

Sensitive and Fast Determination of Endothall in Water Samples by IC-MS/MS

Leo Wang and Bill Schnute,
Thermo Fisher Scientific, Sunnyvale, CA, USA

Introduction

Endothall is a widely used herbicide for both terrestrial and aquatic weeds. Major uses of endothall include defoliation of cotton, the control of aquatic weeds and algae, and as a desiccating agent for lucerne and potatoes. Human exposure to endothall in excess of the maximum contamination level (MCL) may cause gastrointestinal problems. Endothall is regulated by the United States Environmental Protection Agency (EPA) with an MCL at 0.1 mg/L or 100 ppb for drinking water;¹ and the California EPA developed the Public Health Goal (PHG) of 0.58 mg/L or 580 ppb for this compound.²

Current analytical methods for quantitation of endothall in water samples are described in EPA method 548.1³ as gas chromatography with mass spectrometry or flame ionization detection (GC-MS or GC-FID). These methods involve ion-exchange solid phase extraction, sample enrichment, and dimethyl ester derivatization, followed by a 20 min GC separation and MS or FID detection.

This study describes the direct analysis of trace-level endothall in water samples by ion chromatography mass spectrometry (IC-MS). Water samples were directly injected without labor-intensive sample preparation and chromatographic separation was achieved in 10 min, thus significantly improving method throughput. An MS/MS instrument was operated in selected reaction monitoring (SRM) mode, requiring minimal sample cleanup and ensuring highly sensitive (low ppb) and selective quantitation. Isotope labeled glutaric acid (glutaric acid- d_6) was used as an internal standard to ensure quantitation accuracy. This method has been successfully used for quantification of endothall in various water matrices including fresh creek water, lake water, and high salt content lake water. Method performance parameters such as linearity, calibration range, precision, accuracy, detection limits, and recovery were evaluated and will be presented in this application note



Equipment

- Thermo Scientific™ Dionex™ ICS-2000 or ICS-2100 Reagent-Free™ Ion Chromatography (RFIC™) system
- Thermo Scientific™ TSQ Quantum Access™ Triple Stage Quadrupole Mass Spectrometer
- Thermo Scientific™ XCalibur™ 2.1 software with integrated Thermo Scientific™ Dionex™ DCMSLink™ Software Plug-In 2.8 for instrument control, data acquisition and processing, and report generation
- Two Thermo Scientific Dionex AXP-MS auxiliary pumps

Conditions																	
Chromatographic																	
Columns:	Thermo Scientific™ Dionex™ IonPac™ AG16/AS16 hydroxide selective anion-exchange set (2 mm)																
Eluent Source:	Thermo Scientific Dionex EGC II KOH Eluent Generator Cartridge with Thermo Scientific Dionex CR-ATC Continuously-Regenerated Anion Trap Column, 2 mm																
Eluent:	Hydroxide gradient <table border="1"> <thead> <tr> <th>Min</th> <th>Concn</th> </tr> </thead> <tbody> <tr> <td>-4.0</td> <td>15 Mm</td> </tr> <tr> <td>0.0</td> <td>15</td> </tr> <tr> <td>5.0</td> <td>15</td> </tr> <tr> <td>6.0</td> <td>80</td> </tr> <tr> <td>9.0</td> <td>80</td> </tr> <tr> <td>9.5</td> <td>15</td> </tr> <tr> <td>10.0</td> <td>5</td> </tr> </tbody> </table>	Min	Concn	-4.0	15 Mm	0.0	15	5.0	15	6.0	80	9.0	80	9.5	15	10.0	5
Min	Concn																
-4.0	15 Mm																
0.0	15																
5.0	15																
6.0	80																
9.0	80																
9.5	15																
10.0	5																
Flow Rate:	400 µL/min																
Temperature:	30 °C																
Solvent:	200 µL/min acetonitrile delivered by a Dionex AXP-MS pump																
Detector 1:	Suppressed Conductivity with Thermo Scientific™ Dionex™ ASRS™ 300 Anion Self-Regenerating Suppressor,™ 2 mm (external water at 0.5 mL/min delivered by a Dionex AXP-MS pump)																
Detector 2:	TSQ Quantum Access Mass Spectrometer																

Mass Spectrometric			
Interface:	Negative Electrospray Ionization (ESI)		
Spray Voltage:	3500 V		
Sheath Gas:	50 Arbitrary units		
Auxiliary Gas:	30 Arbitrary units		
Capillary Temp.:	350 °C		
Collision Gas:	Argon at 1.5 mTorr		
SRM Acquisition:	4.2 to 6 min		
Operating Mode:	SRM		
Analyte	Parent Ion (<i>m/z</i>)	Product Ion (<i>m/z</i>)	Collision Energy (V)
Endothall-1	185	141	17
Endothall-2	185	123	19
Glutaric Acid-d ₆	137	74	21

These conditions apply to Figure 3.

Reagents and Standards

Endothall standard solution, 1 mg/mL in methanol (AccuStandard)

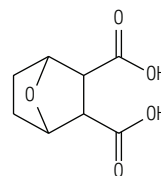
Isotope labeled internal standard (IStd) glutaric acid-d₆ (C/D/N Isotopes)

Seven-anion standard solution with various concentrations from 20 ppm (fluoride) to 150 ppm (sulfate and phosphate)

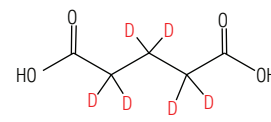
Acetonitrile (HPLC grade, Burdick & Jackson)

Deionized (DI) water, 18.2 MΩ-cm resistance or better (Millipore Corporation)

The chemical structures of endothall and isotope labeled internal standard glutaric acid-d₆ are shown in Figure 1.



Endothall
 Chemical Formula: C₈H₁₀O₅
 Exact Mass: 186.05
 Molecular Weight: 186.16



Glutaric Acid-d₆
 Chemical Formula: C₅H₂D₆O₄
 Exact Mass: 138.08
 Molecular Weight: 138.15

Figure 1. Chemical structures of studied compounds.

Standard Preparation

Dilute endothall standard in DI water to 10 ppm and 1 ppm as the two working standards used for preparation of calibration standards. Dilute the internal standard stock solution (1000 ppm) to 10 ppm for calibration standard preparation and for spiking unknown samples.

Prepare calibration standards by serial dilution from the working standards at 10 concentrations: 1 ppb, 2 ppb, 5 ppb, 10 ppb, 20 ppb, 50 ppb, 100 ppb, 200 ppb, 500 ppb, and 1000 ppb with the IStd spiked at 100 ppb for each concentration with the presence of diluted (100× dilution) seven common anions (fluoride 0.2 ppm, chloride 0.3 ppm, nitrite 1 ppm, bromide 1 ppm, nitrate 1 ppm, phosphate 1.5 ppm, and sulfate 1.5 ppm).

Sample Preparation

Ground water samples were collected from a local creek and the Lake Tahoe west shore. A lake water sample with a high salt content was kindly provided by a customer. These samples were spiked with IStd at 100 ppb and injected directly for IC-MS/MS quantification. (For samples with observed particulates, filtration is required prior to injection.) The salty water sample was diluted with DI water at a 1 to 10 ratio (v/v) due to the observed high concentration of ionic species.

Results and Discussion

Chromatography

As shown in Figure 1, endothall is a dicarboxylic acid. IC is the preferred separation technique for ionic or ionizable compounds; therefore, it was selected in this study for the separation of endothall from interferences. For IC-MS analysis, the major interferences are inorganic matrix ions such as chloride and sulfate, and other ionizable organics at high concentration. The goals of chromatographic separation were to achieve sufficient chromatographic resolution for endothall from common anions, and to have a short run time to improve method throughput.

In this study, several high-capacity anion-exchange columns were evaluated for their selectivity, namely the Dionex IonPac AS20, AS19, and AS16 columns. The Dionex IonPac AS16 column exhibited the best performance by: 1) providing a wide elution window for endothall between chloride and sulfate, and 2) requiring substantially less chromatographic run time than the Dionex IonPac AS19 column. Note that the Dionex IonPac AS19 column would be a better choice for simultaneous quantification of endothall and seven-anion profiling, as it provided complete chromatographic resolution for all analytes (Figure 2).

Mass Spectrometry

A TSQ Quantum Access triple quadrupole mass spectrometer was selected for sensitivity and selectivity, which allowed minimum sample preparation, such as cleanup and enrichment. A matrix diversion valve was used to divert high concentrations of inorganic anions to waste to prevent MS entrance fouling and to maintain long-term system stability. It is important to adjust the diversion window for specific matrices to ensure the analyte signal is collected by the mass spectrometer and inorganic salts are diverted to waste.

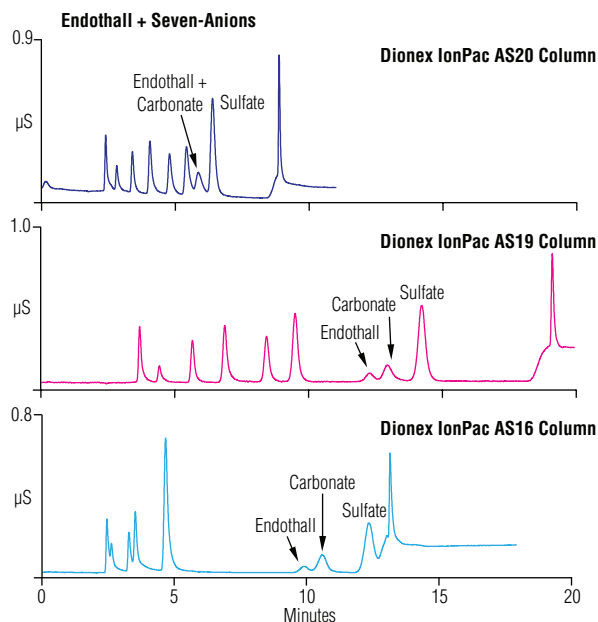


Figure 2. Column selection for endothall separation.

