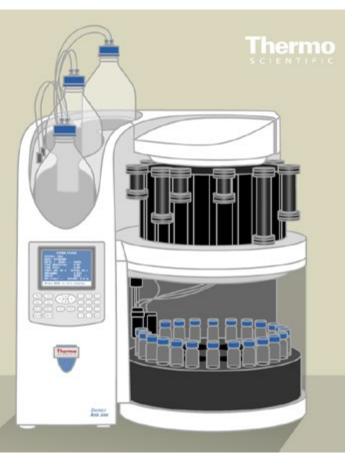


# Accelerated Solvent Extraction Environmental Applications Summary

Pesticides • Chlorinated Compounds • Persistent Organic Pollutants (POPs)



## Thermo Scientific Dionex ASE Accelerated Solvent Extractor Systems



Watch the video to learn more about the Thermo Scientific<sup>™</sup> ASE<sup>™</sup> Accelerated Solvent Extraction System.

Additional product information available at www.thermoscientific.com/samplepreparation

## Contents

<b>Introduction</b> The Accelerated Solvent Extraction System
Dioxins and Furans
Extraction of Polychlorinated Dibenzo- <i>p</i> -dioxins (PCDDs) and Polychlorinated Dibenzofurons (PCDFs) from Environmental Samples Using Accelerated Solvent Extraction
Herbicides
Extraction of Chlorinated Herbicides Using Accelerated Solvent Extraction
Organochlorinated Pesticides (OCPs) Extraction of Chlorinated Pesticides Using Accelerated Solvent Extraction
Organophosphorus Pesticides (OPPs) Extraction of Organophosphorus Pesticides Using Accelerated Solvent Extraction
Persistent Organic Pollutants (POPs)
Rapid Determination of POPs Using Accelerated Solvent Extraction
<b>Polybrominated Diphenyl Ethers (PBDEs)</b> Analyzing PBDEs in House Dust Samples with the Thermo Scientific TSQ Quantum XLS Ultra GC-MS/MS in El-SRM Mode
Polychlorinated Biphenyls (PCBs)
Accelerated Solvent Extraction of Polychlorinated Biphenyls from Polyurethane Foam Adsorbent Cartridges
Extraction of Polychlorinated Biphenyls from Environmental Samples Using Accelerated Solvent Extraction
References
Accelerated Solvent Extraction References

### **The Accelerated Solvent Extraction System**



#### Better extractions in less time using less solvent

Accelerated solvent extraction is an established technique used for fast, accurate sample preparation of air filters, soils, solids, and wastes. Coupled with rapid solvent reduction systems, the often labor-intensive steps of sample preparation can be automated in your laboratory. Accelerated solvent extraction eliminates many of the manual steps involved in preparing samples for analysis, which helps ensure increased reproducibility and accelerates the process significantly.

The Thermo Scientific<sup>™</sup> Dionex<sup>™</sup> ASE<sup>™</sup> Accelerated Solvent Extractor systems use a combination of elevated temperature and pressure with common solvents to increase the efficiency of the extraction process. The result is faster extraction times and a significant reduction in solvent use. Many of the organic solvents used in extractions boil at relatively low temperatures. This is a limitation to techniques such as Soxhlet or automated Soxhlet as the highest temperatures at

which extractions take place in these techniques will be the solvent's boiling point. If sufficient pressure is exerted on the solvent during the extraction, temperatures above the boiling point can be used. Therefore all of the advantages of working at elevated temperature can be realized even with solvents with relatively low boiling points. Operating at elevated pressures also accelerates the extraction process. Pumping solvent through a packed bed is easier at elevated pressures; pressurized solvent is forced into the pores of the sample matrix. Hence, the combination of elevated temperatures and pressures allows extractions to occur rapidly and completely.

When extractions are achieved at elevated temperatures, several factors contribute to improved speed, efficiency, and reduced solvent use: 1) solvent strength is higher, 2) diffusion rates are faster, 3) solvent viscosity is decreased, and 4) solute-matrix interactions (dipole attractions, Van der Waals forces, hydrogen bonding, etc.) are more easily disrupted allowing the analytes to be removed from the matrix.

For environmental applications, accelerated solvent extraction is proven to produce equivalent to or better than traditional methods. Accelerated solvent extraction is accepted for use in EPA Method SW-846 3545 for the extraction of pesticides and herbicides, PAHs and semivolatile compounds, PCBs, dioxins and furans, and explosive compounds. Accelerated solvent extraction is also accepted for use in EPA Method SW-846 6860 for the determination of perchlorate and in CLP OLM 04.2A for semivolatiles and pesticides. The Application Briefs compiled here show conditions for extracting pesticides, chlorinated compounds, and POPs from a variety of environmental samples. To view the complete Application Notes, visit our website at www.thermoscientific.com/ samplepreparation.



Thermo Scientific Dionex ASE 150 and ASE 350 Accelerated Solvent Extractor systems.

Extraction of Polychlorinated Dibenzo-*p*-dioxins (PCDDs) and Polychlorinated Dibenzofurans (PCDFs) from Environmental Samples Using Accelerated Solvent Extraction



#### Introduction

Regulatory agencies are concerned with the high toxicity of PCDDs and PCDFs; great efforts are expended to monitor their presence in environmental samples. The low concentrations at which these compounds are regulated and the strength of their binding to certain matrices pose special challenges to analytical chemists. Accelerated solvent extraction has been applied to the extraction of PCDD and PCDFs from ground chimney brick, urban dust, fly ash and sediment samples. Accelerated solvent extraction complies with U.S. EPA Method 3545 for these compounds.

U.S. EPA Method 3545A: Dioxins and furans from ground chimney brick, urban dust, and fly ash

#### Equipment

Dionex ASE 200 Accelerated Solvent Extractor\*

Gas Chromatograph (GC) with Mass Spectrometer (MS) (low resolution or high resolution MS-MS)

Low pressure LC system for sample clean-up Preparative HPLC for sample clean-up

#### **Solvents**

Toluene, acetic acid, hydrochloric acid. All solvents are pesticidegrade or equivalent and are available from Fisher Scientific.

\*Dionex ASE 150 and 350 can be used for equivalent results.

Solvent:	Toluene or toluene/acetic acid
	(5%, v/v) if HCl pretreatment currently used
Temperature:	175–200 °C
Pressure:	1500 psi**
Static Time:	5–15 min
Static Cycles:	2 or 3
Flush Volume:	60–70%
Purge Time:	60–100 s

\*\*Pressure studies show that 1500 psi is the optimum extraction pressure for all accelerated solvent extraction applications.

#### Analysis

GC-MS

#### Results

Left table: Average values (ng/kg) from ground chimney brick—comparison of Soxhlet vs. ASE. Right table: Average values (ng/kg) from urban dust—comparison of Soxhlet vs. ASE.

Group Totals	Soxhlet (n=1)	ASE <sup>ª</sup> (n=2)
Total T <sub>4</sub> CDD	182	325
Total P <sub>5</sub> CDD	175	221
Total H <sub>6</sub> CDD	86.7	81.7
Total H <sub>7</sub> CDD	221	217
Total O <sub>8</sub> CDD	445	314
Total T <sub>4</sub> CDF	333	419
Total P <sub>5</sub> CDF	146	179
Total H <sub>6</sub> CDF	65.9	122
Total H <sub>7</sub> CDF	13.2	29.4
Total 0 <sub>8</sub> CDF	n.d. (10)	n.d. (10)
Congeners	Soxhlet (n=1)	ASE <sup>ª</sup> (n=2)
2,3,7,8-T <sub>4</sub> CDD	3.3	3.2
1,2,3,7,8-P <sub>5</sub> CDD	11.8	13.1
1,2,3,4,7,8-H <sub>6</sub> CDD	9.8	8.0
1,2,3,6,7,8-H <sub>6</sub> CDD	11.5	9.5
1,2,3,7,8,9-H <sub>6</sub> CDD	n.d. (8)	n.d. (8)
1,2,3,4,6,7,8-H <sub>7</sub> CDD	113	107
2,3,7,8-T <sub>4</sub> CDF	12.5	18.6
1,2,3,7,8(+1,2,3,4,8)-P <sub>s</sub> CDF	9.9	12.0
2,3,4,7,8-P <sub>5</sub> CDF	13.9	18.1
1,2,3,4,7,8(+1,2,3,4,7,9)-H <sub>6</sub> CDF	18.7	23.7
1,2,3,6,7,8-H <sub>6</sub> CDF	10.7	15.8
2,3,4,6,7,8-H <sub>6</sub> CDF	3.3	8.7
1,2,3,7,8,9-H <sub>6</sub> CDF	n.d. (2)	n.d. (2)
1,2,3,4,6,7,8-H <sub>7</sub> CDF	13.2	29.4
1,2,3,4,7,8,9-H <sub>7</sub> CDF	n.d. (3)	n.d. (3)
Toxicity equivalent (NATO)	25.0	28.7
Toxicity equivalent (BgVV)	24.2	29.2

Group Totals	Soxhlet (n=1)	ASE <sup>ª</sup> (n=2)
Total T <sub>4</sub> CDD	440	530
Total P <sub>5</sub> CDD	900	940
Total H <sub>6</sub> CDD	1800	2000
Total H <sub>7</sub> CDD	2000	2100
Total O <sub>8</sub> CDD	2900	2600
Total T <sub>4</sub> CDF	2300	2600
Total P <sub>5</sub> CDF	4100	4300
Total H <sub>6</sub> CDF	4700	4700
Total H <sub>7</sub> CDF	2800	2600
Total O <sub>8</sub> CDF	2000	2000
Congeners	Soxhlet (n=1)	ASE <sup>ª</sup> (n=2)
2,3,7,8-T <sub>4</sub> CDD	6.0	6.0
1,2,3,7,8-P <sub>s</sub> CDD	52	57
1,2,3,4,7,8-H <sub>6</sub> CDD	46	52
1,2,3,6,7,8-H <sub>6</sub> CDD	120	130
1,2,3,7,8,9-H <sub>6</sub> CDD	97	1000
1,2,3,4,6,7,8-H <sub>7</sub> CDD	1000	820
2,3,7,8-T <sub>4</sub> CDF	160	180
1,2,3,7,8(+1,2,3,4,8)-P <sub>5</sub> CDF	430	470
2,3,4,7,8-P <sub>5</sub> CDF	390	390
1,2,3,4,7,8(+1,2,3,4,7,9)-H <sub>6</sub> CDF	1100	1100
1,2,3,6,7,8-H <sub>6</sub> CDF	540	570
2,3,4,6,7,8-H <sub>6</sub> CDF	400	360
1,2,3,7,8,9-H <sub>6</sub> CDF	42	42
1,2,3,4,6,7,8-H <sub>7</sub> CDF	2100	2000
1,2,3,4,7,8,9-H <sub>7</sub> CDF	140	120
Toxicity equivalent (NATO)	540	540
Toxicity equivalent (BgVV)	490	510

Values are corrected for recovery of <sup>13</sup>C-labeled surrogates.

n.d.=not detected. Detection limit, in ppt, given in parentheses. "Sum of two extractions of each sample. Values are corrected for recovery of <sup>13</sup>C-labeled surrogates. <sup>a</sup>Sum of two extractions of each sample.

