



Accelerated Solvent Extraction Applications Summary

Fat Determination • Food Safety

Thermo
SCIENTIFIC

Contents

Introduction

The Accelerated Solvent Extraction (ASE) System.....	4
--	---

Fats

Determination of Unbound Fat in Various Food Matrices Using Accelerated Solvent Extraction (ASE)	5
Determination of Total Fat in Powdered Infant Formula Using Accelerated Solvent Extraction (ASE)	6
Rapid Determination of Fat in Meat Using Accelerated Solvent Extraction (ASE).....	7
Determination of Fat in Dried Milk Products Using Accelerated Solvent Extraction (ASE)	8
Extraction of Fat from Chocolate Using Accelerated Solvent Extraction (ASE)	9
Extraction of Fat from Dairy Products (Cheese, Butter, and Liquid Milks) using Accelerated Solvent Extraction (ASE).....	10
Extraction of Total Fat from Food Samples After Acid Hydrolysis Using Accelerated Solvent Extraction (ASE) with GC-MS Analysis	11

Food Safety

Selective Extraction of PCBs from Fish Tissue Using Accelerated Solvent Extraction (ASE).....	12
Accelerated Solvent Extraction (ASE) of Pesticide Residues in Food Products	13
Extraction of Lipids and Polychlorinated Biphenyls from Fish Tissue in a Single Run Using Accelerated Solvent Extraction (ASE)	14
Determination of PCBs in Large-Volume Fish Tissue Samples Using Accelerated Solvent Extraction (ASE).....	15
Determination of Pesticides in Large-Volume Food Samples Using Accelerated Solvent Extraction (ASE).....	16
Rapid Determination of Organochlorine Pesticides in Animal Feed Using Accelerated Solvent Extraction (ASE).....	17
Extraction of Zearalenone from Wheat and Corn by Accelerated Solvent Extraction (ASE).....	18
Rapid Determination of Persistent Organic Pollutants (POPs) Using Accelerated Solvent Extraction (ASE)	19
Rapid Extraction and Determination of Arsenicals in Fish Tissue and Plant Material Using Accelerated Solvent Extraction (ASE).....	20
Determination of Perchlorate in Vegetation Samples Using Accelerated Solvent Extraction (ASE) and Ion Chromatography.....	21
Extraction and Cleanup of Acrylamide in Complex Matrices Using Accelerated Solvent Extraction (ASE) Followed by Liquid Chromatography Tandem Mass Spectrometry (LC-MS/MS)	22
Extraction of Contaminants, Pollutants, and Poisons from Animal Tissue Using Accelerated Solvent Extraction (ASE).....	23

References

Accelerated Solvent Extraction (ASE) References.....	24
--	----

The Accelerated Solvent Extraction (ASE) System



Better extractions in less time with less solvent

Accelerated Solvent Extraction (ASE) eliminates many of the manual steps involved in preparing food samples for analysis, which helps ensure increased reproducibility and accelerates the process significantly. ASE is an automated extraction technique that uses elevated temperatures and pressures to achieve extractions in very short periods of time (a 10-g sample can be extracted in less than 15 minutes using less than 15 mL solvent).

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 the highest temperatures at which extractions take place in these techniques will be the boiling point of the solvent.

If sufficient pressure is exerted on the solvent during the extractions,

temperatures above the boiling point can be used. This means that 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 helps the extraction process to happen more quickly. 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 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. The net result means performing extractions at elevated temperatures uses less time and with less solvent.

In the food industry, time can be critical. Quality control checks need to be completed quickly and accurately. Whether a laboratory is analyzing the fat content of food for labeling or analyzing food for environmental contaminants, ASE is a powerful tool for preparing these samples in the fastest time possible while ensuring accurate results.

The Application Briefs compiled here show conditions for extracting fats and environmental contaminants from a variety of food and beverage samples. To view the complete Application Notes, visit our website at www.thermoscientific.com/dionex.



Thermo Scientific Dionex ASE 150 and ASE 350 Accelerated Solvent Extraction instruments.

Determination of Unbound Fat in Various Food Matrices Using Accelerated Solvent Extraction (ASE)



Introduction

Accelerated Solvent Extraction (ASE) is an extraction method that significantly streamlines sample preparation. A commonly used solvent is pumped into an extraction cell containing the sample, which is then brought to an elevated temperature and pressure. Minutes later, the extract is transferred from the heated cell to a standard collection vial for cleanup or analysis. The entire extraction process is fully automated and performed in minutes for fast and easy extraction with low solvent consumption. Up to 24 samples can be loaded and extracted sequentially without requiring operator intervention. Recently, the requirements for accurate labeling of fats in foods were revised by the U.S. Food and Drug Administration (U.S. FDA) and the U.S. Department of Agriculture (U.S. DA). This occurred as a result of the Nutrition Labeling and Education Act, which requires the labeling of total saturated and unsaturated fats contained in foods. Though these laws do not directly affect food sold outside of the U.S., there seems to be increased awareness worldwide of fat content in foods. In addition, food manufacturers require a method for the consistent determination of fat content for quality-control purposes.

Accelerated solvent extraction of unsaturated fats in popular snack foods

Equipment

Dionex ASE 200 Accelerated Solvent Extractor* equipped with either 11 or 22 mL stainless steel extraction cells

Analytical balance

Collection vials, 40 mL (P/N 048783) and 60 mL (P/N 048784)

Cellulose filters (P/N 049458)

Solvents

Petroleum ether

Chloroform

Hexane

Isopropanol

Ethanol

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

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

Analysis

Gravimetric

Extraction Conditions

Oven Temperature: 125 °C

Pressure: 1500 psi

Oven Heatup Time: 6 min

Static Time: 5–25 min

Flush Volume: 60%

Purge Time: 60 s

Solvent: Petroleum ether, chloroform, hexane, or hexane/isopropanol (3:2), chloroform/ethanol (1:1), depending on application

Static Cycles: 1 to 3

Results

Extraction of fat from snack crackers: comparison of results by soxhlet and ASE.

Sample	Method	Avg % Fat (wt.%)	Std. Dev.	RSD (%)
Cracker 1	Soxhlet*	15.4	N/A	N/A
Cracker 1	ASE**, n = 3	14.6	0.09	0.65
Cracker 2	Soxhlet*	28–30	N/A	N/A
Cracker 2	ASE**, n = 3	28.1	0.20	0.70

* After acid hydrolysis

** Conditions: 5 g sample, 125 °C, 1500 psi, 6 min heatup, 25 min static, 60% flush, 60 s purge, 1 static cycle, hexane/isopropanol (3:2).

Determination of Total Fat in Powdered Infant Formula Using Accelerated Solvent Extraction (ASE)



Introduction

Accurate determination of fat in certain foods is difficult due to the binding or entrapment of the fat by the matrix. Most methods used to determine fat in these difficult matrices include a pretreatment step to denature or destroy the physical structure of the matrix and allow greater accessibility to the fat.

Common fat determination methods used by the dairy industry are the Roesse-Gottlieb and Modified Mojonnier methods (AOAC Intl. Methods 905.02 and 989.05, respectively). These methods specify the use of ammonium hydroxide to dissolve casein and liberate the fat. Here, Accelerated Solvent Extraction is compared to the Mojonnier method for extraction of fat from infant formula.

Saturated and unsaturated fats

Equipment

Dionex ASE 200 Accelerated Solvent Extractor* equipped with 11-mL stainless steel extraction cells
Cellulose filters (P/N 049458)
Analytical balance (0.001 g or better)
Mortar and pestle (Fisher Scientific or equivalent)
Solvent evaporator
Forced air oven
Ottawa Sand Standard (P/N S23-3)

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

Reagent

ASE Prep DE (diatomaceous earth) (P/N 062819)

Solvents

Acetone
Hexane
Water

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

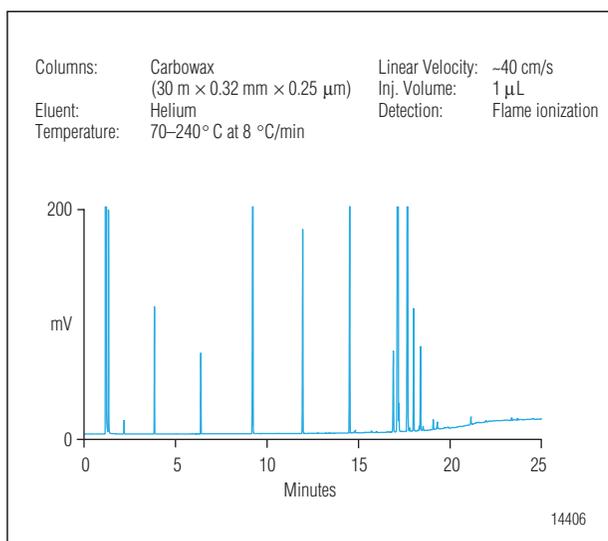
Analysis

GC-FID

Extraction Conditions

Extraction Solvent: Hexane acetone, 4:1 volume
Temperature: 125 °C
Pressure: 1500 psi
Heat Time: 6 min
Static Time: 5 min
Flush Volume: 100%
Purge Time: 60 s
Static Cycles: 3
Total Extraction Time: 24 min per sample

Results



Chromatogram of FAMES from ASE extraction of infant formula.

Rapid Determination of Fat in Meat Using Accelerated Solvent Extraction (ASE)



Introduction

Soxhlet extraction is an accepted technique for extracting fat from meat samples. Though it is simple and robust, there are drawbacks to Soxhlet extraction, such as long drying and extraction times, lack of automation, and the amount of solvent used per sample. ASE is a technique that was developed to replace Soxhlet and other extraction techniques for many samples. The automation and rapid extraction time of ASE overcome the shortcomings of Soxhlet extraction.

Fresh and processed meats

Equipment

ASE 200 Accelerated Solvent Extractor,* with 11 or 22 mL stainless steel extraction cells
 Cellulose Filters (P/N 049458)
 Collection Vials, 40 mL (P/N 048783)
 Ottawa Sand Standard (Fisher Scientific)
 Analytical balance (to read to nearest 0.0001 g or better)
 Mortar and pestle (Fisher Scientific or equivalent)
 Solvent evaporator
 Forced air oven
 Microwave oven (800 W) with carousel
 *ASE 150 and 350 can be used for equivalent results.

Reagent

ASE Prep DE (diatomaceous earth) (P/N 062819)

Solvents

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

Analysis

Gravimetric

Extraction Conditions

Solvent: Petroleum ether or hexane* extraction solvent
 Temperature: 125 °C
 Pressure: 1500 psi
 Heatup Time: 6 min
 Static Time: 1 or 2 min**
 Flush Volume: 60%
 Purge Time: 60 s
 Cycles: 2
 Total Time: 12 min
 Total Solvent: 20 mL

*Petroleum ether and hexane were found to be equivalent as extraction solvents for fat in meat.

**When extracting more than 1 g of a high-fat sample, a 2 min static time may be beneficial.

Results

Percent fat in low-fat processed meat samples (ASE vs soxhlet).

Sample	ASE Run #1	ASE Run #2	ASE Run #3	ASE Average	Standard Deviation	Soxhlet
Beef	2.82	2.90	2.83	2.85	0.046	2.81
Chicken	0.84	0.84	0.79	0.82	0.025	0.75
Ham	1.85	1.74	1.87	1.82	0.069	1.72
Franks	1.90	1.97	1.97	1.94	0.041	1.54
Turkey	1.04	1.00	1.04	1.02	0.026	0.94

Determination of Fat in Dried Milk Products Using Accelerated Solvent Extraction (ASE)



Introduction

Many extraction techniques for the determination of fat in food are labor-intensive or require long extraction times. The Roese-Gottlieb (RG) method requires alkaline pretreatment of the sample before a labor intensive liquid-liquid extraction. The Schmidt-Bondzynski-Ratzlaff (SBR) method calls for acid digestion before liquid-liquid extraction of the sample. Some Soxhlet methods require extraction times from 4 to 24 hours in duration.

