

Determination of Oil Content in Biodiesel Feedstock by Accelerated Solvent Extraction

Prapisala Thepsithar, Yi Zhang, Zhenpeng Zhang, Rong Yan
Institute of Environmental Science and Engineering, Nanyang Technological University, Singapore

Key Words

Accelerated Solvent Extraction, ASE, Jatropha Oil, Jatropha seeds, British standard (BS EN ISO 659:1999), biodiesel feedstocks

Goal

To produce a method that reduces the time required to extract oil from Jatropha seeds when compared to the Soxhlet technique.

Introduction

The determination of oil content in biodiesel feedstock can be performed by several methods, including mechanical press, solvent extraction, and Nuclear Magnetic Resonance (NMR). For feedstock quality control in terms of oil content, it is important that the applied method is universally accepted so that the results can be compared with those from different sources. Although the European standard has specified two methods for the determination of oil content in oilseeds, conventional Soxhlet extraction and NMR imaging, these methods have disadvantages, including excessive time consumption, labor-intensive input, need for more highly skilled labor, a requirement for significant number of samples, high cost, and being harmful to the environment.

Table 1. British Standard ISO 659:1999

Sample size	10 g ground seeds
Solvent	n-hexane
Total volume solvent	~150 mL
Extraction 1	4 h, heat and grind
Extraction 2	2 h, heat and grind
Extraction 3	2 h
Total extraction time	10 h



The accelerated solvent extraction method developed by Dionex, now part of Thermo Fisher Scientific, has great potential in overcoming these constraints. Research by the Institute of Environmental Science and Engineering in Singapore evaluated this potential. The results demonstrate that the accelerated solvent extraction method can be used efficiently to extract oil from Jatropha seeds with comparable results, in terms of accuracy and repeatability, to the results obtained from the European standard. The oil extraction using the accelerated solvent extraction method requires less than 1 h compared to the 10 h consumed by the Soxhlet extraction method (Table 1). The accelerated solvent extraction method can also be used effectively with smaller sample sizes.

Equipment

- Thermo Scientific™ Dionex™ ASE™ Accelerated Solvent Extractor equipped with 10 or 34 mL cells
- Thermo Scientific Dionex vials for extraction collection (60 mL)
- Analytical balance
- Cellulose filters for 1, 5, 10 or 22 mL cells, Pkg of 100, Dionex ASE 150/350 (P/N 068093)
- Cellulose filters for 34, 66 or 100 mL Cells, Pkg of 100, Dionex ASE 150/350 (P/N 056780)
- Diatomaceous Earth Dispersant for Dionex ASE, 1 kg Bottle (P/N 062819)

Solvent

Analytical grade n-Hexane (or light petroleum ether especially composed of hydrocarbons with 6 carbon atoms).

Accelerated Solvent Extraction Conditions

Temperature	80 °C
Time	10 min
Static Time	10 min
Flush Volume	60%
Purge Time	60 s
Solvent	n-Hexane
Static Cycle	3

Sample Preparation

The moisture content of all samples should be less than 10% (w/w). If greater than 10%, the moisture content should be reduced by drying the sample in an oven with a constant temperature of <80 °C. It is also important that the particle size of all samples be in the range of 2–5 mm. The samples should be ground using a mechanical mill without heating or changing moisture and oil content. It is essential that the oil extractions be carried out within 30 min of grinding, especially if the free fatty acid content of the extracted oil is to be determined.

Procedure

Grind the sample into a particle size, specified above, using a mechanical grinder. Place a cellulose filter at the bottom of the extraction cell (10 mL or 34 mL cell). Fill the cell with 5–10 g of the sample. Put Thermo Scientific™ Dionex™ ASE™ Prep DE on top of the sample and leave the head space of the cell around 1/5 of the cell length. Place another cellulose disk on top of the Dionex ASE Prep DE and assemble the extraction cell. Place the extraction cell into the Dionex ASE system. Set the method conditions on the Dionex ASE system and start the extraction. Upon completion of the extraction, transfer the extracts to a round-bottom flask and assemble the evaporation apparatus. Weigh the extracted oil and calculate the percentage of oil content using the following equation:

$$\% \text{ Oil Content} = \frac{\text{Weight of Extracted Oil (g)}}{\text{Weight of Sample (g)}} \times 100$$

Results and Discussions

Accuracy and Reproducibility

The extraction in Table 1 was conducted on Jatropha seeds using both the British standard (BS EN ISO 659:1999) and the accelerated solvent extraction method. The percentage of oil in the Jatropha seeds was determined using a conventional Soxhlet extraction apparatus according to BS EN ISO 659:1999 and the result was $47.2 \pm 0.7\%$ weight oil / weight kernel ($n = 4$) with a total extraction time of 10 h. For accelerated solvent extraction method, the percentage of oil extracted was approximately $48.5 \pm 1.3\%$ weight oil / weight kernel ($n = 3$) when extracted at 80 °C and around $50.6 \pm 0.5\%$ weight oil / weight kernel ($n = 3$) when extracted at 100 °C. It should be noted that the sizes of sample portion for BS EN ISO 659:1999 and ASE were 10 g and 5 g, respectively. It can be concluded that the results obtained from the accelerated solvent extraction method are consistent with those obtained from the standard method (Figure 1).