# Extraction of Chlorinated Herbicides Using Accelerated Solvent Extraction

新教育計

#### Introduction

The use of accelerated solvent extraction in the extraction of chlorinated herbicides from solid wastes is more convenient, faster, and less solvent-intensive than previous methods. The procedures described in this Application Note meets the requirements for sample extraction as prescribed by U.S. EPA Method 3545. This method is applicable to the extraction of water-insoluble or slightly water-soluble chlorinated herbicides in preparation for gas chromatographic measurements. U.S. EPA Method 8150A with 3545A: Chlorinated herbicides from soil, sludge, and sediments

#### Equipment

Dionex ASE 200 Accelerated Solvent Extractor\* equipped with 11 mL or larger stainless steel extraction cells

Gas Chromatograph (GC) with Electron Capture Detector (ECD)

Vials for collection of extracts (40 mL, P/N 049465; 60 mL, P/N 049466)

\*Dionex ASE 150 and 350 can be used for equivalent results.

#### **Solvents and Reagents**

Diatomaceous Earth (DE) Dispersant for ASE<sup>™</sup>, 1 kg Bottle (P/N 062819)

Dichloromethane

Acetone

Phosphoric acid

All solvents are pesticide-grade or equivalent and are available from Fisher Scientific.

Average recovery of chlorinated herbicides from three soil types\*—ASE compared to wrist-shaking.

Chlorinated Herbicide Target Compound	Average Recovery (% of Shaking Method)
2,4-D	116.2
2,4-DB	112.9
2,4,5-T	106.6
2,4,5-TP	117.4
Dalapon	101.8
Dicamba	108.1
Dichlorprop	107.7
Dinoseb	118.4

<sup>a</sup>Averages from extraction of sand, loam, and clay soils.

#### **Extraction Conditions**

Solvent:	Dichloromethane/acetone (1:2, v/v), with 4% (v/v) $H_3PO_4/H_2O$ (1:1)
Temperature:	100 °C
Pressure:	1500 psi**
Static Time:	5 min
Static Cycles:	1
Flush Volume:	60% of extraction cell volume
Purge Time:	60 s

\*\*Pressure studies show that 1500 psi is the optimum extraction pressure for all accelerated solvent extraction applications.

#### Analysis

GC-ECD

#### Results

See tables below.

Average recovery and precision for extraction of chlorinated herbicides from
three soil types by ASE method.

Matrix <sup>a</sup>	ASE (% of Spike)	ASE <sup>b</sup> (RSD,%)	Shaking (% of Spike)	Shaking <sup>b</sup> (RSD,%)	ASE as % of Shaking
Clay (low)	36.1	54.5	42.2	25.2	89.8
Clay (high)	71.1	15.0	61.6	21.6	112.8
Loam (low)	56.7	11.2	36.9	78.7	126.6
Loam (high)	59.9	14.1	43.9	14.7	132.8
Sand (low)	51.1	12.6	49.7	13.3	111.2
Sand (high)	69.2	39.5	66.3	35.5	104.4

 $^{a}\text{Low}$  spiking levels ranged from 50 to 500  $\mu\text{g/kg}.$  High spiking levels ranged from 500 to 5000  $\mu\text{g/kg}.$ 

<sup>b</sup> Each precision (RSD,%) value is the average of seven replicate measurements for each compound, then averaged for all compounds.

### Extraction of Chlorinated Pesticides Using Accelerated Solvent Extraction



#### Introduction

Accelerated solvent extraction provides a more convenient, faster, and less solvent intensive method than previously available for the extraction of chlorinated pesticides from environmentally important samples. Recoveries of these analytes by accelerated solvent extraction are equivalent to or better than other more solvent intense methods such as Soxhlet. Accelerated solvent extraction also avoids the problem of multiple washing procedures associated with sonication.

The procedure described in this Application Note meets the requirement for sample extraction as prescribed by EPA Method 3545. This method is applicable to the extraction of water-insoluble or slightly water-soluble volatiles and semivolatiles in preparation for gas chromatographic or GC/MS measurement. The method is applicable to the extraction of chlorinated pesticides from soils, clays, wastes, and sediments containing from 5 to 250 µg/kg of the target compounds.

U.S. EPA Method 3545A: Organochlorinated Pesticides (OCPs) from soil, clay, sludge, and sediments

#### Equipment

Dionex ASE 200 Accelerated Solvent Extractor\* equipped with 11 mL or larger stainless steel extraction cells

Gas Chromatograph (GC) with Mass Spectrometer (MS) or GC Vials for collection of extracts (40 mL, P/N 049465; 60 mL, P/N 049466)

\*Dionex ASE 150 and 350 can be used for equivalent results.

#### Solvents and Reagents

Diatomaceous Earth (DE) Dispersant for ASE, 1 kg Bottlet (P/N 062819)

#### Hexane

#### Acetone

All solvents are pesticide-grade or equivalent and are available from Fisher Scientific.

#### **Extraction Conditions**

Solvent:	Acetone/hexane (1:1 v/v)
Temperature:	100 °C
Pressure:	1500 psi**
Static Time:	5 min
Static Cycles:	1
Flush Volume:	60%
Purge Time:	60 s

\*\*Pressure studies show that 1500 psi is the optimum extraction pressure for all accelerated solvent extraction applications.

#### Analysis

#### GC-MS

#### Results

Average recovery of pesticides from three soil types\*—ASE compared to automated Soxhlet.

Pesticide	Average Recovery (% of Soxhlet)
Alpha BHC	93.3
Gamma BHC-Lindane	95.6
Beta BHC	98.6
Heptachlor	88.0
Delta BHC	99.5
Aldrin	94.9
Heptachlor Epoxide	100.7
Gamma Chlordane	99.5
Alpha Chlordane	102.0
Endosulfan 1	100.3
p,p'-DDE	98.6
Dieldrin	101.2
Endrin	97.2
p,p-DDD	104.6
Endosulfan II	105.6
p,p'-DDT	74.9
Endrin Aldehyde	104.0
Endosulfan Sulfate	105.2
Methocychlor	79.6
Endrin Ketone	102.9

<sup>a</sup> Averages from extraction of sand, loam, and clay soils.

## Extraction of Organophosphorus Pesticides (OPPs) Using Accelerated Solvent Extraction



#### Introduction

The use of accelerated solvent extraction in the extraction of OPP compounds from solid wastes is a more convenient, faster, and less solvent-intensive method than previously available. OPP recoveries by accelerated solvent extraction are equivalent to recoveries from Soxhlet and other methods. Accelerated solvent extraction also avoids the problem of multiple washing procedures associated with sonication.

The procedures described in this Application Note meet the requirements for sample extraction as prescribed by U.S. EPA Method 3545A. This method is applicable to the extraction of water-insoluble or slightly water-soluble OPPs in preparation for gas chromatographic measurements. This method is applicable to soils, solid wastes, and sediments containing 250–2500 µg/kg of OPPs.

#### U.S. EPA Method 3545A: OPPs from soil, sludge, and sediments

#### Equipment

Dionex ASE 200 Accelerated Solvent Extractor\* equipped with 11 mL or larger stainless steel extraction cells

Gas Chromatograph (GC) with Nitrogen Phosphorous Detector (NPD)

Vials for collection of extracts (40 mL, P/N 049465, 60 mL, P/N 049466)

\*Dionex ASE 150 and 350 can be used for equivalent results.

#### Solvents and Reagents

Diatomaceous Earth (DE) Dispersant for ASE, 1 kg Bottle (P/N 062819)

Dichloromethane

Acetone

All solvents are pesticide-grade or equivalent and are available from Fisher Scientific.

#### **Extraction Conditions**

Solvent:	Dichloromethane/acetone (1:1 v/v)
Temperature:	100 °C
Pressure:	1500 psi**
Static Time:	5 min
Static Cycles:	1
Flush Volume:	60% of extraction cell volume
Purge Time:	60 s

\*\*Pressure studies show that 1500 psi is the optimum extraction pressure for all accelerated solvent extraction applications.

#### Analysis

GC-NPD

#### Results

Average recovery and precision for OPP extractions from three soil types.

Matrix <sup>a</sup>	ASE (% of Spike)	ASE⁵ (RSD,%)	Shaking (% of Spike)	Shaking <sup>b</sup> (RSD,%)	ASE as % of Shaking
Clay (low)	55.0	6.2	56.4	7.6	98.8
Clay (high)	69.2	5.2	72.3	16.3	96.3
Loam (low)	61.3	11.6	60.4	8.6	103.6
Loam (high)	61.4	7.8	64.2	6.3	96.5
Sand (low)	59.0	13.0	63.3	6.7	95.0
Sand (high)	64.1	11.9	63.2	4.8	101.2

 $^{\rm a}\text{Low}$  spiking level is approximately 250 µg/kg. High spiking level is approximately 2500 µg/kg.

<sup>b</sup> Each recovery and precision (%RSD) value is the average of seven replicate measurements for each compound, then averaged for all compounds.

# Rapid Determination of Persistent Organic Pollutants (POPs) Using Accelerated Solvent Extraction



#### Introduction

The United Nations Environmental Program (UNEP) has been implemented in an effort to combat the release of selected POPs. POPs are found in environmental samples such as soils, sludges, solid and semisolid waste, and sediments. POPs are also found in biological samples such as human breast milk, and fish tissue. UNEP is interested in eliminating POPs from the environment because these compounds are considered toxic, carcinogenic, and mutagenic, and degrade slowly in the environment, posing a threat to the global environment.

Accelerated solvent extraction is equivalent to U.S. EPA Methods 3540, 3541, 3550, and 8150 for the extraction of organochlorine pesticides (OCPs), organophosphorous pesticides (OPPs), semivolatiles or base neutral acids (BNAs), chlorinated herbicides, polycyclic aromatic hydrocarbons (PAHs), and polychlorinated biphenyls (PCBs). Accelerated solvent extraction complies with U.S. EPA Method 3545A for these compounds.

## U.S. EPA Methods 3540, 3541, 3550, and 8150: POPs from various environmental samples

#### Equipment

Dionex ASE 200 Accelerated Solvent Extractor\* with Solvent Controller (P/N 048765) Use either:

> 22 mL stainless steel extraction cells (P/N 048764) 11 mL stainless steel extraction cells ( P/N 048765) 33 mL stainless steel extraction cells (P/N 048766)

Cellulose filters (P/N 0469458) Vials for collection of extracts (40 mL, P/N 049465; 60 mL, P/N 049466)

Analytical balance (to read to nearest 0.0001 g or better)

\*Dionex ASE 150 and 350 can be used for equivalent results.

#### Solvents and Reagents

Diatomaceous Earth (DE) Dispersant for ASE, 1 kg Bottle) (P/N 062819) Hexane

Dichloromethane

Acetone

Toluene

All solvents are pesticide-grade or equivalent and are available from Fisher Scientific.

#### **Extraction Conditions**

#### Pesticides and PCBs (8081/8082)

Solvent:Hexane/acetone (1:1), (v/v)Temperature:100 °CPressure:1500 psi\*\*Static Time:5 minStatic Cycles:1-2Flush Volume:60%Purge Time:60-120s

#### Hexachlorobenzene (8270)

Solvent:Dichloromethane/acetone (1:1), (v/v)Temperature:100 °CPressure:1500 psi\*\*Static Time:5 minStatic Cycles:1-2Flush Volume:60%Purge Time:60-120 s

#### Dioxins (PCDD and PCDF) (8290)

Solvent: Toluene (100%) or toluene/acetic acid (5%, v/v) if HCl pretreatment currently used

Temperature:	75–200 °C
Pressure:	1500 psi**
Static Time:	5–15 min
Static Cycles:	2–3
Flush Volume:	60-70%
Purge Time:	60–120 s

\*\*Pressure studies show that 1500 psi is the optimum extraction pressure for all accelerated solvent extraction applications.

#### Analysis

GC-MS

GC-MS/MS

#### Results

Average recovery of pesticides from three soil types<sup>a</sup>—ASE compared to automated Soxhlet.

Pesticide	Average Recovery (% of Soxhlet)
Heptachlor	88.0
Aldrin	94.9
Gamma Chlordane	99.5
Alpha Chlordane	102.0
Dieldrin	101.2
Endrin	97.2
p,p'-DDT	74.9

<sup>a</sup> Averages from extraction of sand, loan, and clay soils.

## Analyzing PBDEs in House Dust Samples with the Thermo Scientific TSQ Quantum XLS Ultra GC-MS/MS in EI-SRM Mode



#### Introduction

Polybrominated diphenyl ethers (PBDEs) were introduced in the 1960s as flame retardants and are used today in a wide variety of household apparatuses, consumer electronics, furniture and more. Environmental levels of PBDEs have been continuously increasing due to their special persistence. Certain congeners have been banned completely and are currently in the list of the Stockholm convention's POPs. Sources of human intake are typically via ingestion (nutritional) and the inhalation of indoor and house dust.

Due to growing concerns over the health risks from constant exposure to this class of compounds and the accumulation effects in the food chain, suitable analytical methods are required to quantify flame retardant compounds at low levels in a variety of matrices.

#### U.S. EPA Method 1614: PBDEs from House Dust

#### Equipment

Dionex ASE 350 Accelerated Solvent Extractor equipped with 22 mL or larger stainless steel extraction cells

Gas Chromatograph (GC) with Mass Spectrometer (MS)

#### Solvents

Methylene chloride

Silica gel

All solvents are pesticide-grade or equivalent and are available from Fisher Scientific.

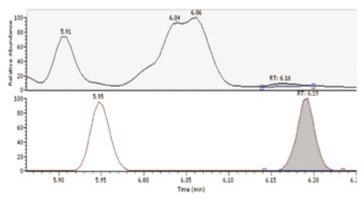
#### **Extraction Conditions**

Solvent:	Methylene chloride (100%)
Temperature:	100 °C
Pressure:	1500 psi*
Static Time:	5 min
Static Cycles:	3
Flush Volume:	60%
Purge Time:	120 s

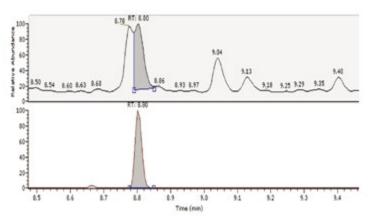
\*Pressure studies show that 1500 psi is the optimum extraction pressure for all accelerated solvent extraction applications.

#### Analysis GC-MS

#### Results



House dust sample in NCI (top) and EI-SRM (bottom) showing BDE47 at a level of 0.07 ppb.



House dust sample in NCI (top) and EI-SRM (bottom) showing BDE 183 at a level of 0.17 ppb.

# Extraction of Polybrominated Biphenyls (PBDEs) from Sediments Using Accelerated Solvent Extraction



#### Introduction

Polybrominated diphenyl ethers (PBDEs) were developed in the early 1970s and are used today as flame retardants for various consumer products, including clothing, furniture, and plastics. Many of these consumer products are disposed in municipal landfills where the PBDEs then leach into groundwater and accumulate in certain biosystems. Recent studies indicate that PBDE concentrations in these biosystems are on the rise. In the 1980s, PBDEs were discovered in European waterways, which led the European Union to ban their production and use. Although the toxicity of PBDEs is still under investigation, evidence suggests that PBDEs may compromise endocrine or hepatic functions.

In this Application Note, PBDEs were extracted from human breast milk (freeze-dried), sediments, fish tissues, and polymers.

#### U.S. EPA Method 1614: PBDEs from sediments

#### Equipment

Dionex ASE 200 Accelerated Solvent Extractor\* with Solvent Controller (P/N 048765)\* 22 mL stainless steel extraction

cells (P/N 048764)

Gas Chromatograph (GC) with Mass Spectrometer (MS) Cellulose filters (P/N 0469458) Vials for collection of extracts, 60 mL (P/N 049466)

\*Dionex ASE 150 and 350 can be used for equivalent results.

#### Solvents and Reagents

Diatomaceous Earth (DE) Dispersant for ASE, 1 kg Bottle (P/N 062819)

Methylene chloride

All solvents are pesticide-grade or equivalent and are available from Fisher Scientific.

#### **Extraction Conditions**

Sediments	
Solvent:	Methylene chloride (100%)
Temperature:	100 °C
Pressure:	1500 psi**
Static Time:	5 min
Static Cycles:	2
Flush Volume:	60%
Purge Time:	120 s

\*\*Pressure studies show that 1500 psi is the optimum extraction pressure for all accelerated solvent extraction applications.

#### Analysis GC-MS

#### Results

With a detection limit of  $0.5 \mu g/kg$ , the GC/MS analysis found BDE-47 in 22% of the 332 sediment sample extracts. BDE-47 is one of two major constituents of Penta, which is a commercial product used to flame retard polyurethane foam.

## Accelerated Solvent Extraction of Polychlorinated Biphenyls from Polyurethane Foam Absorbent Cartridges

Appl



#### Introduction

U.S. EPA Method TO-41 has been used for a number of years for the determination of polychlorinated biphenyls (PCBs) in air. The traditional extraction method for PCBs has been Soxhlet, as typified by U.S. EPA Method 3540. This method requires 12-14 hours of extraction time and 350 mL of solvent per sample to perform. The size and volume requirements of the PUF cartridges have made alternative extraction methods difficult to employ. Accelerated solvent extraction has been found to be a successful alternative method with several advantages over traditional methods.

#### U.S. EPA Method TO-41: PCBs from air sampling PUF cartridges

#### Equipment

Dionex ASE 200 Accelerated Solvent Extractor\* equipped with either:

22 mL stainless steel extraction cells (P/N 048764)

11 mL stainless steel extraction cells (P/N 048765)

Vials for collection of extracts (40 mL, P/N 049465; 60 mL, P/N 049466)

Gas Chromatograph (GC) Kimwipes<sup>®</sup> (Kimberly-Clark

Corporation)

2.5 inch diam. × 3.0 inch PUF cartridge

\*Dionex ASE 150 and 350 can be used for equivalent results.

#### Solvents

Hexane

All solvents are pesticide-grade or equivalent and are available from Fisher Scientific.

