Application Note: 50723

# Rapid Analysis of Trans Fat Content Using a Fourier Transform Infrared Spectrometer

Joni Miller, Mike Garry, Mike Bradley, Thermo Fisher Scientific, Madison, WI, USA

# **Key Words**

- ATR
- FT-IR
- GasChromatography
- Infrared
- Oil Analysis
- Trans Fat

## Introduction

Evidence continues to mount on the adverse affects of human consumption of *trans* fatty acids. A report by the US Food and Drug Administration (FDA) concluded that consumption of *trans* fat contributes to increased LDL ("bad") cholesterol levels, which increase the risk of coronary heart disease. The FDA estimates that publicizing the *trans* fat content of processed food products would prevent up to 1,200 cases of heart disease and up to 500 deaths each year.<sup>1</sup>

*Trans* fats are found in many common food products in the form of hydrogenated oils. The use of hydrogenated oils is widespread because they improve the consistency and shelf-life of many processed food products, especially baked goods.

Nutrition Facts Serving Size 1 cup (228g) Servings Per Container 2						
Amount Per Serving Calories 260			Calories from Fat 120			
				%Da	ily Value*	
Total Fat 13g			20%			
Saturate				25%		
Trans Fat 2g						
Cholester			10%			
Sodium 660mg					28%	
Total Carbohydrate 31 109						
Dietary Fiber 0g					0%	
Sugars 5g						
Protein 5g						
Vitamin A	A 4%	•	Vi	tamir	n C 2%	
Calcium 15% •			Iron 4%			
Percent Daily Values are based on a 2,000 calorie diet. Your Daily Values me be higher or lower depending on your calorie needs:						
	Calories		2,000	1	2,500	
Total Fat	Less tha		65g		80g	
Sat Fat	Less tha		20g		25g	
Cholesterol	Less tha		300g		300g	
Sodium	Less tha	ın	2,400	mg	2,400 mg	
Total Carbohydrate					375g	
Dietary Fiber 2					30g	
Calories per gr Fat 9 *		ohydrate	4		Protein 4	

Figure 1: Revised label, from the FDA, which includes information on *trans* fat

Countries around the world, led by Denmark, the United States, and Canada, have begun to enact strict guidelines requiring food manufacturers to list trans fatty acid content on the packaging, as shown in Figure 1. The importance of this labeling task places strict requirements on the analysis - speed, reliability, and robustness are essential. Certified methods by the American Oil Chemist Society (AOCS) or the Association of Official Analytical Chemists (AOAC)) entail infrared (IR) spectroscopic or gas chromatographic (GC) analysis.

Most present analyses require extraction to isolate the fats. GC methods are commonly used, due to the availability of GC equipment and its sensitivity. GC methods (AOCS Ce 1f-96 and AOAC 996.06) have good sensitivity down to 0.5%, but the triglycerides must be broken down to release the fatty acids, which must then be converted to fatty acid methyl esters (FAME) for injection. In addition, overlap of *cis* and *trans* peaks causes fat values to be underestimated, so fractionation of the *cis* and *trans* isomers using TLC or HPLC must be done, adding significant time and effort.

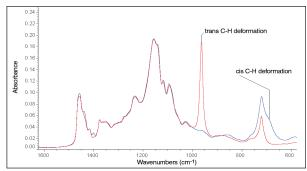


Figure 2

The infrared features of the *cis* and *trans* molecular configurations occur in different spectral regions, so no interference occurs. This can be seen in Figure 2, which shows overlaid infrared spectra of trielaidin (100% trans) and triolein (100% cis). However, the traditional infrared analysis method, Cd 14-95, also uses FAME derivatives, and requires carbon disulfide (CS<sub>2</sub>) as a dilution solvent. This is cumbersome, limited in sensitivity, and the use of  $CS_2$  is objectionable due to its odor. Although the method specifies applicability down to 0.5%, the *trans* peak is a shoulder on a large peak in the fat spectrum, making accurate measurement difficult, so the method may only work to 5%.

A newer infrared method was developed based on a heated horizontal attenuated total reflectance (ATR) accessory and is specified in methods AOCS Cd14d-99 and AOAC 2000.10. The ATR method is easy, rapid, and reproducible. Direct analysis of the *trans* isomer in the fat sample can be completed without weighing or preparation of FAME derivatives, and no smelly solvents are needed. The small sampling area of modern single bounce ATR accessories only require sample volumes of 50 µL or less, allowing for a reasonable sample size when extracting fat from food.

One key requirement for the ATR method is the use of *trans*-free reference fat for background correction, to eliminate the sloping baseline and shoulder peak measurement seen in the Cd 14-95 method. The method states it can be applied down to *trans* levels of 1%. However, the reference must approximate the fatty acid profile of the fat sample or it can adversely affect the methods accuracy, particularly near the limit of quantification. This means the user must select a *trans*-free fat that is similar to the fat sample being analyzed.



#### **Materials and Methods**

Data was collected on a Thermo Scientific Nicolet™ 380 FT-IR spectrometer. Samples were run on a heated diamond ATR accessory (100 scans, 4 cm<sup>-1</sup> resolution, at 65 °C). Trielaidin (100% trans) and triolein (0% trans) standards supplied by Nu-Chek Prep were used to prepare the calibration standards per the official method AOCS Cd 14d-99. The area of the peak at 966 cm<sup>-1</sup> due to the outof-plane C-H deformation about trans double bonds, was used to quantify the trans content, using a linear regression generated in the TQ Analyst™ software. Standards from 1 to 50% trans are shown overlaid in Figure 3 along with a 100% trielaidin, and the resultant calibration is shown in Figure 4.

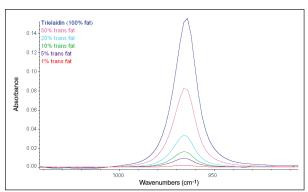


Figure 3

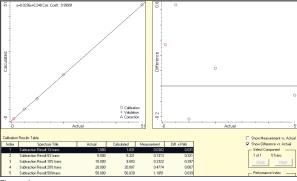


Figure 4

### Conclusion

Food manufacturers can use the infrared ATR technique for rapid determination of the trans fat content of the fats and oils used in the manufacture of food products. This analysis is instrumental in helping them comply with the food labeling requirements set by various countries throughout the world to help promote healthy eating habits. To help manufacturers meet these requirements, instrumentation companies have developed systems that can be used to quantify the trans fat content of edible fats and oils using an FT-IR spectrometer.

The high quality of the data, and the resulting calibration curve, show the excellent dependability of the FT-IR method. The speed and sensitivity of modern FT-IR spectrometers coupled with the ease of sampling of horizontal ATR accessories has moved the technique back into the mainstream. For the analysis of lipid material, FT-IR provides unique information difficult to obtain using other techniques. The implementation of diamond as an ATR crystal material has further enhanced the technique by providing a rugged sampling interface.

#### Resources

1. FDA Web site: http://www.fda.gov/oc/initiatives/transfat/

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

Austria +43 1 333 50340 Belgium Canada China India Italy +39 02 950 591 Spain

Australia

**Japan** +81 45 453 9100 **Latin America** +1 608 276 5659

**Netherlands** +31 76 587 98 88 South Africa

+27 11 570 1840 4 91 657 4930

Sweden/Norway/ **Finland** +46 8 556 468 00

**Switzerland** +41 61 487<u>84 00</u> UK +44 1442 233555

+1 800 532 4752

www.thermo.com



no Electron Scientific ments LLC, Madison, WI

are the property of Thermo Fisher Scientific Inc. and its

pricing are subject to change. Not all products are available

AN50723\_E 05/07M