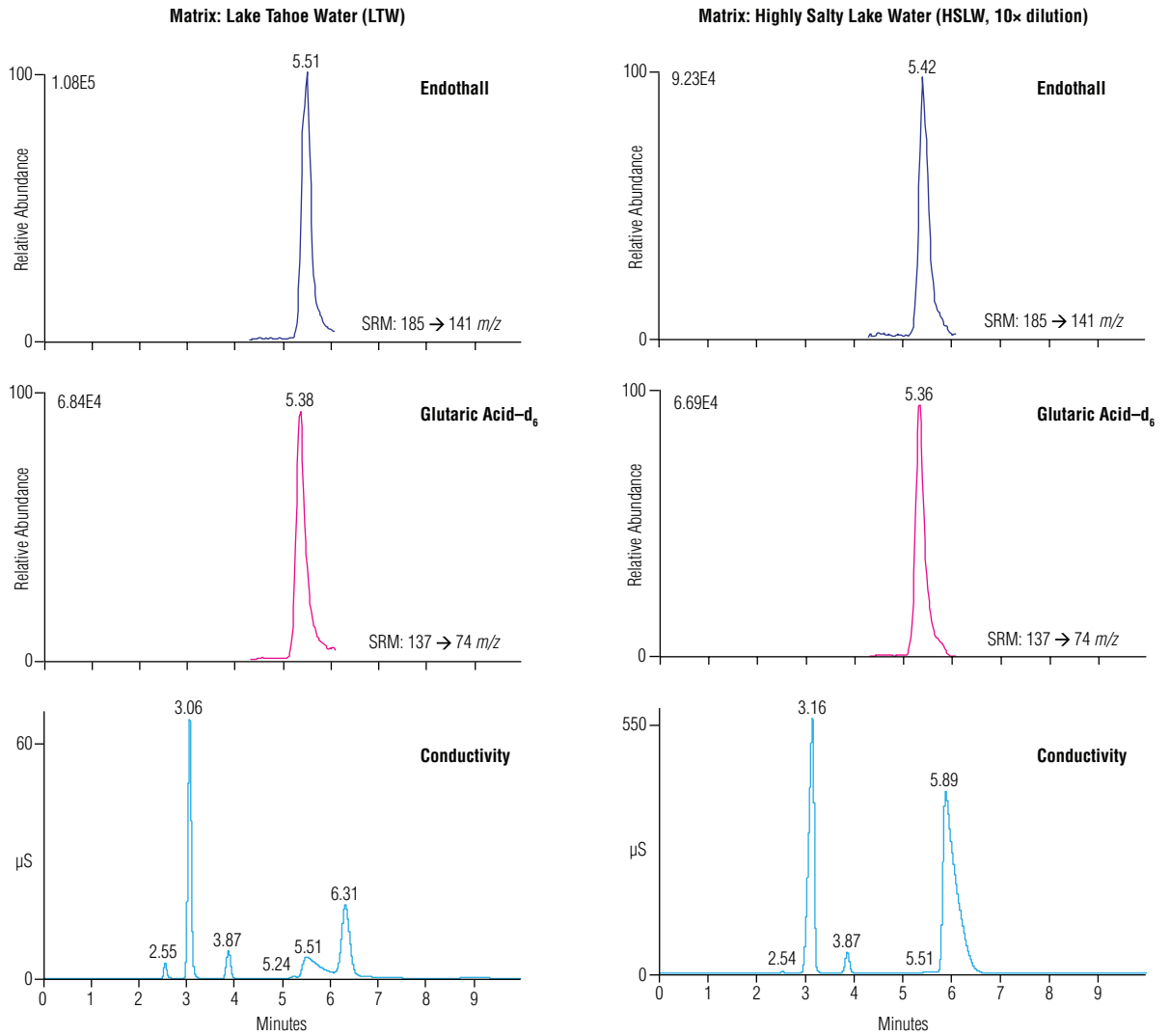


Figure 3. Conductivity and SRM chromatograms of endothall spiked in different samples: left, 5 ppb endothall spiked in LTW; right, 50 ppb endothall in salty lake water with 10-fold dilution.

As seen in Figure 3, two different matrices, LTW and HSLW, were spiked with endothall at 5 ppb and analyzed by this method. The HSLW is ~1000 ppm sulfate, (estimated by conductivity peak area), and the retention time of endothall in this 10× diluted matrix was slightly shifted (0.09 min earlier than in the LTW matrix). In addition, the retention time for sulfate shifted significantly earlier by 0.42 min; thus, the diversion window had to be adjusted accordingly to avoid MS entrance contamination. In rare situations where dilution of the original sample is not applicable, such as when endothall is present in very low concentration in a high-salt matrix, the gradient program will have to be adjusted, i.e., decrease the gradient slope from 5 to 9 min in order to increase the resolution of endothall from sulfate (as seen in Figure 2).

The quantitation of endothall was carried out in SRM mode, the precursor ion was observed as the deprotonated molecular ion $[M-H]^-$ at 185 m/z , and the predominant product ion was observed as $[M-H-CO_2]^-$ at 141 m/z and used as the quantitative SRM transition. Another characteristic product ion was also observed as $[M-H-CO_2-H_2O]^-$ at 123 m/z with less intensity and was used as a confirmative SRM transition. Isotope labeled glutaric acid- d_6 was used as the internal standard due to its similarity in chemical structure and chromatographic retention to endothall.

Because the IC eluent is 100% aqueous postsuppression, acetonitrile was introduced to assist desolvation/ionization and was delivered by an auxiliary pump at 0.2 mL/min.

Method Performance

Typical conductivity and SRM chromatograms are shown in Figure 3. Sulfate, chloride, and nitrate were observed as the main anionic species in tested ground water, and were chromatographically separated from the target analyte, endothall, and the internal standard, glutarate- d_6 . The MS/MS instrument operated in SRM mode provided sensitive and selective detection: each target analyte was observed as a single peak in each SRM channel with excellent intensity.

This method was used to determine endothall in various water samples mentioned in the Sample Preparation section, and no quantifiable level of target analyte was observed in any of the tested samples. Thus, these three matrices were used to evaluate the recovery from matrices.