#### **Extraction Conditions**

Solvent:	Hexane
Temperature:	100 °C
Pressure:	1500 psi**
Static Time:	5 min
Static Cycles:	1
Flush Volume:	60%
Purge Time:	100 s

\*Pressure studies show that 1500 psi is the optimum extraction pressure for all accelerated solvent extraction applications.

Method detection limit study for ASE and Soxhlet.

#### Analysis

GC

#### Results

See tables below.

Initial demonstration of laboratory performance for ASE and Soxhlet extraction of PUF cartridges\*.

Sample #	ASE (µg)	Soxhlet (µg)	US EPA Method 608 Acceptance Criteria
LCS 1	4.63	4.87	
LCS 2	4.44	4.93	
LCS 3	4.72	5.16	
LCS 4	4.10	4.26	
Average (µg)	4.47	4.80	3.42 - 8.40 µg
Average Recovery	90%	96%	38 - 158%
Standard Deviation	0.274	0.384	< 1.23

\*Recoveries are based on a spiked value of 5.00 µg Aroclor 1248.

Sample #	ASE (µg)	Soxhlet (µg)
MDL 1	0.92	0.79
MDL 2	0.97	0.90
MDL 3	0.91	0.95
MDL 4	0.95	0.78
MDL 5	0.80	0.79
MDL 6	1.11	0.88
MDL 7	0.95	0.90
MDL 8	1.21	0.82
Average	0.98	0.85
Standard Deviation	0.13	0.06

0.38

0.19

\*Recoveries are based on a spiked value of 1.00 µg Aroclor 1248.

**MDL (**3δ)

## Extraction of PCBs from Environmental Samples Using Accelerated Solvent Extraction



#### Introduction

Polychlorinated biphenyls (PCBs) are found in many solid waste materials worldwide. This Application Note describes the application of accelerated solvent extraction to the extraction of PCBs from sewage sludge, river sediments, marine sediments, and marine tissue (oyster). The procedures described in this Application Note meet the requirements for sample extraction as determined by U.S. EPA Method 3545 for solid samples.

Accelerated solvent extraction is a direct replacement for solvent-intensive techniques such as Soxhlet and sonication. For the preparation of solid waste samples, containing PCBs, accelerated solvent extraction provides more convenient, faster extractions with significantly less solvent usage than these other methods.

#### U.S. EPA Method 3545A: PCBs from soil, sludge, and sediments

#### Equipment

Dionex ASE 200 Accelerated Solvent Extractor\* equipped with 11 mL or larger stainless steel extraction cells

Gas Chromatograph (GC) with Electron Capture Detector (ECD) Vials for collection of extracts (40 mL, P/N 049465; 60 mL, P/N 049466)

\*Dionex ASE 150 and 350 can be used for equivalent results.

#### **Solvents and Reagents**

Diatomaceous Earth (DE) Dispersant for ASE, 1 kg Bottle (P/N 062819) Acetone

#### Hexane

All solvents are pesticide-grade or equivalent and are available from Fisher Scientific.

#### **Extraction Conditions**

Solvent:	Hexane/acetone (1:1), (v/v)
Temperature:	100 °C
Pressure:	1500** psi
Static Time:	5 min
Static Cycles:	1
Flush Volume:	60% of extraction cell volume
Purge Time:	60 s

\*\*Pressure studies show that 1500 psi is the optimum extraction pressure for all accelerated solvent extraction applications.

#### Analysis

GC-ECD

#### Results

PCB recoveries from sewage sludge\*.

PCB Congener	Average Recovery, n = 6 (as % of Soxhlet)	RSD (%)
PCB 28	118.1	2.5
PCB 52	114.0	4.7
PCB 101	142.9	7.4
PCB 153	109.5	5.8
PCB 138	109.6	3.9
PCB 180	160.4	7.5

\*Analyte concentration range: 160-200 µg/kg per component

#### PCB recovery from river sediment (SRM 1939)\*.

PCB Congener	Average, n = 6 (as % of Soxhlet)	RSD (%)
PCB 101	89.2	3.7
PCB 153	62.3	4.1
PCB 138	122.1	2.3
PCB 180	111.5	5.9

\*Analyte concentration range: 170-800 µg/kg per component

#### Recovery of Aroclor 1254 from soil (CRM911-050).

Run Number Aroclor Found (µg/kg)		
1	1290.0	
2	1365.8	
3 1283.4		
4	1368.6	
Average	1327.0 (99.0%)	
RSD	3.51%	

### Literature References

J. L. Ezzell, B. E. Richter, W. D. Felix, S. R. Black, J. E. Meikle. A comparison of ASE with conventional solvent extraction for organophosphorus pesticides and herbicides. *LC/GC*. **1995**: 13, 390–398.

J. L. Ezzell, B. E. Richter. Automated sample preparation for environmental laboratories using ASE. *American Environmental Laboratory*. **1996**, 16–18.

B. E. Richter, B. A. Jones, J. L. Ezzell, N. L. Porter, N. Avdalovic, C. Pohl. ASE: A technique for sample preparation. *Analytical Chemistry*. **1996**: 68, 1033–1039.

B. E. Richter, J. L. Ezzell, D. E. Knowles, F. Höfler, A. K. R. Mattulat, M. Scheutwinkel, D. S. Waddell, T. Gobran, V. Khurana. Extraction of PCDDs and PCDFs from environmental samples using ASE. *Chemosphere*. **1997**: 34, 5–7, 975–987.

P. Popp, P. Keil, M. Möder, A. Paschke, U. Thuss. Application of ASE followed by GC, HPLC and GC-MS for the determination of polycyclic aromatic hydrocarbons, chlorinated pesticides and PCDDs and PCDFs in solid wastes. Journal of Chromatography A. **1997**: 774, 203-211.

H. Bautz, J. Polzer, L. Stieglitz. Comparison of pressurized liquid extraction with Soxhlet extraction for the analysis of PCDDs and PCDFs from fly ash and environmental matrices. *Journal of Chromatography A*. **1998**: 815, 231-241.

S. P. Frost, J. R. Dean, K. P. Evans, K. Harradine, C. Cary, M. H. I. Comber. Extraction of hexaconazole from weathered soils: a comparison between Soxhlet extraction, microwave-assisted extraction SFE and ASE. *Analyst.* **1997**: 122, 895–898.

E. Conte, R. Milani, G. Morali, F. Abballe. Comparison between ASE and traditional extraction methods for the analysis of the herbicide difluenican in soil. *Journal of Chromatography A*. **1997**: 765, 121–125.

S. R. Sumpter, B. A. Peterson. Determination of pyrithiobac sodium in soil using subcritical water extraction followed by HPLC/UV analysis in support of a Staple® herbicide small-scale prospective groundwater study. 1997 Pittsburgh Conference. **1997**: 16–21 March, Paper 1264, .

L. Guzzella, F. Pozzoni. ASE of herbicides in agricultural soil samples. *International Journal of Environmental Analytical Chemistry.* **1998**: 71, 3–4, 1–9.

J. R. Donnelly, A. H. Grange, N. R. Herron, G. R. Gregory, R. Nichol, J. L. Jeter, R. J. White, W. C. Brumley, J. Van Emon. Modular methodology for determination of PCBs in soil as aroclors and individual congeners. *Journal of AOAC International.* **1996**: 79, 953–961.

O. Zuloaga, N. Etxebarria, L. A. Fernández, J. M. Madariaga. Comparison of ASE with microwave-assisted extraction and Soxhlet for the extraction of PCBs in soil samples. *Trends in Analytical Chemistry.* **1998:** 17, 10, 642–647.

W. C. Brumley, E. Latorre, V. Kelliher, A. Marcus, D. E. Knowles. Determination of chlordane in soil by LC/GC/ECD and LC/GC/EC NIMS with comparison of ASE, SFE, and Soxhlet extraction. *Journal of Liquid Chromatography and Related Technology*. **1998**: 21, 1199–216. B. A. Tomkins, G. A. Sega, S. J. Macnaughton. The quantitation of sulfur mustard by-products, sulfur containing herbicides, and organo-phosphonates in soil and concrete. *Analytical Letters*. **1998**: 31, 1603–1622.

S. M. Pyle, A. B. Marcus. Rapid and sensitive determination of pesticides in environmental samples by accelerated solvent extraction and tandem mass spectrometry. *Journal of Mass Spectrometry*. **1997:** 32, 897–898.

S. Heise, N. Litz. Extraction of surfactants out of solid matrices. *Tenside, Surfactants, Detergents.* **1999:** 36, 185–191.

T. S. Reighard, S. V. Olesik. Bridging the gap between supercritical fluid extraction and liquid extraction techniques: alternative approaches to the extraction of solid and liquid environmental matrices. Critical Reviews in Analytical Chemistry. 1996: 26, 2 and 3, 61–99.

C. Bandh, E. Bjorklund, L. Mathiasson, C, Naf, Y. Zebuhr. Determination of PCBs in baltic sea sediments using accelerated solvent extraction. *Organonbalogen Compounds*. **1998**: 35, 17–19.

G. S. Chen, K. W. Schramm, B. Henkelmann, Y. Xu, Y. Y. Zhang, T. Wottgen, A. Kettrup. Comparative evaluation of "Soxhlet-" and "Accelerated Solvent-" Extraction in the Determination of PCDD/F in Sediment. *Organohalogen Compounds*. **1997**: 31, 114–118.

H. Wagenaar, N. Pronk, H. Olthof. Evaluation of accelerated solvent extraction of PCDDs and PCDFs from native contaminated samples. *Organohalogen Compounds*. **1996**: 27, 265–268.

J. C. Johnson, J. M. Van Emon. Re-evaluation study of polychlorinated biphenyl contaminated gencorp lawrence soil using quantitative immunoassay. *Book of Abstracts, 214th ACS National Meeting, Las Vegas, NV.* **1997, Sept.** 7–11:029, .

W. Knoth. Comparison of pressurized fluid and Soxhlet extraction of PCDD/F from soil. 4th Symposium on Extraction for Sample Preparation SFE-()SE-SPME. 1999: 22–23 Sept., 71–74.

J. Gan, S. K. Papiernik, W. C. Koskinen, S. R. Yates. Evaluation of accelerated solvent extraction (ASE) for analysis of pesticide residues in soil. *Environmental Science and Technology*. **1999**: 33, 3249–3253.

D. E. Kimkbrough, R. Chin, J. Wakakuwa. Wide-spread and systematic errors in the analysis of soils for polychlorinated biphenyls: Part 2. Comparison of extraction systems. *Analyst.* **1994**: 119, 1283–1292.

E. Bjorklund. Determination of persistent organic pollutants in solid environmental sample using ASE and SFE. *US DOE Facsimile Report*. **1999:** 25-Feb, 1–52.

M. Krappe, F. Hoefler. A fully automated system for extraction, clean-up and analysis of solid matrices using ASE 200 combined with an automated SPE sampler. 4th Symposium on Extraction for Sample Preparation SFE-()SE-SPME. **1999**: 22–23 Sept., 80–81.

A. Hubert, K. D. Wenzelm, M. Manz, L. Weissflog, W. Engewald, G. Schuurmann. Improvement of extraction efficiency for POPs in real contaminated soil samples using ASE. 4th Symposium on Extraction for Sample Preparation SFE-()SE-SPME. **1999**: 22–23 Sept, Poster. J. Gan, S. K. Papiernik, W. C. Koskinen, S. R. Yates. Evaluation of accelerated solvent extraction (ASE) for analysis of pesticide residues in soil. *Environmental Science and Technology*. **1999**: 33, 18, 3249–3253.

D. D. McCant, L. S. Inouye, V. A. McFarland. A one-step ASE extraction method for TCDD TEQ determination. *Bulletin of Environmental Contamination and Toxicology*. **1999:** 63, 282–288.

L. J. Fitzpatrick, J. R. Dean, M. H. I. Comber, K. Harradine, K. P. Evans. Extraction of DDT and its metabolites DDE and DDD from aged contaminated soil. *Journal of Chromatography A*. 2000: 874, 257–264.

B. Henkelmann, T. Wottgen, G. Chen, K. W. Schramm, A. Kettrup. Accelerated solvent extraction (ASE) of different matrices in the analysis of polychlorinated dibenzo-*p*-dioxins and dibenzofurans: Method development and comparison to soxhlet extraction. *Organohalogen Compounds.* **1999**: 40, 133–136.

A. Hubert, K. D. Wenzel, M. Manz, L. Weissflog, W. Engewald, G. Schuurmann. High extraction efficiency for POPs in real contaminated soil samples using accelerated solvent extraction. *Analytical Chemistry.* 2000: 72, 1294–1300.

M. D. David, S. Campell, Q. X. Li. Pressurized fluid extraction of nonpolar pesticides and polar herbicides using in situ derivation. *Analytical Chemistry*. **2000**: 72,15, 3665–3670.

D. L. Poster, M. M. Schantz, S. A. Wise, M. G. Vangel. Analysis of urban particulate standard reference materials for the determination of chlorinated organic contaminants and additional chemical and physical properties. *Fresenius Journal of Analytical Chemistry.* **1999:** 363, 380–390.

I. Windal, D. J. Miller, E. De Pauw, S. B. Hawthorne. Supercritical fluid extraction and accelerated solvent extraction of dioxins from high-and low-carbon fly ash. *Analytical Chemistry*. **2000**: 72, 3916–3921.

C. M. Reddy, L. J. Heraty, B. D. Holt, N. C. Sturchio, T. I. Eglinton, N. J. Drenzek, L. Xu, J. L. Lake, K. A. Maruya. Stable chlorine isotopic compositions of aroclors and aroclor-contaminated sediments. Environmental Science and Technology. 2000: 34, 13, 2866–2870.

B. E. Richter, L. Covino. New environmental applications of accelerated solvent extraction. *LCGC*. **2000**: 18, 10, 1068–1073.

L.J. Fitzpatrick, O. Zuloaga, N. Etxebarria, J. R. Dean. Environmental applications of pressurized fluid extraction. Reviews in Analytical Chemistry. 2000: 19, 75-122.

J. S. Yang, S. K. Lee, Y. H. Park, D. W. Lee. Analytical method for dioxin and organo-chlorinated compounds: (II) Comparison of extraction methods of dioxins from XAD-2 adsorbent. *Bull. Korean Chem. Soc.* **1999:** 20, 689–695.

Y. Zhu, K. Yanagihara, F. Guo, Q. X. Li. Pressurized fluid extraction for quantitative recovery of chloroacetanilide and nitrogen heterocyclic herbicides in soil. *Journal of Agricultural and Food Chemistry*. 2000: 48, 4097–4102.

T. Bergloef, W. C. Koskinen, H. Kylin, T. B. Moorman. Characterization of triadimefon sorption in soils using supercritical fluid extraction (SFE) and accelerated solvent extraction (ASE) techniques. *Pest Management Science.* **2000**: 56, 10, 927–931.

H. Miyamoto, K. Ohtsuka, Y. Fukuda, Y. Ishibashi. Rapid extraction of dioxins from soil, fly ash and XAD-2 resin using ASE and hot extraction. *Organobalogen Compounds*. **1999:4** 0, 215–218.

A. Rubel, R. Bierl. Routine analysis of vinicultural relevant fungicides, insecticides and herbicides in soil samples using enhanced solvent extraction (ESE). *Fresenius J Anal Chem.* **1999:** 364, 648–650.

S. G. Reddy, K. E. Keeler, M. W. Coyle. A sample preparation method for PCB's in PUFs by an accelerated solvent extraction system: A pollution prevention and waste minimization initiative. *American Environmental Labs.* **1999.** 

J. L. Ezzell. Pressurized fluid extraction (PFE) in organic analysis. Extraction Methods in Organic Analysis: Book edited by Alan J handley, Sheffield CRS Press Boca Raton Fl 33431 **1999.** 147–165.

O. Zuloaga, N. Etxebarria, L. A. Fernandez, J. M. Madariaga. Optimization and comparison of MAE, ASE and Soxhlet extraction for the determination of HCH isomers in soil samples. *Fresenius J analytical Chemistry.* **2000**: 367, 733–737.

C. Bandh, E. Bjorklund, L. Mathiasson, C. Naf, Y. Zebuhr. Comparison of accelerated solvent extraction and soxhlet extraction for the determination of PCB's in baltic sea sediments. *Environmental Science and Technology.* **2000**: 34, 23, 4995–5000.

Y. Abrha, D. Raghavan. Polychlorinated biphenyl recovery from spiked organic matrix using accelerated solvent extraction and soxhlet extraction. Journal of Hazardous Materials. 2000: B80, 147-157.

M. Weichbrodt, W. Vetter, B. Luckas. Microwave-assisted extraction and ASE with ethyl acetate-cyclohexane before determination of organochlorine in fish tissue by gas chromatography with electron capture detector. *Environmental Analysis*. 2000: 83, 1334–1343.

W-H. Ding, J. C. H. Fann. Determination of linear alkylbenzenesulfonates in sediments using pressurized fluid extraction and ion-pair derivitization gas chromatography-mass spectrometry. *Analitica Chimica Acta.* **2000**: 408, 408, 291–297.

E. Bjorkland, T. Nilsson, S. Bowardt. Pressurized liquid extraction of persistent organic pollutants in environmental analysis. *Trends in Analytical Chemistry.* 2000: 19, 434–445.

A. A. Mackay, P. M. Gschwend. Enhanced concentrations of PAHs in groundwater at a coal tar site. *Environmental Science Technology.* 2001: 35, 1320–1328.

S. Marchese, D. Perret, A. Gentili, R. Curini, A. Marino. Development of a method based on accelerated solvent extraction and liquid chromatography/mass spectrometry for the determination of arylphenoxypropionic herbicides in soil. *Rapid Communications in Mass Spectrometry.* 2001: 15, 393–400.

S. Raccanelli, V. Bonamin, M. Favotto, V. Di Marco, W. Tirler. Comparative PCDD/Fs analysis in different matrices (fish harbor sediment, Industrial sludge, MSWI ash) extracted by pressurized fluid extraction and by traditional soxhlet. Preliminary results. *Oganohalogen Compounds.* **1999:** 40, 239–242.

E. Braekevelt, G. T. Tomy, G. A. Stern. Comparison of individual congener standard and a technical mixture for the quantitation of toxaphene in environmental matrixes by HRGC/ENCI-HRMS. *Environmental Science and Technology.* **2001**: 35 ,17, 3515–3518.

R. E. Clement, P. W. Young. Environmental Analysis: 2001 application review. *Analytical Chemistry*. **2001**: 73, 2761–2790.

G. LeBlanc. A review of EPA sample preparation techniques for organic compound analysis of liquid and solid samples. *LCGC*. **2001**: November 19, 11, 1120–1130.

S. Campbell, Q.X. Li. NA+ EDTA assisted insitu derivitization pressurized fluid extraction of polar herbicides in soil. *Analytica Chimica Acta*. **2001**: 434, 283–289.

S. Valsechchi, S. Polesello, S. Cavalli. Recovery of 4-nonylphenol and 4-nonylphenol ethoxylates from river sediment by pressurized fluid extraction. *Journal of Chromatography A*. **2001**: 925, 297–301.

R. C. Hale, M. J. La Guardia, E. P. Harvey, T. Matteson Mainor, W. H. Duff, M. O. Gaylor. Polybrominated diphenyl ether flame retardant in Virginia freshwater fishes. *Environmental Science and Technology*. 2001: 35, 23, 4585–4591.