Low- and high-fat samples

Equipment

Dionex ASE 200 Accelerated Solvent Extractor* equipped with 11-mL stainless steel extraction cells
Cellulose filters (P/N 049458)
Analytical balance (0.0001 g or better)
Solvent evaporator
Forced air oven

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

Solvents

Hexane
Dichloromethane
Methanol
Petroleum ether (40-60 °C boiling range)
All solvents are pesticide-grade or equivalent and available from Fisher Scientific.

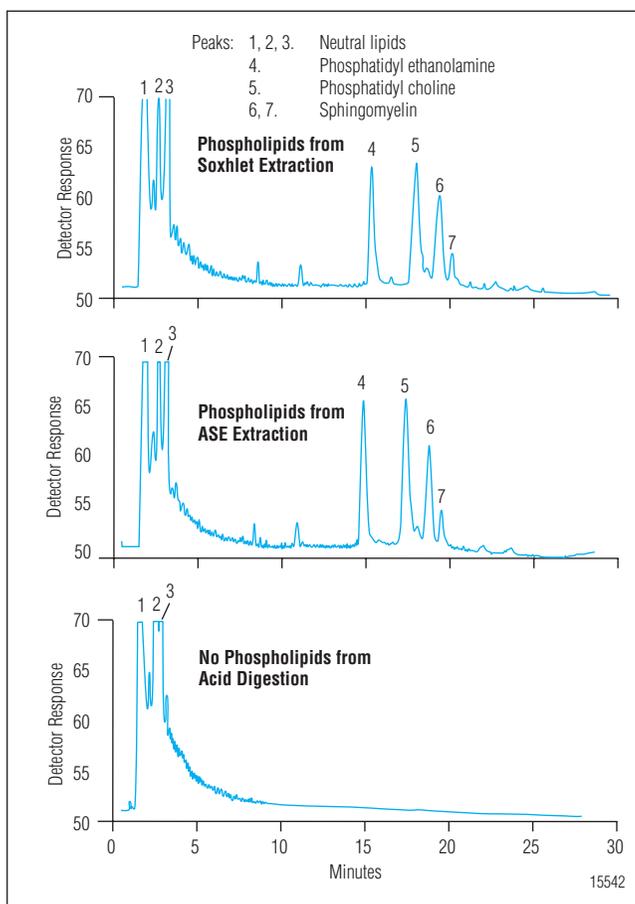
Analysis

HPLC

Extraction Conditions

Temperature: 80 °C
Pressure: 1500 psi
Heatup Time: 5 min
Static Time: 1 min
Flush Volume: 100%
Purge Time: 40 s
Static Cycles: 3
Total Time: 11 min
Total Solvent: <30 mL
Solvent: Mixtures of hexane, dichloromethane, and methanol in various volume ratios

Results



HPLC with ELSD traces of lipids extracted from whey protein concentrate.

Extraction of Fat from Chocolate Using Accelerated Solvent Extraction (ASE)



Introduction

Accelerated Solvent Extraction (ASE) is a new way to speed up gravimetric fat determination of chocolate products and greatly reduce the amount of solvent used. The ASE system uses a combination of elevated temperature and pressure to increase the extraction kinetics, thus decreasing time and solvent consumption. Current methods for determining the fat content in chocolate are labor-intensive and require large amounts of solvent and time. For example, the Mojonnier ether extraction method takes 2–3 h and over 110 mL of solvent and requires the laboratory technician to be present for most of the extraction. Using ASE, extraction time is reduced to 18 min and solvent use to 20 mL. ASE has been shown to produce comparable if not better results than the current methods. Furthermore, the ASE process is fully automated, making it possible to extract up to 24 samples unattended.

Fats in solid and powdered chocolate samples

Equipment

Dionex ASE 200 Accelerated Solvent Extractor* with 11-mL stainless steel extraction cells
Cellulose Filters (P/N 049458)
Collection Vials, 40 mL (P/N 048783)
Analytical balance (to read to the nearest 0.0001 g or better)
Mortar and pestle (Fisher Scientific or equivalent)
Solvent evaporator
Forced air oven

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

Reagent and Solvent

ASE Prep DE (diatomaceous earth) (P/N 062819)
Petroleum ether (pesticide grade or equivalent; Fisher Scientific)

Analysis

Gravimetric

Extraction Conditions

Solvent: Petroleum ether 100%
Temperature: 125 °C
Pressure: 1500 psi
Heatup Time: 6 min
Static Time: 3 min
Flush Volume: 60%
Purge Time: 60 s
Cycles: 3
Total Time: 18 min
Total Solvent: 20 mL

Results

Baking chocolate (top) and milk chocolate (bottom) % fat* recovery: ASE vs mojonnier method (n = 3).

	ASE	Mojonnier
Average	52.80	51.69
SD	0.35	0.26
RSD	0.67	0.50

*%fat = (residue/sample wt.) × 100

	ASE	Mojonnier
Average	31.80	32.34
SD	0.32	0.33
RSD	1.02	1.02

*%fat = (residue/sample wt.) × 100

Extraction of Fat from Dairy Products (Cheese, Butter, and Liquid Milks) Using Accelerated Solvent Extraction (ASE)



Introduction

The current methods for determining fat in dairy products, though acceptable, have several drawbacks. Many dairy-based products require a pretreatment prior to extraction. This denatures the casein, allowing greater exposure of the fat to the solvents. For example, a 1-g sample of cheese must be pretreated with ammonium hydroxide followed by hydrochloric acid and boiled for several minutes. The standard fat extraction methods, including the pretreatment, are very time consuming. Large amounts of solvents are required to remove the fat from each sample matrix, which can be costly. For example, manual extraction of pretreated cheese usually requires 2–3 h and more than 110 mL of solvent per sample. Therefore, the standard fat extraction methods are not time- or cost-efficient and can expose laboratory technicians to potentially dangerous solvents.

Extraction of dairy sample matrices made simple!

Equipment

ASE 200 Accelerated Solvent Extractor* equipped with 11- and 33-mL stainless steel extraction cells

Cellulose Filters (P/N 049458)

Collection Vials, 40 mL (P/N 048783) and 60 mL (P/N 048784)

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

Mortar and pestle (Fisher Scientific or equivalent)

Solvent evaporator

Heated block

Forced air oven

Ottawa Sand Standard (P/N S23-3)

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

Reagent

ASE Prep DE (diatomaceous earth) (P/N 062819)

Solvents

Petroleum ether

Acetone I

Isopropanol

Hexane

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

Analysis

Gravimetric

Extraction Conditions

Cheese

Solvent: Hexane: isopropanol (3:2)

Temperature: 110 °C

Pressure: 1500 psi

Cell Heatup Time: 6 min

Static Time: 2 min

Flush Volume: 100%

Purge Time: 60 s

Cycles: 3

Total Time: 10 min

Total Solvent: <30 mL

Butter

Solvent: Petroleum ether: acetone (3:2)

Temperature: 100 °C

Pressure: 1500 psi

Heatup Time: 5 min

Static Time: 2 min

Flush Volume: 60%

Purge Time: 60 s

Cycles: 1

Total Time: 8 min

Total Solvent: <30 mL

Milk and Cream

Sample: Cream (40%)

Solvent: Petroleum ether:acetone:isopropanol (3:2:1)

Sample: Whole milk (4–6%)

Solvent: Petroleum ether:isopropanol (2:1)

Sample: Homogenized/UHT milk (3%)

Solvent: Petroleum ether:isopropanol (3:2)

Sample: Skim milk (0.1%)

Solvent: Petroleum ether:isopropanol (3:2)

Temperature: 120 °C

Pressure: 1500 psi

Cell Heatup Time: 6 min

Static Time: 1 min

Flush Volume: 100%

Purge Time: 60 s

Cycles: 3

Total Time: 10 min

Total Solvent: <30 mL

Results

Milk and cream, % fat recovery: ASE vs. RG method.

Sample	ASE Mean ± SD (n)	RG Method
Cream	40.62 ± 0.06 (3)	40.58
Whole Milk	4.42 ± 0.02 (4)	4.50
Homogenized Milk	3.39 ± 0.03 (6)	3.39
Skim Milk	0.053 ± 0.010 (7)	0.053

Extraction of Total Fat from Food Samples After Acid Hydrolysis Using Accelerated Solvent Extraction (ASE) with GC-MS Analysis



Introduction

Sample preparation—specifically, solvent extraction—is an important step in the analytical process. For many years, analysts have used an array of solvent extraction techniques including Soxhlet, shaking, sonication, and blending.

ASE technology provides a flow-thru solvent extraction system that increases productivity while decreasing cost and providing a platform for automation. Complex matrices such as food typically require acid hydrolysis or pretreatment prior to solvent extraction.

Pretreatment or hydrolysis of these matrices is often necessary to facilitate complete extraction of lipids from the sample. Time-consuming and labor-intensive liquid extraction techniques such as Soxhlet, automated Soxhlet, and Mojonnier extraction are typically used to extract fatty acids after acid hydrolysis.

Turnkey extraction and analysis of snack and processed foods

Equipment

Dionex ASE 150 or 350 with pH-hardened pathway (P/N 066401 or 066230)
 Dionium extraction cells (100 mL) (P/N 068103)
 Glass fiber filters (P/N 056781)
 Collection bottles (250 mL) (P/N 056284)
 Collection vials (40 mL) (P/N 048783)
 GC-MS
 Carbowax capillary GC column
 Pressure tubes (ACE Glass Inc.)

Solvents and Reagents

Chloroform
 Pyrogallol
 Alcohol; reagent-grade
 Hexane
 Ethyl ethern
 ASE Prep DE P/N 062819)
 ASE Prep CR P/N 080024)
 8.3 M HCL
 Toluene
 12% BF3 in MeOH
 Na₂SO₄

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

Analysis

GC-MS

Extraction Conditions

Pressure: 1500 psi*
 Temperature: 100 °C
 Solvent: Hexane
 Static Time: 5 minutes
 Static Cycles: 3
 Flush Volume: 70%
 Purge Time: 120 sec

*Pressure studies show that 1500 psi is the optimum extraction pressure for all ASE extraction.

Results

Extraction results using Mojonnier techniques and ASE (n = 3).

Mayonnaise	Average	RSD	%RSD
Mojonnier	75.1	0.89	1.18
ASE	74.2	0.43	0.575
Corn Chips			
Mojonnier	30.41	0.37	1.21
ASE	29.85	0.33	1.10
Parmesan Cheese			
Mojonnier	26.41	0.284	1.08
ASE	26.27	0.220	0.839
Baked Shortbread			
Mojonnier	13.95	0.033	0.238
ASE	14.07	0.451	3.20
Bologna			
Mojonnier	25.58	0.275	0.968
ASE	28.60	0.375	1.31

Selective Extraction of PCBs from Fish Tissue Using Accelerated Solvent Extraction (ASE)



Introduction

Accelerated Solvent Extraction (ASE) is a new extraction method that significantly streamlines sample preparation. A solvent is delivered into an extraction cell containing the sample, which is then brought to an elevated temperature and pressure. Minutes later, the extract is transferred from the heated cell to a standard collection vial for cleanup or analysis. The entire extraction process is fully automated and performed in minutes for fast and easy extraction with low solvent consumption.

