union permitted levels. Overall, the method performed very well for the parameters investigated. Trueness (recovery) ranged from 80% to 94%, precision (intra-laboratory reproducibility), ranged from 5% to 22%, and LOQs were from 0.9 to 4.8 pg on column. Overall method uncertainty of 38% was estimated for OA, using certified reference material and a top-down approach considering contributions arising from the trueness and precision studies. The authors concluded that the Orbitrap mass spectrometer-based method enables full-scan acquisition with good sensitivity and better selectivity than other approaches, and can help avoid false positives when confirmation criteria are used.

- Exceed EU-specified levels for accuracy, linearity, LOQ, and specificity
- Enjoy full-scan data acquisition with enhanced sensitivity and selectivity
- Avoid false positives using confirmation criteria

16. Domènech, A.; Cortés-Francisco, N.; Palacios, O.; Franco, J. M.; Riobó, P.; Llerena, J. J., Vichi, S.; Caixach, J. Determination of lipophilic marine toxins in mussels. Quantification and confirmation criteria using high resolution mass spectrometry. J. Chromatogr. A. 2014, Feb 7;1328:16-25.



Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

White wines can contain dyes added to produce a richer color and these dyes can cause reactions in susceptible people. **Jia et al**. developed a method combining QuEChERS sample preparation and extraction with Q Exactive mass spectrometer detection for accurate and sensitive screening of 69 dyes in wine.¹⁷ After optimization of the QuEChERS procedure, the maximum predicted recovery was 99.48% for canacert indigo carmine. Recovery rates of the other 68 compounds ranged from 87.2 to 107.4%, with a coefficient of variation < 6.4%. The mass accuracy obtained was routinely better than 1.6 ppm with once-perweek calibration. The LODs ranged from 1 to 1000 µg/kg. The authors concluded that the method is very useful for fast screening of dyes in commercial wines.

Screen rapidly, accurately, and with high sensitivity

17. Jia, W.; Chu, X.; Ling, Y.; Huang, J.; Lin, Y.; Chang, J. Simultaneous determination of dyes in wines by HPLC coupled to quadrupole orbitrap mass spectrometry. J. Sep. Sci. 2014, Apr;37(7):782-91.



Dyes in wine

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

The active ingredients in herbal medicines and dietary supplements can augment or antagonize the actions of other prescription and non-prescription drugs. In the US, antidiabetics are prescription-only medicines designed to be taken only with physician supervision. Guo et al. developed a Q Exactive mass spectrometer-based method for rapid screening, confirmation, and quantitation of 11 illegal antidiabetic adulterants in herbal medicines and dietary supplements.¹⁸ Sixty-three batches of herbal medicine, and 34 batches of dietary supplement samples were tested. The mass spectrometer was operated in the full MS/ddMS² mode. Full-scan data was used for identification and provided superior accuracy, precision, and sensitivity for quantitation. The data-dependent MS/MS scans produced product ion spectra for unambiguous compound confirmation. Quantitation was performed using matrix-matched standard calibration curves with phenacetin as the internal standard. Method validation parameters were selectivity, sensitivity, calibration curve, accuracy and precision, recovery, matrix effects, and stability.

Herbal medicines and dietary supplements

Response was linear over wide analyte concentration ranges (e.g., 0.0004 to 1 µg/g for metformin) with coefficients of correlation $r^2 > 0.9991$. The LODs ranged from 0.05 to 0.5 ng/g. Recoveries were higher than 74.3%. Accuracy ranged from -6.75 to 3.85% and the intraand inter-day precision ranged from 0.048 to 11.5%. The adulterants metformin, phenformin, and glibenclamide were detected in seven of the dietary supplements tested, but in none of the herbal medicines. The authors concluded that the method demonstrated very good performance for the identification, confirmation, and quantitation of antidiabetics in herbal medicines and dietary supplements.

- Use full-scan data for compound ID as well as accurate quantitation
- Confirm compounds confidently using data-dependent MS/MS

18. Guo, C.; Shi, F.; Jiang, S.; Gong, L.; Zhao, Y.; Zhang, J.; Zeng, S. Simultaneous identification, confirmation and quantitation of illegal adulterated antidiabetics in herbal medicines and dietary supplements using high-resolution benchtop quadrupole-Orbitrap mass spectrometry. J. Chromatogr. B. Analyt. Technol. Biomed. Life. Sci. 2014, Sep 15;967:174-82.



Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Herbal medicines and dietary supplements

Vaclavik et al. developed and validated a Q Exactive mass spectrometer method for the simultaneous determination of 96 pharmaceuticals, plant toxins, and other plant metabolites in herbal dietary supplements.¹⁹ Target analytes were extracted using the QuEChERS method. The mass spectrometer was operated in full MS/ddMS² acquisition mode, which provided high-resolution full-scan data for quantitation and high-resolution MS/MS data for confirmation in a single analytical run. The method provided excellent selectivity in both full MS and ddMS² modes. Confirmation of analytes was with a high degree of confidence. Method LODs and LOQs differed significantly depending on the sample matrix tested. Across the five different matrices, the LODs \leq 10 µg/kg and LOQs \leq 50 µg/kg were obtained for 48 to 81% of the target compounds. With the exception of the highly polar analytes, the QuEChERS extraction provided acceptable recoveries in the range of 70% to 120%. The precision of the method, defined as the relative standard deviation (RSD, n = 5), was \leq 25% and \leq 18% at spiked concentrations of 50 µg/kg and 500 µg/kg, respectively. Because of sample-to-sample variation in matrix effects of the extracts, the method of standard additions and an approach based on dilution of matrix components followed by quantitation using solvent standards were applied to the quantitative analysis.

Quantitate and confirm in a single run

19. Vaclavik, L.; Krynitsky, A. J.; Radar, J. Targeted analysis of multiple pharmaceuticals, plant toxins and other secondary metabolites in herbal dietary supplements by ultra-high performance liquid chromatography-quadrupole-orbital ion trap mass spectrometry. Anal. Chim. Acta. 2014, Jan 31;810:45-60.



Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Environmental contaminants analysis

Targeted screening and quantitation of environmental contaminants: workflow

Contaminants in drinking and wastewater

Contaminants in surface water

60





software.

