

Infrared Microspectroscopy in Forensic Science, Part 1 – Hair Fiber Analysis

Pamela A. Martoglio, Ph.D., Spectra-Tech, Inc., Shelton, CT, USA

KEYWORDS

FT-IR/forensic science/hair analysis/microscope/microspectroscopy/ATR objective

INTRODUCTION

Forensic science has benefited greatly from the many advantages that are inherent with infrared microspectroscopy. First of all, microspectroscopic techniques require only a small amount (down to a picogram or below) of sample for analysis. In addition, infrared microspectroscopy is a non-destructive technique. Both of these advantages make infrared microspectroscopy ideal for any sample which is available in a limited quantity, such as evidence in forensic investigations. This application note will discuss the infrared microspectroscopic analysis of one particular type of forensic sample: hair fibers.

PROBLEM

Hair fibers are a very common evidence type found at crime scenes. Light microscopy is routinely used in forensics to determine if an unknown hair sample could have originated from a known source. If chemical information from the hair can also be included, the ability to match an unknown hair with a known source greatly increases; thus, infrared microspectroscopy is quickly becoming a necessary tool in most forensic laboratories.

By using infrared spectroscopy, much more information about the sample can be obtained. For example, if a fiber has been degraded by burial or by time, its surface features which lead to its identification may no longer be present. In this case, infrared microspectroscopy could identify the fiber as proteinaceous, cellulosic, or synthetic. Additionally, infrared techniques can detect any chemical treatments done on the hair (such as bleaching or permanent-waving). Infrared microspectroscopy can be used to detect the chemical damage along the length of individual hair fibers, which could be useful in determining the extent of

natural weathering of the hair or the frequency of chemical treatments. Finally, ATR microscopy enables discrete surface areas on individual hair fibers to be analyzed to detect any residues left behind from styling aids, such as hair spray and conditioner.

All of these techniques will supply the investigator with even more information to aid in the identification and the determination of a possible source of the unknown fibers.

SOLUTION

A single hair fiber was cut to a length of approximately 100 microns and flattened with a roller knife on a clean glass slide. The flattened fiber was picked up with a tungsten probe and transferred to the bottom KBr salt plate in a micro-compression cell, along with a small crystal of KBr. A second KBr salt plate was placed on top of the bottom plate, and the micro-compression cell was tightened until optical contact was made between the fiber and the salt plates. The cell was placed on the microscope stage of a Nic-Plan™ microscope, interfaced to a Magna-IR® FT-IR spectrometer. A background spectrum was obtained through the KBr crystal, and the sample spectrum was obtained through the fiber. Both spectra were collected with 4 cm⁻¹ spectral resolution, and 64 scans were co-added for each. The sample size was 50 x 100 microns.

Figure 1 shows the infrared spectra obtained from a normal, untreated hair and a hair which had been bleached and permanent-waved. The increase in absorbance of the bands at 1175 and 1040 cm⁻¹ is indicative of disulphide oxidation of cystine in keratin. This oxidation can be caused by treatment with alkaline hydrogen

peroxide, or bleaching. The 1040 cm⁻¹ band is due to the symmetric S=O stretch in cysteic acid, and the 1175 cm⁻¹ band is due to the asymmetric S=O stretch. Table 1 shows the band positions indicative of different chemical treatments on hair fibers. Clearly, this information can be very useful in determining possible sources of unknown hair fibers.

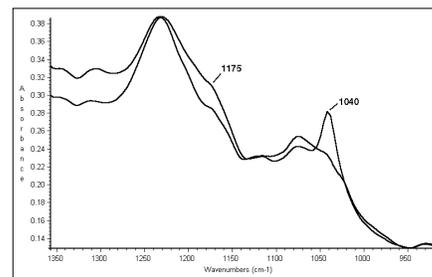


Figure 1: Infrared spectra of a normal, untreated hair fiber and a chemically damaged hair fiber