The data presented in this application brief demonstrate that selective extractions can be performed using ASE with the proper choice of solvent and sorbent in the extraction cell. Results are given for the recovery of PCBs from contaminated fish tissue showing that extracts can be obtained using ASE that do not require further cleanup prior to analysis by gas chromatography.

Environmental contaminants in fish

Equipment

Dionex ASE 200 Accelerated Solvent Extractor* equipped with 11-, 22-, or 33-mL stainless steel extraction cells

Analytical balance

Collection vials, 40 mL (P/N 048783) and 60 mL (P/N 048784)

Cellulose filter (P/N 049458)

Gas chromatograph (GC) with electron capture detector (ECD)

*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.

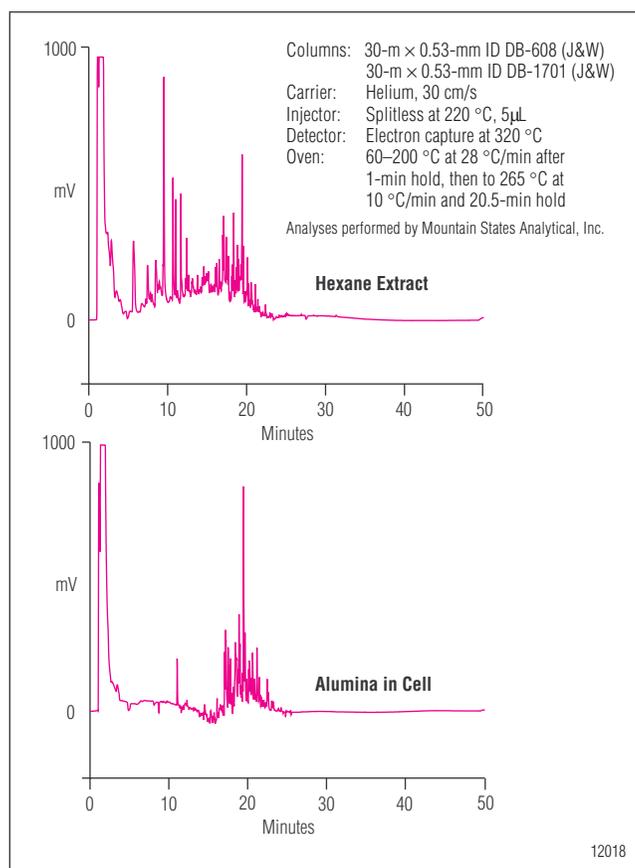
Analysis

GC-ECD

Extraction Conditions

Extraction Solvent:	Hexane
Temperature:	100 °C
Pressure:	1500 psi
Heat Time:	5 min
Static Time:	5 min
Flush Volume:	60%
Purge Time:	90 s
Static Cycles:	2
Total Extraction Time:	17 min per sample

Results



Chromatograms obtained from the nonselective ASE extraction of the fish tissue (top) and from the selective ASE extraction of a portion of the same sample (bottom).

Accelerated Solvent Extraction (ASE) of Pesticide Residues in Food Products



Introduction

Residue analysis in crops and food products is routinely performed in regulatory and industrial laboratories around the world. Many of the traditional procedures used to perform these extractions are time-consuming and solvent-intensive. Accelerated Solvent Extraction (ASE) is an extraction technique that speeds the extraction process and reduces the total amount of solvent used. The system uses conventional liquid solvents at elevated temperatures and pressures, which results in increased extraction kinetics. Extraction of samples ranging from 1 to 30 g typically requires 12–17 min and 15–50 mL of solvent.

Organochlorine residues in bananas and potatoes

Equipment

Dionex ASE 200 Accelerated Solvent Extractor* equipped with 11, 22, or 33 mL stainless steel extraction cells

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

Cellulose filter (P/N 049458)

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

Reagents

Acetonitrile, Optima grade (Fisher Scientific)

ASE Prep DE (diatomaceous earth) (P/N 062819)

Solvents

Acetone

Acetonitrile

Hexane

Sodium sulfate, anhydrous (added after extraction)

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

Analysis

GC-ECD

Extraction Conditions

Temperature:	100 °C
Pressure:	1500 psi
Heatup Time:	5 min
Static Time:	5 min
Flush Volume:	60%
Purge Time:	100 s
Static Cycles:	1–2
Total Extraction Time:	14–18 min per sample
Total Solvent Used:	15–45 mL per sample

Results

Recovery of organochlorine pesticides spiked onto raw banana at the 100 ppm level.*

Compound	Av. Recovery (%)	SD (µg/kg)	RSD (%)
α-BHC	100.3	2.3	2.3
α-BHC	102.2	2.3	2.3
α-BHC	98.9	3.2	3.2
Heptachlor	89.2	7.6	8.5
Aldrin	89.4	2.2	2.5
Heptachlor Epoxide	93.5	2.1	2.2
Dieldrin	93.7	1.6	1.7
4,4'-DDE	92.1	1.8	1.9
2,4'-DDD	95.4	2.5	2.6
Endrin	94.4	2.7	3.0
4,4'-DDD	88.0	2.7	3.0
4,4'-DDT	89.6	5.8	6.4

*N = 3

Extraction of Lipids and Polychlorinated Biphenyls from Fish Tissue in a Single Run Using Accelerated Solvent Extraction (ASE)



Introduction

The presence of polychlorinated biphenyls (PCBs) in fish and other marine organisms is of immediate environmental and regulatory concern. To determine the concentrations of PCBs indicative of contaminant exposure, and lipids, which also characterize physiological conditions of fish, the analytes must first be extracted. Traditionally, this has been done using Soxhlet methods that are both time consuming and use large volumes of solvent. This application note summarizes the use of ASE to quickly and efficiently extract lipids and PCBs in a single 20 min extraction using only 40 mL of solvent.

Fats and contaminants in fish

Equipment

ASE 200 Accelerated Solvent Extractor* equipped with 33 mL stainless steel extraction cells
 ASE Solvent Controller (optional)
 Gas Chromatograph
 Methylsiloxane column
 Microwave oven (800 W) with carousel
 Analytical balance
 Graduated Conc. Vial (P/N 055442)
 Analytical Evaporator
 Screw-on Stainless Steel Funnel (P/N 049288)
 Cellulose Filter Insertion Tool (P/N 049495)
 Cellulose Filters (P/N 049458)
 *ASE 150 and 350 can be used for equivalent results.

Solvents

Hexane
 PCB Aroclor Standard
 Sulfuric Acid, ACS grade or equivalent

Reagent

ASE Prep DE (diatomaceous earth) (P/N 062819)

Analysis

GC-ECD

Extraction Conditions

Accelerated Solvent Extractor

Solvent: Hexane
 System Pressure: 1500 psi*
 Oven Temperature: 25 °C
 Sample Size: 10 g
 Heatup Time: 6 min
 Static Time: 5 min
 Static Cycles: 2
 Flush Volume: 60% of extraction cell volume
 Nitrogen Purge: 1 MPa (150 psi) for 60 s
 Total Extract Volume: 40 mL
 Total Extraction Time: 20 min

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

Gas Chromatograph

Column: Methylsiloxane
 Carrier Gas: Helium (16 psi)
 Detector: Electron Capture Detector
 Temperature: 300 °C
 Temperature Program: 100 °C (hold for 2 min), 100–160 °C at 15 °C/min, followed by 160–270 °C at 5 °C/min
 Injector Temperature: 225 °C

Results

PCBs in high-fat fish tissue: comparison of soxhlet and ASE results.

Method	Solvent	Avg. PCBs (µg/g)	Std. Dev. (µg/g)	RSD (%)
Soxhlet	Hexane/Acetone (1:1)	0.19	NA	NA
ASE, N = 3	Hexane	0.21	0.01	4.8

Determination of PCBs in Large-Volume Fish Tissue Samples Using Accelerated Solvent Extraction (ASE)



Introduction

Extraction and analysis of PCBs from fish and other marine tissues continues to be a necessary step in the monitoring of the aquatic food supply. Due to the large number of samples requiring analysis, automated extraction systems have proven useful in this area. ASE technology has been shown to produce good recoveries of naturally occurring PCBs from fish tissue samples, and is approved for use in U.S. EPA SW-846 Method 3545 for the extraction of PCBs, OCPs, BNAs, OPPs, herbicides, and dioxins and furans. ASE was designed to replace time-consuming and solvent-intensive methods such as Soxhlet and sonication in the environmental area. ASE operates at temperatures higher than those possible in traditional techniques, thus increasing the efficiency of the extraction process.

Large-volume capacity yields increased sample throughput

Equipment

Dionex ASE 300 Accelerated Solvent Extractor* equipped with 100-mL stainless steel extraction cells

Collection vials, 250 mL (P/N 056785)

Cellulose filters (P/N 056780)

Gas Chromatograph equipped with electron capture detector (ECD)

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

Analysis

GC-ECD

Reagents and Standards

Methylene chloride (Optima Grade, Fisher Scientific)
ASE Prep DE (diatomaceous earth) (P/N 062819)
Alumina (basic, Brockman activity I, Fisher Scientific)
PCB standards (ULTRA Scientific Inc.)

Extraction Conditions

Extraction Solvent: Methylene chloride
Temperature: 125 °C
Pressure: 1500 psi
Heatup Time: 5 min
Static Time: 3 min
Flush Volume: 60%
Purge Time: 120 s
Static Cycles: 3
Total Extraction Time: 18 min per sample

Results

Recovery of spiked PCB congeners from 30-g fish tissue samples using selective ASE extraction conditions.

Cogener	BZ #	Spike (µg)	% Recovery	% RSD
2-Chlorobiphenyl	1	2.5	99.8	3.0
2,3-Dichlorobiphenyl	5	2.5	103.8	8.8
2,4,5-Trichlorobiphenyl	29	2.5	107.1	3.1
2,2',4,6-Tetrachlorobiphenyl	50	5	98.4	2.4
2,2',3,4,5'-Pentachlorobiphenyl	87	5	92.3	7.9
2,2',4,4',5,6'-Hexachlorobiphenyl	154	5	89.0	5.9
2,2',3,4',5,6,6'-Heptachlorobiphenyl	186	7.5	91.1	8.5
2,2',3,3',4,5',6,6'-Octachlorobiphenyl	200	7.5	96.0	6.5
Decachlorobiphenyl	209	12.5	94.2	8.7

Determination of Pesticides in Large-Volume Food Samples Using Accelerated Solvent Extraction (ASE)



Introduction

Pesticide residue analysis in crops and food products is performed in regulatory and industrial laboratories around the world. Many of the traditional procedures used to perform the extractions for these analyses are time consuming and solvent intensive. Accelerated Solvent Extraction (ASE) is an extraction technique that speeds the extraction process and reduces the total amount of solvent used. The system uses conventional solvents at elevated temperatures and pressures, which results in improved extraction kinetics. The extraction of samples ranging from 1 to 30 g typically requires 12–17 min and 15–50 mL of solvent.

Extraction of samples up to 30 g have been reported using the Dionex ASE 200 extractor with an upper limit sample cell size of 33 mL. However, for many pesticide residue analyses, this volume is insufficient. Food samples such as fruit and vegetables have very high water contents and must be mixed with desiccants such as sodium sulfate to achieve quantitative pesticide recovery. In this case, the actual weight of the sample extracted will be much less than 30 g. The Dionex ASE 300 has the capability to extract samples with volumes as large as 100 mL. This capability allows the direct extraction of food and vegetable samples with weights in the 30 to 50-g range.

High-volume capacity for ultralow detection levels of pesticide residues

Equipment

Dionex ASE 300 Accelerated Solvent Extractor* with 34-, 66-, or 100 mL stainless steel extraction cells

Dionex vials (250 mL) for collection of extracts (P/N 056785)

Cellulose filters (P/N 056780)

*ASE 150 and 350 can be used for equivalent results

Analysis

GC

Solvents and Reagents

Acetone, Optima grade (Fisher Scientific)

Methylene chloride, Optima grade (Fisher Scientific)

Ethyl acetate, Optima grade (Fisher Scientific)

Hexane, Optima grade (Fisher Scientific)

Cyclohexane, Optima grade (Fisher Scientific)

ASE Prep DE (diatomaceous earth) (P/N 062819)

Sodium sulfate, anhydrous (Fisher Scientific)

Extraction Conditions

Temperature: 100 °C

Pressure: 1500 psi

Solvent: Ethyl acetate/cyclohexane or MeCl₂/acetone (1:1, v/v)

Heatup Time: 5 min

Static Time: 5 min

Flush Vol.: 60%

Purge Time: 180 s

Static Cycles: 1–2

Total Extraction Time: 14–20 min per sample

Total Solvent: 135–145 mL per sample

Results

Percent recovery of organophosphorus pesticides from apple puree fortified at 50 ppb.

Apples	1	2	3	4	5	6	7	8	9	10	11	12	Mean (%)	SD	RSD (%)
Dichlorvos/Naled	76	80	93	82	80	97	102	95	97	90	67	92	87	10	12
Mevinphos	91	96	105	93	90	108	115	111	110	104	71	106	100	12	12
TEPP	117	141	124	120	96	126	144	137	150	107	79	115	121	20	16
Demeton-O	64	78	47	44	64	67	83	71	75	51	64	77	65	12	19
Ethoprophos (Ethoprop)	84	86	105	91	87	106	110	106	103	91	70	97	95	11	12
Sulfotep	94	100	101	95	88	102	105	101	108	87	72	90	95	10	10
Phorate	80	84	85	77	83	88	100	93	95	83	71	89	86	8	9
Demeton-S	60	68	45	46	72	55	65	55	72	41	63	73	59	11	18
Dimethoate	128	125	146	133	106	141	148	140	149	115	81	121	128	19	15
Diazinon	87	92	101	86	86	99	107	101	104	91	73	93	93	9	10
Disulfoton	59	73	46	44	66	59	78	63	75	52	66	80	63	11	18
Parathion-methyl	91	95	104	89	88	103	108	104	101	94	70	95	95	10	10
Fenclorophos	82	89	101	86	85	100	103	99	96	91	71	93	91	9	10
Malathion	87	96	106	97	84	106	116	104	105	82	62	89	94	14	15
Fenthion	82	89	87	79	83	91	98	93	94	82	71	90	86	7	8
Chlorpyrifos	89	97	94	82	84	101	99	100	95	89	70	87	91	9	10
Parathion-ethyl	100	104	105	99	87	104	109	106	118	92	75	91	99	11	11
Trichloronat	80	91	98	83	82	99	98	95	96	90	68	90	89	9	10
Tetrachlorvinphos	87	90	100	87	85	95	100	97	98	91	71	94	91	8	9
Prothiofos	76	87	93	78	78	93	93	91	97	85	64	90	85	9	11
Merphos	78	79	91	76	74	88	89	88	96	82	63	83	82	9	10
Fensulfothion	91	92	113	93	90	106	110	106	105	100	71	95	98	11	11
Sulprofos	70	85	80	70	76	86	92	82	88	76	64	85	80	8	10
EPN	96	103	100	97	88	105	111	103	111	89	70	93	97	11	11
Azinphos-methyl	95	99	106	87	84	105	111	104	111	99	75	104	98	11	11
Coumaphos	95	98	102	92	89	102	110	102	106	101	80	96	98	8	8

Rapid Determination of Organochlorine Pesticides in Animal Feed Using Accelerated Solvent Extraction (ASE)



Introduction

Animal feed contaminated with organochlorine pesticides (OCPs) has begun to attract worldwide attention. When ingested, the OCPs from animal feed tend to accumulate in certain animal products, especially those rich in fat, such as meat, milk, and butter. Because these types of animal products are widely consumed by humans, methods are needed that quickly extract and determine OCPs in the feeds of animals used to produce products for human consumption. Traditional methods used to extract OCPs from animal feed require large amounts of organic solvents and take from one to several hours per extraction. Also, many of the traditional methods are very labor intensive and require constant analyst attention.

Environmental contaminants in feedstocks

Equipment

Dionex ASE 200 Accelerated Extractor* with Solvent Controller
11-mL stainless steel extraction cells (P/N 055422)
Dionex cellulose filters (P/N 049458)
Collection vials, 40 mL (P/N 048783)
Analytical balance (accurate to the nearest 0.0001 g or better)
Laboratory grinder
Ottawa Sand Standard (P/N S23-3)
Dichloromethane silica gel, 0.063–0.200 mm, water content 2.62% (Merck, Darmstadt, Germany)
S-X3 Bio-Beads® (Bio Rad Laboratories)

*ASE and 350 can be used for equivalent results.

Analysis

GC

Reagents

For reagents, use either:
Bulk Isolute® Sorbent (International Sorbent Technology Ltd., UK)
ASE Prep DE (diatomaceous earth) (P/N 062819)

Solvent

Petroleum ether (All solvents are pesticide-grade or equivalent and available from Fisher Scientific.)

Extraction Conditions

Solvent: Petroleum ether 100%
Temperature: 125 °C
Pressure: 1500 psi
Heatup Time: 6 min
Static Time: 3 min
Flush Volume: 60%
Purge Time: 60 s
Cycles: 3
Total Time: 18 min
Total Solvent: 20 mL

Results

Concentration values (ng g⁻¹) and RSD (%) for the extraction of CRM BCR 115.

Compounds	Certified Value		ASE (n = 3)	
	C (ng g ⁻¹)	RSD (%)	C (ng g ⁻¹)	RSD (%)
α-HCH	*	*	21.5 ± 0.5	2.5
HCB	19.4 ± 1.4	7.2	20.6 ± 0.4	1.8
β-HCH	23 ± 3	13.0	26.0 ± 2.3	8.7
γ-HCH	21.8 ± 2	9.2	27.1 ± 1.4	5.3
Heptachlor	19 ± 1.5	7.9	20.0 ± 0.5	2.7
Aldrin	*	*	56.0 ± 3.1	5.5
p,p'-DDE	47 ± 4	8.5	54.6 ± 2.6	4.7
Dieldrin	18 ± 3	16.7	22.0 ± 0.6	2.6
Endrin	46 ± 6	13.0	52.1 ± 1.9	3.6
p,p'-DDD	*	*	91.8 ± 2.6	2.8
o,p'-DDT	46 ± 5	10.9	49.8 ± 0.5	1.1
p,p'-DDT	*	*	59.4 ± 1.8	3.1

* Present but not certified.

Extraction of Zearalenone from Wheat and Corn by Accelerated Solvent Extraction (ASE)



Introduction

Zearalenone (ZON) is a mycotoxin produced by the Fusarium fungus. ZON can be found in a wide variety of plants and soils, and can have negative health effects on animal husbandry and humans. Traditional methods for extracting ZON from soils or animal feed include wrist shaking or blending. These methods normally take 30–60 min per sample with constant lab technician attendance. Because of the time-consuming nature of these traditional extraction techniques, many sample prep labs experience large bottlenecks that hinder the flow of samples to the analytical lab.

Mycotoxins in grains

Equipment

Dionex ASE 200 Accelerated Solvent Extractor* equipped with 22 mL stainless steel extraction cells (P/N 048764)

Cellulose Filters (P/N 049458)

Collection Vials, 60 mL (P/N 048784)

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

Ottawa Sand Standard (P/N S23-3)

Laboratory grinder or blender (Fisher Scientific)

Tyler Sieve 0.5 mm (Fisher Scientific)

PTFE Syringe Filter 0.45 µm (Fisher Scientific)

**ASE 150 and 350 can be used for equivalent results*

Analysis

LC-MS

Reagents

ASE Prep DE (diatomaceous earth) (P/N 062819)

Solvents

Methanol Acetonitrile (All solvents are pesticide-grade or equivalent and available from Fisher Scientific.)

Extraction Conditions

Solvent: 50% methanol, 50% acetonitrile

Temperature: 80 °C

Pressure: 1500 psi

Heatup Time: 5 min

Static Time: 5 min

Static Cycles: 2

Flush Volume: 75%

Purge Time: 100 s

Total Extraction Time: 15 min

Volume of Solvent Use: 25–35 mL

Results

Results of extraction of ZON from wheat and corn using ASE.

Sample	Target Value (ng/g)	Average Recovery (ng/g) n=3	Percent Recovery	Percent RSD
Wheat	112	132	118	5.2
Corn	285	305	107	2.2

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



Introduction

The United Nations Environmental Program (UNEP) has been implemented in an effort to combat the release of selected persistent organic pollutants (POPs). POPs are found in environmental samples such as soils, sludges, solid and semi-solid 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. The following compounds are listed by UNEP to be POPs:

- Pesticides: Aldrin, Chlordane, DDT, Dieldrin, Endrin, Heptachlor, Mirex, and Toxaphene
- Industrial chemicals: Hexachlorobenzene, and PCB (polychlorinated biphenyl)
- Chemical by-products (Dioxins): Polychlorinated dibenzo-p-dioxins and polychlorinated dibenzofurans (PCDD and PCDF)

Persistent organic pollutants in environmental samples

Equipment

Dionex ASE 200 Accelerated Extractor* with Solvent Controller

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 049458)

Collection Vials 60 mL (P/N 048784) or 40 mL (P/N 048783)

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

*ASE 150 and 350 can be used for equivalent results

Solvents

Hexane

Dichloromethane

Acetone

Toluene

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

Reagents and Standards

Methylene chloride (Optima Grade, Fisher Scientific)

ASE Prep DE (diatomaceous earth) (P/N 062819)

Alumina (basic, Brockman activity I, Fisher Scientific)

PCB standards (ULTRA Scientific)

Analysis

GC-ECD

Extraction Conditions

Pesticides and PCBs (8081/8082)

Solvent: Hexane/acetone (1:1) (v/v)

Temperature: 100 °C

Pressure: 1500 psi*

Static Time: 5 min

Static Cycles: 1–2

Flush Vol.: 60%

Purge Time: 60–120s

Hexachlorobenzene (8270)

Solvent: Dichloromethane/acetone (1:1), (v/v)

Temperature: 100 °C

Pressure: 1500 psi

Static Time: 5 min

Static Cycles: 1–2

Flush Vol.: 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: 175–200 °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 ASE applications.