Environmental contaminants analysis

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

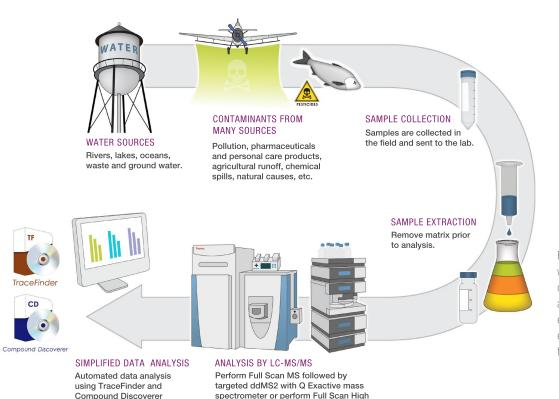
Targeted screening and quantitation of environmental contaminants: workflow

Ultimately the goal of any environmental analysis is the quantitative determination of contaminants. Water, soil, and air have regulated levels of contaminants established to protect human health. These levels are based on the degree of toxicity along with prevalence and cost of cleanup. Monitoring for and removing toxic compounds from the environment, especially drinking water, also increases costs. Therefore, quantitation of exact concentration is crucial to both preserving health and minimizing costs. Also important are investigations of non-regulated, emerging contaminants to assess if and when a future regulatory determination is to be made.

Ensuring the safety of drinking water is one area of interest. Because lakes, rivers, streams, and oceans are impacted by human and natural activities, monitoring these resources continues to be of interest. HRAM analysis is ideally suited to rapid screening and quantitation of water samples for targeted contaminants such as pesticides, herbicides, and other pollutants. Orbitrap mass spectrometer-based LC-MS analyses of water are performed using variations of the workflow shown in Figure 8.

> DidyOUknow? Screening, identification, and quantitation of environmental contaminants is faster and more certain with Thermo Scientific[™] High-Resolution Accurate-Mass MS/MS

> > Spectral Libraries. To learn more: click here



Speed Screening with Exactive Plus.

Figure 8. Q Exactive mass spectrometer workflow for targeted screening and quantitation of contaminants in water. Samples are processed and contaminants are extracted using various extraction techniques, from online solid-phase extraction to offline extraction, prior to analysis on the mass spectrometer.





Environmental contaminants analysis

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

1,2 N-nitrosamines (NAs) are receiving increased attention due to their high carcinogenic and mutagenic potential at low concentrations. Though a countrywide maximum level in drinking water is not yet established in North America, California and Ontario have set limits of 10 ng/L and 9 ng/L, respectively. N-Nitrosamine compounds are produced by industrial activity and resist removal during wastewater treatment. In addition, N-Nitrosamines such as NDMA are by-products of disinfection by chlorination and chloramination. Because of their hydrophobicity and polarity, extraction and detection of NAs at low levels is a challenge. **Ngongang et al.** developed a selective and robust Q Exactive mass spectrometer-based SPE-UHPLC-MS method for the analysis of nine N-nitrosamines in drinking and wastewater matrices.²⁰ SPE was employed as a cost-effective method that helps in achieving low detection limits and high sample throughput.

Contaminants in drinking and wastewater

The authors validated the method in HPLC grade water, drinking water and wastewater matrices and determined that the method sensitivity is comparable with that of GC-MS and LC-MS/MS methods. HRAM data produced by the Q Exactive mass spectrometer helped the authors identify and quantify the target NAs unambiguously, and the selectivity of the HRAM method eliminated matrix interferences. The authors concluded that although GC-MS can provide better sensitivity for N-nitrosamines, LC-MS provides significant time savings because of the longer retention times of GC-MS. Furthermore, LC-MS enabled detection of both GC-detectable and GC-undetectable NAs such as NDPhA.

• Identify and quantify with confidence using HRAM data

• Eliminate matrix interferences with high selectivity

20. Ngongang, A.D.; Vo Duy, S.; Sauve', S. Analysis of nine N-nitrosamines using liquid chromatography-accurate mass high resolution mass spectrometry on a Q-Exactive instrument. Anal. Methods. 2015.





Environmental contaminants analysis

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

New ways to correlate the amounts of drug residues in wastewater with drug consumption are driving growing interest in their quantitation. Due to this interest, **Fedorova et al**. compared the quantitative performance of hybrid quadrupole-Orbitrap-based mass spectrometry with that of QQQ MS for LC-MS analysis of drug residues in wastewater.²¹ The Q-Exactive mass spectrometer was operated at resolution of 70,000 (FWHM) at *m/z* 200 in full-scan (HRFS) mode and 17,500 (FWHM) at *m/z* 200 in product ion scan (HRPIS) mode. For the SRM analyses, the first and third quadrupoles of the QQQ mass spectrometer were set to 0.7 FWHM. A mass-extracted window of 5 ppm around the theoretical *m/z* of each target analyte was used to construct mass chromatograms.

While all three methods showed good linearity and repeatability, the HRPIS mode delivered better sensitivity and selectivity for certain compounds. The selectivity and sensitivity advantages of the Q Exactive mass spectrometer result from quadrupole precursor ion selection combined with HRAM Orbitrap mass-analyzer detection. The LOQs ranged from 0.46 to 20 ng/L. Both the QQQ and the HRPIS Orbitrap mass spectrometer-based MS/MS methods showed good selectivity. The HRFS-based method suffered from more interferences and

Contaminants in drinking and wastewater

showed some false positive results when confronted with co-extracted matrix interferences. The authors concluded that the Q Exactive mass spectrometer is suited for trace-level detection and quantitation of most of the drugs tested in complex wastewater matrices and offers potentially better selectivity than QQQ MS instruments for certain compounds.

When exposed to chemical disinfectants used in wastewater treatment, many pharmaceuticals degrade into by-products that can be more toxic than the parent compound itself. To address this problem, **Negregaria et al**. investigated the stability of a widely used cytostatic etoposide in chlorinated water using the HRAM capability of the Q Exactive mass spectrometer.²² The authors identified two new etoposide oxidation by-products and were able to measure the time course of etoposide degradation into its by-products at different pH values and free chlorine concentrations, and in different water matrices.

- Detect and quantify trace-level compounds in wastewater matrices
- Experience selectivity and sensitivity surpassing QQQ instruments

21. Fedorova, G.; Randak, T.; Lindberg, R.H.; Grabic, R. Comparison of the quantitative performance of a Q-Exactive high-resolution mass spectrometer with that of a triple quadrupole tandem mass spectrometer for the analysis of illicit drugs in wastewater. Rapid. Commun. Mass Spectrom. 2013, Aug 15;27(15):1751-62.





22. Negreira, N.; López de Alda, M.; Barceló, D. Degradation of the cytostatic etoposide in chlorinated water by liquid chromatography coupled to quadrupole-Orbitrap mass spectrometry: identification and quantification of by-products in real water samples. Sci. Total Environ. 2015, Feb 15;506-507:36-45.



Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Moschet et al. developed a Q Exactive mass spectrometer-based LC-MS method that relies on exact mass rather than reference standards. to assess rarely-investigated pesticides and their transformation products (TPs) in 76 surface water samples.²³ One hundred eighty five watersoluble insecticides, fungicides and their major TPs were analyzed. A SPE LC-MS method was developed using 45 known, persistent, and highsales-volume pesticides. Seventy percent of these targets had LOQs < 5 ng/L. This compound set was then used to develop and optimize a screening method using only exact mass as a priori information. This method resulted in a 70% success rate. False negatives were mainly due to low intensity peaks. The authors then applied the method to the remaining 140 compounds. Nineteen additional substances were detected including two TPs that had never been found in the environment before. The authors concluded that the screening approach was fast, successful, and easily expanded to other micropollutant classes for which reference standards are not accessible.

Contaminants in surface water

Due to the large number of potential contaminants, comprehensive assessment of pesticides in surface waters is challenging. Most scientific studies and routine monitoring programs include only 15 to 40 pesticides. Routine, comprehensive screening has not been feasible due to labor-intensive analysis with QQQ mass spectrometers. For this reason, Moschet et al. developed a comprehensive LC-MS screening method relying on the HRAM capability of the Q Exactive mass spectrometer.²⁴ The method covered 86% of all polar organic pesticides sold in Switzerland that are applied to agricultural or urban land (249 compounds) and 134 TPs. Between March and July 2012, the authors regularly drew samples from five medium-sized rivers containing large areas of diverse crops and urban settlements within their catchments. The authors detected over 100 parent compounds and 40 TPs and the sum of the pesticide concentrations was above 1000 ng/L in 78% of samples. The chronic environmental quality standard was exceeded for 19 single substances and the "exceedances" occurred over the entire measurement period in all rivers.

- Screen for surface water contaminants routinely and comprehensively
- Avoid the labor-intensive method set-up of QQQ instruments

23. Moschet, C.; Piazzoli, A.; Singer, H.; Hollender, J. Alleviating the reference standard dilemma using a systematic exact mass suspect screening approach with liquid chromatography-high resolution mass spectrometry. Anal. Chem. 2013, Nov 5;85(21):10312-20.



24. Moschet, C.; Wittmer, I.; Simovic, J.; Junghans, M.; Piazzoli, A.; Singer, H.; Stamm, C.; Leu, C.; Hollender, J. How a complete pesticide screening changes the assessment of surface water quality. Environ. Sci. Technol. 2014, May 20;48(10):5423-32.





Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Moschet et al. also applied the Q Exactive mass spectrometer to a large field study designed to evaluate the in-situ calibration of a passive sampler (styrene divinylbenzene (SDB) covered by a polyether sulfone (PES) membrane for 322 polar organic micropollutants.²⁵ As before, 5 rivers with different agricultural and urban influences were sampled between March and July 2012 using two methods: i) two-week time-proportional composite water samples, and ii) two-week passive sampler deployment. All compounds of different compound classes (logKow -3 to 5, and neutral, anionic, cationic, and zwitterionic species) were analyzed using the Q Exactive mass spectrometer. Because the number of detected substances was similar, LOQs were comparable (median: 1.3 ng/L vs. 1.6 ng/L), and handling was fast and easy. The authors concluded the SDB passive samplers are well-suited for qualitative screening of polar micropollutants.

Contaminants in surface water

Of concern because it is poorly biodegradable, iodinated contrast media (ICM) is used to aid in visualization of human tissue, organs, and the cardiovascular system during radiographic medical procedures. **Zonja et al**. studied the applicability of Orbitrap mass spectrometer-based LC-MS analysis for screening and quantitation of six ICMs and their photo TPs in surface waters.²⁶ The authors began by performing a photodegradation study of ICMs using a sunlight lab-scale simulator. Differential analysis was used to determine the exact masses of molecular ions and retention times of the TPs. Hits were manually filtered to produce a list of 108 suspected TPs.

Next, solid-phase extraction of real surface water samples followed by LC-MS analysis was used to screen for the compounds detected in the photodegradation study. The eleven TPs detected in more than 50% of the samples were selected for structural elucidation using LC-MS and NMR. The median concentrations of parent ICMs ranged from 110 ng/L to 6 μ g/L. TPs were detected at concentrations of 8 ng/L to 0.4 μ g/L. The authors concluded that the LC-MS method facilitates characterization of ICM degradation products and detection of TPs without standards.

Detect, characterize, and quantitate without standards

25. Moschet, C.; Vermeirssen, E. L.; Singer, H.; Stamm, C.; Hollender, J. Evaluation of in-situ calibration of Chemcatcher passive samplers for 322 micropollutants in agricultural and urban affected rivers. Water Res. 2015, Mar 15;71:306-17.



26. Zonja, B.; Delgado, A.; Pérez, S.; Barceló, D. LC-HRMS suspect screening for detection-based prioritization of iodinated contrast media photodegradates in surface waters. Environ. Sci Technol. 2015, Mar 17;49(6):3464-72.



Environmental contaminants analysis

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Anatoxin-a (ANA-a) is highly neurotoxic cyanotoxin produced by cyanobacteria that is a threat to humans and livestock when ingested. **Roy-Lachapelle et al**. developed a rapid method that relies on HRAM Orbitrap mass spectrometer-based detection to distinguish ANA-a from the natural amino acid phenylalanine (PHE).²⁷ The Q Exactive mass spectrometer was equipped with laser diode thermal desorption-atmospheric pressure chemical ionization (LDTD-APCI). Full-scan and targeted ion fragmentation modes were compared to determine which achieve highest selectivity and sensitivity. The method was then applied to eight lake water samples that showed signs of cyanobacterial blooms.

Even though the resolving power of the Q Exactive mass spectrometer was sufficient to distinguish the ANA-a and PHE, the targeted ion fragmentation mode greatly increased selectivity. Internal calibration with standard addition was validated using isotopically labeled phenylalanine (PHE-D5) as the internal standard. The method was validated and determined to be linear with correlation coefficients (r^2) above 0.999. The method yielded better signal-to-noise and thus the lower detection and quantification limits of 0.2 and 0.6 µg/L, respectively, for real samples

than in a previous study of a QQQ-based method. Accuracy, inter-day, and intra-day relative standard deviations were below 15%, and signal recovery in the targeted ion fragmentation mode showed no significant matrix effects with values ranging from 96 to 108%. Using the lock mass feature, mass accuracy was below 1 ppm with low variation. In sum, the LDTD-APCI Q Exactive mass spectrometer system enabled ultra fast screening—in less than 15 seconds per sample–of anatoxin-a at concentrations below established guidelines (3.7 μ g/L) in water matrices, with simplified sample preparation and high selectivity.

Contaminants in surface water

- Screen at concentrations below established guidelines, rapidly, with simplified sample preparation and high selectivity.
- Achieve lower detection and quantification limits than QQQ instruments

27. Roy-Lachapelle, A.; Solliec, M.; Sinotte, M.; Deblois, C.; Sauvé, S. High resolution/accurate mass (HRMS) detection of anatoxin-a in lake water using LDTD-APCI coupled to a Q-Exactive mass spectrometer. Talanta. 2015, Jan;132:836-44.





Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes



Clinical research



Metabolism and Drug Monitoring Research



Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Because it offers the required analytical sensitivity and selectivity, QQQ LC-MS/MS analysis in the SRM mode has been routinely employed for targeted analyte quantitation in a range of forensic toxicology and clinical research applications. However, the SRM duty cycle makes monitoring and quantitating of large numbers of analytes very difficult, if not impossible. For clinical research and forensic toxicology applications where quantitation of multiple targeted and untargeted analytes are required, Orbitrap mass spectrometers offer an outstanding orthogonal solution with QQQ-like analytical selectivity, sensitivity, and dynamic range, which are critical for performing successful quantitative analysis of both endogenous and exogenous compounds in plasma, serum, whole blood, urine, cerebrospinal fluid, and oral fluid. Accurate mass measurements help eliminate the risk of over- and underestimations, reduce the rate of false results, and improve sensitivity by minimizing background noise. Moreover, the ability to analyze and guantitate in the same analytical run, together with the potential for retrospective data analysis, makes Orbitrap mass spectrometers a must-have for clinical research, including drug monitoring research, forensic toxicology, and sports anti-doping (particularly for emerging drugs of abuse) applications.

Clinical research

Clinical research overview

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



Clinical research

Clinical research overview

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Clinical Research

Being able to accurately quantify both endogenous and exogenous compounds in biological samples is of critical importance to clinical researchers. High-resolution MS technologies offer high identification power and easy development of quantitative research methods. Specifically, the advantages of high-resolution MS compared to QQQ MS for clinical research are:

- Comparable sensitivity can be achieved with easier research method development because MS optimization for each analyte is not needed.
- Identification of analytes using isotopic patterns and high-resolution product ion spectra is possible.
- Particularly with Orbitrap mass spectrometer technology, much higher analytical selectivity can be obtained with the ability to determine inaccurate detection of an analyte that is not present and is caused by interferences.

- High-resolution full-scan quantitation methods can monitor a much larger number of compounds and data can be re-evaluated later to address analytical questions that arise.
- Often a smaller amount of sample and a shorter sample preparation time can be used.

Though both Orbitrap and TOF mass analyzers provide high to medium resolving power, higher mass spectral noise can be a major limiting factor in the detection threshold and dynamic range of TOFs instruments.

Didyouknow?

Screening, identification, and quantitation of endogenous and exogenous compounds in plasma, serum, whole blood and other bodily fluids is faster and more certain with Thermo Scientific[™] High-Resolution Accurate-Mass MS/MS Spectral Libraries. To learn more: <u>click here</u>

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





Clinical research

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Though drug metabolism is thought to impact drug toxicity, determination of drug metabolites is not fully explored due to the challenges faced in metabolite profiling. Streamlining analysis of *in vivo* drug metabolism and pharmacovigilance studies could offer a more comprehensive understanding of drug biotransformation. To understand if the analytical challenges can be addressed, **Dahmane et al**. compared the quantification of the pharmaceutical drug tamoxifen, and three of its metabolites in plasma using high-resolution Orbitrap and QQQ mass spectrometer technologies.²⁸ Of particular interest was the ability of the Orbitrap mass spectrometer to collect full-scan mass spectral data with comparable analytical selectivity and sensitivity to the SRM data collected using a QQQ mass spectrometer.

Using simultaneous Quan/Qual capability, the Orbitrap mass spectrometerbased research method enabled identification and relative quantitation of 37 more tamoxifen metabolites than the QQQ method. Using multivariate analysis, the researchers associated metabolite patterns and ratios with the administration of tamoxifen and the CYP2D6 genotype,

Metabolism and drug monitoring research

an enzyme important in drug metabolism. Two hydroxylated metabolites were identified as putative CYP2D6 substrates. Relative quantitation was demonstrated to be reasonably precise at < 20% and suggested that the metabolites found were in consequential amounts.

Mueller et al. presented a method to quantify the five main components of the therapeutic drug teicoplanin.²⁹ The method coupled turbulent-flow-based online extraction with LC-MS analysis performed using a Q Exactive mass spectrometer.⁴ The results of the LC-MS method were compared to the commercially available immunoassay QMS[®] teicoplanin. Results obtained using the Q Exactive mass spectrometer-based LC-MS method correlated well with the results obtained using immunoassay, demonstrating within- and between-day precision and accuracy suitable for quantitative analysis of therapeutic drugs for clinical research.

Detect, characterize and quantitate without standards

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

 \bigcirc

28. Dahmane, E.; Boccard, J.; Csajka, C.; Rudaz, S.; Décosterd, L.; Genin, E.; Duretz, B.; Bromirski, M.; Zaman, K.; Testa, B.; Rochat, B. Quantitative monitoring of tamoxifen in human plasma extended to 40 metabolites using liquid-chromatography high-resolution mass spectrometry: new investigation capabilities for clinical pharmacology. Anal. Bioanal. Chem. 2014, Apr;406(11):2627-40



29. Mueller, D. M; von Eckardstein, A.; Saleh, L. Quantification of teicoplanin in plasma by LC-MS with online sample clean-up and comparison with QMS® assay. Clin. Chem. Lab. Med. 2014, Jun:52(6):879-87.



Clinical research

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Henry et al. compared the full-scan HRAM performance of the Q Exactive mass spectrometer with that of the Thermo ScientificTM TSQ Quantum DiscoveryTM and Thermo ScientificTM Quantum UltraTM triple quadrupole mass spectrometers operating in the SRM mode for LC-MS quantitation of drugs in plasma for clinical research.³⁰ The Q Exactive mass spectrometer was set to mass resolution of 50,000 (FWHM) at *m/z* 200 and a mass extracted window of 5 ppm around the theoretical *m/z* of each analyte was used to construct chromatograms for quantitation. Seventeen drugs, including eight antifungal agents (anidulafungin, caspofungin, fluconazole, itraconazole, hydroxyitraconazole posaconazole, voriconazole, and voriconazole-N-oxide), four immunosuppressants (ciclosporine, everolimus, sirolimus, and tacrolimus) and five protein kinase inhibitors (dasatinib, imatinib, nilotinib, sorafenib, and sunitinib), were analyzed.

