

The Determination of Organotins in Water Using Triple Quadrupole GC-MS/MS

Inge de Dobbeleer, Thermo Fisher Scientific, Breda, Netherlands
Hans-Joachim Huebschmann, Thermo Fisher Scientific, Bremen, Germany
Anton Mayer, Joachim Gummersbach, Thermo Fisher Scientific, Dreieich, Germany

Key Words

- Comparison of EI-SIM and EI-SRM
- Detection Limits
- Organotin
- PTV
- Timed-SRM

Introduction

The European Union Water Framework Directive seeks to lower detection and reporting limits for many contaminants.¹ One class of these is organotins. These include mono-, di-, tri- and tetrabutyl and triphenyl tin compounds. Tributyltin compounds are considered the most hazardous, and several studies have shown the effect on shell malformation of oysters, imposex of marine snails and reduced resistance to infections.^{2,3} Exposure to these compounds has been reported to cause acute kidney and central nervous disorders in humans.

Organotins are used widely in industry for a variety of applications, including as antifouling agents on underwater structures for both tributyl and triphenyltin; and triphenyltin as a fungicide in crop protection. Mono- and dibutyltin have uses as stabilizers in plastics and catalysts in soft foam production.^{1,2}

Concentration levels of organotin compounds in seawater are in the ppb to ppt range. Higher concentrations are found in sediments and biological samples. Organotin compounds are lipophilic and get absorbed into adipose tissue. They can also be adsorbed onto particulate matter. The toxicity of these compounds at low concentrations drives the requirement for accurate and sensitive analytical methods for their detection, quantitation and research for less-toxic replacements.^{4,5,6}

The method described in this paper uses the Thermo Scientific TSQ Quantum XLS triple quadrupole GC-MS/MS system to provide detection and quantitation limits that go below regulatory requirements with the use of timed selected reaction monitoring (t-SRM).

Sample Preparation

Derivatization and extraction of organotin compounds from water samples:^{7,8,9} From the water sample, 400 mL was taken and the pH was adjusted to 5 using a 1 M acetic acid/sodium acetate buffer prior to derivatization. For the ethylation of the organotin compounds, a 2% w/v sodium tetraethyl borate solution in 0.1 M NaOH was added.

Extraction was performed by adding pentane, followed by shaking for at least ten minutes. The organic phase was transferred and slowly evaporated to 400 µL. A 3 µL injection was made on the GC.

Method

All sample analyses were carried out using the TSQ Quantum XLS™ GC-MS/MS system, equipped with a Thermo Scientific TRACE GC Ultra gas chromatograph. The TRACE GC Ultra™ was configured with a programmable temperature vaporizer (PTV) injector. Sample introduction was performed using the Thermo Scientific TriPlus AS liquid autosampler. The capillary column was a Thermo Scientific TraceGOLD TG-5MS column (5% phenyl film) of 30 m length, 0.25 mm inner diameter and 0.25 µm film thickness. An uncoated 2 m DMTPS deactivated 0.53 mm ID pre-column was used as a guard column. Selected instrument parameters are shown in Table 1.

TRACE GC Ultra

| | |
|-------------------|--|
| Injection Volume: | 3 µL injection |
| Liner: | Thermo Scientific SurfaSil treated glass liner p/n PTV straight liner: 45352054 p/n SurfaSil™: TS-42800 (*) |
| Carrier Gas: | He, constant flow, 1.4 mL/min |
| Column Type: | TraceGOLD™ TG-5MS column (5% phenyl film) of 30 m length, 0.25 mm inner diameter and 0.25 µm film thickness. p/n: 26099-1420 |
| Oven Temperature: | Initial 45 °C, Hold 2 min, Ramp 55.0 °C/min – 175 °C, Ramp 35.0 °C/min – 300 °C Hold 2.0 min |
| Transfer Line: | 300 °C |

TRACE GC Ultra PTV Injector

| | |
|-----------------------|--|
| Injector Temperature: | 50 °C, hold 0.1 min, splitless injection |
| PTV Transfer: | 8 °C/sec to 280 °C, hold 1 min |
| PTV Cleaning Step: | 350 °C, 11 min, clean flow 50 mL/min |

TSQ Quantum XLS Mass Spectrometer

| | |
|---------------------|------------------------------|
| Source Temperature: | 250 °C, closed EI ion volume |
| Ionization: | EI, 70 eV |
| Emission Current: | 50 A |
| Resolution: | 0.7 Da Q1, Q3 |
| Collision Gas: | Argon, 1.0 mTorr |