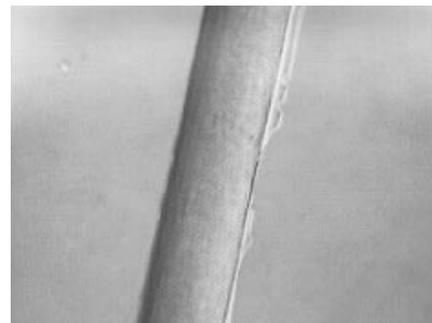


Figure 2: Video image capture of a hair fiber with hair spray visible on the surface

The second example involves the detection of cosmetic treatments found on hair fiber surfaces. Figure 2 shows a video image capture of a hair fiber with hair spray on the surface. If this sample were to be analyzed by a transmission method, as in the first example, the spectrum would predominantly show protein. Hair is a very strong infrared absorber, and detecting something on its surface is very difficult by traditional transmission techniques. However, by using ATR microscopy, a spectrum of

Hair Treatment	Oxidation Products
Alkaline hydrogen peroxide	Cysteic acid (1040 cm ⁻¹ and 1175 cm ⁻¹)
Metabisulfite treatment	S-sulfonate (Bunte-salt) (1022 cm ⁻¹)
Natural weathering	Cystine monoxide (1071 cm ⁻¹), cysteic acid (1040 cm ⁻¹ and 1175 cm ⁻¹), Bunte-salt (1022 cm ⁻¹)

TABLE 1: Oxidation products resulting from different hair treatments¹

predominantly the hair spray can easily be obtained. For this example, a hair fiber was mounted on a glass slide and held in place at both ends with double-sided sticky tape. By placing the sample on a glass slide, the illumination from below the sample can be used to aid in viewing the sample with the survey mode of the ATR objective. A ZnSe crystal ($n = 2.4$) was used in the ATR objective, which yields a sampling diameter of $42 \mu\text{m}$ when using the 2.5 mm diameter upper aperture ($100/2.4 = 42 \mu\text{m}$). The background spectrum was obtained through air (the crystal in contact with nothing), and 64 sample scans were co-added and ratioed against 64 background scans to obtain the infrared spectra at a resolution of 8 cm^{-1} .

Figure 3 shows the infrared spectra obtained from a clean hair and a hair spray coated hair. Clearly, major differences exist between the spectra. By performing a spectral subtraction of the clean hair from the hair sprayed hair, the difference spectrum in Figure 4 results. The main resin in the hair spray, poly(vinylacetate) is easily identified.

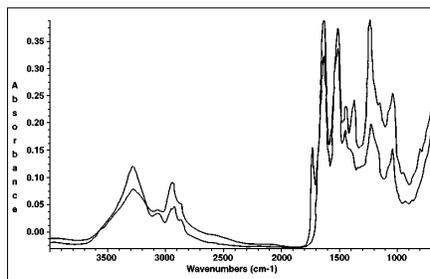


Figure 3: Infrared spectra of a clean hair fiber and a hair spray coated hair fiber

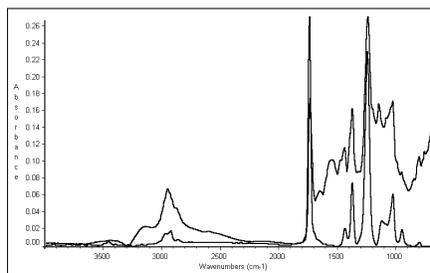


Figure 4: Infrared difference spectrum of the clean hair from the hair sprayed hair, and a reference spectrum of poly(vinylacetate)

CONCLUSION

Infrared microspectroscopy is clearly a very important tool for forensic science. Individual hair fibers can easily be analyzed, and differences due to chemical damage, natural weathering, and cosmetic treatments are readily apparent. Clearly, the information supplied by infrared microspectroscopy will greatly aid forensic scientists in their investigations.

REFERENCES

1. M. Joy and D. M. Lewis, *Int. J. Cosmet. Sci.*, **13**, 1991.
Magna-IR is a registered trademark and *Nic-Plan* is a trademark of Nicolet Instrument Corporation.