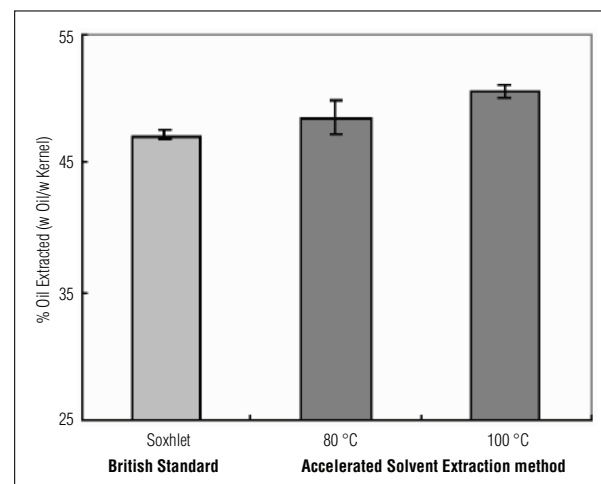


Figure 1. Percentage of oil extracted using the British standard and Accelerated Solvent Extraction method.

In addition, the repeatability of the test results was very good. The accelerated solvent extraction method can also be effectively used when the sample portion is only 3 g. The percentage of extracted oil is around $49.5 \pm 2.6\%$ weight oil / weight kernel ($n = 3$). Although the standard deviation is slightly higher compared to the results obtained from BS EN ISO 659:1999 and accelerated solvent extraction when using 5 g of sample, the results from 3 g are consistent with the 5 g results (Figure 2).

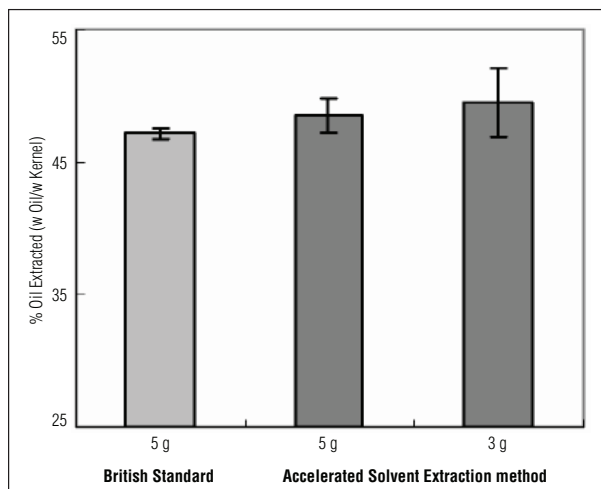


Figure 2. Percentage of oil extracted using different amounts of sample.

Conclusions

Accelerated solvent extraction method is a very effective method for oil extraction in the determination of oil content in *Jatropha* seeds. The test results are very consistent with the results obtained from the British standard. The results obtained from the accelerated solvent extraction method have high accuracy and precision. In addition, the accelerated solvent extraction method can be applied in the determination of oil content when a limited amount of sample is available. It should also be emphasized that oil extraction using the accelerated solvent extraction method requires less than 1 h compared to the 10 h consumed by Soxhlet extraction.

www.thermoscientific.com/samplepreparation

©2014 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific 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 Fisher Scientific,
Sunnyvale, CA USA is
ISO 9001:2008 Certified.

Africa +43 1 333 50 34 0
Australia +61 3 9757 4300
Austria +43 810 282 206
Belgium +32 53 73 42 41
Brazil +55 11 3731 5140
Canada +1 800 530 8447
China 800 810 5118 (free call domestic)
 400 650 5118

Denmark +45 70 23 62 60
Europe-Other +43 1 333 50 34 0
Finland +358 9 3291 0200
France +33 1 60 92 48 00
Germany +49 6103 408 1014
India +91 22 6742 9494
Italy +39 02 950 591

Japan +81 6 6885 1213
Korea +82 2 3420 8600
Latin America +1 561 688 8700
Middle East +43 1 333 50 34 0
Netherlands +31 76 579 55 55
New Zealand +64 9 980 6700
Norway +46 8 556 468 00

Russia/CIS +43 1 333 50 34 0
Singapore +65 6289 1190
Sweden +46 8 556 468 00
Switzerland +41 61 716 77 00
Taiwan +886 2 8751 6655
UK/Ireland +44 1442 233555
USA +1 800 532 4752

CAN-71162-EN-0614S

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand