

Determination of Lipophilic Marine Biotoxins in Molluscs by LC-MS/MS using Offline Extraction

Francesca G. Bellagambi, Department of Chemistry and Industrial Chemistry, University of Pisa, Italy

Cláudia P. B. Martins, Thermo Fisher Scientific, Barcelona, Spain

Michal Godula, Thermo Fisher Scientific, Prague, Czech Republic

Alessio Ceccarini, Department of Chemistry and Industrial Chemistry, University of Pisa, Italy

Key Words

TSQ Quantum Ultra, Food Safety, Marine Biotoxins

Introduction

In recent years many countries have had to deal with the consequences of toxic microalgal blooms in both marine and fresh water, such as the deaths of wild animals and domestic livestock.

Several cases of poisoning in humans have been associated with the direct consumption of shellfish, fish, or water contaminated by algal toxins. People may also come into contact with toxins during recreational activities along sea coasts affected by episodes of algal blooms. Depending on the type of toxin involved, there are forms of mild and usually self-limiting symptoms, characterized by gastrointestinal disorders or allergic-like episodes. Much more severe symptoms of the neurological type can lead to death.

The foods most frequently involved in episodes of human poisoning are represented by bivalve molluscs. These shellfish can accumulate and concentrate any biotoxins present in the plankton they ingest through filtering large quantities of water for trophical reasons. It is not possible to evaluate shellfish edibility by an organoleptic examination alone. While human ingestion of contaminated food with biotoxins can lead to the onset of different clinical symptoms, in shellfish it usually has only marginal effects. An important risk factor lies in the thermostability of such molecules which are not completely inactivated by common physical treatments for fish products (cooking, smoking, salting, freezing, housing), but remain virtually unchanged in the finished product.



There are a series of regulations issued by the European Union (EU) that relate to marine biotoxins. One is Regulation (EC) No 853/2004 which concerns the control of lipophilic toxins, establishing maximum levels for lipophilic toxins in bivalve molluscs destined for human consumption:¹

- For okadaic acid, dinophysistoxins, and pectenotoxins together – 160 micrograms of okadaic acid equivalents per kilogram
- For yessotoxin – 1 milligram of yessotoxin equivalent per kilogram
- For azaspiracids – 160 micrograms of azaspiracid equivalents per kilogram

In the past, aside from bioassays on mice, most analytical techniques developed for the determination of marine biotoxins in bivalve molluscs have been based on offline methodologies. These include methods involving solid phase extraction (SPE) or liquid-liquid extraction (LLE) followed by high pressure liquid chromatography (HPLC) with fluorimetric or UV-diode array detection, as well as detection by liquid chromatography coupled with mass spectrometry (LC-MS).

The EU Commission Regulation (EC) No 15/2011, amending Regulation (EC) No 2074/2005 about the testing methods for detecting marine biotoxins in bivalve molluscs, describes an LC-MS/MS procedure as the reference method for the quantification of lipophilic marine biotoxins – namely okadaic acid, pectenotoxin 2, azaspiracid 1, and yessotoxin.^{2,3} Moreover, dinophysistoxin 1 (DTX-1) and dinophysistoxin 2 (DTX-2) can be quantified by the calibration curve of okadaic acid, pectenotoxin 1 by calibration of pectenotoxin 2, azaspiracid 2 and 3 by calibration of azaspiracid 1 and 45-OH-, and 45-homo-OH-yessotoxin by the calibration of yessotoxin.

In accordance with current European regulations, we propose a quick, selective, sensitive, and accurate analytical method for the determination of lipophilic marine biotoxins in bivalve molluscs using an LC-MS/MS method.

Goal

Our goal is to validate analytical procedures proposed in “EU-Harmonised Standard Operating Procedure for determination of Lipophilic marine biotoxins in molluscs by LC-MS/MS – Version 3” by LC-MS/MS using offline extraction.⁴

Experimental

Sample Preparation

About 1 kg of bivalve molluscs (*Mytilus Galloprovincialis*) were cleaned with water and put in a solution of NaCl (3.5 g/L). After opening, the molluscs were washed with fresh water, their flesh was removed and placed on a stainless steel net, and they were washed again with deionized water. The whole collected raw tissue, not less than 150 g, was chopped and blended by a mixer.

Extraction procedure

9 mL of 100% methanol (gradient quality) were added to 2.00 ± 0.05 g of blended tissue, put into a centrifuge tube, and mixed by vortex for 3 minutes at maximum speed. After centrifugation at 4000 rpm for 10 minutes, the supernatant solution was transferred into a vial.