Table 1: Selected instrumental conditions for the TRACE GC Ultra and TSQ Quantum XLS mass spectrometer

* Treating liners with SurfaSil is described in instructions on this website www.separatedbyexperience.com/literature/pierce_instructions.pdf

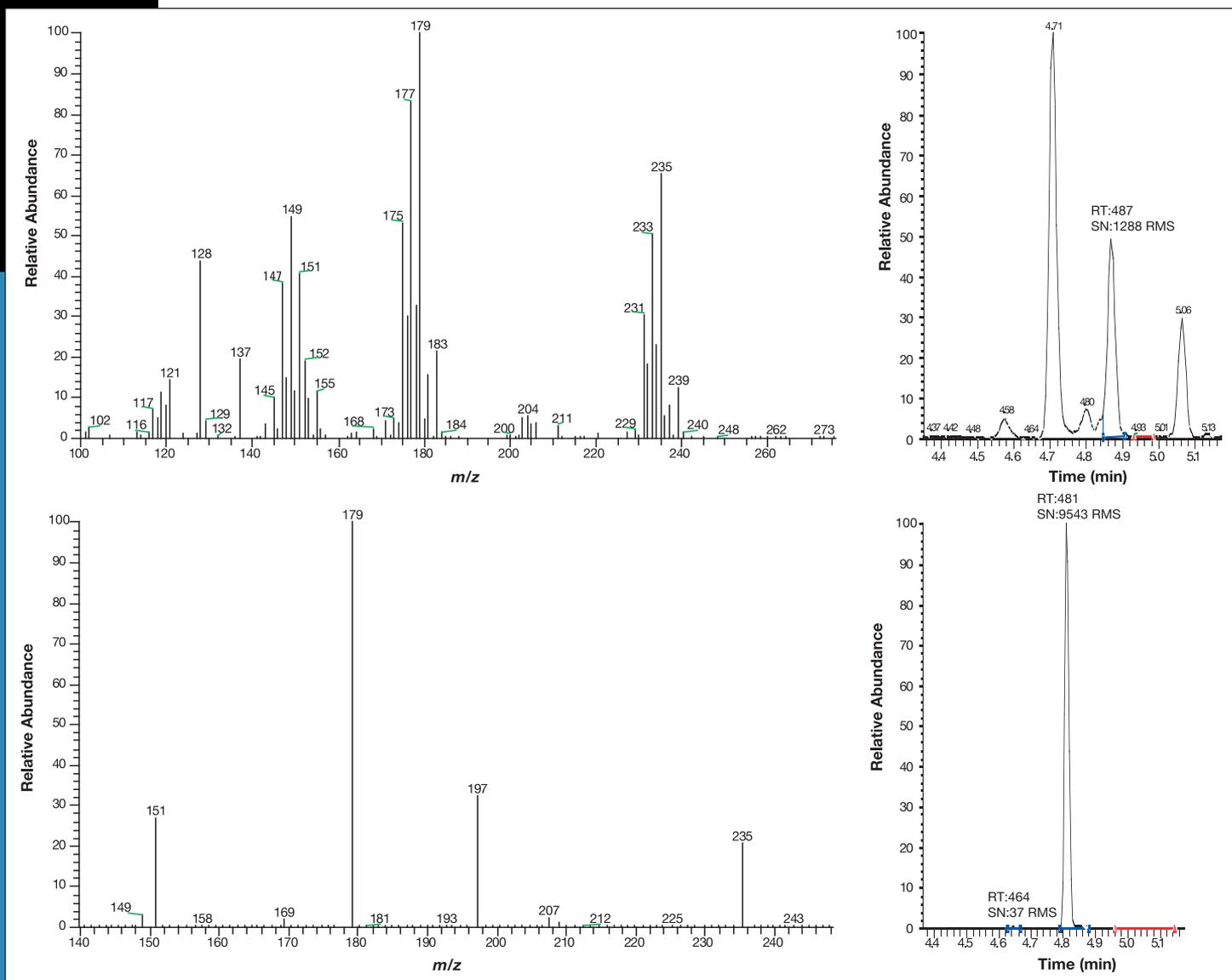


Figure 1: Clockwise: full scan spectrum of monobutyltin; S/N of m/z 235.08 in full scan mode; S/N of 235.08 > 178.95 in SRM mode; product spectrum of monobutyltin

EI-SRM Method Development

A higher-level standard was used for optimizing the transitions on the TSQ Quantum XLS operated in electron ionization (EI)-SRM mode. An example of one SRM optimization is given in Figure 1. The same procedure was followed for all organotin compounds, resulting in at least two transitions each.