Results

PCB recovery from oyster tissue^a

PCB Congener	Average Recovery, n = 6 (as % of Soxhlet)	RSD (%)
PCB 28	90.0	7.8
PCB 52	86.9	4.0
PCB 101	83.3	1.5
PCB 153	84.5	3.5
PCB 138	76.9	3.0
PCB 180	87.0	4.3

^a Analyte concentration range: 50–150 µg/kg per component

Rapid Extraction and Determination of Arsenicals in Fish Tissue and Plant Material Using Accelerated Solvent Extraction (ASE)



Introduction

The toxicity of arsenic is species dependent. Inorganic arsenic species such as arsenite (As[III]) and arsenate (As[V]) have been classified as carcinogens. Methylated forms such as monomethylarsonic acid (MMA) and dimethylarsinic acid (DMA) have recently been labeled as cancer promoters. Arsenobetaine (AsB), arsenocholine (AsC), and arseno sugars have been found to be relatively nontoxic. Two major pathways for toxic arsenic exposure include drinking water and diet. Seafood (including fish and seaweed) accounts for the majority of ingested arsenic, most of which is nontoxic, however, fruits and vegetables grown in contaminated soils and sediments contribute another significant source. Due to the variable levels of toxicity associated with arsenic species in foods, total arsenic determination is not sufficient to assess potential harmful contamination. Determination of individual arsenic species is necessary. This has increased the need to improve separation and detection methods for organo-metallic speciation.

Arsenic species extracted from a variety food sample matrices

Equipment

Dionex ASE 200 Extractor*
equipped with 11-mL stainless steel extraction cells
Cellulose Glass-fiber Filters (P/N 049458 or 047017)
Collection Vials (40 or 60 mL) (P/N 048783 or 048784)
Analytical Balance (to read to the nearest 0.0001 g or better)
Solvent Evaporator
**ASE 150 and 350 can be used for equivalent results*

Reagents

Methanol (HPLC grade)
HPLC water
Ottawa Sand Standard (P/N S23-3)
ASE Prep DE (diatomaceous earth) (P/N 062819)

Solvents

Hexane
Dichloromethane
Acetone
Toluene
(All solvents are pesticide-grade or equivalent and available from Fisher Scientific.)

Analysis

LC-ICP-MS

Extraction conditions

Fish Tissue

Solvent: Methanol 100%
Temperature: 100 °C
Pressure: 1500 psi Cell
Heatup Time: 5 min
Static Time: 2 min
Flush Volume: 60%
Purge Time: 60 s
Cycles: 5
Total Time: 7 min
Total Solvent: <30 mL

Ribbon Kelp

Solvent: 30/70 (w/w) Methanol/ H₂O
Temperature: Ambient
Pressure: 1500 psi
Heatup Time: N/A
Static Time: 1 min
Flush Volume: 90%
Purge Time: 120 s
Cycles: 3
Total Time: 7 min
Total Solvent: <30 mL

Carrots

Solvent: Water
Temperature: 100 °C
Pressure: 1500 psi
Cell Heat-up Time: 5 min
Static Time: 1 min
Flush Volume: 100%
Purge Time: 90 s
Cycles: 3
Total Time: 18 min
Total Solvent: <30 mL

Results

Results of ASE extraction of fish tissue CRMs (n=6). Data obtained for AsB in two certified reference materials and a candidate reference material* extracted with ASE.

	Measured Value	Certified Value
DORM-2 (dogfish muscle)	16.3 ± 0.9 (±1s)	16.4 ± 1.1 (±95% C.I.)
BCR 627 (tuna fish)	3.69 ± 0.21 (±1s)	3.90 ± 0.22 (±95% C.I.)
BCR 710** (oyster tissue)	31.8 ± 1.1 (±1s)	32.7 ± 5.1 (±1s)

* Expressed as mg/kg As, unless otherwise stated.

** Concentration as species. The data shown for this material is based on the consensus mean of the final certification round after the removal of statistical outliers.

Determination of Perchlorate in Vegetation Samples Using Accelerated Solvent Extraction (ASE) and Ion Chromatography



Introduction

Perchlorate (ClO_4^-) is an environmental contaminant that has been found in drinking, ground, and surface waters in several states within the United States. Most of the contaminated sites have been traceable to sources near military installations or manufacturing sites where perchlorate salts are used to manufacture rocket propellant, munitions, or fireworks. The solubility, mobility, and persistence of perchlorate have resulted in the contamination of drinking water, soil, and vegetation in several areas.

Environmental contaminants in a variety of vegetable matrices

Equipment

Dionex ASE 200 or ASE 300* system

Collection vials, 60 mL (P/N 048784)

Collection bottles, 250 mL (P/N 056284)

Glass fiber filters (P/N 047017 for ASE 200, P/N 056781 for ASE 300)

OnGuard II Sample Pretreatment Cartridges

Ag (P/N 057089)

Ba (P/N 057093)

H (P/N 057085)

RP (P/N 057083)

ASE Prep DE (P/N 062819)

Analytical balance with 0.1 mg resolution

Dionex ICS-2500 chromatography system consisting of:

GP50 Gradient Pump with vacuum degas option

EG50 Eluent Generator with EluGen EGC II

NaOH cartridge (P/N 058908)

AS40 Autosampler

LC30 Chromatography Oven

CD25 Conductivity Detector with conductivity cell

*ASE 150 and 350 can be used for equivalent results

Analysis

IC

Reagents and Standards

Deionized water ($\text{DI H}_2\text{O}$), Type I reagent grade, 18 Ω -cm resistance or better

Sodium perchlorate, 98% ACS reagent grade or better (Fisher Scientific)

ACS reagent grade sodium salts (Mallinckrodt, Fisher)

Sodium Hydroxide (NaOH) 50% w/w (Fisher Scientific)

ASE Prep DE (diatomaceous earth) (P/N 062819)

Extraction Conditions

Extraction Solvent: Water

Pressure: 1,500 psi

Temperature: 80 °C

Equilibration Time: 5 min

Extraction Time: 5 min (static)

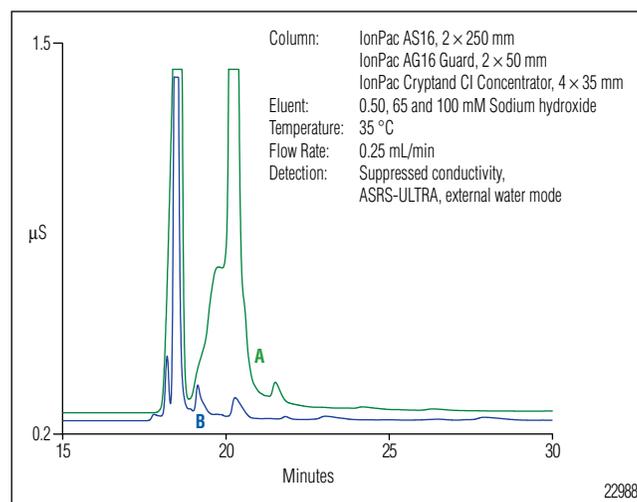
Solvent Flush: 30% (of cell volume)

Nitrogen Purge: 1, 20 s (after extraction)

Extraction Cycles: 3

Cell Sizes: 33 mL and 100 mL

Results



Alfalfa extracts obtained using (A) no in-line cleanup and (B) OnGuard resins combined with basic alumina in the ASE extraction cell.

Extraction and Cleanup of Acrylamide in Complex Matrices Using Accelerated Solvent Extraction(ASE) Followed by Liquid Chromatography-Tandem Mass Spectrometry (LC-MS/MS)



Introduction

Acrylamide is formed during the cooking process of certain plant-based foods which are rich in carbohydrates and low in protein. Specifically, it forms when asparagines reacts with sugars such as glucose at high temperatures. Acrylamide was detected in fried foods by the Swedish National Food Authority in 2002. Since then, many food laboratories have successfully performed determinations for this compound on a variety of different food matrices. Acrylamide is a known carcinogen in animals.

Contaminants in coffee and chocolate

Equipment

Dionex ASE 200* equipped with 33-mL stainless steel extraction cells
 Cellulose filters (P/N 049458)
 Collection vials, 60 mL (P/N 048784)
 Dionex SE 500 Solvent Evaporator (P/N 063221, 120 v) (P/N 063218, 240 v)
 Standard laboratory tissue homogenizer
 Standard laboratory centrifuge (rated to 10,000 rpm or greater)
 Centrifuge tubes (40–50 mL)

*ASE 150 and 350 can be used for equivalent results

Analysis

LC-MS/MS

Chemicals and Reagents

Acrylamide, purity 99% (Fisher Scientific)
d3-Acrylamide (2,3,3-*d3*-2-propenamide) (Cambridge, Isotope Laboratories USA)
 Florisil, 60–100 mesh (Fisher Scientific)
 Potassium hexacyanoferrate (II) trihydrate (Carrez I) (Fisher Scientific)
 Zinc sulfate heptahydrate (Carrez II) (Fisher Scientific)
 ASE Prep DE (diatomaceous earth) (P/N 062819)
 Termamyl® 120 L (Type L thermostable amyloglucosidase enzyme) (Novozymes, Denmark)
 Ethyl acetate (Fisher Scientific, HPLC Grade)
 Dichloromethane (Fisher Scientific, HPLC Grade)
 Methanol (Fisher Scientific, HPLC Grade)

Extraction Conditions

Solvent: Ethyl acetate (100%)
 Temperature: Ambient
 Pressure: 1500 psi
 Static Time: 3 min
 Static Cycles: 3
 Flush Volume: 100%
 Purge Time: 60 s

Results

Comparison of manual extraction versus ASE for quantification of acrylamide spiked samples in soluble chocolate powder (n = 6).

Spiking Levels	Manual Extraction		ASE Extraction	
	Recovery %	%RSD	Recovery %	%RSD
12.7 µg/kg	103.7	17.2	94.6	4.3
304.7 µg/kg	108.0	6.3	102.2	7.0
2504 µg/kg	104.3	5.3	101.5	2.4

ASE of roast ground coffee, soluble coffee, coffee surrogate, and cocoa.

Materials	Acrylamide Level (µg/kg)			
	Spiked at 150 µg/kg			
	Incurred ^a	Expected ^b	Measured ^a	CV%
R&G coffee	136	286	298	3.1
Soluble coffee powder	299	449	435	2.9
Coffee surrogate	632	782	782	1.0
Cocoa powder	192	342	343	1.1

Extraction of Contaminants, Pollutants, and Poisons from Animal Tissue Using Accelerated Solvent Extraction (ASE)



Introduction

ASE also be used also to extract organic materials from matrices such as milk, foodstuffs, plant material, plasma, serum, and tissue. This application brief details procedures for extracting the following contaminants from animal tissues:

- Dioxins/Furans
- Polybrominated Flame Retardants (PBDE)
- PCBs
- Pesticides
- PAHs
- Organotin

Environmental contaminants in fish and egg samples

Equipment

Dionex ASE 200 Accelerated Solvent Extractor* with ASE Solvent Controller
 Choose either 11 mL stainless steel extraction cells (P/N 049560) or 22 mL stainless steel extraction cells (P/N 049561) or 33 mL stainless steel extraction cells (P/N 049562)
 Cellulose filters (P/N 049458)
 Collection vials, 40 mL (P/N 048783) or 60 mL (P/N 048784)
 Dionex SE 500 Solvent Evaporation system (P/N 063221)
 Analytical balance (to read to nearest 0.0001 g or better)
 Tissue homogenizer (Buchi B-400 or equivalent)
 Freeze drier (for PCB extraction)
 Centrifuge (for organotin extraction)
 Mechanical shaker (for organotin extraction)
**ASE 150 and 350 can be used for equivalent results*

Analysis

GC
 GC-MS
 GC-ECD
 HPLC

Extraction Conditions

Pressure: 1500 psi
 Temperature: 175 °C
 Solvent: 100% Toluene
 Static Time: 10 min
 Static Cycles: 2
 Flush Volume: 60%
 Purge Time: 60 sec
 Static Time: 1 or 2 min**
 Flush Volume: 60%
 Purge Time: 60 sec
 Cycles: 2
 Total Time: 12 min
 Total Solvent: 20 mL

** Petroleum ether and hexane were found to be equivalent as extraction solvents for fat in meat.*

***When extracting more than 1 g of a high-fat sample, a 2 min static time may be beneficial.*

Results

PCDDs/PCDFs in fish tissue samples (ng/kg or ppt)using ASE.