A second aliquot of 9 mL of 100% methanol was further added to the residual tissue pellet and homogenized for 1 minute by Ultra-turrax® (IKA®, USA) at 12,000 rpm and the mixture was centrifuged at 4000 rpm for 10 minutes. Then the supernatant solution was transferred and

combined with the first extract and made up to 20 mL with 100% methanol. When not immediately analyzed, the solution was stored at -20 °C.

Spikes of toxin standard solutions can be added to the blended tissue before the extraction procedure.

Purification Procedure

The organic extract was purified by being passed through a C18 SPE cartridge preliminarily conditioned with 1 mL of 100% methanol. A 0.45 µm syringe filter was placed at the end of the cartridge to improve purification.

LC Conditions for the Thermo Scientific Hypersil GOLD Column

| | |
|-------------------|--|
| System | Thermo Scientific Accela UHPLC |
| Solvent A | 100% water with 2 mM ammonium formate and 50 mM formic acid |
| Solvent B | 95% acetonitrile + 5% water with 2 mM ammonium formate and 50 mM formic acid |
| Flow Rate | 200 µL/min |
| Gradient | The mixture started at 30% solvent B (8.0 min) followed by a linear gradient up to 90% solvent B in 3.0 min. It went up to 30% of solvent B in 0.5 min. This composition was maintained for 5.5 min. |
| Analytical Column | Hypersil GOLD™; 50 × 2.1 mm, particle size 1.9 µm, part number 25002-052130 |

H-ESI II Source Conditions

| Ion Source Polarity | Positive Ion Mode | Negative Ion Mode |
|--|-------------------|-------------------|
| Spray Voltage | 3000 V | 2700 V |
| Capillary Temperature | 270 °C | 270 °C |
| Vaporizer Temperature | 240 °C | 240 °C |
| Sheath Gas Pressure (N ₂) | 15 units | 15 units |
| Auxiliary Gas Pressure (N ₂) | 5 units | 5 units |

MS/MS Setup

MS analysis was carried out on a Thermo Scientific TSQ Quantum Ultra triple quadrupole mass spectrometer equipped with a heated electrospray ionization probe (H-ESI II).

| Collision Gas (Ar) | 1.5 mTorr |
|-----------------------|------------------------------|
| Q1/Q3 Peak Resolution | 0.7 u (unit mass resolution) |
| Scan Time | 0.100 s |
| Scan Width | 0.500 m/z |
| Data Acquisition Mode | SRM |

| Analyte | ESI Mode | Parent Mass | Product Mass | Collision Energy | Tube Lens |
|---------|----------|-------------|--------------|------------------|-----------|
| AZA-1 | ESI+ | 842.3 | 806.1 | 51 | 207 |
| | | | 824.2 | 42 | 207 |
| AZA-2 | ESI+ | 856.3 | 838.1 | 42 | 214 |
| | | | 820.2 | 49 | 214 |
| AZA-3 | ESI+ | 828.3 | 792.2 | 48 | 192 |
| | | | 810.0 | 40 | 192 |
| PTX-2 | ESI+ | 876.3 | 841.3 | 35 | 205 |
| | | | 823.0 | 40 | 205 |
| | | | 805.3 | 41 | 205 |
| DTX-1 | ESI- | 817.0 | 255.0 | 69 | 197 |
| | | | 113.1 | 67 | 197 |
| DTX-2 | ESI- | 803.15 | 255.3 | 61 | 207 |
| | | | 113.1 | 50 | 207 |
| YTX | ESI- | 1141.5 | 1061.7 | 50 | 240 |
| | | 570.2 | 467.3 | 40 | 240 |
| OA | ESI- | 803.3 | 254.9 | 68 | 216 |
| | | | 113.1 | 50 | 216 |

Table 1: Selected ion transitions (m/z) of the studied compounds and optimized collision energy and tube lens value for the TSQ Quantum Ultra triple quadrupole mass spectrometer

The optimization of selective reaction monitoring (SRM) parameters was performed by direct infusion of standards. Collision-induced dissociation (CID) data were recorded for each analyte including optimum collision energies for the selected ion transitions.

Table 1 summarizes all the mass transitions found for each analyte and its relative collision energy (CE) and tube lens values.

Results and Discussion

To ensure thorough validation of the method, neat standard solutions, standard addition on purified extracts, and spiked blank tissue extracts were prepared and compared.

