Using Accelerated Solvent Extraction in Alternative Fuel Research

Richard Carlson, David Knowles, and Bruce Richter Thermo Fisher Scientific, Salt Lake City, UT, USA

Introduction

As global interest in alternate fuel sources increases, many laboratories are researching effective ways to test, develop, and produce fuel from renewable energy sources. One of these is the production of alcohol from biomass, called bioalochol. Alcohol produced from biomass has several benefits over fossil fuels because a) it is produced from crops, a renewable energy source, and b) when burned, it causes little environmental pollution. As bioalcohol is produced from biomass via sugar fermentation, it is necessary to research which types of plants produce the best yields of usable sugars.

Thermo Scientific[™] Dionex[™] ASE[™] Accelerated Solvent Extractors provide a fast and efficient way to extract various biomass samples for sugar analysis to determine sample viability for alcohol production. The Dionex ASE 150 and 350 Accelerated Solvent Extractors with the pH-hardened pathways provide additional benefits to the method by allowing pre-hydrolyzed biomass samples or those with acidified solvents to be extracted. This allows samples to be hydrolyzed in the extraction cell.¹ This application note describes methods for the extraction of sugars from biomass samples using acidified solvents (in-cell hydrolysis) and extracting pre-hydrolyzed samples.²

Equipment

- Dionex ASE 350 Accelerated Solvent Extractor (P/N 057794)
- Thermo Scientific[™] Dionium[™] Extraction Cells 66 mL (P/N 068102)
- Collection Bottles 250 mL (P/N 056284)
- Glass Fiber Filters (P/N 056781)
- Ottawa Sand (Fisher Scientific)
- Standard Laboratory Sample Evaporation System
- Standard Laboratory Grinder or Mill

Reagents

- HPLC Grade Water (Fisher Scientific)
- H₂SO₄ 8 M (Sigma-Aldrich)

Sample Prep

Grind raw sample material using a standard laboratory grinder or mill.

Sample Prep for Acidified Solvent Extraction (In-Cell Hydrolysis):

Weigh approximately 1.0–10.0 g of ground sample into a tared 66 mL Dionium extraction cell containing a glass fiber filter. It is important not to compact the sample into the cell as a tightly compacted sample will not hydrolyze well. Ottawa Sand can be added to the cell to take up any dead volume. This will help reduce the amount of solvent used per extraction. Weigh the appropriate number of 250 mL collection vials and place them onto the Dionex ASE Accelerated Solvent Extractor.

Sample Prep for Pre-Hydrolyzed Samples

Weigh approximately 2 g of sample and hydrolyze using 0.5% H₂SO₄ at 200 °C. Transfer the acidified sample mixture to a 66 mL Dionium extraction cell containing a glass fiber filter. Weigh the appropriate number of 250 mL collection bottles and place them onto the Dionex ASE Accelerated Solvent Extractor.

Conditions

Accelerated Solvent Extraction Conditions Using Acidified Solvent		
Pressure:	1500 psi*	
Temperature:	140–150 °Cª	
Solvent:	0.1 M $\rm H_2SO_4$ solution with HPLC water	
Static Time:	5 min	
Static Cycles:	3	
Flush:	10%	
Purge:	120 s	

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

^aTests concluded that an extraction temperature of no higher than 150 °C was optimal for the corn stover samples. Other biomass sample matrices may vary and could require an extraction temperature lower than 150 °C.



Accelerated Solvent Extraction Conditions for Pre-Hydrolyzed Samples²

Pressure:	1500 psi
Temperature:	100 °C
Solvent:	HPLC water
Static Time:	7 min
Static Cycles:	3
Flush:	50%
Purge:	120 s

Extraction

Place the extraction cells containing the samples onto the Dionex ASE Accelerated Solvent Extractor cell tray. Create an extraction method based on the desired method listed above. Once the extractions are completed, prepare the extracts according to the analytical method needed. For gravimetric determination, evaporate the solvent to dryness using a standard laboratory sample evaporation system and weigh the collection vials to determine % residue.

(Weight of evaporated vial – tared weight of vial)/ sample weight \times 100.

If sample is to be analyzed chromatographically for sugar content, then the appropriate volume adjustment should be made before injection into the chromatographic system.

Results and Discussion

Acidified Solvent Extraction (In-Cell Hydrolysis)

Corn stover^b samples were extracted for the quantifiable determination of sugars used for ethanol production. Two different extraction temperatures were examined for determining the best method (Table 1). Results showed that the samples extracted at 190 °C produced more total extractables. However, HPLC analysis showed that this higher temperature caused a breakdown of the sugars. The higher temperature was too aggressive and caused an extraction of unwanted coextractables resulting in the increased final weight. Analysis of the extracts at 150 °C showed no breakdown of sugars and no increase in unwanted coextractables.

^bCorn stover consists of the leaves and stalks of maize plants left in a field after harvest and includes the residue stalk, the leaf, husk, and cob.

Table 1. Corn stover extracted with acidified solvent.

Extraction Temperature	% Total Extractables
150 °C	48.7
190 °C	87.3

Pre-Hydrolyzed Samples

Two different concentrations of acidic solvents were examined for hydrolysis of switchgrass samples to determine the optimal concentration for recovery of glucose (Table 2). The HPLC analysis showed that for samples pretreated with $0.2\% H_2SO_4$, xylose was the predominant sugar present (Figure 1). Samples pretreated with $0.5\% H_2SO_4$ had a significant increase in glucose levels (Figure 2). Table 3 lists the HPLC conditions.

Table 2. Pre-hydrolyzed switchgrass samples extracted with HPLC-grade water.

Acidic Solvent Concentration	% Total Extractables
0.2% H ₂ SO ₄	76.76
0.5% H ₂ SO ₄	72.39



Figure 1. HPLC characterization of switchgrass pretreated with 0.2% H₂SO₄.



Figure 2. HPLC characterization of switchgrass pretreated with 0.5% $\rm H_{9}SO_{4}.$

Table 3. HPLC conditions for Figures 1 and 2.

Column Type	Aminex [®] HPX-87H
Flow Rate	0.6 mL/min
Temperature	65 °C
Mobile Phase	0.0005 N H ₂ SO ₄
Detector	RI
Run Time	26 min
Post Run Time	10 min

Conclusions

The new pH-hardened pathways of the Dionex ASE 350 and 150 Accelerated Solvent Extractors allow the extraction of samples that have been pretreated with acid, and also allows in-cell hydrolysis with acidified solvents. This new feature, along with the rapid extraction time and minimal solvent usage, provides users with a way to rapidly quantify biomass samples for biofuel research.

Acknowledgements

The accelerated solvent extraction method for extraction of corn stover samples was developed by NREL in Golden, Colorado.

The authors would like to thank Dr. Doug Raynie from South Dakota State University (SDSU) Dept. of Chemistry and Biochemistry for the data on switchgrass samples.

List of Manufacturers

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

Sigma-Aldrich Chemical Company, 3050 Spruce St., St. Louis, MO 63103 USA, Tel: 800-325-3010, www. sigmaaldrich.com.

References

- 1. Accelerated Solvent Extraction ASE Systems Brochure, LPN 2012-02, 2008. Dionex Corporation (now part of Thermo Scientific), Sunnyvale, CA.
- 2. Minnesota Chromatography Forum, Poster Presentation (SDSU) May 7, 2008.

©2013 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. Aminex is a registered trademark of Bio-Rad Laboratories, Inc. 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. and its subsidiaries. This information is presented as in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

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

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



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