The quantitative results obtained from the Q Exactive mass spectrometer demonstrated detection selectivity, assay precision, accuracy, linearity, and sensitivity comparable and with good correlation to that obtained using the QQQ MS instruments in the SRM mode. The Q Exactive mass spectrometer method was compatible with sample preparation approaches used for QQQ-based SRM analyses. The authors commented that the Q Exactive mass spectrometer method development with no optimization of SRM parameters, ability to rapidly transfer the analysis from one mass

Metabolism and drug monitoring research

spectrometer to another, more complete full-scan qualitative information, and easier troubleshooting. Considering the advantages, the authors suggested that there should be a shift in how routine quantitative analyses of small molecules are performed for clinical research.

Dafachronic acid (DA) is a steroid hormone required for the reproductive development of *Caenorhabditis elegans* into fertile adults. **Li et al.** developed a simple, sensitive research method for the absolute quantification of DA that employs derivatization with 2-picolylamine and the high-resolution SIM capability of the Q Exactive mass spectrometer.³¹ Due to the research method sensitivity, only relatively small amounts of worms were needed for the analysis. Orbitrap mass spectrometer-based high-resolution SIM analyses outperformed targeted MS/MS analysis (on the same instrument) and QQQ-based SRM. The LOQ was determined to be as low as 1 pg of DA, enabling absolute quantification of endogenous DA during the worms' reproductive development. The DA levels at different developmental stages and in different Daf-c mutants showed that during the L2 larval stage, DA is highly elevated to ensure complete reproductive development of *C. elegans*.

• Obtain high-resolution SIM results surpassing QQQ-based SRM

• Reduce the amount of sample needed for high-sensitivity analysis

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

30. Henry, H.; Sobhi, H. R.; Scheibner, O.; Bromirski, M.; Nimkar, S. B.; Rochat, B. Comparison between a high-resolution single-stage Orbitrap and a triple quadrupole mass spectrometer for quantitative analyses of drugs. Rapid Commun. Mass Spectrom. 2012, Mar 15;26(5):499-509.



31. Li, T. M.; Chen, J.; Li, X.; Ding, X.J.; Wu, Y.; Zhao, L.F.; Chen, S.; Lei, X.; Dong, M.Q. Absolute quantification of a steroid hormone that regulates development in Caenorhabditis elegans. Anal. Chem. 2013, Oct 1;85(19):9281-7.

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Forensic toxicology and sports doping

Quantitation of drugs and metabolites in urine: workflow

- Equine doping
- Cannabis detection
- Designer drugs in urine





Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

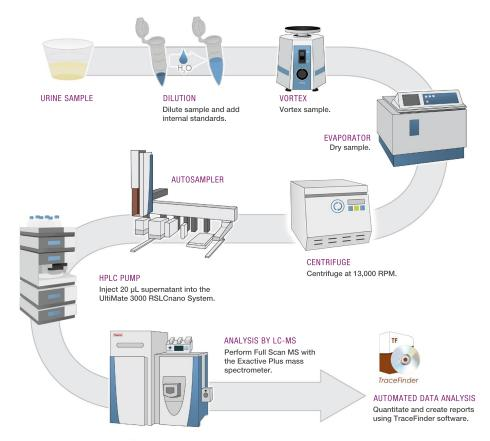
Pharmaceutical Discovery

Conclusion

Application Notes

Quantitation of drugs and metabolites in urine: workflow

In the fields of forensic toxicology and doping, quantitation of drugs in body fluids is challenging due to varying concentrations and substantially different chemical and physical properties of drugs, interfering matrices, occasionally small volumes of sample to test, and the presence of many similar compounds. Further, the constant evolution of drugs and their analogs makes it harder to identify and quantitate them. Orbitrap mass spectrometer-based LC-MS analyses of body fluids are performed using variations of the workflow shown in Figure 9.



Didyouknow?

Screening, identification, and quantitation of toxins and drugs of abuse is faster and more certain with Thermo Scientific[™] High-Resolution Accurate-Mass MS/MS Spectral Libraries. To learn more: <u>click here</u>

For Forensic Use Only.

Figure 9. Q Exactive mass spectrometer-based workflow for quantitation of drugs and metabolites in urine. Currently, urine is a matrix of choice for quantitative analysis in equine drug testing. Many drugs and their metabolites are present in conjugated form in urine and include isomers that need to be identified after separation using chromatography. The analytical workflow must handle a wide range of polarities, many similar molecules, matrix interferences, wide dynamic ranges and sample-to-sample differences caused by the varying nature of the matrix.





Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

The cobra venom toxin α -cobratoxin (α -Cbtx) has analgesic potency greater than morphine. After being found in the facilities of a thoroughbred trainer, the lack of a detection method for the protein became a serious problem in horseracing. To address the problem, **Bailly-Chouriberry et al.** developed the first method, a method that relies on the Q Exactive mass spectrometer, to detect and quantitate α -Cbtx in equine plasma.³² Prior to LC-MS/MS analysis, 3 mL of equine plasma sample was treated with ammonium sulphate precipitation, methanol precipitation, SPE extraction, concentration via filtration, and then digestion with trypsin. LC-MS/MS analysis of the product ions of the doubly-charged precursor of the target peptide was performed at 70,000 resolution. The method was validated and proved to be sufficiently specific, robust and sensitive to enable confirmation of the presence of α -Cbtx in 18 different equine plasma samples spiked at 5 g/L (640 pmol/L), thus meeting the Association of Official Racing Chemists (AORC) requirements. The LOD was determined to be 1 µg/L (130 pmol/L). The method makes it possible to confirm the presence of α -Cbtx in horse plasma from 30 minutes up to 24 hours after administration with an upper limit of 48 hours.

For Forensic Use Only.

Equine doping

Confirm compounds at regulated levels with sensitivity, specificity, and robust operation

32. Bailly-Chouriberry, L.; Cormant, F.; Garcia, P.; Kind, A.; Popot, M. A.; Bonnaire, Y. Identification of -cobratoxin in equine plasma by LC-MS/MS for doping control. Anal. Chem. 2013, May 21:85(10):5219-25.





Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Due to advancements in analytical technology and because it is possible to obtain a specimen quickly and easily, oral fluid is gaining acceptance as an alternative matrix for forensic toxicology applications. Though Δ 9-tetrahydrocannabinol (THC) is the primary target for detecting cannabis use in oral fluid, THC carboxylic acid (THCA) has been demonstrated more reliable because its presence in oral fluid does not occur from passive exposure. However, THCA quantitation is more difficult because it is found in very low concentrations in oral fluid. LC–MS/MS methods provide the requisite sensitivity, but can involve complicated sample preparation procedures. For this reason **He et al.** developed a sensitive LC–MS/MS forensic method for the simultaneous quantitation of THC and THCA in oral fluid using online sample extraction low-flow LC coupled to a Q Exactive mass spectrometer.³³

The method proved simple, robust, and efficient with a total runtime of 12.5 minutes. The HRAM capability of the Q Exactive mass spectrometer enabled the method to achieve high specificity and low-pg/mL sensitivity. Quantitative results were linear from 7.5 to 300 pg/mL for THCA, with a lower limit of 7.5 pg/mL. Intra- and inter-batch precision of ranged from 3.3% to 9.3% for both THC and THCA.