Table 2 lists the correlation coefficients (r^2) indicating the linearity of the calibration curves for the three types of samples analyzed; five concentrations of the sample solution are considered (2, 5, 10, 20, and 50 $\mu\text{g}/\text{kg}$ or similar).

To assess the inter-day repeatability of the method, ten replicates of spiked samples were analyzed between days. Solutions were prepared containing all the toxins in the five different concentrations used to perform the calibration curves (2, 5, 10, 20, and 50 $\mu\text{g}/\text{kg}$ or similar).

The repeatability of the method expressed as the coefficient of variation percentage (CV %) has been rated less than 20% as shown in Table 3.

| Analyte | Neat Solution | Spiked Purified Extract | Spiked Extract |
|---------|---------------|-------------------------|----------------|
| AZA-1 | 0.9932 | 0.9965 | 0.9970 |
| AZA-2 | 0.9973 | 0.9964 | 0.9901 |
| AZA-3 | 0.9972 | 0.9958 | 0.9993 |
| DTX-1 | 0.9964 | 0.9995 | 0.9953 |
| DTX-2 | 0.9973 | 0.9966 | 0.9965 |
| YTX | 0.9999 | 0.9923 | 0.9988 |
| PTX-2 | 1.0000 | 0.9977 | 0.9927 |
| OA | 0.9955 | 0.9924 | 0.9927 |

Table 2: Correlation Coefficient (r^2) of the calibration curves for the three types of samples analyzed in the concentration range of 2–50 $\mu\text{g}/\text{kg}$

| Analyte | Standard Deviation | Repeatability | CV% |
|---------|--------------------|---------------|-----|
| AZA-1 | 0.25 | 0.79 | 20 |
| AZA-2 | 0.29 | 0.90 | 18 |
| AZA-3 | 0.43 | 1.37 | 17 |
| DTX-1 | 0.18 | 0.55 | 2 |
| DTX-2 | 0.22 | 0.68 | 9 |
| YTX | 0.12 | 0.40 | 12 |
| PTX-2 | 0.39 | 1.36 | 20 |
| OA | 0.16 | 0.48 | 4 |

Table 3: Values of CV% obtained for the repeatability of the lower concentrated curve point (2 $\mu\text{g}/\text{kg}$)

The calculations, of limit of detection (LOD) and limit of quantification (LOQ) were made in accordance with the *UNICHIM Manual N. 179/0* where the calculation of the limit of detection is made through the calibration curve of the instrument used for analysis.⁵

To estimate the LOD and LOQ of the method (Table 4), ten samples were prepared by adding standard solution to 500 mg of homogenized mussel flesh and repeating the extraction procedure according to the method in “*EU-Harmonised Standard Operating Procedure for determination of Lipophilic marine biotoxins in molluscs by LC-MS/MS – Version 3*”. LOD and LOQ are expressed in µg/Kg. Recoveries are shown in Table 5.

| Hypersil GOLD | | | |
|---------------|-------------|-------------|-----------------------|
| Analyte | LOD (µg/Kg) | LOQ (µg/Kg) | Outliers (Huber Test) |
| AZA-1 | 0.56 ± 0.18 | 1.11 | NO |
| AZA-2 | 0.93 ± 0.31 | 1.86 | NO |
| AZA-3 | 1.28 ± 0.43 | 2.57 | NO |
| DTX-1 | 5.66 ± 1.02 | 10.45 | NO |
| DTX-2 | 0.71 ± 0.23 | 1.42 | NO |
| YTX | 1.67 ± 0.56 | 3.33 | NO |
| PTX-2 | 1.40 ± 0.46 | 2.79 | NO |
| OA | 3.95 ± 1.32 | 7.91 | NO |

Table 4: LOD and LOQ of the method

| Analyte | Spiked Purified Extract | Spiked Extract |
|---------|-------------------------|----------------|
| AZA-1 | 96 ± 11 | 97 ± 11 |
| AZA-2 | 101 ± 9 | 94 ± 14 |
| AZA-3 | 104 ± 10 | 99 ± 6 |
| DTX-1 | 101 ± 6 | 101 ± 7 |
| DTX-2 | 101 ± 6 | 108 ± 42 |
| YTX | 99 ± 15 | 102 ± 17 |
| PTX-2 | 103 ± 13 | 102 ± 20 |
| OA | 95 ± 7 | 93 ± 18 |

Table 5: Recovery values, where
 $R\% = \left[\frac{(\mu\text{g/Kg})_{\text{CALCULATED}}}{(\mu\text{g/Kg})_{\text{THEORETICAL}}} \right] * 100$

