

Measurement of Samples by Transmission Spectroscopy with the Thermo Scientific Antaris FT-NIR Analyzer

William J. McCarthy, Thermo Fisher Scientific, Madison, WI, USA

Abstract

Near-infrared spectroscopy has significant fundamental advantages that allow the use of transmission for solid samples such as polymers and pharmaceutical formulations in the near-infrared.

Introduction

Near-infrared (identified in 1800 by Sir William Herschel) was the second region of electromagnetic spectrum discovered after the visible, yet little analytical work was performed in the near infrared until the 1950's when Karl Norris at the United States Department of Agriculture (USDA) began investigating agricultural products. Finding transmission difficult, he worked primarily with diffuse reflectance. When transmission is used in NIR, it is more commonly transfectance, in which the beam is reflected from a mirror-like substrate. Yet, the major advantages of Fourier transform near-infrared (FT-NIR) spectroscopy readily lend themselves to true transmission spectroscopy. The intensity of near-infrared bands is approximately an order of magnitude weaker than the fundamental mid-infrared features allowing for relatively long pathlengths. Glass does not absorb strongly in the near infrared, allowing the use of inexpensive culture tubes and glass vials. Cross contamination in the analysis is eliminated since the tubes can be thrown away after use. Glass sampling also makes it possible to easily obtain high quality spectra by heating samples with low melting points so that the spectrum can be obtained in liquid state. Thick samples, particularly packaging materials (typically from thickness of ~0.125 to 3 mm), do not have to be pressed or dissolved and cast as films but can be run as processed.

Experimental

All spectra were acquired on a Thermo Scientific Antaris™ FT-NIR analyzer with a 20-second measurement time at a resolution of 8 cm⁻¹. The background spectra were obtained automatically as part of the collect using the internal background position. The three-position (non-heated) cardholder (Figure 1) was used for the polymer film studies. The polymer film was a commercially purchased protective sheet holder approximately 0.120 mm thick. The cast films (0.190-0.225 mm thick) were polymer alloys of polystyrene and polyethylene. A quantitative model for the cast films was built using Classical Least Squares (CLS) in Thermo Scientific TQ Analyst™ software.

To study the effect of temperature, a proprietary waxy pharmaceutical

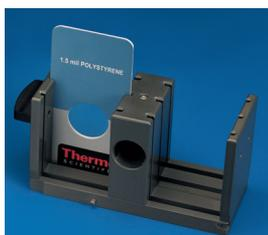


Figure 1: Three-position cardholder

formulation was examined in transmission. The waxy formulation was broken into small (< 1 mm³) pieces and placed in a 6 mm-diameter disposable glass vial.

The tube was placed in the sample position of the heated cuvette/culture tube holder in the Antaris analyzer (Figure 2). The temperature was raised from 28 °C to 60 °C at 2-degree increments. (Figure 3) Prior to collecting the spectrum of the sample, the temperature was allowed to equilibrate and then a background was collected using the computer controlled sample shuttle.



Figure 2: Heated three-position cuvette/culture tube holder

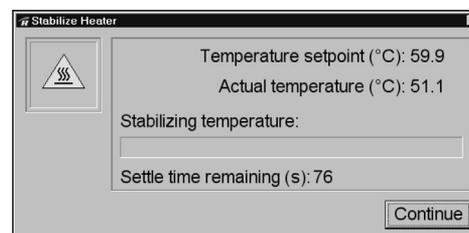


Figure 3: Temperature control in Thermo Scientific RESULT™ software

Results and Discussion

Transmission of Polymers

The near-infrared transmission of the polymer sheet demonstrates that excellent spectra can be obtained from commercial samples run as received. All of the absorbance peaks are in the linear range (absorbances are directly proportional to concentration) and can be readily observed and quantified (Figure 4). The absorbance peaks in the mid-infrared spectra of the same sheet are either in the non-linear range, which would make quantitation difficult or are completely absorbing (Figure 5).

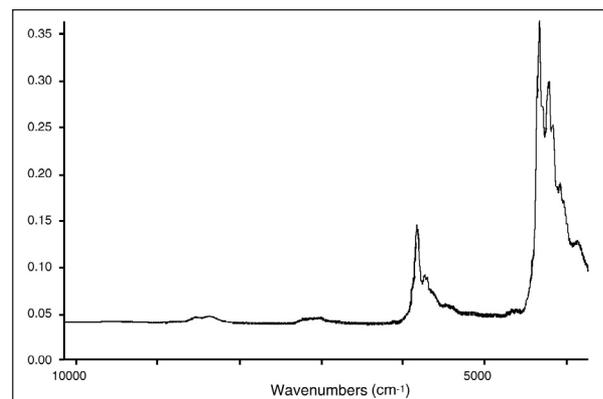


Figure 4: FT-NIR spectrum of plastic sheet

Key Words

- FT-NIR
- Pharmaceutical
- Polymers
- Transmission

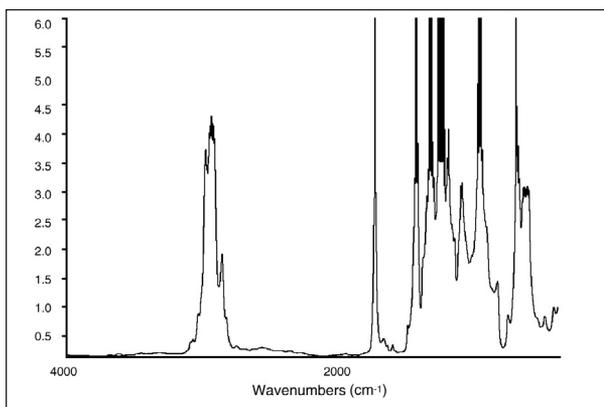


Figure 5: Mid-infrared plastic sheet

Quantification of Polymers

The cast films (0.190-0.225 mm thick) used for quantitative analysis were polymer alloys of polystyrene and polyethylene. Each component (polystyrene and polyethylene) was varied from 0 to 100%. Ten samples were selected as calibration standards and an additional standard was used as a validation standard. The standards were collected using the Collect Standards feature of RESULT Operation software. The CLS method was created using the Build TQ Analyst feature of RESULT Integration software. Three regions were used. As can be seen in Figure 6, the quality of the fit is excellent. For the polystyrene component, the correlation coefficient R is 0.99970 and the (root mean square) error of calibration (RMSEC) is 0.813%. For the polyethylene, R=0.99977 and RMSEC=0.712%.

