

# Automated Solid-Phase Extraction (SPE) of Endocrine Disruptors in Water

## Introduction

Endocrine disrupting compounds (EDCs) include a wide range of compounds that have the potential to interfere with the body's endocrine system by disrupting the function of natural or endogenous hormones. Examples of potential EDCs include pesticides, such as DDT, dioxins and dioxin-like compounds, polychlorinated biphenyls, flame retardants, and plastic additives. Drinking water supplies need to be tested for the presence of EDCs because of their effect on hormones. The solid-phase extraction (SPE) method described in this Application Brief provides a quick and simple means of concentrating the components using controlled flow rates and volumes that will consistently produce good recoveries.

## Instrumentation Used for Sample Preparation

Thermo Scientific™ Dionex™ AutoTrace™ 280 Solid-Phase Extraction instrument.

## Sample Preparation

- A 500 mL sample of water is taken for the analysis.
- Screen for the presence of octylphenol, nonylphenol, octachlorostyrene, butylbenzylphthalate, diazinon, BPA, BHA, and BHT.
- Thermo Scientific™ Dionex™ SolEx™ SPE C18, 3 mL Cartridge, with 0.5 g of packing (P/N 074412).
- Dionex SolEx SPE C8, 3 mL Cartridge, with 0.5 g of packing (P/N 074413).
- The solvents methyltertbutylethylene (MTBE) and dichloromethane are used as eluents.
- The eluent can be dried using an in-line anhydrous sodium sulphate cartridge.

## Solvents

- Isopropyl alcohol (pesticide quality or equivalent)
- Cyclohexane (pesticide quality or equivalent)
- Extraction solvent: 2.5% (v/v) cyclohexane in isopropyl alcohol (IPA)

## Dionex AutoTrace 280 SPE Method and Conclusion

All solvent lines are purged and primed with solvent first. A maximum of five solvents can be used to run a wide range of methods. Independent lines separate the aqueous and organic waste solvents. The instrument sample rack has six positions that can take volumes from 10 mL to 2000 mL and six sample collection positions for the eluent. This method offers an automated rugged and reproducible solution for cleaning up the samples to concentrate and remove interferences.

No.	Method: Estimated time 1 h 44 min
1	Process six samples using the following method steps:
2	Condition the column with 5.0 mL of dichloromethane into solvent waste.
3	Condition column with 5.0 mL of MTBE into solvent waste.
4	Condition column with 5.0 mL of CH <sub>3</sub> OH into solvent waste.
5	Condition column with 5.0 mL of DI water into aqueous waste.
6	Load 500.0 mL of sample into column.
7	Rinse column with 5.0 mL of DI water into aqueous waste.
8	Wash syringe with 5.0 mL of CH <sub>3</sub> OH/MTBE 10:90.
9	Dry column with gas for 15.0 min.
10	Collect 5.0 mL fraction into sample tube using CH <sub>3</sub> OH/MTBE 10:90.
11	Collect 5.0 mL fraction into sample tube using dichloromethane.
12	Wash syringe with 5.0 mL of CH <sub>3</sub> OH.
13	End.

<b>Parameters</b>	
<b>Flow Rates</b>	
Cond Flow:	15.0 mL/min
Load Flow:	10.0 mL/min
Rinse Flow:	20.0 mL/min
Elute Flow:	5.0 mL/min
Cond Air Push:	15.0 mL/min
Rinse Air Push:	20.0 mL/min
Elute Air Push:	5.0 mL/min
<b>SPE Parameters</b>	
Push Delay:	5 sec
Air Factor:	1.0
Autowash Vol.:	1.00 mL
<b>Instrument Parameters</b>	
Max. Elution Vol.:	12.0 mL
Exhaust Fan On:	Yes
Beeper On:	Yes

### [www.thermoscientific.com/autotrace](http://www.thermoscientific.com/autotrace)

©2013 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. 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.

**Australia** +61 3 9757 4486  
**Austria** +43 1 333 50 34 0  
**Belgium** +32 53 73 42 41  
**Brazil** +55 11 3731 5140  
**China** +852 2428 3282

**Denmark** +45 70 23 62 60  
**France** +33 1 60 92 48 00  
**Germany** +49 6126 991 0  
**India** +91 22 6742 9494  
**Italy** +39 02 51 62 1267

**Japan** +81 6 6885 1213  
**Korea** +82 2 3420 8600  
**Netherlands** +31 76 579 55 55  
**Singapore** +65 6289 1190  
**Sweden** +46 8 473 3380



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

**Switzerland** +41 62 205 9966  
**Taiwan** +886 2 8751 6655  
**UK/Ireland** +44 1442 233555  
**USA and Canada** +847 295 7500

**Thermo**  
 SCIENTIFIC

Part of Thermo Fisher Scientific