The transitions that were found in the method development process were added to the t-SRM table, shown in Table 2. The start and stop times show partial overlap. This function of t-SRM allows the instrument to monitor SRM transitions more efficiently by only monitoring at the specific elution times for the compounds analyzed.

Results and Discussion

A single-point calibration was used at a level of 10 ng/L. The concentration is calculated on the actual levels in the water sample, not in neat solutions. The result shown in Figure 2 demonstrates the instrument's capability to reach 0.05 ng/L of organotins in the water sample and below. This is four times lower than the annual average stated in the Water Framework Directive. The actual amount injected on column was 0.2 pg for each organotin compound. Calculations were performed using tripropyltin as an internal standard. The ratio deviation was calculated and is shown in Table 3.

Peak identification is a key element in mass spectrometry. In SRM mode, at least two product ion transitions for each compound are needed, with peaks positively identified if the ion ratios in standards and samples are the same. The deviation for the calculated ion ratios must therefore be established by running their respective standards. Demonstration of ion ratios is shown in Table 3.

| Name | Parent Ion | Product Ion | Collision Energy (V) | Start Time (min) | Stop Time (min) |
|----------------------------------|------------|-------------|----------------------|------------------|-----------------|
| Monobutyltin | 235.08 | 150.98 | 6 | 7.44 | 8.44 |
| Monobutyltin | 233.08 | 176.95 | 6 | 7.44 | 8.44 |
| Monobutyltin | 235.08 | 178.95 | 6 | 7.44 | 8.44 |
| Tripropyltin internal standard | 249.08 | 164.91 | 8 | 8.06 | 9.71 |
| Tripropyltin internal standard | 245.08 | 160.91 | 8 | 8.06 | 9.71 |
| Tripropyltin internal standard | 247.08 | 162.91 | 8 | 8.06 | 9.71 |
| Tetrapropyltin internal standard | 249.08 | 164.91 | 8 | 8.06 | 9.71 |
| Tetrapropyltin internal standard | 245.08 | 160.91 | 8 | 8.06 | 9.71 |
| Tetrapropyltin internal standard | 247.08 | 162.91 | 8 | 8.06 | 9.71 |
| Dibutyltin | 261.03 | 205.03 | 8 | 8.92 | 9.92 |
| Dibutyltin | 263.03 | 150.98 | 8 | 8.92 | 9.92 |
| Dibutyltin | 263.03 | 207.03 | 8 | 8.92 | 9.92 |
| Monoheptyltin internal standard | 178.95 | 150.98 | 8 | 10.05 | 11.05 |
| Monoheptyltin internal standard | 275.07 | 176.95 | 8 | 10.05 | 11.05 |
| Monoheptyltin internal standard | 277.07 | 178.95 | 8 | 10.05 | 11.05 |
| Tributyltin | 287.09 | 174.94 | 8 | 10.17 | 12.24 |
| Tributyltin | 291.08 | 235.08 | 8 | 10.17 | 12.24 |
| Tributyltin | 289.09 | 176.95 | 8 | 10.17 | 12.24 |
| Monooctyltin | 287.09 | 174.94 | 8 | 10.17 | 12.24 |
| Monooctyltin | 289.09 | 176.95 | 8 | 10.17 | 12.24 |
| Monooctyltin | 291.08 | 178.95 | 8 | 10.17 | 12.24 |
| Tetrabutyltin | 287.09 | 174.94 | 8 | 10.17 | 12.24 |
| Tetrabutyltin | 289.09 | 176.95 | 8 | 10.17 | 12.24 |
| Tetrabutyltin | 291.08 | 235.08 | 8 | 11.24 | 12.24 |
| Diheptyltin internal standard | 275.07 | 176.95 | 8 | 13.21 | 14.21 |
| Diheptyltin internal standard | 277.07 | 178.95 | 8 | 13.21 | 14.21 |
| Diheptyltin internal standard | 245.08 | 145.98 | 8 | 13.21 | 14.21 |
| Dioctyltin | 261.03 | 148.98 | 8 | 14.44 | 15.44 |
| Dioctyltin | 263.03 | 150.98 | 8 | 14.44 | 15.44 |
| Dioctyltin | 375.17 | 263.03 | 8 | 14.44 | 15.44 |
| Triphenyltin | 196.94 | 119.9 | 18 | 15.71 | 16.71 |
| Triphenyltin | 349.15 | 194.98 | 22 | 15.71 | 16.71 |
| Triphenyltin | 351.15 | 196.94 | 22 | 15.71 | 16.71 |
| Tricyclohexyltin | 233.08 | 150.98 | 8 | 15.78 | 16.78 |
| Tricyclohexyltin | 287.09 | 205.03 | 8 | 15.78 | 16.78 |
| Tricyclohexyltin | 315.04 | 150.98 | 8 | 15.78 | 16.78 |