Though Δ (9)-Tetrahydrocannabinol (THC) is the primary target in oral fluid for detecting cannabis use, additional biomarkers such as 11-nor-9-carboxy-THC (THCCOOH), cannabidiol (CBD), and cannabinol (CBN) are needed to address the possibility of passive inhalation. **Concheiro et al.** developed and validated a Q Exactive mass spectrometer-based microflow LC-MS method for the simultaneous quantitation of THC, THCCOOH, CBD, and CBN in oral fluid.³⁴ Authentic oral fluid specimens were collected using the Oral-Eze([®]) and Quantisal[™] collection systems. Proteins were precipitated and the supernatant extracted using CEREX[™] Polycrom[™] THC SPE. Target compounds were identified and quantified in targeted MS/MS experiments.

The forensic method was linear from 0.5 to 50 ng/mL for THC, CBD and CBN, and from 15 to 500 pg/mL for THCCOOH. Intra- and inter-day and imprecision were < 10.8% CV, and bias was from 86.5 to 104.9%. Extraction efficiency ranged from 52.4 to 109.2%, process efficiency from 12.2 to 88.9%, and matrix effect from -86 to -6.9%. The authors concluded that the method provides rapid simultaneous quantitation of THCCOOH, THC, CBD, and CBN with good selectivity and sensitivity, and thus presents an opportunity to improve interpretation of the results obtained from cannabinoid analyses of oral fluids.

For Forensic Use Only.

Quantify challenging low-level compounds, with high selectivity and sensitivity

33. He, X.; Marta Kozak M.; Nimkar, S. Ultra-Sensitive Measurements of 11-Nor-∆9-Tetrahydrocannabinol-9-Carboxylic Acid in Oral Fluid by Microflow Liquid Chromatography–Tandem Mass Spectrometry Using a Benchtop Quadrupole/Orbitrap Mass Spectrometer. Anal. Chem. 2012, 84 (18), pp 7643–7647.



34. Concheiro, M.; Lee D.; Lendoiro, E.; Huestis, M. A. Simultaneous quantification of ∆(9)-tetrahydrocannabinol, 11-nor-9-carboxy-tetrahydrocannabinol, cannabidiol and cannabinol in oral fluid by microflow-liquid chromatography-high resolution mass spectrometry. J. Chromatogr. A. 2013, Jul 5;1297:123-30.

Cannabis detection



Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

New designer drugs and the various drug analogues developed to circumvent legislation have varying effects and potencies, which can complicate interpretation of forensic toxicology cases. Synthetic cathinones are a group of designer stimulants with amphetamine or cocaine-like effects. Comprehensive multi-analyte confirmation methods are needed due to the wide spectrum of synthetic cathinones available. **Concheiro et al.** developed an Orbitrap mass spectrometer-based method for simultaneous quantitation of 28 synthetic cathinones and four metabolites in urine.³⁵ The target cathinones included cathinone, methcathinone, and synthetic cathinones position-3'-substituted, N-alkyl-substituted. Extraction was via solid phase cation exchange extraction (SOLA SCX) followed by reverse-phase LC. Target compounds were identified and quantified using targeted MS/MS experiments. The method was then applied to the urine specimens containing synthetic cathinones.

The authors determined that the method was linear from 0.5-1 to 100 µg/L, with LODs of 0.25-1 µg/L. Imprecision (n = 20) was < 15.9% and accuracy was (n = 20) 85.2 to 118.1%. Extraction efficiency was 78.9 to 116.7% (CV 1.4 to 16.7%, n = 5), process efficiency was 57.7 to 104.9%, and matrix effects were from -29.5% to 1.5% (CV 1.9 to 13.1%, n = 10). The confirmation method proved comprehensive for the 28 synthetic cathinones, with good analytical specificity. The authors concluded that the research would help in the interpretation of test results in forensic toxicology cases, and in the evaluation of the toxicity of designer cathinone drugs.

For Forensic Use Only.

Designer drugs in urine

Confirm and quantitate of a wide range of compounds, simultaneously and with high specificity

35. Concheiro, M.; Anizan, S.; Ellefsen, K.; Huestis, M. A. Simultaneous quantification of 28 synthetic cathinones and metabolites in urine by liquid chromatography-high resolution mass spectrometry. Anal. Bioanal. Chem. 2013, Nov;405(29):9437-48.



Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Pharmaceutical discovery assays

Drug discovery process: workflow

- Bioanalysis
- Discovery-stage screening
- ADME screening and DMPK research





Drug discovery process: workflow

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

As potential leads for successful drug candidates move through the drug discovery process, there is a need for quantitative *in vitro* and *in vivo* analysis at each step. Though the requirement for quantitative LC-MS assays that provide sensitivity, ruggedness, and linear response has remained relatively constant over time, other considerations are emerging as MS technology improves. Ease of use, simplified method development, and troubleshooting tools are growing in importance when choosing the most appropriate MS technology for a particular assay and its corresponding method development. Orbitrap mass spectrometer-based LC-MS analyses of plasma are performed using variations of the workflow shown in Figure 10.

SAMPLE PREPARATION Cleaning up the often Biological samples from in vitro or in vivo studies complex matrixes of typical UHPLC samples is an important step SPE, protein precipitation, With analysis for in vivo mass filtration and centrifugation balance and profiling studies are common spanning days or months, the need for robust, reproducible UHPLC performance is key. MASS SPECTROMETRY SEPARATION By combining robust accurate mass, The need to selectively retain and extreme resolution, and MSⁿ fragmentaseparate a wide array of pharmaceutical tion with multiple techniques. Orbitrap compounds requires a range of high mass spectrometers provide superior performing stationary phases. structural information CD METABOLITE DETECTION LC CHROMATOGRAM By combining accurate mass, isotope pattern, elemental composition, and Compound Discove fragmentation searching, metabolites can DATA ANALYSIS be easily detected in complex samples. In addition to detecting the metabolites within a sample assigning a structure is aided through fragmentation

> techniques within Compound Discoverer

TIME

Quantitation using HRAM Orbitrap mass analyzer technology produces reliable analytical results at extremely low concentration levels and in complex matrices. The performance, in particular sensitivity and linear response, and the HRAM results produced by Orbitrap mass spectrometers are comparable to and in some cases exceed, that of QQQ mass spectrometers. Additionally, HRAM capability reduces the time and effort needed to develop quantitative methods. Full-scan and SIM modes of operation enable direct analysis and monitoring of target analytes without the need to remove interferences through fragmentation or MRM. Data collection in both full-scan and SIM modes also provides additional information about possible matrix interferences near the target analyte m/z that can facilitate quantitative method development and troubleshooting.