V. Camel. Recent extraction techniques for solid matrixes-supercritical fluid extraction, pressurized fluid extraction and microwave-assisted extraction: their potential and pitfalls. *The Analyst.* 2001: 126, 1182–1193.

A. Hubert, K-D Wenzel, W. Engelwald, G. Schuurman. ASE- More efficient extraction of POPs and PAHs from real contaminated plant and soils. *Reviews in Analytical Chemistry.* **2001**, 101–143.

C. Schroder, P. Castle. Life history parameters and PCB in long finned pilot whales from New Zealand. A report to WWF. **1998**.

M. Papagiannopolus, B. Zimmerman, A. Mellinthin, M. Krappe, G. Maio, R. Galensa, ASE-ASPEC-HPLC kopplung - automatic bestimmung von proanthocycidien in malz. *GIT Labor Fachzeitschrift*. 2001.

E. Bjorklund, S. Bowadt, T. Nilsson, L. Mathiasson. Pressurized fluid extraction of polychlorinated biphenyls in solid environmental samples. *Journal of Chromatography A.* **1999**: 836, 285–293.

I. Ferrer, E. T. Furlong. ASE followed by on line solid phase extraction coupled to Ion trap LC/MS/MS for analysis of benzalkonium chlorides in sediment samples. *Analytical Chemistry.* **2002**: 74, 1275–1280.

D. Martens, M. Gfrerer, T. Wenzel, A Zhang, B.M. Gawlik, K.-W. Schramm, E. Lankmayr, A. Kettrup. Comparison of different extraction techniques for the determination of polychlorinated organic compounds in sediment. *Analytical Bioanal Chem.* **2002**: 372, 562–568.

E.M. Golet, A. Strehler, A.C. Alder, W. Giger. Determination of flouroquinolone antibacterial agents in sewage sludge and sludge-treated soil using accelerated solvent extraction followed by solid-phase extraction. *Analytical Chemistry.* **2002**: 74, 5455–5462.

I. C. Ryu, J.h. Uom, Y. G. Lee, S. W. Eom, J. y. Shin. Comparison of the extraction efficiency of polychlorinated dibenzo-p-dioxins and polychlorinated dibenzofurans from soils using ASE and Soxhlet. *Organohalogen Compounds.* 2000: 45, 78–80.

H. Eun, N. Seike, K. Baba, R. Uegaki, Y. Ishii, M. Kuwahara, M. Ueji. Comparative efficiency of PCDD/FS with a variety of extraction methods. *Organobalogen Compounds*. **2002**: 55, 163–165.

Y. Kemmochi, K. Tsutsumi. Rapid PCDD/PCDF screening method for fly ash with ion trap MS/MS. *Chemosphere*. **2001**:43, 433–437.

H. Preud'homme, M. Potin-Gautier. Optimization of accelerated solvent extraction for polyhalogenated dibenzo-*p*-dioxins and benzo-*p*-furans in mineral and environmental matrixes using experimental designs. *Analytical Chemistry.* **2003**: 75, 22, 6109–6118.

K.Saito, M. Takekuma, M. Ogawa, S.Kobayashi, Y. Sugawara, M. Ishizuka, H. Nakazawa, Y. Matsuki. Extraction and cleanup methods of dioxins in house dust from two cities in Japan using accelerated solvent extraction and a disposable multi-layer silica-gel cartridge. *Chemosphere*. 2003: 53, 137–142.

R. Wahlen, C. Wolff-Briche. Comparison of GC-IP-MS and HPLC-ICP-MS for species-specific isotope dilution analysis of tributyltin in sediment after accelerated solvent extraction. Anal. Bioanal. Chem. 2003: 377, 140-148.

Y-h. Lang, X. Jiang, D. Martens, L. Sun, J-w. Yang. Accelerated solvent extraction and GC-MS analysis of chlorophenols in the sediment. *Journal of Analytical Science*. **2003**: 19, 228–230.

C. H. Marvin, S. Painter, G. T. Tomy, G. A. Stern, E. Braekevelt, D. C. G. Muir. Spatial and temporal trends in short-chain chlorinated paraffins in Lake Ontario sediments. *Environmental Science and Technology.* 2003: 37, 20, 4561–4568.

Y. Cheng, S-M. Li. Analytical method development of long-chain ketones in PM 2.5 aerosols using Accelerated Solvent Extraction. Intern. J. Environ. *Anal. Chem.* **2004**: 84, 367–378.

T. Yarita, A. Nakama, M. Numata, Y. Aoyagi, M. Yamazaki, A. Takatsu. Determination of polychlorinated biphenyls in sediment by isotope-dilution gas chromatography/mass spectrometry with pressurized fluid extraction. *Japan Society for Analytical Chemistry.* **2003**: *52*, 1011–1017.

G. MacInnis, G. Tomy, M. Alaee, C. Marvin. Analysis of hexabromocyclododecane (HBCD) in sediments using negative ESI LC/MS/MS. Poster.

E. Kremer, M. Rompa, B. Zygmunt. Extraction of acid herbicides from soil by means of accelerated solvent extraction. *Chromatographia Supplement.* **2004**: 60, S169–S174.

W. Wilcke, W. Amelung. Persistent organic pollutants in native grassland soils along a climosequence in North America. *Soil Sci. Soc. Am. J.* 2000: 64, 2140–2148.

J. S. Yang, S. K. Lee, Y. H. Park, D. W. Lee. Analytical method for dioxin and organo-chlorinated compounds: (II) comparison of extraction methods of dioxins from XAD-2 adsorbent. *Korean Chem. Soc.* **1999:** 20, 689–695.

C. Robinson, P. Blow, F. Dorman. Rapid dioxin analysis using accelerated solvent extraction (ASE), multi-column sample cleanup and RTx-Dioxin2 gas chromatography column. Organohalogen Compounds. 2004: 66, 101-106.

A. Sjodin, O. Papke, E. McGahee III, R. Jones, J-F Focant, T. Pless-Mulloli, L-M Toms, R. Wang, Y. Zhang, L. Needham, T. Herrmann, D. Patterson Jr. Concentration of polybrominated diphenyl ethers (PBDEs) in house hold dust from various countries-inhalation a potential route of human exposure. *Organohalogen Compounds*. 2004: 66, 3770–3775.

J. R. Kucklick, K. J.S. Tuerk, S. S. Vander Pol, M. M. Schantz, B. J. Porter, S. A. Wise. Concentration of polybrominated diphenyl ether congeners and toxaphene in selected marine standard reference materials. *Organohalogen Compounds*. **2003**: 60, 126–129.

W. W. Brubaker Jr, M. M. Schantz, S. A. Wise. Determination of non-ortho polychlorinated biphenyls in environmental standard reference materials. *Fresenius J Anal Chem.* **2000**: 367, 401–406.

S. Tao, F. L. Xu, X. j. Wang, W. X. Liu, Z. M. Gong, J. Y. Fang, L. Z. Zhu, Y. M. Luo. Organochlorine pesticides in agricultural soil and vegetables from Tianjin, China. *Environ. Sci. Technol.* **2005**: 39, 2494–2499.

Holscher, A. Maulshagen, H. Shirkhan, G. Lieck, P. A. Behnisch. Automated rapid analysis for dioxins and PCBs in food, feedingstuff and environmental matrices. *Organohalogen Compounds*. 2004: 66, 116–124.

M. R. Burkhardt, R. C. ReVello, S. G. Smith, S. D. Zaugg. Pressurized liquid extraction using water/isopropanol coupled with solid-phase extraction cleanup for industrial and anthropogenic waste-indicator compounds in sediment. *Analytica Chimica Acta*. 2005: 534, 89–100.

H. Stapleton, M. Schantz, S. Wise. Measurement of polybrominated diphenyl ethers in environmental matrix standard reference materials. *Organohalogen Compounds*. **2004**: 66, 3696–3699.

D. R. Oros, D. Hoover, F. Rodigari, D. Crane, J. Sericano. Levels and distribution of polybrominated diphenyl ethers in water, surface sediments and bivalves from the San Francisco estuary. *Environmental Science and Technology*. 2005: 39, 33–41.

H. M Stapleton, N. G. Dodder, J. H. Offenberg, M. M. Schantz, S. A. Wise. Polybrominated diphenyl ethers in house dust and clothes dryer lint. *Environmental Science and Technology*. **2005**:3 9, 925–931.

J. R. Kucklick, K. J. S. Tuerk, S. S. Vander Pol, M. M. Schantz, S. A. Wise. Polybrominated diphenyl ether congeners and toxaphene in selected marine standard reference. *Anal Bioanal Chem.* **2004:** 378, 1147–1151.

L. Turrio-Baldassarri, S. Alivernini, C. L. Battisteli, S. Carasi, M. Casella, I. Fochi, N. Iacovella, A. L. Iamiceli, A. Indelicato, C. La Rocca, A. Mariani, C. Scarcella. A study on PCB, PCDD/PCDF industrial contamination in an urban/agricultural area. Part I: soil. *Organohalogen Compounds.* **2004**: 66, 1346–1350.

H. Stapleton, N. Dodder, M. Schantz, S. Wise. Measurements of the flame retardants polybrominated diphenyl ethers (PBDEs) and hexabromocyclododecane (HBSDD) in house dust. *Organobalogen Compounds*. **2004:** 66, 3691–3695.

G. Larsen, Z. Fan, F. Casey, H. Hakk. Sorption. Mobility and fate of 1,4,7,8-tetrachlorodibenzo-p-dioxin in soils. *Organohalogen Compounds*. 2004: 66, 2292–2296.

E. Eljarrat, A. de la Cal, D. Raldua, D. Barcelo. Fate and behaviour of brominated flame retardants (PBDEs and HBCD) in Ebro River basin. Department of Environmental Chemistry, Spain. 74–77.

E. M. Krummel, I. Gregory-Eaves, R. W. MacDonald, L. E. Kimpe, M. J. Demers, J. P. Smol, B. Finney, J. M. Blais. Concentrations and fluxes of salmon-derives polychlorinated biphenyls (PCBs) in lake sediments. *Environmental Science Technology*. **2005**: 39, 18, 7020–7026.

S. Sporring, S. Bowadt, B. Svensmark, E. Bjorklund. Comprehensive comparison of classis soxhlet extraction with soxtec extraction, ultrasonication extraction, supercritical fluid extraction, microwave assisted extraction and accelerated solvent extraction for the determination of polychlorinated biphenyls. *Journal of Chromatography A.* 2005: 1090, 1–9.

Y. Ashizuka, R. Nakagawa, K. Tobiishi, T. Hori, T. Iida. Determination of polybrominayed diphenyl ethers and polybrominated dibenzo-pdioxins/dibenzofurans in marine products. *Journal of Agricultural and Food Chemistry.* **2005**: 53, 3807–3813.

K. Kawata, T. Asada, K. Oikawa. Determination of pesticides in compost by pressurized liquid extraction and gas chromatography-mass spectrometry. *Journal of Chromatography A*. 2005: 1090, 10–15.

M. I. H. Helaleh, A. Al-Omair, N. Ahmed, B. Gevao. Quantitive determination of organochlorine pesticides in sewage sludges using soxtex, soxhlet and pressurized liquid extractions and ion trap mass-mass spectrometric detection. *Anal Bioanal Chem.* **2005** :382, 1127–1134.

S. Campbell, L. Chen, J. Yu, Q. X. Li. Adsorption and analysis of the insecticides thiamethoxam and indoxacarb in Hawaiian soils. *Journal of Agricultural and Food Chemistry.* 2005: 53, 5373–5376.

T. Dagnac, S. Bristeau, R. Jeannot, C. Mouvet, N. Baran. Determination of chloroacetanilides, triazines and phenylureas and some of their metabolites in soils by pressurized liquid extraction, GC-MS/MS, LC-MS and LC-MS/MS. *Journal of Chromatography A*. 2005: 1067, 225–233.

X. Pan, B. Zhang, S. B. Cox, T. A. Anderson, G. P. Cobb. Determination of N-nitroso derivatives of hexahydro-1,3,5-trinitro-1,3,5-triazine (RDX) in soils by pressurized liquid extraction and liquid chromatography-electrospray ionization mass spectrometry. *Journal of Chromatography A*. 2006: 1107, 2–8.

A. Hussen, R. Westborn, N. Megersa, L. Mathiasson, E. Bjorklund. Development of a pressurized liquid extraction and clean-up procedure for the determination of a-endosulfan, B-endosulfan and endosulfan sulfate in aged contaminated Ethiopian soil. *Journal of Chromatography A.* **2006**: 1103, 202–210.

E. Eljarrat, A. de la Cal, D. Barcelo. Potential chlorinated and brominated interferences on the polybrominated diphenyl ether determinations by gas chromatography-mass spectrometry. *Journal of Chromatography A.* **2003**: 1008, 181–192.

A. Marklund, B. Andersson, P. Haglund. Organophosphorus flame retardants and plasticizers in Swedish sewage treatment plants. *Environmental Science and Technology*. **2005**: 39, 19, 7423–7429.

S. Josefsson, R. Westbom, L. Mathiasson, E. Bjorklund. Evaluation of PLE exhaustiveness for the extraction of PCBs from sediments and the influence of sediment characteristics. *Analytica Chimica Acta*. 2006: 560, 94–102.

O. Kiguchi, T. Kobayashi, K. Saitoh, N. Ogawa. Pressurized liquid extraction of polychlorinated dibenzo-p-dioxins, polychlorinated dibenzofurans and coplanar polychlorinated biphenyls from contaminated soil. *Journal of Chromatography A.* **2006**: 1108, 176–182.

R. C. Hale, M. J. La Guardia, E. Harvey, T. M. Mainor. Potential role of fire retardant-treated polyurethane foam as a source of brominated diphenyl ethers to the US environment. *Chemosphere*. 2002: 46, 729–735.

E. Eljarrat, A. de la Cal, D. Barcelo. Determination of decabromodiphenyl ether in sediments using selective pressurized liquid extraction followed by GC-NCI-MS. *Anal Bioanal Chem.* **2004**: 378, 610–614.

A. de la Cal, E. Eljarrat, D. Barcelo. Determination of 39 polybrominated diphenyl ether congeners in sediment samples using fast selective pressurized liquid extraction and purification. *Journal of Chromatography A.* **2003**: 1021, 165–173.

E. Eljarrat, A. del la Cal, D. Raldua, C. Duran, D. Barcelo. Brominated flame retardants in Alburnus alburnus from Cinca River Basin (Spain). *Environmental Pollution.* **2005**: 133, 501–508.

G. T. Tomy, W. Budakowski, T. Halldorson, D. M. Whittle, M. J. Keir, C. Marven, G. MacInnis, M. Alaee. Biomagnification of ol- and y-hexabromocyclododcane isomers in a Lake Ontario food web. *Environmental Science and Technology.* **2004**:3 8, 8, 2298–2303.

S. Polati, M. Roz, S. Angioi, V. Gianotti, F. Gosetti, E. Marengo, C. Rinaudo, M.C. Gennaro. Statistical evaluation of recovery of 3,4-dichloroaniline in soil as function of particle size and analyte concentration. *Talanta*. **2005**: 68, 93–98.

R. Rodil, P. Popp. Development of pressurized subcritical water extraction combined with stir bar sorptive extraction for the analysis of organochlorine pesticides and chlorobenzenes in soil. *Journal of Chromatography A.* 2006:1 124, 82–90.

S. Harrad, S. Hazrati, C. Ibarra. Concentrations of polychlorinated biphenyls in indoor air and polybrominated diphenyl ethers in indoor air and dust in Birmingham, United Kingdom: Implications for human exposure. *Environmental Science and Technology.* **2006**: 40, 4633–4638.

M. S. Diaz-Cruz, D. Barcelo. Highly selective sample preparation and gas chromatographic-mass spectrometric analysis of chlorpyrifos, diazinon and their major metabolites in sludge and sludge-fertilized agricultural soils. *Journal of Chromatography A*. 2006: 1132, 21–27.

R. Koppen, R. Becker, C. Jung. Investigation of extraction procedure and HPLC-DAD/MS for the determination of the brominated flame retardant tetrabromobisphenol of the brominated flame retardant tetrabromobisphenol A bis(2,3-dibromopropylether) in environmental samples. *Anal Bioanal Chem.* **2006**: 384, 1485–1492.

R. C. Brandli, T. D. Bucheli, T. Kupper, F. X. Stadelmann, J. Tarradellas. Optimized accelerated solvent extraction of PCBs and PAHs from compost. *International Journal of Environmental Analytical Chemistry.* **2006:** 86, 7, 505–525.

M. Nording, M. Nichkova, E. Spinnel, Y. Persson, S. J. Gee, B. D. Hammock, P. Haglund. Rapid screening of dioxin-contaminated soil by accelerated solvent extraction/purification followed by immunochemical detection. *Anal Bioanal Chem.* **2006**: 385, 357–366.

M. M. Schantz. Pressurized liquid extraction in environmental analysis. *Anal Bioanal Chem.* **2006**: 386, 1043–1047.

H. M. Stapleton, N. G. Dodder, J. R. Kucklick, C. M. Reddy, M. M. Schantz, P. R. Becker, F. Gulland, B.J. Porter, S. A. Wise. Determination of HBCD, PBDEs and MeO-BDEs in California sea lions (Zalophus californianus) stranded between 1993 and 2003. *Marine Pollution Bulletin.* 2006: *52*, *522–531*.

I. Johansson, K. Heas-Moisan, N. Guiot, C. Munschy, J. Tronczynski. Polybrominated diphenyl ethers (PBDEs) in mussels from selected French coastal sites: 1981–2003. *Chemosphere*. **2006**: 64, 296–305.

D. Chen, B. Mai, J. Song, Q. Sun, Y. Luo, X. Luo, E. Y. Zeng, R. C. Hale. Polybrominated diphenyl ethers in birds or prey from Northern China. *Environmental Science and Technology*. 2007: 41, 1828–1833.

Ahmed Hussen, Rikard Westbom, Negussie Megersa, Negussie Retta, Lennart Mathiasson, Erland Bjorklund. Optimisation of pressurised liquid extraction for the determination of p,p'-DDT and p,p'-DDE in aged contaminated Ethiopian soils. *Anal. Bional Chem.* **2006**: 386, 1525–1533.

Ahmed Hussen, Rikard Westbom, Negussie Megersa, Lennart Mathiasson, Erland Bjorklund. Selective pressurized liquid extraction for multi-residue analysis of organochlorine pesticides in soil. *Journal of Chromatography A.* 2007: 1152, 247–253.

Alain Hildebrandt, Silvia Lacorte, Damia Barcelo. Assessment of priority pesticides, degradation products, and pesticide adjuvants in groundwaters and top soils from agricultural areas of the ebro river basin. Anal Bioanal Chem. 2007: 387, 1459-1469.

Masahiko Numata, Takashi Yarita, Yoshie Aoyagi, Yoko Tsuda, Misako Yamazaki, Akiko Takatsu, Keiichiro Ishikawa, Koichi Chica, Kensaku Okamaoto. Sediment certified reference materials for the determination of polychlorinated biphenyls and organochlorine pesticides from the National Metrology Institute of Japan. *Anal Bioanal Chem.* **2007**: 387, 2313–2323.

Osamu Kiguchi, Katsumi Saitoh, Nobuaki Ogawa. Simultaneous extraction of polychlorinated dibenzo-p-dioxins. Polychlorinated dibenzofurans and coplanar polychlorinated biphenyls from contaminated soil using pressurized liquid extraction. *Science Direct.* 2007: 1144, 262–268.

Pedro Antunes, Paula Viana, Tereza Vinhas, J.L. Capelo, J. Rivera, Elvira M.S. Gaspar.Optimization of pressurized liquid extraction (PLE) of dioxin-furans and dioxin-like PCB's from environmental samples. *Science Direct*. 2007: 75, 916–925. Wentao Wang, Bingjun Meng, Xiaoxia Lu, Yu Liu, Shu Tao. Extraction of polycyclic aromatic carbons and oranochlorine pesticides from soils: A comparison between Sochlet extraction, microwave-assisted extraction and accelerated solvent extraction techniques. *Science Direct.* **2007**: 602, 211–221.

Clara Coscolla, Vincent Yusa, Pedro Marti, Augustin Pastor Analysis of currently used pesticides in fine airborne particulate matter (PM 2.5) by pressurized liquid extraction and liquid chromatography-tandem mass spectrometry. *Journal of Chromatography A*. 2008: 1200, 100–107.

Rikard Westbom, Sune Sporring, Louise Cederberg, Lars-Ola Linderoth, Erland Bjorklund Selective pressurized liquid extraction of polychlorinated biphenyls in sediment. *Analytical Sciences*. **2008**: 24, 531–533.

Robert C. Hale, Stacy L Kim, Ellen Harvey, Mark J. La Guardia, T. Matt Mainor, Elizabeth O. Bush, Elizabeth M Jacobs Anarctic reasearch bases: Local sources of polybrominated diphenyl ether (PBDE) flame retardants. *Environmental Science and Technology*. 2007.

Adrain Covaci, Stefan Voorspoels, Lourdes Ramos, Hugo Neels, Ronny Blust. Recent developments in the analysis of brominated flame retardants and brominated natural compounds. *Journal of Chromatography A*. 2007: 1153,145–171.

Heather Stapleton, Nathan Dodder, Michele Schantz, Stephen Wise. Measurement of the Flame Retardants polybrominated diphenyl ehters (PBDEs) and hexabromocyldodecane (HBCCD) in house dust. *Organohalogen Compounds.* **2004**: 66, 3691–3695.

Augustina de la Cal, Ethel Eljarrat, Damia Barcelo. Determination of 39 polybrominated diphenyl ether congeners in sediment samples using fast selective pressurized liquid extraction and purification. *Journal of Chromatography A*. 2003: 1021,165–173.

Hanna Giergielewicz-Mozajska, Lukasz Da browski, and Jacek Namiesnik. Accelerated Solvent Extraction (ASE) in the analysis of environmental solid samples- some aspects of theory and practice. *Critical Reviews in Analytical Chemistry.* **2001**: 31(3), 149–165.

E. Kremer, B. Zygmunt. Accelerated solvent extraction-gas chromatographic determination of acidic herbicides in soil. *Chem. Anal.* (*Warsaw*). 2007: 52, 673–687.

E. Concha-Grana, V. Fernandez-Gonzalez, M. I. Turnes-Carou, S. Muniategui-Lorenzo, P. Lopez-Mahia. Pressurized liquid extraction of organochlorine pesticides from certified solid materials. *Journal of Chromatographic Science*. 2007: 45, 369–374.

Dongli Wang, Xiusheng Miao, Qing X. Li. Analysis of Organochlorine pesticides in coral (Porites evermanni) Samples using accelerated solvent extraction and gas chromatography/Ion trap mass spectrometry. *Arch Environ Contam Toxicol.* 2008: 54, 211–218.

S. P. J. van Leeuwen, J. de Boer. Advances in the gas chromatographic determination of persistent organic pollutants in the aquatic environment. *Journal of Chromatography A*. 2008: 1186, 161–182.

C. Lesueur, M. Gartner, A. Mentler, M. Fuerhacker. Comparison of four extraction methods for the analysis of 24 pesticides in soil samples with gas chromatography-mass spectometry and liquid chromatography-ion trap-mass spectrometry. *Talanta*. 2008: 75, 284–293.

Dongli Wang, Shannon Atkinson, Anne Hoover-Miller, Seilin L. Shelver, Qing X. Li. Simultaneous use of gas chromatography/ion trap mass spectometry- electron capture detection to improve the analysis of bromodiphenyl ethers in biological and environmental samples. *Rapid Communications in Mass Spectrometry.* 2008: 22, 647–656.

Ralf Bernhard Schaefer, Ralf Mueller, Werner Brack, Kalus-Dieter Wenzel, Georg Streck, Wolfgang Ruck, Matthias Liess. Determination of 10 particle-associated multiclass polar and semi-polar pesticides from small streams using accelerated solvent extraction. *Chemosphere*. 2008: 70 (11), 1952–1960. Anthony F. Lagalante, Timothy D. Oswald. Analysis of polybrominated diphenyl ethers (PBDE's) by liquid chromatography with negative-ion atmospheric pressure photoionization tandem mass spectrometry (LC/ NI-APPI/MS/MS): application to house dust. *Analytical and Bioanalytical Chemistry*. 2008: 391 (6), 2249–2256.

Jean-Baptiste Baugros, Cecile Cren-Olive, Barbara Giroud, Jean-Yves Gauvrit, Pierre Lanteri, Marie-Florence Grenier-Loustalot. Optimization of pressurized liquid extraction by experimental design for quantification of pesticides and alkyl phenols in sludge, suspended materials and atmospheric fallout by liquid chromatography-tandem mass spectrometry. *Journal of Separation Science.* **2009**: 32 (9), 1383–1389.

E. Schreck, F. Geret, L. Gontier, M. Treilhou. Development and validation of a rapid multiresidue method for pesticide determination using gas chromatography-mass spectrometry: A realistic case in vineyard soils. *Food Science and Technology*. **2009:** 42 (1) 233–238.

Jing You, Dongli Wang, Michael J. Lydy. Determination of pyrethroid insecticides in sediment by gas chromatography-Ion trap tandem mass spectrometry. *Talanta*. **2010**: 81 (1, 2) 0039–9140.

Pu Wang, Qinghua Zhang, Yawei Wang, Thanh Wang, Xiaomin Li, Lei Ding, Guibin Jiang. Evaluation of Soxhlet extraction, accelerated solvent extraction and microwave-assisted extraction for the determination of polychlorinated biphenyls and polybrominated diphenyl ethers in soil and fish samples. *Analytica Chimica Acta*. **2010**: 663 (1) 43–48.

T. L. Do, S. Lundstedt, P. Haglund. Selective pressurized liquid extraction of PCDD/Fs in soil: Optimization using experimental design. *Organohalogen Compounds*. **2010**: (72) 530–533.

G. Mascolo, V. Locaputo, G. Mininni. New perspective on the determination of flame retardants in sewage sludge by using ultrahigh pressure liquid chromatography-tandem mass spectrometry with different ion sources. *Journal of Chromatography, A.* **2010**: (27) 4601–4611.

Rahel C. Brandli, Thomas D. Bucheli, Thomas Kupper, Franz X. Stadelmann, Joseph Tarradellas. Optimised accelerated solvent extraction of PCBs and PAHs from compost. *International Journal of Environmental Analytical Chemistry.* **2006**: 86 (7) 505–525

Amal Al-Rashdan, Murad I.H. Helaleh, A. Nisar, Z. Al-Ballam. Simultaneous determination of polycyclic aromatic hydrocarbons (PAHs) and organochlorinated pesticides (OCPs) in sewage sludge using gas chromatography tandem mass spectrometry. *Current Analytical Chemistry.* **2010**: 6 (2) 177–183.

Donald C. Hilton, Richard S. Jones, Andreas Sjoedin. A method for rapid, non-targeted screening for environmental contaminants in household dust. *Journal of Chromatography*, A. **2010**: 1217 (44) 6851–6856.

Zulin Zhang, Murali Shanmugam, Stewart M. Rhind. PLE and GC-MS determination of Polybrominated Diphenyl Ethers in soils. *Chromatographia.* **2010**: 72 (7,8) 777.

Dongli Wang, Jing You, Michael J. Lydy. Sediment matrix effects in analysis of pyrethroid insecticides using gas chromatography-mass spectrometry. *Archives of Envirnomental Contamination and Toxicology*. **2010:** 59 (3) 382–392.

P. Lopez, S. A. Brandsma, P. E. G. Leonards, J. de Boer. Optimization and development of analytical methods for the determination of new brominated flame retardants and polybrominated diphenyl ethers in sediments and suspended particulate matter. *Analytical and Bioanalytical Chemistry*. **2011:** 400 (3) 871–883.

J. Malavia, F. J. Santos, M. T. Galceran. Simultaneous pressurized liquid extraction and clean-up for the analysis of polybrominated biphenyls by gas chromatography-tandem mass spectrometry. *Talanta*. **2011**: 84 (4) 1155–1162.

Kresinova, Milan Murzikar, Jana Olsovska, Tomas Cajthaml. Determination of 15 isomers of chlorobenzoic acid in soil samples using accelerated solvent extraction followed by liquid chromatography. *Talanta.* **2011:** 84 (4) 1141–1147.

## Additional product information available at www.thermoscientific.com/samplepreparation

©2012 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 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.

 Australia
 +61
 3
 9757
 4486

 Austria
 +43
 1
 616
 51
 25

 Benelux
 +31
 20
 683
 9768

 +32
 3
 353
 42
 94

 Brazil
 +55
 11
 3731
 5140

Ireland +353 1 644 0064 Italy +39 02 51 62 1267 Japan +81 6 6885 1213 Korea +82 2 3420 8600 Singapore +65 6289 1190



Sweden +46 8 473 3380 Switzerland +41 62 205 9966 Taiwan +886 2 8751 6655 UK +44 1276 691722 USA and Canada +847 295 7500

