

US EPA Method 525.2 — Determination of Semivolatile Organics in Water by Solid-Phase Extraction and GC/MS Detection

INTRODUCTION

Method 525.2, determination of organic compounds in drinking water by SPE and capillary column gas chromatography/mass Spectrometry, is a drinking water method for the determination of a wide variety of semi-volatile organic compounds including PAHs, pesticides, and PCBs.

Method 525.2 is performed by loading a one-liter sample of water onto a conditioned C-18 SPE cartridge. The cartridge is dried and eluted with a combination of ethyl acetate (EtOAc) and methylene chloride (DCM). The extract is dried by passing through 5–7 g of fired sodium sulfate and concentrated under nitrogen to 1 mL. For this work, we used the optional cartridge SPE (6 mL syringe body, 1000 mg of C18 packing), and a Varian Saturn Ion Trap GC/MS.

AUTOTRACE EXTRACTION PROCEDURES

Loading Procedure (Method One):

- Step 1: Process 6 samples using the following procedure
- Step 2: Wash syringe with 2 mL of CH₃OH
- Step 3: Rinse column with 5 mL of EtOAc into solvent waste
- Step 4: Rinse column with 5 mL of DCM into solvent waste
- Step 5: Condition column with 10 mL of CH₃OH into solvent waste
- Step 6: Condition column with 10 mL of water into aqueous waste
- Step 7: Load 1000 mL of sample onto column
- Step 8: Dry column with gas for 10 minutes

Flow Rate:

Cond. Flow: 40 mL/min
Load Flow: 20 mL/min
Rinse Flow: 40 mL/min
Elute Flow: 5 mL/min
Cond. Air Push: 15 mL/min
Rinse Air Push: 20 mL/min
Elute Air Push: 5 mL/min

Eluting Procedure (Method Two):

Step 1: Process 6 samples using the following procedure

Step 2: Manually rinse the sample container with 7 mL to collect. Use Ethyl acetate to rinse the sample container. See Elution details section.

Step 3: Manually rinse sample container with 10 mL to collect. Use Dichloromethane to rinse the sample container. See Elution details section.

Step 4: Elute to soak 3 mL fraction using EtOAc

Step 5: Elute to collect 2 mL fraction into sample tube using EtOAc

Step 6: Elute to soak 3 mL fraction using DCM

Step 7: Elute to collect 2 mL fraction into sample tube using DCM

Flow Rate:

Cond. Flow: 40 mL/min
Load Flow: 5 mL/min
Rinse Flow: 40 mL/min
Elute Flow: 5 mL/min
Cond. Air Push: 15 mL/min
Rinse Air Push: 20 mL/min

Manual Rinse Details:

In Step 2, add 5 mL of EtOAc to sample container, rinsing the walls, and reposition tube, and press COND. Again in Step 3, add 5 mL of DCM to sample container, rinsing the walls, and reposition tube, and press COND. The extra volume pulled allows air to follow each rinsing. Note that the "Load Flow" parameter is different for the the eluting procedure (Method Two) and should be set at 5mL/min.

VARIAN SATURN ION TRAP CONDITIONS

Column: Restek Rtx[®]-5MS 0.25 mm ID by 30 m with 0.25 µm df.

Injection Temp.: 245 °C, splitless for 1 min.

Temp. Prog.: 70 °C (1 min), 50 °C/min, 130 °C (0.8 min), 12 °C/min, 180°, 7 °C/min, 240 °C, 12 °C/min, 320 C° (1.61 min) for total of 24 min.

Pressure Temp: 20 psig (1 min), -3.6 psig/min, 13.1 psig, 0.6 psig/min, 15.7 psig, 0.4 psig/min, 18.7 psig, 0.5 psig/min, 22.7 psig (1.2 min) for total of 24 min.

Table 1. Analytical Results of Spiked Samples

Compound	%REC.	%RSD	Compound	%REC.	%RSD
Chlordane	110%	7.8%	Methoxychlor	111%	4.5%
Toxaphene	92%	7.4%	Heptachlor	109%	3.4%
Aroclor 1248	104%	3.6%	Heptachlor epoxide	106%	2.8%
Atrazine	119%	12.8%	Hexachlorobenzene	120%	5.2%
Alachlor	109%	4.8%	Hexachlorocyclopentadiene	85%	7.6%
Simazine	142%	5.8%	Aldrin	104%	3.1%
Metolachlor	128%	2.8%	Dieldrin	106%	4.0%
Propachlor	147%	5.4%	DDT	91%	4.6%
Butachlor	102%	14.3%	Benzo(a)pyrene	71%	3.3%
Metribuzin	126%	7.2%	bis(2-ethylhexyl)adipate	120%	3.4%
Lindane	119%	4.1%	bis(2-ethylhexyl)phthalate	79%	2.7%
Endrin	111%	4.3%			

SUMMARY

Using AutoTrace® system effectively automates sample extraction using a cartridge.

Solid Phase Extraction provides reduced solvent consumption and sample handling when compared to traditional liquid-liquid extraction techniques. The AutoTrace instrument automates the conditioning loading rinse and eluting steps of Manual SPE that normally require extensive involvement by an operator. AutoTrace provides both Automation of the SPE steps and positive pressure flow of sample and solvent through the cartridges. Positive pressure flow of the sample and solvent ensure an uniform extraction and elution. These features contribute greatly to increased laboratory productivity and reliable results for determination of organic pollutants in drinking water.

ACKNOWLEDGEMENTS

Data submitted by David Larabee-Zierath of the University Hygienic Laboratory, Iowa's Environmental and Public Health Laboratory, Iowa City, IA.

AutoTrace is a registered trademark of Dionex Corporation.
Restek Rtx is a registered trademark of Restek.

Passion. Power. Productivity.

**Dionex Corporation**

1228 Titan Way
P.O. Box 3603
Sunnyvale, CA
94088-3603
(408) 737-0700

North America

U.S. (847) 295-7500
Canada (905) 844-9650

South America

Brazil (55) 11 3731 5140

Europe

Austria (43) 1 616 51 25 Benelux (31) 20 683 9768 (32) 3 353 4294
Denmark (45) 36 36 90 90 France (33) 1 39 30 01 10 Germany (49) 6126 991 0
Ireland (353) 1 644 0064 Italy (39) 02 51 62 1267 Sweden (46) 8 473 3380
Switzerland (41) 62 205 9966 United Kingdom (44) 1276 691722

Asia Pacific

Australia (61) 2 9420 5233 China (852) 2428 3282 India (91) 22 2764 2735
Japan (81) 6 6885 1213 Korea (82) 2 2653 2580 Singapore (65) 6289 1190
Taiwan (886) 2 8751 6655

www.dionex.com



LPN 2162 PDF 01/09
©2009 Dionex Corporation