Didyouknow?

Screening, identification, and quantitation of pharmaceuticals is faster and more certain with Thermo Scientific[™] High-Resolution Accurate-Mass MS/MS Spectral Libraries. To learn more: click here



Figure 10. Orbitrap mass spectrometer-based workflow for plasma-protein binding (PPB), an early ADME in vitro study that indicates the likelihood of a test compound binding to proteins in blood plasma. The sample matrices analyzed require solvent addition and centrifugation to remove the proteins prior to injection to the LC-MS system. Use of on-line sample preparation such as Thermo Scientific[™] Turboflow[™] Technology allows direct injection of the plasma matrix, thereby increasing sample throughput and reducing cost.



Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Bioanalytical scientists strive to achieve lower limits of quantitation. The reasons range from smaller sample volumes available for analysis, to more potent analytes and the growth of biologics in drug development. As a result, scientists are exploring new LC-MS techniques, involving both high- and low-flow rate LC devices and HRAM. Because most biological samples are matrices comprised of plasma, serum, whole blood, urine, and various tissues, desirable new analytical technology should provide greater sensitivity and robustness. **Wang et al.** assessed application of a Q Exactive mass spectrometer combined with microspray ionization and microflow separation.³⁶ A standard flow rate sprayer needle was compared to a microspray needle and Q Exactive mass spectrometer performance was compared to QQQ MS. Specifically, the sensitivity and noise level of four ionization modes were evaluated: QQQ MS SRM, HRAM full scan, HRAM SIM, and HRAM MS/MS.

Pharmaceutical discovery assays

Bioanalysis

The authors determined that microflow LC-MS showed less matrix effects than UHPLC-MS, but the extent of suppression for both was compound-dependent and the amount of reduction was not proportional for each analyte. Because the precision was improved by using microflow LC-MS, relative quantitation might be more reliable. Microflow LC-MS did provide a substantial increase in sensitivity for some compounds without additional sample preparation or chromatographic method development. Carryover of high-concentration samples was a disadvantage of microflow LC-MS. When using SIM, the Q Exactive mass spectrometer was found to be as or more sensitive than the Thermo Scientific TSQ Vantage Triple Stage Quadrupole Mass Spectrometer, but was also compounddependent. The authors also concluded that the HRAM full-scan capability of the Q Exactive mass spectrometer allowed observation of co-eluting compounds during method development, which in turn enabled reduction or elimination of undesirable matrix effects.

- Obtain sensitivity equal to or exceeding QQQ instruments
- Observe co-eluting compounds, reduce matrix effects, using full-scan HRAM data

36. Wang, H.; Bennett, P. Performance assessment of microflow LC combined with high-resolution MS in bioanalysis. Bioanalysis. 2013, May;5(10):1249-67.





Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

MS/MS with CID enables QQQ and ion trap mass spectrometers to quantify biological analytes in complex matrices with high sensitivity and selectivity. However, the molecular ions produced from cyclic and large disulfide-containing peptides, sterols, and fatty acids are not amenable to CID, making their analysis using these technologies difficult. In particular, cyclic peptides are of growing interest in drug discovery because of their stability in blood and potential for oral dosing. As a solution, **Ciccimaro et al.** demonstrated a Q Exactive mass spectrometer method where the target ion is selectively isolated for quantitation while interfering matrix components undergo MS/MS fragmentation by CID (HCD)—an approach the authors termed HRAM survivor-SIM.³⁷ Fundamentally different than

the traditional MS/MS with CID where a target analyte's unique fragments are monitored following CID (HCD) fragmentation, the new approach significantly enhanced selectivity by removing isobaric interferences. Comparative QQQ MS experiments were performed using a Triple Quad 6500 LC/MS/MS System (AB Sciex). Because molecular ion detection is more sensitive than fragment ion detection, the authors observed a fiveto ten-fold improvement in LLOQs compared to traditional approaches for CID-resistant peptides in plasma extracts. The results demonstrated that the Q Exactive mass spectrometer is an ideal orthogonal platform for quantitation in situations where compounds are not amenable to MS/MS with HCD/CID fragmentation.

- Quantitate using HRAM survivor-SIM when compounds are not amenable to HCD/CID fragmentation
- Eliminate isobaric interferences and enhance selectivity

37. Ciccimaro, E;. Ranasinghe, A.; D'Arienzo, C.; Xu, C.; Onorato, J.; Drexler, D. M.; Josephs, J. L.; Poss, M.; Olah, T. Strategy to improve the quantitative LC-MS analysis of molecular ions resistant to gas-phase collision induced dissociation: application to disulfide-rich cyclic peptides. Anal. Chem. 2014, Dec 2;86(23):11523-7.





Bioanalysis



Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Though high-throughput screening using LC-MS is routinely implemented throughout the drug-discovery process, using it earlier in the lead discovery stage can provide significant benefits. For example, avoiding labeling agents can reduce costly sample preparation, and LC separation of analytes of interest from interferences and can increase selectivity and sensitivity. For this reason **Murphy et al.** applied high-resolution MS-LC multiplexing to a screening assay of phosphorylated peptides.³⁸ Phosphorylated peptide standards were prepared using common enzyme buffers and these were plated into a 96-well plate format prior to LC-MS analysis using a Q Exactive mass spectrometer.

Laboratories supporting the lead discovery stage must analyze several thousand samples per day, generally with complex matrixes that vary from project to project. Therefore, assay methods must be fast and robust. The authors concluded that Q Exactive mass spectrometer coupled with LC multiplexing provides robust, high-quality results at rapid sampling rates (up to 18 seconds per sample). Samples analyzed in both simple and complex sample matrices demonstrated an LOQ of 5 nM with linear response across the working range of the assay.

Discovery-stage screening

Achieve robust, high-throughput screening of complex sample matrices

38. Murphy, K.; Bennett, P. K.; Duczak, N. Jr. High-throughput quantitation of large molecules using multiplexed chromatography and high-resolution/accurate mass LC-MS. Bioanalysis. 2012, May;4(9):1013-24.





Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Application of high-resolution MS to absorption, distribution, metabolism and excretion (ADME), and drug metabolism, and pharmacokinetics (DMPK) studies has generated considerable interest within pharmacokinetics and pharmacology laboratories. **Zhang et al**. explored the benefits of high-resolution MS for quantitative bioanalysis using full-scan data acquisition.³⁹ Of particular interest is the lack of compound-specific MS method development, simultaneous data collection for both targeted and non-targeted components, and suitability for simultaneous quantitation of an unlimited number of compounds, which is not possible using QQQ SRM methods. In addition, the information obtained from HRAM at the ADME phase can be transferred to reduce method development time during downstream routine quantitation analysis using QQQ MS. ADME screening and DMPK research

The authors developed an *in vitro* ADME workflow involving cassette incubation of as many as 32 compounds, followed by quantitative analysis using an Orbitrap mass spectrometer in full-scan mode. The workflow was evaluated for serum protein-binding and parallel artificial membrane permeability (PAMPA) assays. The workflow was found to have acceptable sensitivity, selectivity, and linearity for all compounds, and the biological results obtained were similar to those obtained from discrete incubation and analysis, demonstrating the feasibility of the workflow.

