

# Extraction and Cleanup of Acrylamide in Complex Matrices Using Accelerated Solvent Extraction 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 asparagine 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.

Accelerated solvent extraction is an excellent technique for extraction of acrylamide from various fried food products; until recently, however, extraction of this compound from matrices such as coffee and chocolate has proven difficult. Traditional extraction techniques are time consuming, and may cause bottlenecks in sample preparation. This application note describes a new accelerated solvent extraction method that combines the extraction of low-levels of acrylamide from coffee and chocolate with an in-cell, solid-phase cleanup step.

## Equipment

---

Thermo Scientific Dionex ASE 200 accelerated solvent extraction system with 33 mL stainless steel extraction cells (P/N 048763)

---

Thermo Scientific Dionex cellulose filters (P/N 049458)

---

Thermo Scientific Dionex collection vials 60 mL (P/N 048784)

---

Thermo Scientific 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)

---

\*Thermo Scientific Dionex ASE 150 and 350 accelerated solvent extraction systems can be used for equivalent results.

## Chemicals and Reagents

---

Acrylamide, purity 99% (Sigma-Aldrich®, St. Louis, USA)

---

*d3*-Acrylamide (2,3,3-*d3*-2-propenamamide)

---

Florisil®, 60–100 mesh (U.S. Silica Co., Frederick, USA)

---

Potassium hexacyanoferrate (II) trihydrate (Carrez I)

---

Zinc sulfate heptahydrate (Carrez II)

---

Thermo Scientific Dionex ASE Prep DE (P/N 062819)

---

TermamyI® 120 L (Type L thermostable amyloglucosidase enzyme) (Novozymes, Bagsvaerd, Denmark)

---

Fisher Scientific HPLC grade ethyl acetate

---

Fisher Scientific HPLC grade dichloromethane

---

Fisher Scientific HPLC grade methanol

---

## Food Samples

The coffee and chocolate samples were purchased from a local grocery store and stored at room temperature.

## Reagent Solutions

0.68 M potassium hexacyanoferrate (II) trihydrate (Carrez I) solution

Dissolve 28.722 g of  $K_4Fe(CN)_6 \cdot 3H_2O$  in 100 mL of water.

2 M zinc sulfate heptahydrate (Carrez II) solution

Dissolve 57.512 g of  $ZnSO_4 \cdot 7H_2O$  in 100 mL of water.

## Standard Solutions

Prepare aqueous stock solutions of acrylamide and *d3*-acrylamide at concentrations of 50 and 5 µg/mL, respectively.

To make 50 µg/mL acrylamide, add 5 mg to 100 mL water.

To make 5 µg/mL *d3*-acrylamide, add 0.5 mg to 100 mL water.

Dilute the acrylamide solutions in water to obtain the following matrix-equivalent levels: 0, 10, 50, 200, 500, and 2500 µg/kg. The matrix-equivalent concentration of *d3*-acrylamide should be 250 µg/kg.

## Sample Preparation

### Hydrolysis

Weigh 2.0 g of sample into a centrifuge tube and add 10 mL of water (heated to 60 °C) then add 50 µL of Termamyl. Place the tube in a water bath at 90°C for 45 min. Homogenize the mixture for 1 min. To precipitate the proteins, add 1 mL of the 0.68 M potassium hexacyanoferrate (II) trihydrate solution and 1 mL of the 2 M zinc sulfate heptahydrate solution to the centrifuge tube, swirling constantly for 1–2 min. Add 5 mL dichloromethane and swirl for an additional minute. Centrifuge for 15 min at 10,000 rpm.

### Preparing the Accelerated Solvent Extraction Cell

Prepare the 33-mL extraction cell by successively inserting: (1) a cellulose filter, (2) 6 g Florisil deactivated with 3% deionized water, (3) a second cellulose filter, and (4) at least 8 g Dionex ASE™ Prep DE so that there is approximately 0.5 cm of empty space at the top of the cell.

Transfer 6 mL of the supernatant from the centrifuge tube and drip onto the Dionex ASE Prep DE layer in the prepared extraction cell. Fill the extraction cell to the top with additional Dionex ASE Prep DE and cover with a third cellulose filter. Carefully place the extraction cell cap on the cell and tighten by hand.

## Extraction Conditions

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

\*Although the authors of this application note used ambient temperature for the extraction, this is not a typical extraction temperature for accelerated solvent extraction. Normal extraction temperatures range from 40–200 °C. If recoveries are lower than expected using ambient temperature, increasing temperature may improve results.