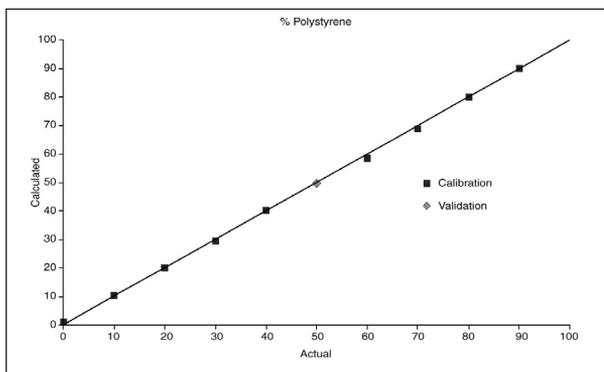


Figure 6: Calibration Plot for Polystyrene

Pharmaceutical Formulation

Diffuse reflection is the most common method for analyzing pharmaceutical compounds. The reflectance spectrum of the waxy pharmaceutical formulation acquired with the integrating sphere (Figure 7) shows that this is not an appropriate analysis technique for this type of sample. The surface of the formulation does not provide a diffuse medium. Consequently the signal-to-noise level is low as

evidenced by the noise (small, sharp peaks) throughout the spectrum. The peaks themselves are broad and ill-defined. Transmission at ambient temperature did not give acceptable results as can be seen in the top spectrum of Figure 8. At ambient temperature, the material has a diffuse, white appearance due to scattering. Because of the short wavelength in the near-infrared, small particles reflect much of the near-infrared light away from the detector. Consequently, the baseline is high (> 2.5 absorbance units) and the signal-to-noise is low. As the material softens, it becomes clearer in appearance, the baseline drops, peaks sharpen and signal-to-noise rises. At 60 °C, the sample spectrum will produce good results. The spectra are presented in common scale with no baseline offset.

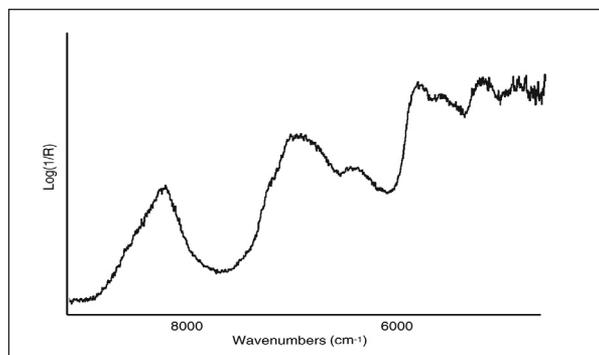


Figure 7: Reflectance spectrum of waxy formulation

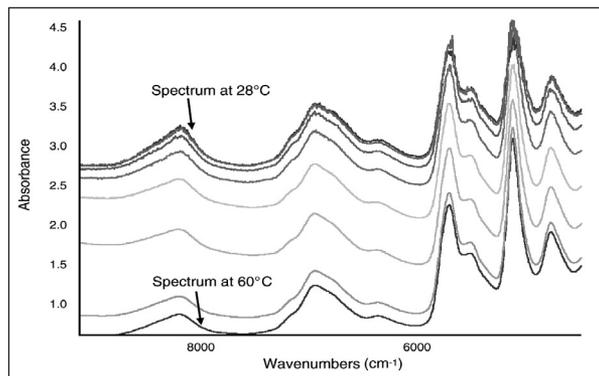


Figure 8: Spectra of waxy formulation acquired at different temperatures displayed on common scale

Conclusion

The Antaris FT-NIR analyzer transmission module provides reproducible quantifiable spectra from thick plastics, liquids and solid samples with a low melting point. The transmission module clearly demonstrates that commercially produced plastic films can be measured without sample preparation. The ability to measure thick samples in a controlled temperature environment with automatic background collection makes transmission spectroscopy with the Antaris FT-NIR analyzer a high-precision and easy-to-implement analytical technique.

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

Africa
+43 1 333 5034 127

Australia
+61 2 8844 9500

Austria
+43 1 333 50340

Belgium
+32 2 482 30 30

Canada
+1 800 530 8447

China
+86 10 8419 3588

Denmark
+45 70 23 62 60

Europe-Other
+43 1 333 5034 127

France
+33 1 60 92 48 00

Germany
+49 6103 408 1014

India
+91 22 6742 9434

Italy
+39 02 950 591

Japan
+81 45 453 9100

Latin America
+1 608 276 5659

Middle East
+43 1 333 5034 127

Netherlands
+31 76 579 55 55

South Africa
+27 11 570 1840

Spain
+34 914 845 965

Sweden/Norway/Finland
+46 8 556 468 00

Switzerland
+41 61 48784 00

UK
+44 1442 233555

USA
+1 800 532 4752

www.thermo.com



Thermo Electron Scientific Instruments LLC, Madison, WI
USA is ISO Certified.

AN51668_E 05/08M

©2001, 2008 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.