- Achieve a dynamic range exceeding Q-TOF instruments
- Attain scan speed equal to that of Q-TOF systems
- Conserve time, space, and budget with Quan/Qual capability
- Quantify with the instrument for the lab of the future

39. Zhang, J.; Maloney, J.; Drexler, D. M.; Cai, X.; Stewart, J.; Mayer, C.; Herbst, J.; Weller, H.; Shou, W. Z. Cassette incubation followed by bioanalysis using high-resolution MS for in vitro ADME screening assays. Bioanalysis. 2012, Mar;4(5):581-93.





Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

The ability to acquire quantitative data, along with qualitative data for parallel or retrospective data analysis—the quan/qual approach—is powerful in conserving time, laboratory space, and budgets. For this reason, **King et al**. described their evaluation and implementation of Q Exactive mass spectrometer-based Quan/Qual analyses at UCB's research DMPK department.⁴⁰ The authors compared the quantitative performance of Q-TOF and Q Exactive instruments. Both types of instruments performed equally in terms of mass accuracy, sensitivity, robustness, and scan speed. The key difference between the Q-TOF and Orbitrap instruments was the linear quantitative dynamic range achieved, four and greater than four orders of magnitude, respectively, for the compounds tested, a feature considered important in measuring all time points at all doses.

ADME screening and DMPK research

The authors concluded by summarizing the benefits of adopting the Q Exactive mass spectrometer-based Quan/Qual workflow: the additional information obtained produced a more integrated understanding of bioanalytical processes and the DMPK properties of compounds; a large amount of time is saved by not having to develop targeted MS/MS methods for individual compounds; and the ability to obtain both Quan/Qual data from a single analysis has reduced the need for *in vivo* and *ex vivo* studies involving animals, sample preparation and processing, consumables, and laboratory space. The authors expect that over time the development of less expensive high-resolution MS systems will lead to the demise of the use of QQQ mass spectrometers in routine bioanalytical applications.

40. King, L.; Kotian, A.; Jairaj, M. Introduction of a routine quan/qual approach into research DMPK: experiences and evolving strategies. Bioanalysis. 2014, 6(24):3337-48.



Conclusion

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Quantitate for the future, close your uncertainty gap

Quadrupole-Orbitrap mass spectrometer technology has been successfully used for the quantitation of both small and large molecule analytes in a range of clinical research, forensic toxicology, pharmaceutical discovery, environmental analysis, and food testing applications. As observed from numerous peer-reviewed publications, the popularity of performing quantitation using Orbitrap mass spectrometer technology is growing rapidly due to its unique ability to provide HRAM Quan/Qual MS and MS/MS data in a single analytical run. The quality of data produced provides ultimate confidence in the results obtained in a wide variety of qualitative and quantitative applications. High resolving power increases analytical selectivity for compounds in complex matrices and thus reduces the uncertainty associated with the chance of false detection of the analyte of interest. The unique ability to capture all relevant data in using full scan allows retrospective data analysis, reducing the need for additional sample injections. In addition, quadrupole-Orbitrap mass spectrometers are enabling clinical research, forensic toxicology, food safety, and environmental laboratories to raise their productivity to new levels. Compared to QQQ MS-based quantitation methods that involve time consuming optimization of hundreds of SIM or SRM transitions in numerous time windows, quadrupole-Orbitrap mass spectrometer methods for the quantitation of virtually unlimited numbers of compounds are fast and easy to set up. Calibrations are easy to perform and remain stable for days. Fast scanning and spectral multiplexing capabilities make the instrument fully compatible with UHPLC and fast chromatography techniques. With several operating modes such as full-scan MS, AIF, MS/MS, SIM, HCD, PRM, and rapid polarity switching, the Q Exactive and Q Exactive Focus mass spectrometers provide unmatched analytical versatility to meet the needs of laboratories today and into the future.



Thermo Scientific Application Notes

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Detection of Stanozolol Glucuronides in Human Sports Drug Testing by Means of High-Resolution, Accurate-Mass Mass Spectrometry

Quick and Sensitive Analysis of Multiclass Veterinary Drug Residues in Meat, Plasma, and Milk on a Q Exactive Focus LC-MS System

Quantitation of Opiates to Low ng/mL Levels in Urine for Forensic Use Using an Affordable, High-Resolution, Accurate-Mass Mass Spectrometer

Quantitative and Qualitative Confirmation of Pesticides in Beet Extract Using a Hybrid Quadrupole-Orbitrap Mass Spectrometer

Demonstrating High-Performance Quantitative Analysis of Benzodiazepines using Multiplexed SIM with High-Resolution, Accurate Mass Detection on the Q Exactive LC/MS

HRAM Targeted Peptide Quantitation on a Q Exactive MS: A Unique Combination of High Selectivity, Sensitivity and Throughput

Low pg/mL Detection of rHuEPOs in Horse Plasma Employing High-Resolution MS

Full-Scan Fragmentation Options for the Detection of Food Contaminants by an Affordable LC-Q-Orbitrap MS

Variable Data-Independent Acquisition (vDIA) Delivers High Selectivity and Sensitivity in Combined Targeted and Untargeted Analyses for Small Molecules



Thermo Scientific Application Notes

Table of Contents

Introduction

Q Exactive Series Mass Spectrometer

Quantitative Capability

Growing Body of Publications

Food Safety Testing

Environmental Contaminants Analysis

Clinical Research

Forensic Toxicology

Pharmaceutical Discovery

Conclusion

Application Notes

Analysis of Five Barbiturates in Urine Using an Affordable High-Resolution Mass Spectrometer

Quantitation of 25-Hydroxyvitamin D3 and 25-Hydroxyvitamin D2 in Plasma for Clinical Research Using an Affordable **High-Resolution Mass Spectrometer**

Quantification of Steroids in Human Plasma by Liquid Chromatography Coupled with the Affordable Q Exactive Focus HRMS

Bulk Substance Analysis Using DART Source Coupled to LTQ XL Ion Trap and Q Exactive Focus Mass Spectrometers

Analysis of Immunosupressant Drugs in Whole Blood Using a High-Efficiency Method on an Affordable High-Resolution **Mass Spectrometer**