Compound	Soxhlet	ASE	Certified
2,3,7,8-TCDD	7.6	7.6	6.6
1,2,3,4,8-PCDD	4.3	4.3	4.4
1,2,3,4,7,8-HCDD	1.4	1.4	1.9
2,3,4,7,8-TCDF	13.4	12.6	11.9
1,2,3,7,8-PCDF	5.4	5.1	5.0
1,2,3,4,7,8-HCDF	12.5	12.2	12.2
OCDD	12.4	6.4	6.3
Total TEQ	21.4	21.1	21.0

References

- H. Obana, K. Kikuchi, M. Okihashi, and S. Hori. Determination of Organophosphorus Pesticides in Foods Using an ASE System. *Analyst* **1997**, *122*, 217–220.
- J. W. McKiernan, J. T. Creed, C. A. Brockhoff, J. A. Caruso, and R. M. Lorenzana. A Comparison of Automated and Traditional Methods for the Extraction of Arsenicals from Fish. *Journal of Analytical Atomic Spectrometry* **1999**, *14*, 607–613.
- M. Okihashi, H. Obana, and S. Hori. Determination of *N*-methylcarbamate Pesticides in Foods Using an ASE with a Mini-column Cleanup. *The Analyst* **1998**, *123*, 711–714.
- M. Davis, J. Paper, H. Shafei, and J. Stanley. Comparison and Evaluation of a Method for the Extraction of Polychlorinated Dibenzo-*p*-dioxins and Polychlorinated Dibenzofurans from Chicken Eggs Using Accelerated Solvent Extraction. *Organohalogen Compounds* **1999**, *35*, 105–108.
- E. Francis, B. Joyce. Accelerated Solvent Extraction (ASE): A Low-volume Solvent Extraction Technique. *AOAC International* **1999**, July, 18–20.
- J. Toth, J. Ezzell, L. Covino, and B. Richter. The Use of ASE for Analysis of Trace Contaminants in Foods. Poster.
- Accelerated Solvent Extraction for Rapid Extraction of Pesticide Residues. *LaborPraxis* **2000**, March, 44–46.
- A. Franchi, L. Bensi, S. Viti, and F. Abballe. Use of Accelerated Solvent Extraction in Multiresidual Analysis of Pesticides in Vegetable Food. *Bollettino dei Chemical Igienisti, Parte Scientifica* **1998**, *49*, 119–125.
- J. Ezzell. Pressurized Fluid Extraction: Non-environmental Applications. *Encyclopedia of Separation Science* **2000**, 3993–3999.
- M. S. S. Curren and J. W. King. Accelerated Solvent Extraction of Triazine Pesticides from Beef Kidney Using Subcritical Water. Pittcon **2001**.
- B. E. Richter, N. Rothe, J. L. Ezzell, E. S. Francis, L. Covino, and J. Toth. Expanding the Capabilities of Automated Sample Preparation. *American Laboratory* **2001**, March, 56–60.
- B. Richter, F. Hoefler, and M. Linkerhaegner. Determining Organophosphorus Pesticides in Foods Using Accelerated Solvent Extraction with Large Sample Sizes. *LCGC* **2001**, *19*, 410–412.
- N. P. Vela, D. T. Heitkemper, and K. R. Stewart. Arsenic Extraction and Speciation in Carrots Using Accelerated Solvent Extraction, Liquid Chromatography, and Plasma Mass Spectrometry. *The Analyst* **2001**, *126*, 1011–1017.
- M. Papagianopoulos, B. Zimmermann, A. Mellenthin, M. Krappe, G. Maio, and R. Galensa. Online Coupling of Pressurized Liquid Extraction, Solid-phase Extraction, and HPLC for Automated Analysis of Proanthocyanidins in Malt. *Journal of Chromatography A* **2002**, *958*, 9–16.
- M. S. S. Curren and J. W. King. Ethanol-modified Subcritical Water Extraction Combined with Solid Phase Micro-extraction for Determining Atrazine in Beef kidney. *Journal of Agriculture Food Chemistry* **2001**, *49*, 2175–2180.
- K.-O. Honikel. Method for Determination of Nitrogen in Meat and Meat Products. Federal Centre for Meat Research **2002**.
- R. M. Alonso-Salces, E. Korta, A. Barranco, L. A. Berrueta, B. Gallo, and F. Vicente. Pressurized Liquid Extraction for the Determination of Polyphenols in Apple. *Journal of Chromatography A* **2001**, *933*, 37–43.
- M. Papagiannopoulos and A. Mellenthin. ASE in the Investigation of Polyphenols in the Brewing Process. *Special publication-Royal Society of Chemistry* **2001**, *269 Biological*, 199–201.
- J. C. Chuang, K. Hart, J. S. Chang, L. E. Boman, J. M. Van Emon, and A. W. Reed. Evaluation of Analytical Methods for Determining Pesticides in Baby Food and Adult Duplicate-diet Samples. *Analytica Chimica Acta* **2001**, *444*, 87–95.
- T. Pihlstrom, G. Isaac, M. Waldeback, B.-G. Osterdahl, and K. E. Markides. Pressurized Fluid Extraction (PLE) as an Alternative General Method for the Determination of Pesticide Residues in Rape Seed. *The Analyst* **2002**, *127*, 554–559.
- B. Kim, D.-H. Kim, E. Hwang, and H. Y. So. Comparison of Sample Extraction Methods for the Determination of Bis(*p*-chlorophenyl) Dichloroethylene (*p,p*-dde) in Rice Flour Using Isotope Dilution Mass Spectrometry. *Bulletin Korean Chem Society* **2002**, *23*, 935–936.
- J. Ezzell. Accelerated Solvent Extraction (ASE) of Mycotoxins Contaminants in Food Matrixes. Pittcon Presentation **2001**, 1099.
- J. F. Lawrence, B. Niedzwiedek, and P. M. Scott. Effects of Temperature and Solvent Composition on Extraction of Fumonisin B1 and B2 from Corn Products. *Journal of AOAC* **2000**, *83*, 604–611.
- R.A. Moreau, A. Nunez, and V. Singh. Diferuloylputrescine and *p*-coumaroyl-feruloylputrescine, Abundant Polyamine Conjugates in Lipid Extracts of Maize Kernels. *AOCS Press Lipids* **2001**, *36*, 839–844.
- L. Pallaroni and C. von Holst. Determination of Zearalenone from Wheat and Corn by Pressurized Liquid Extraction and Liquid Chromatography-electrospray Mass Spectrometry. *Journal Chromatography A* **2003**, *993*, 39–45.
- I. Pecorelli, R. Galarini, R. Bibi, Al. Floridi, E. Casciarri, and A. Floridi. Simultaneous Determination of 13 Quinolones from Feeds Using Accelerated Solvent Extraction and Liquid Chromatography. *Analytica Chimica Acta* **2003**, *483*, 81–89.
- O. Ezquerro, A. Garrido-Lopez, and M. T. Tena. Determination of 2,4,6-trichloroanisole and Guaiacol in Cork Stoppers by Pressurized Fluid Extraction and Gas Chromatography-mass Spectrometry. *Journal of Chromatography A* **2006**, *1102*, 18–24.
- C. Pirard, J.-F. Focant and E. De Pauw. An Improved Clean-up Strategy for Simultaneous Analysis of Polychlorinated Dibenzo-*p*-dioxins (PCDD), Polychlorinated Dibenzofurans (PCDF), and Polychlorinated Biphenyls (PCB) in Fatty Food Samples. *Anal Bioanal Chem* **2002**, *372*, 373–81.