Method performance was evaluated against quality parameters such as linearity, reproducibility, precision, accuracy, detection limits, and recovery from matrix. Calibration was performed by analyzing calibration standards at 10 concentrations from 1 to 1000 ppb using internal calibration with IStd at 100 ppb in each standard. Excellent linearity was achieved through three orders of magnitude with the coefficient of determination (r^2) at 0.9996. The calibration curve is shown in Figure 4, and the insert shows the calibration curve at lower concentrations

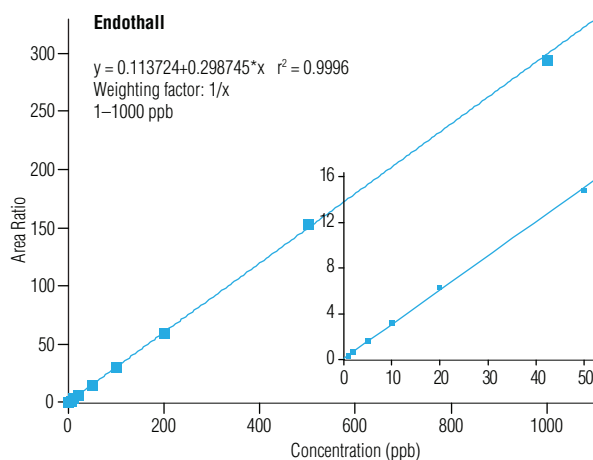


Figure 4. Calibration curve of endothall using glutaric acid- d_6 as an internal standard.

Table 1. Recovery of endothall from three matrices.

Sample	5 ppb			500 ppb		
	Mean	Standard Deviation	% Recovery	Mean	Standard Deviation	% Recovery
Creek Water	5.00	0.15	100	551	31.6	110
LTW	5.20	0.08	104	540	7.32	108
HSLW	4.76	0.06	95.1	535	50.3	107

Precision and accuracy were evaluated by seven replicate assays of a standard at 5 ppb with the calculated mean at 4.89 ppb (97.8% accuracy) and the standard deviation at 0.18 (3.63% RSD). The method detection limit (MDL) was calculated following the equation $MDL = s \times t$, where s is the standard deviation and t is the Student's t at 99% confidence interval. The calculated MDL is 0.56 ppb in DI water ($n = 7$).

Recovery was evaluated by spiking three samples—creek water, LTW, and a HSLW (10-fold dilution)—at two levels: 5 ppb and 500 ppb with IStd spiked at 100 ppb. The results are summarized in Table 1.

Conclusion

A fast and sensitive IC-MS/MS method was presented for the quantitative determination of trace level endothall in environmental water samples. The detection limit was estimated at 0.56 ppb, and linear response was observed from 1 ppb to 1000 ppb. The significantly improved sensitivity enables direct analysis of water samples without labor-intensive sample enrichment and derivatization. The total chromatographic run time was halved from the 20 min GC run in EPA Method 548.1.

References

1. United States Environmental Protection Agency, *Basic Information about Endothall in Drinking Water*. <http://water.epa.gov/drink/contaminants/basicinformation/endothall.cfm> (accessed September, 2010).
2. Pesticide and Environmental Toxicology Section, Office of Environmental Health Hazard Assessment, California Environmental Protection Agency, *Public Health Goal for ENDOTHALL in Drinking Water*, December 1997.
3. United States Environmental Protection Agency, *Method 548.1: Determination of Endothall in Drinking Water by Ion-Exchange Extraction, Acidic Methanol Methylation and Gas Chromatography/Mass Spectrometry*. Available from http://www.epa.gov/waterscience/methods/method/files/548_1.pdf (accessed January 31, 2011).

www.thermoscientific.com/dionex

©2014 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. Burdick & Jackson is a registered trademark of Honeywell International Inc. 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 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.



Thermo Fisher Scientific, Sunnyvale, CA
USA is ISO 9001:2008 Certified.

Africa +43 1 333 50 34 0
Australia +61 3 9757 4300
Austria +43 810 282 206
Belgium +32 53 73 42 41
Brazil +55 11 3731 5140
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 9 3291 0200
France +33 1 60 92 48 00
Germany +49 6103 408 1014
India +91 22 6742 9494
Italy +39 02 950 591

Japan +81 6 6885 1213
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
Sweden +46 8 556 468 00
Switzerland +41 61 716 77 00
Taiwan +886 2 8751 6655
UK/Ireland +44 1442 233555
USA +1 800 532 4752
AN70945_E 01/14S

Thermo
SCIENTIFIC

Part of Thermo Fisher Scientific