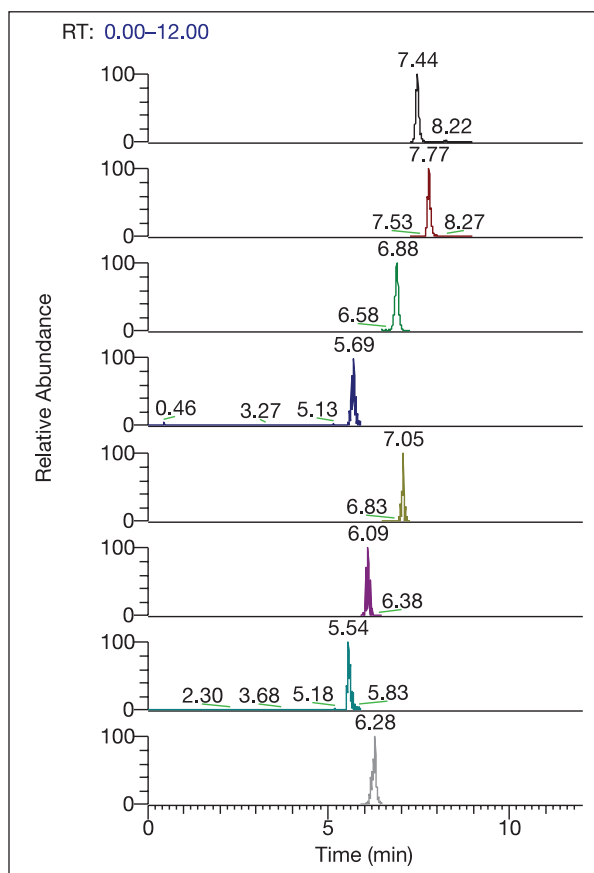


Figure 1: Chromatogram of sample containing 40 ppb of toxins (Retention Time: 7.44 min – AZA-1; 7.77 min – AZA-2; 6.88 min – AZA-3; 5.69 min – OA; 7.05 min – DTX-1; 6.06 min – DTX-2; 5.54 min – YTX; 6.28 min – PTX-2)

Conclusion

This method proved to be selective, sensitive, accurate, and reproducible. It can be successfully applied for the quantitative determination of several classes of lipophilic marine biotoxins in bivalve mollusc samples.

References

1. Regulation (EC) No 853/2004 of the European Parliament and of the Council of 29 April 2004. *Official Journal of the European Communities* L139, 55-255.
2. Commission Regulation (EU) No 15/2011 of 10 January 2011. *Official Journal of the European Union*. L6, 3-9.
3. Commission Regulation (EC) No 2074/2005 of 5 December 2005. *Official Journal of the European Communities*, L388, 27-59.
4. European Union Reference Laboratory for Marine Biotoxins (EU-RL-MB), Agencia Española de Seguridad Alimentaria y Nutrición (AESAN). *EU-Harmonised Standard Operating Procedure for determination of Lipophilic marine biotoxins in molluscs by LC-MS/MS – Version 3*. 2011.
5. UNICHIM, Manual N. 179/0. *Linee guida per la convalida di metodi analitici nei laboratori chimici - criteri generali*. Milano. 1999.

www.thermoscientific.com

©2012 Thermo Fisher Scientific Inc. All rights reserved. ULTRA-TURRAX and IKA are registered trademarks of IKA-Werke GmbH & Co. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific Inc. products. It is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Africa-Other +27 11 570 1840
Australia +61 3 9757 4300
Austria +43 1 333 50 34 0
Belgium +32 53 73 42 41
Canada +1 800 530 8447
China +86 10 8419 3588
Denmark +45 70 23 62 60
Europe-Other +43 1 333 50 34 0
Finland/Norway/Sweden +46 8 556 468 00
France +33 1 60 92 48 00
Germany +49 6103 408 1014
India +91 22 6742 9434
Italy +39 02 950 591

Japan +81 45 453 9100
Latin America +1 561 688 8700
Middle East +43 1 333 50 34 0
Netherlands +31 76 579 55 55
New Zealand +64 9 980 6700
Russia/CIS +43 1 333 50 34 0
South Africa +27 11 570 1840

Spain +34 914 845 965
Switzerland +41 61 716 77 00
UK +44 1442 233555
USA +1 800 532 4752

AN63552_E 07/12M

Thermo
 SCIENTIFIC
 Part of Thermo Fisher Scientific