- M. Biedermann, S. Biedermann-Brem, A. Noti, K. Grob, P. Egli, and M. Veritas. Two GC-MS Methods for the Analysis of Acrylamide in Foodstuffs. *Mitteilungen aus Lebensmitteluntersuchung und Hygiene*, 1–12.
- M. Schlummer, G. Wolz, L. Gruber, and A. Moller. Accelerated Dioxin Sample Preparation. Paper by the Fraunhofer Institute for Process Engineering and Packaging.
- J. Focant and H. Shirkhan. Power-prep Automated Extraction and Clean-up System for PCDDs, PCDFs, and PCBs. Power Point Presentation.
- A. Muller, E. Bjorklund, and C. von Holst. On-line Clean-up of Pressurized Liquid Extracts for the Determination of Polychlorinated Biphenyls in Feedingstuffs and Food Matrices Using Gas Chromatography-mass Spectrometry. *Journal of Chromatography A* 2001, 925, 197–205.
- B. Zimmermann, M. Papagiannopoulos, A. Mellenthin, M. Krappe, G. Maio, and R. Gelensa. Coupling of ASE, ASPEC and HPLC-automated Determination of Proanthocyanidines in Malt. *G.I.T. Laboratory Journal* 2002, 175–177.
- J. Haib, I. Hofer, and J-M. Renaud. Analysis of Multiple Pesticide Residues in Tobacco Using Pressurized Liquid Extraction, Automated Solid-phase Extraction Clean-up and Gas Chromatography-tandem Mass Spectrometry. *Journal of Chromatography A* 2002, 173–187.
- K. Kitamura, Y. Takazawa, S. Hashimoto, J-W. Choi, H. Ito, and M. Morita. Effective Extraction Method for Dioxin Analysis from Lipid-rich Biological Matrices Using a Combination of Pressurized Liquid Extraction and Dimethyl Sulfoxide/Acetonitrile/Hexane Partitioning. *Analytica Chimica Acta* 2004, 512, 27–37.
- M. Papagiannopoulos and A. Mellenthin. Automated Sample Preparation by Pressurized Liquid Extraction-Solid-phase Extraction for the Liquid Chromatographic-Mass Spectrometric Investigation of Polyphenols in the Brewing Process. *Journal of Chromatography A* 2002, 976, 345–348.
- T. Hori, K. Tabiishi, Y. Ashizuka, R. Nakagawa, T. Iida, T. Tsutsumi, and K. Sasaki. Comparison of Accelerated Solvent Extraction and Standard Shaking Extraction for Determination of Dioxins in Foods. *Organohalogen Compounds* 2004, 66, 537–541.
- M. Bonoli, E. Marconi, and M. F. Caboni. Free and Bound Phenolic Compounds in Barley Flours. Evaluation of the extraction capability of different solvent mixtures and pressurized liquid methods by micellar electrokinetic chromatography and spectrophotometry. *Journal of Chromatography A* 2004, 1057, 1–2.
- J. L. Gomez-Ariza, M. A. Caro de la Torre, I Giraldez, and E. Morales. Speciation Analysis of Selenium Compounds in Yeasts Using Pressurized Liquid Extraction and Liquid Chromatography-Microwave-Assisted Digestion-Hydride Generation-Atomic Fluorescence Spectrometry. *Analytica Chimica Acta* 2004, 524, 305–314.
- E. Korta, A. Bakkali, A. Berrueta, B. Gallo, and F. Vicente. Study of an Accelerated Solvent Extraction Procedure for the Determination of Acaricide Residues in Honey by High-performance Liquid Chromatography-Diode Array Detector. *Journal of Food Protection* 2002, 65, 161–166.
- A. Kot-Wasik and A. Wasik. Determination of Eobenedine in Animal Feeds by Liquid Chromatography Coupled with Diode-array Detection and Mass Spectrometry after Accelerated Solvent Extraction. *Analytica Chimica Acta* 2005, 543, 46–51.
- M. Marsin Sanagi, H. H. See, W. Aini Wan Ibrahim, A. and Abu Naim. Determination of Carotene, Tocopherols and Tocotrienols in Residue Oil from Palm Pressed Fiber Using Pressurized Liquid Extraction-Normal Phase Liquid Chromatography. *Analytica Chimica Acta* 2005, 538, 71–76.
- R. Carabias-Martinez, E. Rodriguez-Gonzalo, P. Revilla-Ruiz, and J. Hernandez-Mendez. Pressurized Liquid Extraction in the Analysis of Food and Biological Samples. *Journal of Chromatography A* 2005, 1089, 1–17.
- J. L. Urraca, E. Benito-Pena, C. Perez-Conde, M.C. Moreno-Bondi, and J.J. Pestka. Analysis of Zearalenone in Cereal and Swine Feed Samples Using Automated Flow-through Immunosensor. *Journal of Agricultural and Food Chemistry* 2005, 53, 3338–3344.
- H. Obana, M. Furuta, and Y. Tanaka. Analysis of 2-alkylcyclobutanones with Accelerated Solvent Extraction to Detect Irradiated Meat and Fish. *Journal of Agricultural and Food Chemistry* 2005, 53, 6603–6608.
- J. L. Gomez-Ariza, T. Garcia-Barrera, F. Lorenzo, and A. Gustavo Gonzalez. Optimization of a Pressurized Liquid Extraction Method for Haloanisoles in Cork Stoppers. *Analytica Chimica Acta* 2005, 540, 17–24.
- Y. X. Gong, S. P. Li, Y. T. Wang, P. Li, and F. Q. Yang. Simultaneous Determination of Anthraquinones in Rhubarb by Pressurized Liquid Extraction and Capillary Zone Electrophoresis. *Electrophoresis* 2005, 26, 1778–1782.
- O. Pardo, V. Yusa, N. Leon, and A. Pastor. Determination of Bisphenol Diglycidyl Ether Residues in Canned Foods by Pressurized Liquid Extraction and Liquid Chromatography-Tandem Mass Spectrometry. *Journal of Chromatography A* 2006, 1107, 70–78.
- C. Blasco, G. Font, and Y. Pico. Analysis of Pesticides in Fruits by Pressurized Liquid Extraction and Liquid Chromatography-Ion Trap-Triple Stage Mass Spectrometry. *Journal of Chromatography A* 2005, 1098, 37–43.
- C. Juan, L. Gonzalez, J. M. Soriano, J. C. Molto, and J. Manes. Accelerated Solvent Extraction of Ochratoxin A from Rice Samples. *Journal of Agricultural and Food Chemistry* 2005, 1–4.
- L. Gonzalez Osnaya, J. M. Soriano del Castillo, J. C. Molto Cortes, and J. Manes Vinuesa. Extraction and Analysis of Ochratoxin A in Bread Using Pressurized Liquid Extraction and Liquid Chromatography. *Journal of Chromatography A* 2006, 1113, 32–36.
- Collection of Official Methods under Article 35 of the German Federal Food Act. Analysis of foodstuffs. Modular multiple analytical method for the determination of pesticide residues in foodstuffs. L00.00-34 1999.
- G.-F. Pang, Y.-M. Liu, C.-L. Fan, J.-J. Zhang, Y.-Z. Cao, X.-M. Li, Z.-Y. Li, Y.-P. Wu, and T.-T. Guo. Simultaneous Determination of 405 Pesticide Residues in Grain by Accelerated Solvent Extraction then Gas Chromatography-Mass Spectrometry or Liquid Chromatography-Tandem Mass Spectrometry. *Anal Bioanal Chem* 2006, 384, 1366–1408.
- C. Soler, J. Manes, and Y. Pico. Determination of Carbosulfan and Its Metabolites in Oranges by Liquid Chromatography Ion-trap Triple-stage Mass Spectrometry. *Journal of Chromatography A* 2006, 1109, 228–241.
- L. Gonzalez, C. Juan, J.M. Soriano, J. C. Molto, and J. Manes. Occurrence and Daily Intake of Ochratoxin A of Organic and Non-organic Rice and Rice Products. *International Journal of Food Microbiology* 2006, 107, 223–227.
- S.-H. Nam, J. Cheng, W. R. Mindak, and S. G. Capar. Preliminary Results of Extraction, Separation and Quantitation of Arsenic Species in Food and Dietary Supplements by HPLC-ICP-MS. *Bull. Korean Chem. Soc.* 2006, 27,6, 903–908.
- S. Kleinhenz, W. Jira, and K-H. Schwind. Dioxin and Polychlorinated Biphenyl Analysis: Automation and Improvement of Clean-up Established by Example of Spices. *Mol. Nutr. Food Res* 2006, 50, 362–367.

- H. Zhao, L. Wang, Y. Qiu, Z. Zhou, X. Li, and W. Zhong. Simultaneous Determination of Three Residual Barbiturates in Pork Using Accelerated Solvent Extraction and Gas Chromatography-Mass Spectrometry. *Journal of Chromatography B* 2006, 840, 139–145.
- G. Morlock and W. Schwack. Determination of Isopropylthioxanthone (ITX) in Milk, Yogurt and Fat by HPTLC-FLD, HPTLC-ESI/MS and HPTLC-DART/MS. *Anal Bioanal Chem.* 2006, 385, 586–595.
- A. Cascone, S. Eerola, A. Ritieni, and A. Rizzo. Development of Analytical Procedures to Study Changes in the Composition of Meat Phospholipids Caused by Induced Oxidation. *Journal of Chromatography A* 2006, 1120, 211–220.
- V. Yusa, G. Quintas, O. Pardo, P. Marti, and A. Pastor. Determination of Acrylamide in Foods by Pressurized Fluid Extraction and Liquid Chromatography-Tandem Mass Spectrometry Used for a Survey of Spanish Cereal-based Foods. *Food Additives and Contaminants* 2006, 23,3, 237–244.
- J. K. Houessou, C. Delteil, and V. Camel. Investigation of Sample Treatment Steps for the Analysis of Polycyclic Aromatic Hydrocarbons in Ground Coffee. *Journal of Agricultural and Food Chemistry* 2006, 54, 7413–7421.
- M. M. Delgado-Zamarreno, M. Bustamante-Rangel, M. Garcia-Jimenez, A. Sanchez-Perez, and R. Carabias-Martinez. Off-line Coupling of Pressurized Liquid Extraction and LC/ED for the Determination of Retinyl Acetate and Tocopherols in Infant Formulas. *Talanta* 2006, 70, 1094–1099.
- R. Carabias-Martinez, E. Rodriguez-Gonzalo, and P. Revilla-Ruiz. Determination on Endocrine-disrupting Compounds in Cereals by pressurized Liquid Extraction and Liquid Chromatography-Mass Spectrometry. Study of background contamination. *Journal of Chromatography A* 2006, 1137, 207–215.
- S. Herranz, M. C. Moreno-Bondi, and M. D. Marazuela. Development of a New Sample Pretreatment Procedure Based on Pressurized Liquid Extraction for the Determination of Fluoroquinolone Residues in Table Eggs. *Journal of Chromatography A* 2007, 1140, 63–70.
- K. Wiberg, S. Sporning, P. Haglund, and E. Bjorklund. Selective Pressurized Liquid Extraction of Polychlorinated Dibenzo-p-dioxins, Dibenzofurans and Dioxin-like Polychlorinated Biphenyls from Food and Feed Samples. *Journal of Chromatography A* 2007, 1138, 55–64.
- G. Sagratini, J. Manes, and D. Giardina, Y. Pico. Determination of Isopropyl Thioxanthone (ITX) in Fruit Juices by Pressurized Liquid Extraction and Liquid Chromatography-Mass Spectrometry. *Journal of Agricultural and Food Chemistry* 2006, 54, 7947–7952.
- A.I. Smeds, P. C. Eklund, R. E. Sjöholm, S. M. Willfor, S. Nishibe, T. Deyama, and B. R. Holmbom. Quantification of a Broad Spectrum of Lignans in Cereals, Oilseeds and Nuts. *Journal of Agricultural and Food Chemistry* 2007, 55, 1337–1346.
- M. Bustamante-Rangel, M. M. Delgado-Zamarreno, A. Sanchez-Perez, and R. Carabias-Martinez. Determination of Tocopherols and Tocotrienols in Cereals by Pressurized Liquid Extraction-Liquid Chromatography-Mass Spectrometry. *Analytica Chimica Acta* 2007, 587, 216–221.
- C. Soler, B. Hamilton, A. Furey, K. J. James, J. Manes, and Y. Pico. Liquid Chromatography Quadrupole Time-of-flight Mass Spectrometry Analysis of Carbosulfan, Carbofuran, 3-hydroxycarbofuran, and Other Metabolites in Food. *Analytical Chemistry* 2007, 79,4, 1492–1501.
- O. Pardo, V. Yusa, N. Leon, and A. Pastor. Development of a Pressurized Liquid extraction and Liquid Chromatography with Electrospray Ionization-Tandem Mass Spectrometry Method for the Determination of Domoic Acid in Shellfish. *Journal of Chromatography A* 2007, 1154, 287–294.
- C. Soler, B. Hamilton, A. Furey, K. J. James, J. Manes, and Y. Pico. Liquid Chromatography Quadrupole Time-of-Flight mass Spectrometry Analysis of Carbosulfan, Carbofuran, 3-Hydroxycarbofuran, and other metabolites in food. *Anal. Chem* 2007, 79, 1492–1501.
- J. A. Mendiola, M. Herrero, A.O. Cifuentes, and E. Ibanez. Use of Compressed Fluids for Sample Preparation: Food Applications. *Journal of Chromatography* 2007, 1152, 234–246.
- K. Ridgway, S. P. D Lalljie, and R. M. Smith. Sample Preparation Techniques for the Determination of Trace Residues and Contaminants in Foods. *Journal of Chromatography A* 2007, 1153, 36–53.
- S. Vichi, C. Santini, N. Natali, C. Riponi, E. Lopez-Tamames, and S. Buxaderas. Volatile and Semi-volatile Components of Oak Wood Chips Analysed by Accelerated Solvent Extraction (ASE) Coupled to Gas Chromatography-Mass Spectrometry. *Food Chemistry* 2007, 102, 1260–1269.
- G. Font, A. Juan-Garcia, and Y. Pico. Pressurized Liquid Extraction Combined with Capillary Electrophoresis-Mass Spectrometry as an Improved Methodology for the Determination of Sulfonamide Residues in Meat. *Journal of Chromatography A* 2007, 1159, 233–241.
- J. Wilga, A. Kot-Wasik, and J. Namiesnik. Comparison of Extraction Techniques of Robenidine from Poultry Feed Samples. *Science Direct* 2007, 72, 812–819.
- T. Tanaka, T. Hori, T. Asada, K. Oikawa, and K. Kawata. Simple One-step Extraction and Cleanup by Pressurized Liquid Extraction for Gas Chromatographic-Mass Spectrometric Determination of Pesticides in Leafy Vegetables. *Journal of Chromatography A* 2007, 1175, 181–186.
- P. Arapitsas and C. Turner. Pressurized Solvent Extraction and Monolithic Column-HPLC/DAD Analysis of Anthocyanins in Red Cabbage. *Talanta* 2008, 74, 1218–1223.
- S. Chu and C. D. Metcalfe. Analysis of Paroxetine, Fluoxetine, and Norfluoxetine in Fish Tissues Using Pressurized Liquid Extraction, Mixed Mode Solid Phase Extraction Cleanup and Liquid Chromatography-Tandem MS. *Journal of Chromatography A* 2008, 1163, 112–118.
- M. R. Khan, R. Busquets, F. J. Santos, and L. Puignou. New Method for the Analysis of Heterocyclic Amines in Meat Extracts Using Pressurized Liquid Extraction and Liquid Chromatography-Tandem Mass Spectrometry. *Journal of Chromatography A* 2008, 1194, 155–160.
- S. Lacorte and M. Guillamon. Validation of a Pressurized Solvent Extraction and GC-NCI-MS Method for the Low Level Determination of 40 Polybrominated Diphenyl Ethers in Mothers' Milk. *Chemosphere* 2008, 73, 70–75.
- R. A. Jacques, C. Dariva, J. V. de Oliveira, and E. S. Caramao. Pressurized Liquid Extraction of Mate Tea Leaves. *Analytica Chimica Acta* 2008, 625, 70–76.
- M. Barriada-Pereira, M. J. Gonzalez-Castro, S. Muinatogui-Lorenzo, P.Lopez-Mahia, and D. Prada-Rodriguez. Comparison of Pressurized Liquid Extraction and Microwave Assisted Extraction for the Determination of Organochlorine Pesticides in Vegetables. *Talanta* 2007, 71.
- J. Sokol and E. Matisova. Determination of Tetracycline Antibiotics in Animal Tissues of Food-Producing Animals by High-performance Liquid Chromatography Using Solid-phase Extraction. *Journal of Chromatography A* 1994, 669, 75–80.
- J. Chen, W. Li, B. Yang, X. Guo, F. Sen-Chun Lee, and X. Wang. Determination of Four Major Saponins in the Seeds of *Aesculus Chinensis* Bunge Using Accelerated Solvent Extraction Followed by High-performance Liquid Chromatography and Electrospray-Time of Flight Mass Spectrometry. *Analytica Chimica Acta* 2007, 596, 273–280.

