

# Validated Method for the Analysis of Lipophilic Marine Biotoxins in Bivalves by Liquid Chromatography—High-Resolution Mass Spectrometry

Complete method: Rúbies, A. et al, 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. Chrom. A, 2015 Mar 20;1386:62-73. doi: 10.1016/j.chroma.2015.01.088

## Highlights

- The successfully validated method for the confirmatory, quantitative analysis of lipophilic marine biotoxins meets the requirements of European Union legislation.
- The QuEChERS strategy provides effective and reliable sample preparation.
- UHPLC with HRMS/MS delivers excellent results in the quantitative analysis of MBTXs.
- Calibration with surrogate matrix-matched standards (SMMS) provides reliable quantitative results.
- The lower limit quantitation of the method is 25 µg/kg.

## Introduction

A harmful algal bloom is a serious environmental problem that results when the population of microscopic algae in a water system rapidly increases. The algae have the ability to produce marine biotoxins (MBTXs), which are dangerous to humans and other aquatic life. These biotoxins can accumulate in filter-feeding bivalve mollusks and, in turn, pose significant food safety risks to humans who ingest the shellfish, including gastrointestinal illnesses or neurological disorders.

Adequate testing for biotoxins is necessary to ensure public safety and long-term viability of commercial shellfish markets. Maximum concentration levels for MBTXs in bivalves have been established by many countries. European Union Regulations (EC) No 853/2004<sup>1</sup> and (EU) 786/2013<sup>2</sup> define permitted limits of lipophilic MBTXs in live bivalve mollusks (Table 1). As of 2015, LC-MS/MS is mandatory for official controls.

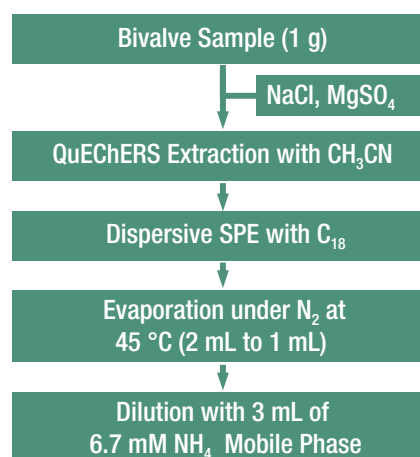
Table 1. Permitted limits for marine biotoxins according to EU regulations.<sup>1,2</sup>

| Marine Biotoxins                                  | Permitted Limit                       |
|---|---------------------------------------|
| Okadaic acid<br>Dinophysistoxins<br>Pectenotoxins | 160 µg/kg of okadaic acid equivalents |
| Yessotoxins                                       | 3.75 mg/kg of yessotoxins equivalents |
| Azaspiracids                                      | 160 µg/kg of azaspiracids equivalents |

## Experimental

### Sample Preparation

A sample to acetonitrile to C<sub>18</sub> proportion of 1:5:0.025 produced the best recovery of analytes. To detect and quantify the total content of the okadaic acid group toxins, an additional hydrolysis step was required to transform the esters of the toxins.



### LC-MS Conditions

UHPLC analysis was performed using the Thermo Scientific™ Accela™ UHPLC system with a C18 column (100 × 2.1 mm, 1.7 µm). Mobile phases were (A) 6.7 mM ammonia aqueous solution and (B) acetonitrile. The injection volume was 5 µL, and the flow rate was 0.25 mL/min. A Thermo Scientific™ Q Exactive™ hybrid quadrupole-Orbitrap mass spectrometer, operating in tandem mass spectrometry mode, was used for analysis.

|                               |   |
|-------------------------------|---|
| Spray voltage:                | 3.5 kV (positive ionization),<br>1.5 kV (negative mode) |
| Sheath gas (N <sub>2</sub> ): | 30 arb. units   |
| Capillary temperature:        | 300 °C  |
| S-lens:                       | 50 arb. units   |
| Heater temperature:           | 200 °C  |
| Normalized collision energy:  | 15% to 34% for the various<br>compounds                 |
| Resolution:                   | 70,000 ( <i>m/z</i> 200, FWHM)                          |
| Scan rate:                    | 2 Hz  |
| Automatic gain control:       | 5 × 104   |
| Maximum injection time:       | 35–100 ms   |

Thermo Scientific™ TraceFinder™ software version 3.1 was used for data processing.

## Data

### Evaluation of QuEChERS Method

The extraction method was tested on canned and fresh samples of various bivalve species. A comparison of the trueness of the spiked samples at 25 µg/kg in the different matrices showed no variations. In addition, the internal standard responses were the same in these samples.

### Validation of Complete QuEChERS UHPLC-HRMS Method

For each analyte, the molecular ion and one or two product ions were acquired with a mass accuracy better than 5 ppm (Figure 1). Quantification was performed using SMMS, with epinomectin as internal standard. A single SMMS calibration curve built from blank mussel samples was shown to be suitable for the analysis of different bivalve species, which is of importance to laboratories that analyze multiple types of shellfish.

The method was validated in mussel muscle. Specificity was evaluated by analyzing blank mussel samples, which were found to be free of endogenous interferences. In addition, blank clam, cockle, and razor clam samples were investigated and, again, no interfering peaks were observed at the retention time of the analytes.

The lower limit of the method was established as 25 µg/kg for the set of analytes. Linearity studies showed excellent results over the range 25–350 µg/kg, with coefficients of regression ( $r^2$ ) greater than or equal to 0.9998. Residuals were below 25% at 25 µg/kg and below 15% for higher concentrations. Recovery ranged between 79% and 114%, while relative standard deviations ranged between 1.5% and 18.7%.

To further test the method, two mussel certified reference materials were analyzed and the data showed good agreement (Table 2).



Working group from Laboratori de l'Agència de Salut Pública de Barcelona in Spain

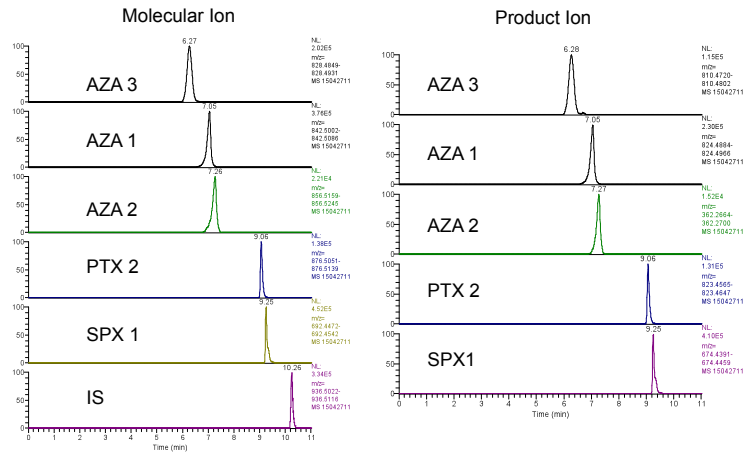


Figure 1. Extracted molecular ion chromatograms and extracted product ion chromatograms for a mixture of MBTXs.

Table 2. Analysis of certified reference materials.

| MBTX              | Certified Value       |                 | Determined Amount     |              |
|-------------------|-----------------------|-----------------|-----------------------|--------------|
|                   | Concentration (mg/kg) | Uncertainty (%) | Concentration (mg/kg) | Trueness (%) |
| Okadaic acid      | 0.361                 | 0.034           | 0.336                 | 93.1         |
| Dinophysistoxin-1 | 0.206                 | 0.019           | 0.189                 | 91.8         |
| Dinophysistoxin-3 | 0.283                 | 0.054           | 0.287                 | 101.4        |
| Azaspiracid-1     | 1.16                  | 0.100           | 1.06                  | 91.4         |
| Azaspiracid-2     | 0.273                 | 0.024           | 0.256                 | 93.8         |
| Azaspiracid-3     | 0.211                 | 0.023           | 0.222                 | 105.2        |

## References

1. Regulation (EC) No 853/2004 of the European Parliament and of the Council of 29 April 2004 Laying Down Specific Hygiene Rules for Food of Animal Origin, OJ L 139, 30.4.2004, p. 55.
2. Commission Regulation (EU) No 786/2013 of 16 August 2013 Amending Annex III to Regulation (EC) No 853/2004 of the European Parliament and of the Council as Regards the Permitted Limits of Yessotoxins in Live Bivalve Molluscs, OJ L 220, 17.8.2013, p. 14.

[www.thermofisher.com](http://www.thermofisher.com)

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific 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** +43 1 333 50 34 0  
**Australia** +61 3 9757 4300  
**Austria** +43 810 282 206  
**Belgium** +32 53 73 42 41  
**Canada** +1 800 530 8447  
**China** 800 810 5118 (free call domestic)  
 400 650 5118

**Denmark** +45 70 23 62 60  
**Europe-Other** +43 1 333 50 34 0  
**Finland** +358 10 3292 200  
**France** +33 1 60 92 48 00  
**Germany** +49 6103 408 1014  
**India** +91 22 6742 9494  
**Italy** +39 02 950 591

**Japan** +81 45 453 9100  
**Korea** +82 2 3420 8600  
**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  
**Norway** +46 8 556 468 00

**Russia/CIS** +43 1 333 50 34 0  
**Singapore** +65 6289 1190  
**Spain** +34 914 845 965  
**Sweden** +46 8 556 468 00  
**Switzerland** +41 61 716 77 00  
**UK** +44 1442 233555  
**USA** +1 800 532 4752

**Thermo**  
 SCIENTIFIC

A Thermo Fisher Scientific Brand