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

## Extraction

Place the prepared extraction cells onto the accelerated solvent extraction carousel. Enter the extraction conditions into the *Method Editor* screen, and save this method with the desired number. Begin the extraction by pushing *Start*. The method can also be set up as a Schedule in the *Schedule Editor* screen. Running the accelerated solvent extraction system under *Schedule* control enables the system to track any errors that may occur throughout an extraction run. This is especially helpful if the system is set up to run unattended overnight. Any problems are logged in the *Error Log* for the user to view the next morning.

When the extraction is complete, evaporate the extracts under vacuum (40 °C, 200 mbar) until only a few droplets remain. Evaporate the residual ethyl acetate under a gentle stream of nitrogen, or transfer the vials to the Dionex SE™ 500 Solvent Evaporation System (P/N 063221~120v; 063218~240v) and evaporate to dryness using standard conditions. Redissolve the extract with 500 µL water.

If necessary, the extract can be filtered through a 0.2-µm cellulose filter. Decant approximately 180 µL of the extract and mix with 90 µL of methanol prior to analysis by LC-MS/MS.

## Results and Discussion

The accelerated solvent extraction method automates the extraction and cleanup steps of extraction of acrylamide from cocoa and coffee. Compared to the manual method, accelerated solvent extraction greatly reduces the time and the amount of sample handling required (Table 1). The addition of Florisil to the extraction cell eliminates the need for an additional cleanup step of the extract. Figure 1 shows coffee extracted with various amounts of Florisil, followed by filtration through SPE cartridges. Method optimization determined that 6 g of Florisil were sufficient to obtain a clear extract, however some samples may require additional filtering before analysis.

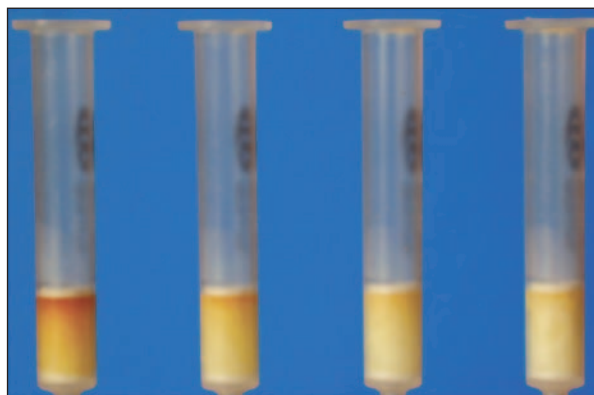


Figure 1. Residual co-extractables from coffee extracts, trapped on an Isolute® (Biotage AB, Uppsala, Sweden) Multimode SPE cartridge. The four extraction cells contained from 0 to 6 g of Florisil. From left to right: (1) no Florisil, (2) 2 g Florisil, (3) 4 g Florisil, and (4) 6 g Florisil.

Table 1. Comparison of the Sample Preparation Steps for the Manual Method and Accelerated Solvent Extraction Method for Extraction of Acrylamide from Coffee and Chocolate

Manual Method	Accelerated Solvent Extraction Method
1. Weigh 2 g of sample, add 100 $\mu$ L of <i>d</i> 3-acrylamide solution (5 $\mu$ g/mL) and 10 mL of water into a centrifuge tube. Homogenize for 1 min.	1. Weigh 2 g of sample, add 100 $\mu$ L of <i>d</i> 3-acrylamide solution (5 $\mu$ g/mL) and 10 mL of water into a centrifuge tube. Homogenize for 1 min.
2. Add 1 mL Carrez I, swirl, add 1 mL Carrez II, swirl. Add 5 mL dichloromethane, shake vigorously for 1 min.	2. Add 1 mL Carrez I, swirl, add 1 mL Carrez II, swirl. Add 5 mL dichloromethane, shake vigorously for 1 min.
3. Centrifuge at 3–5 $^{\circ}$ C, 10,000 rpm for 15 min.	3. Centrifuge at 3–5 $^{\circ}$ C, 10,000 rpm for 15 min.
4. Transfer 6 mL of the supernatant into a centrifuge tube containing 1.8 g of NaCl, swirl to dissolve.	4. Prepare the accelerated solvent extraction cell by successively adding 1 cellulose filter, 6 g of Florisil (deactivated with 3% water), a second cellulose filter and 8 g Dionex ASE Prep DE.
5. Add 13 mL ethyl acetate and shake vigorously for 1 min.	5. Transfer 6 mL of the extract on the Dionex ASE Prep DE, fill the rest of the column with Dionex ASE Prep DE, add a third cellulose filter and close the cell.
6. Centrifuge at 3–5 $^{\circ}$ C, 13800 g for 15 min.	6. Perform the accelerated solvent extraction step.
7. Transfer the organic phase into an amber vial containing 2 mL water. Shake vigorously for 1 min.	7. Evaporate the organic fraction under vacuum (40 $^{\circ}$ C, 200 mbar) to about 500 $\mu$ L and finish the evaporation with a gentle stream of N <sub>2</sub> .
8. Evaporate the organic phase with N <sub>2</sub> at 40 $^{\circ}$ C.	8. Re-dissolve the extract in 500 $\mu$ L of water.
<b>Repeat the ethyl acetate extraction (2<math>\times</math>), steps 5–7</b>	<b>Add 90 <math>\mu</math>L of extract. Proceed with LC-MS/MS analysis (60 <math>\mu</math>L injected).</b>
9. Condition the SPE cartridge with 3 mL methanol, then twice with 3 mL distilled water.	
10. Load the aqueous extract onto the cartridge, elute and rinse with 1 mL water. Collect both fractions.	
11. Reduce the extract volume to approximately 500 $\mu$ L (N <sub>2</sub> , 40 $^{\circ}$ C).	
12. Add 90 $\mu$ L of methanol to 180 $\mu$ L of extract. Proceed with LC-MS/MS analysis (60 $\mu$ L injected).	