- A. Gil-Vergara, C. Blasco, and Y. Pico. Determination of 2-isopropyl thioxanthone and 2-ethylhexyl-4-dimethylaminobenzoate in Milk: Comparison of Gas and Liquid Chromatography with Mass Spectrometry. *Anal Bioanal Chem* 2007, 389, 605–617.
- B. Shao, H. Han, D. Li, Y. Ma, X. Tu, and Y. Wu. Analysis of Alkylphenol and Bisphenol A in Meat by Accelerated Solvent Extraction and Liquid Chromatography with Tandem Mass Spectrometry. *Food Chemistry* 2007, 105, 1236–1241.
- J. L. Martinez Vidal, J. L. Fernandez Moreno, F. J. Arrebola Liebanas, and A. G. Frenich. Application of Low-pressure Gas Chromatography/Tandem Mass Spectrometry to the Determination of Pesticide Residues in Tropical Fruits. *Journal of AOAC International* 2007, 90, 1146–1164.
- I. Garcia, M. Ignacio, A. Moureira, J. Cobias, and N. Carro. Assisted Solvent Extraction an Ion-trap Tandem Mass Spectrometry for the Determination of Polychlorinated Diphenyls in Mussels. Comparison with other extraction techniques. *Analytical and Bioanalytical Chemistry* 2008, 390, 729–737.
- M. Barrida-Pereira, I. Iglesias-Garcia, M. J. Gonzalez-Castro, S. Muniategui-Lorenzo, P. Lopez-Mahia, and D. Prada-Rodriguez. Pressurized Liquid Extraction and Microwave-assisted Extraction in the Determination of Organochlorine Pesticides in Fish Muscle Samples. *Journal of AOAC International* 2008, 91, 174–180.
- R. Gadzala-Kopciuch, A. Filipiak, and B. Buzewski. Isolation, Purification and Determination of 4-n-nonylphenol and 4-tert-octylphenol in Aqueous and Biological Samples. *Talanta* 2008, 74.
- J. Wilga, A. Kot-Wasik, A. Astel, and J. Namiesnik. Statistical Evaluation of the Effectiveness of Different Extraction Techniques for Determining Robenidine Level in Poultry Feed. *Critical Reviews in Analytical Chemistry* 2008, 38, 19–33.
- 1S.-K. Cho, A. M. Abd El-Aty, J.-H. Choi, Y.-M. Jeong, H.-C. Shin, B.-J. Chang, C. Lee, and J.-H. Shim. Effectiveness of Pressurized Liquid Extraction and Solvent Extraction for the Simultaneous Quantification of 14 Pesticide Residues in Green Tea Using GC. *Journal of Separation Science* 2008, 31, 1750–1760.
- F. J. Lara, A. M. Garcia-Campana, F. Ales-Barrero, J. M. Bosque-Sendra. In-line Solid-phase Extraction Preconcentration in Capillary Electrophoresis-Tandem Mass Spectrometry for the Multiresidue Detection of Quinolones in Meat by Pressurized Liquid Extraction. *Electrophoresis* 2008, 29, 2117–2125.
- J.-M. Lee, J.-W. Park, G.-C. Jang, and K.-J. Hwang. Comparative Study of Pesticide Multi-residue Extraction in Tobacco for Gas Chromatography-Triple Quadrupole Mass Spectrometry. *Journal of Chromatography* 2008, 1187, 25–33.
- P. Haglund, P. Korytar, C. Danielsson, J. Diaz, K. Wiberg, P. Leonards, A. T. Udo Brinkman, J. de Boer. GC x GC-ECD: A Promising Method for the Determination of Dioxins and Dioxin-like PCB's in Food and Feed. *Analytical and Bioanalytical Chemistry* 2008, 390, 1815–1827.
- M. Lores Pablo, C. Garcia-Jares, M. Garcia-Chao, T. Dagnac. Simultaneous Extraction and Cleanup Method Based on Pressurized Solvent Extraction for Multiresidue Analysis of Pesticides in Complex Feed Samples. *Journal of Agricultural and Food Chemistry* 2009, 57 (10), 3963–3973.
- S.-K. Cho, A. M. Abd El-Aty, H.-R. Jeon, J.-g.-H. Choi, H.-Ch. Shin, J.-H. Shim. Comparison of Different Extraction Methods for the Simultaneous Determination of Pesticide Residues in Kiwi Fruit Using Gas Chromatography-Mass Spectrometry. *Biomedical Chromatography* 2008, 22.
- W. Jira, K. Ziegenhals, K. Speer. Gas Chromatography-Mass Spectrometry (GC-MS) Method for the Determination of 16 European Priority Polycyclic Aromatic Hydrocarbons in Smoked Meat Products and Edible Oils. *Food Additives and Contaminants* 2008, 25(6), 704–713.
- C. Blasco, A. Di Corcia, and Y. Pico. Determination of Tetracyclines in Multi-specie Animal Tissues by Pressurized Liquid Extraction and Liquid Chromatography-Tandem Mass Spectrometry. *Food Chemistry* 2009, 116 (4), 1005–1012.
- T. Hori, D. Yasutake, K. Tobiishi, Y. Ashizuka, J. Kajiwara, R. Nakagawa, T. Lida, T. Tsutsumi, and K. Sasaki. Comparison of Accelerated Solvent Extraction and Alkaline Digestion-Hexane Shaking Extraction Determination of Dioxins in Animal-origin Food Sample. *Organohalogen Compounds* 2007, 69, 669/1–669/4.
- M. Lund, L. Duedahl-Olesen, and J. H. Christensen. Extraction of Polycyclic Aromatic Hydrocarbons from Smoked Fish Using Pressurized Liquid Extraction with Integrated Fat Removal. *Journal of Pharmaceutical and Biomedical Analysis* 2009, 49 (5), 1172–1178.
- D. Calderon-Preciado, C. Jimenez-Cartagena, G. Penuela, and J. M. Bayona. Development of an Analytical Procedure for the Determination of Emerging and Priority Organic Pollutants in Leafy Vegetables by Pressurized Solvent Extraction Followed by GC-MS Determination. *Talanta* 2009, 79 (1), 10–15.
- T. Otake, Y. Aoyagi, M. Matsuo, N. Itoh, and T. Yarita. Evaluation of Pressurized Liquid Extraction for the Analysis of Four Pesticides in Unpolished Rice. *Journal of Chromatography* 2008, 1203 (1), 81–87.
- H. Berrada, F. Borrull, G. Font, and R. M. Marce. Determination of Macrolide Antibiotics in Meat and Fish Using Pressurized Liquid Extraction and Liquid Chromatography-Mass Spectrometry. *Journal of Environmental Science and Health* 2008, 43 (5), 390–394.
- S. Kinani, S. Bouchonnet, J. Abjean, and C. Campargue. Determination of Polybromodiphenyl Ethers (PBDEs) in Milk Cream by Gas Chromatography-Mass Spectrometry. *Journal of Pharmaceutical and Biomedical Analysis* 2008, 48 (1), 231–235.
- A. R. Mauri-Aucejo, A. Campos-Candel, R. Marin-Saez, and M. Llobat-Estelles. Extraction of Cd, Pb, Ni, Cu, and Zn from Tobacco Samples by Pressurized Fluid Extraction and Determination by Atomic Spectrometry. *Atomic Spectroscopy* 2008, 29 (6), 217–223.
- E. Benito-Pena, J. L. Urraca and M. C. Moreno-Bondi. Quantitative Determination of Penicillin V and Amoxicillin in Feed Samples by Pressurized Liquid Extraction and Liquid Chromatography with Ultraviolet Detection. *Journal of Pharmaceutical and Biomedical Analysis* 2009, 49 (2), 7031–7085.
- O. Pardo, V. Yusa, N. Leon, and A. Pastor. Development of a Method for the Analysis of 7 Banned Azo-dyes in Chilli and Hot Chilli Food Samples by Pressurized Liquid Extraction and Liquid Chromatography with Electrospray Ionization-Tandem Mass Spectrometry. *Talanta* 2009, 78 (1), 178–186.
- K. Verheyden, H. Noppe, J. Vanden Bussche, K. Wille, K. Bekaert, L. De Boever, J. Van Acker, C. R. Janssen, H. F. De Brandbander, and L. Vanhaecke. Characterisation of Steroids in Wooden Crates of Veal Calves by Accelerated Solvent Extraction (ASE) and Ultra-high Performance Liquid Chromatography Coupled to Triple Quadrupole Mass Spectrometry (U-HPLC-QqQ-MS-MS). *Analytical and Bioanalytical Chemistry* 2010.

www.thermoscientific.com/dionex

©2012 Thermo Fisher Scientific Inc. All rights reserved. Arcolor is a trademark of Arco Íris Brasil Industria. Bio-beads is a registered trademark of Bio-Rad Laboratories Inc. Isolute is a registered trademark of International Sorbent Technologies Ltd. Equatherm is a registered trademark of Xaloy Inc. Termamyl is a registered trademark of Novozymes, Denmark. 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.



Thermo Scientific Dionex products are designed, developed, and manufactured under an ISO 9001 Quality System.

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

China +852 2428 3282
Denmark +45 36 36 90 90
France +33 1 39 30 01 10
Germany +49 6126 991 0
India +91 22 2764 2735

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

Thermo
SCIENTIFIC

Part of Thermo Fisher Scientific