

# Monitoring the UV Cure Process of a Polymer Based Ink by FT-IR

## Key Words

- FT-IR/polymers
- Inks
- Macros\Basic
- Paints & Pigments/  
HATR/Quant
- Solid coatings

## Introduction

The degree of cure of a polymer-based ink applied to a Mylar film is readily determined by FT-IR. The ink is screened onto the Mylar film and then exposed to UV light to cure the ink. The determination of percent cure is an important quality control (QC) tool and may also be used to optimize the product manufacturing process.

## Experimental

Spectra were collected using a Thermo Scientific Nicolet™ FT-IR spectrometer and a Smart Multi-Bounce horizontal ATR accessory with a zinc selenide crystal. The ATR sampling technique was chosen for this analysis to enhance the spectral response of the inked surface and to minimize the response of the bulk Mylar film base. The inked side of the Mylar film sample was simply pressed onto the surface of the ZnSe crystal of the ATR accessory. No sample preparation was required. Spectra were collected at 4 cm<sup>-1</sup> with 32 sample scans (40 second sample collection time).

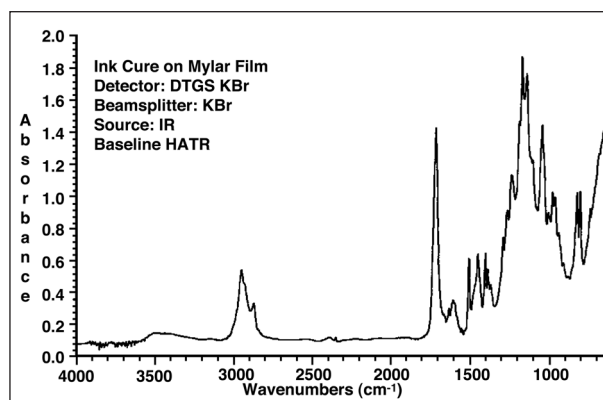


Figure 1

## Results

Figure 1 shows the spectrum of the sample prior to UV irradiation.

For this particular ink, the peak of interest is at 810 cm<sup>-1</sup> and is related to free acrylate monomer. As the ink cures, there will be less of this free monomer, so the intensity of this peak will decrease. In order to quantify the degree of polymerization, an internal standard peak is needed. For this ink, the peak at 830 cm<sup>-1</sup> is unrelated to the cure chemistry and remains unchanged. It is a simple exercise to monitor the ratio of these two peak heights in order to determine the amount of polymerization that has occurred during the cure process.

Figure 2 shows six spectra from the polymerization process with percent of cure ranging from 0 to 87%. Note that for these six spectra autoscaled on the 830 cm<sup>-1</sup> absorbance, the band at 810 cm<sup>-1</sup> varies in intensity. Where the peak heights of the 830 cm<sup>-1</sup> and 810 cm<sup>-1</sup> bands are nearly identical, the ink is uncured. Where the peak height of the 810 cm<sup>-1</sup> band is lowest with respect to the 830 cm<sup>-1</sup> absorbance the ink is 87% cured.

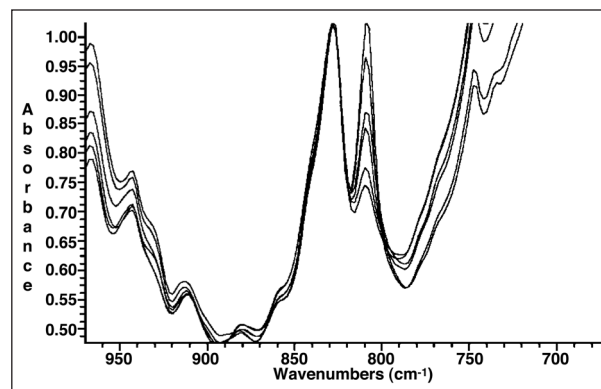


Figure 2

Figure 3 shows a simple macro, written using Thermo Scientific OMNIC™ Macros\Basic™ software, that computes the ratioed peak height of 810 cm<sup>-1</sup>/830 cm<sup>-1</sup> for each sample spectrum using a base-line correction point at 895 cm<sup>-1</sup>.

Start
Peak Height
Store Result
Peak Height
Store Result
Math
Report
End

Figure 3

The plot in Figure 4 shows a linear relationship between percent cure and the ratioed peak area. The physical properties of the cured ink are then related to the percent cure to determine the optimal manufacturing process for the UV cured ink based upon the QC determination of the spectral band ratios.

## Conclusion

The use of FT-IR with ATR sampling provides a fast and easy determination of the quality and state of UV curable polymerization in inks.

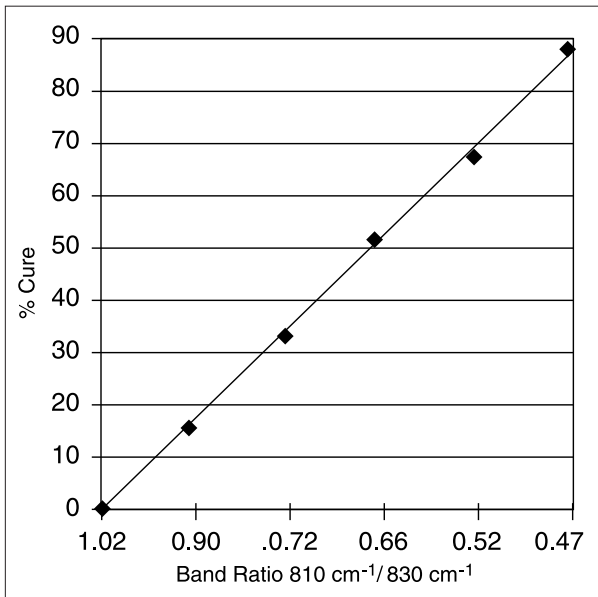


Figure 4

*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 5850 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 76 587 98 88

**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](http://www.thermo.com)

©2007 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.



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

AN50752\_E 11/07M