Table 2 compares the results of manual extraction to accelerated solvent extraction of blank samples spiked with acrylamide standard.

Table 2. Comparison of Manual Extraction and Accelerated Solvent Extraction for Quantification of Acrylamide-Spiked Samples in Soluble Chocolate Powder (n = 6)

Spiking Levels	Manual Extraction		Accelerated Solvent 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

Table 3 shows the results of using accelerated solvent extraction for the extraction of acrylamide from various difficult matrices.

Table 3. Accelerated Solvent Extraction of Roast Ground Coffee, Soluble Coffee, Coffee Surrogate, and Cocoa

Materials	Acrylamide level (µg/kg)			
	Incurred <sup>a</sup>	Spiked at 150 µg/kg		
		Expected <sup>a</sup>	Measured <sup>b</sup>	CV%
R&G coffee	136	286	298	3.1
Soluble coffee powder	299	449	435	2.9
Coffee surrogate	632	762	782	1.0
Cocoa powder	192	342	343	1.1

<sup>a</sup>Mean of two independent determinations. <sup>b</sup>Mean incurred level + spike level.

## Conclusions

Accelerated solvent extraction has consistently proven to be an excellent alternative to the traditional labor-intensive extraction methods used for determination of acrylamide in food. Accelerated solvent extraction allows extraction and cleanup to be performed simultaneously, eliminating the need for a post-extraction cleanup step. Automation of accelerated solvent extraction allows for unattended extraction, and can be set up to run over-night to provide the user with filtered extracts that are ready for analysis in the morning. The advantages of speed and decreased sample handling as compared to manual extraction techniques are clear. Acrylamide recoveries obtained with the accelerated solvent extraction method were comparable to traditional methods, proving that accelerated solvent extraction is an effective tool for the extraction of polar compounds from complex matrices.

## References

- Wenzl, T., de la Calle, B. Anklam, E.; Analytical Methods for the Determination of Acrylamide in Food Products: A Review. *Food Addit. Contam.* 2003, 20, 885–902.
- Delatour, T., Perisset, A., Goldmann, T., Riediker, S., Stadler, R. H.; Improved Sample Preparation to Determine Acrylamide in Difficult Matrixes Such as Chocolate Powder, Cocoa, and Coffee by Liquid Chromatography Tandem Mass Spectroscopy. *J. Agric. Food Chem.* 2004, 52, 4625–4631.
- Cavalli, S., Maurer R., Hoefler F.; Fast Determination of Acrylamide in Food Samples Using Accelerated Solvent Extraction Followed by Ion Chromatography with UV or MS Detection. *LC-GC International* 2003, 16, Applications Book, 9–11.
- Hoefler F., Maurer R., Cavalli S.; Schnelle Analyse von Acrylamid in Lebensmitteln mit ASE und LC/MS, *GIT Labor-Fachzeitschrift*, 2002, 9, 968–970.

## Acknowledgment

Thanks to the authors at the Nestle Research Center Ltd. and the Nestle Product Technology Center in Switzerland for allowing us to use their data for this application note.

## Suppliers

Fisher Scientific, 2000 Park Lane, Pittsburgh, PA 15275-1126 USA, Tel: 800-766-7000, www.fishersci.com.

Novozymes A/S, Krogshoejvej 36, 2880 Bagsvaerd, Denmark, Tel.: +45 88 24 99 99 Fax: +45 88 24 99 98, www.novozymes.com.

[www.thermoscientific.com/dionex](http://www.thermoscientific.com/dionex)

©2012 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. Sigma-Aldrich is a registered trademark of Sigma-Aldrich Co. LLC. Florisil is a trademark of U.S. Silica Co. Termamyl is a trademark of Novozymes. Isolute is a registered trademark of Biotage AB. 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**  
 S C I E N T I F I C

Part of Thermo Fisher Scientific