Table 2: Transitions for EI-SRM

The peaks were all easily detected at the lowest level and confirmed on two transitions. A fast method was used to allow a large number of samples to be processed per day. Figure 2 displays extracted ion chromatograms for all quantifier peaks at the 0.2 pg injected level.

| Injection | Ion Ratio in % |
|----------------|----------------|
| 1 | 72.84 |
| 2 | 66.29 |
| 3 | 70.73 |
| 4 | 70.87 |
| 5 | 70.18 |
| 6 | 72.84 |
| Average | 70.62 |
| Stdev | 2.40 |
| RSD | 3.4% |

Table 3: Ion ratio in percentage for monobutyltin at 10 ng/L; the ratio between 233.08>176.95 and 235.08>178.95 was calculated

Conclusion

This method demonstrates the ability of the TSQ Quantum XLS system to exceed the EU Water Framework Directive requirements for the detection and quantitation of organotins.^{1,10} The t-SRM functionality gives the instrument the ability to automatically determine optimal SRM transition times, even allowing for partially overlapping SRM transitions. This method demonstrates detection and quantitation levels to 0.05 ng/L, exceeding the EU Directive's Annual Allowable average of 0.2 ng/L. The methodology provided here offers the operator an easier implementation of this method in the laboratory.

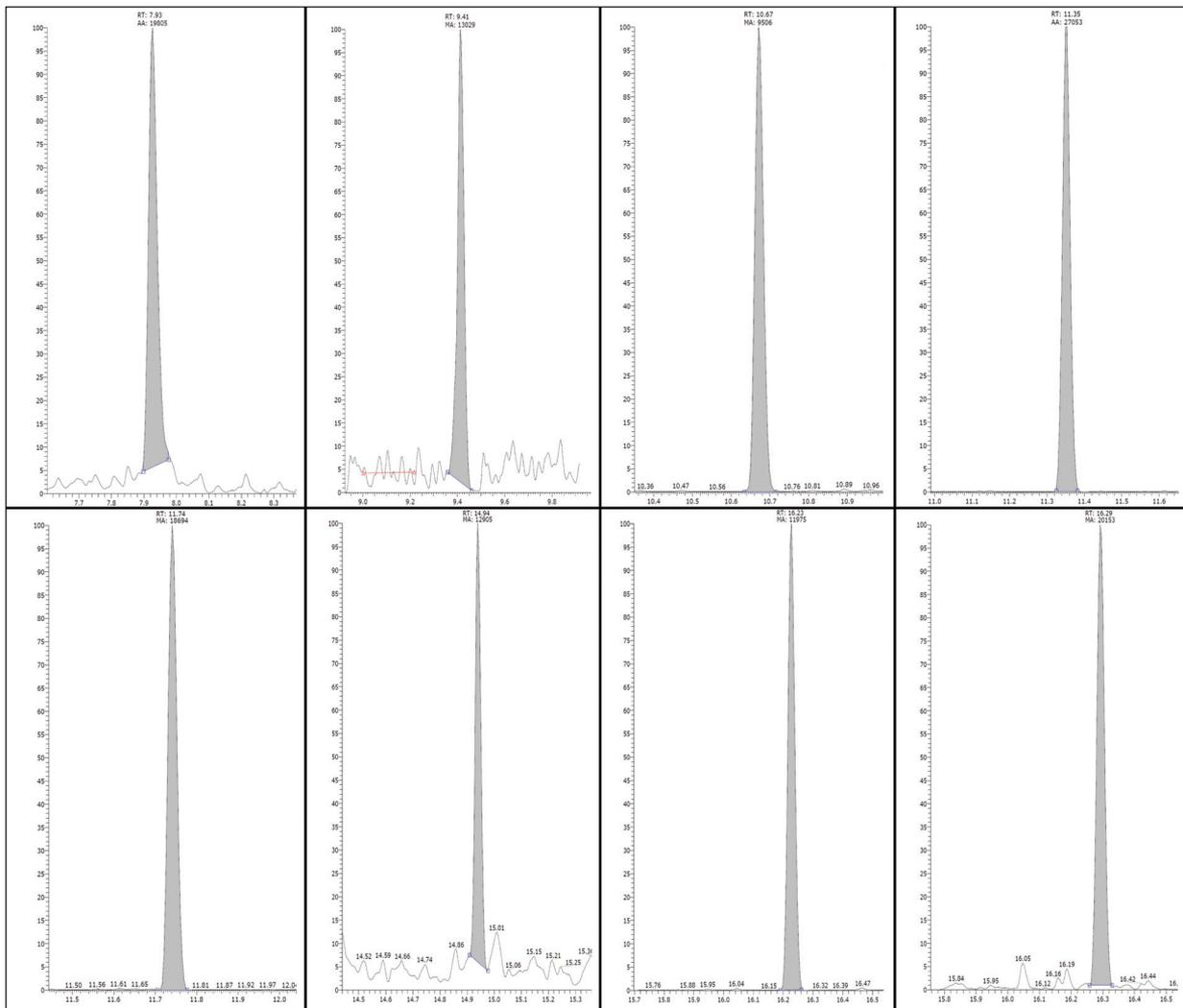


Figure 2: Compound peaks at 0.05 ng/L (0.2 pg on column). The peak names are listed in Table 4. Compounds 1 through 4 are on the top, compounds 5 through 8 are on the bottom (left to right).

| Compound | RT | Transition | Area | S/N |
|--------------------|-------|---------------|-------|------|
| 1 Monobutyltin | 7.93 | 235.08>178.95 | 19805 | 103 |
| 2 Dibutyltin | 9.42 | 263.03>207.03 | 13029 | 47 |
| 3 Tributyltin | 10.67 | 289.09>176.95 | 9506 | 784 |
| 4 Monoctyltin | 11.35 | 289.09>176.95 | 27053 | 2458 |
| 5 Tetrabutyltin | 11.74 | 289.09>176.95 | 18694 | 1651 |
| 6 Dioctyltin | 14.94 | 261.03>148.98 | 12905 | 80 |
| 7 Triphenyltin | 16.23 | 351.15>196.94 | 11975 | 923 |
| 8 Tricyclohexyltin | 16.29 | 233.08>150.98 | 20153 | 187 |

Table 4: Signal-to-noise ratio for all of the organotins at 0.05 ng/L

References

1. Directive 2008/105/EC on Environmental Quality Standards in the Field of Water Policy
2. OSPAR Background document on Organic Tin Compounds
3. Gunnar Nordberg, Handbook on the Toxicology of metals, Third edition p854
4. M. Takeuchi, K. Mizuishi and T. Hobo, *Analytical Sciences*, 2000, 16, 349
5. R. Morabito and Ph. Quevaullier, *Spectroscopy Europe*, 2002, 4, 18
6. Quantification of Organotin Compounds Using S/SL Injection and LVI-GC-MS by R.A. Correa C. and D. Claus
7. G. Centineo, P. Rodriguez-Gonzalez, E. Blanco Gonzalez, J.I. Garcia Alonso, and A. Sanz-Medel, Simultaneous determination of mono-, di- and tributyltin in environmental samples using isotope dilution gas chromatography-mass spectrometry, *J. Mass Spectrom.*, 2004, 39, 485-494
8. Analysis of Organotins by LVI GCMS SIM, *Thermo Fisher Application Note 10305*, by Jessie Butler, Eric Phillips
9. Determination of Organo-Tin Species in sediment extracts using GC coupled with X-series ICP-MS, *Thermo Application Note 40698*, by Dr. Martin Nash, Dr. Eva Krupp
10. Council Directive 96/23/EC concerning the performance of analytical methods and the interpretation of results

www.thermoscientific.com

Legal Notices: ©2011 Thermo Fisher Scientific Inc. All rights reserved. All 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 manners 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.

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

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



Thermo Fisher Scientific, Austin, TX USA is ISO Certified.

AN52099_E 05/11M