

Protein secondary structure elucidation using FTIR spectroscopy

Author

Suja Sukumaran
Thermo Fisher Scientific, USA

Keywords

FTIR, ATR, protein structure elucidation, Biocell calcium fluoride cell, ConcentrateIR2 ATR, transmission

Abstract

Fourier-transform infrared (FTIR) spectroscopy is one of the most versatile analytical tools used across various disciplines. In this study, the Thermo Scientific™ Nicolet™ iS10 and Nicolet iS50 FTIR Spectrometers, equipped with attenuated total reflection (ATR) FTIR and transmission FTIR, were used for the determination of protein secondary structures. Structure calculations based on a protein database as well as spectral deconvolution are discussed. The analyses were quick and easy.

Introduction

Protein secondary structure describes the repetitive conformations of proteins and peptides. There are two major forms of secondary structure, the α -helix and β -sheet, so named for the patterns of hydrogen bonds between amine hydrogen and carbonyl oxygen atoms that create the peptide backbone of a protein.¹ Understanding protein secondary structure is important to gain insight into protein conformation and stability. For example, temperature dependent analysis of the secondary structure is critical in determining storage conditions for maintaining active therapeutic proteins.² Protein secondary structure is also crucial in understanding the structure–function relationship and enzyme kinetics of various proteins.³

FTIR has long been established as a powerful analytical technique to investigate protein secondary structure and local conformational changes.^{1,4} A typical protein infrared (IR) spectrum often contains nine amide bands, with vibrational contributions from both protein backbone and amino acid side chains. Among which, of particular pertinence to protein secondary structure are amide I and amide II bands. The absorptions associated with C=O stretching are denoted as amide I, whereas those associated with N–H bending are amide II. Since both C=O and N–H bonds are involved in the hydrogen bonding between different moieties of secondary structure, the positions of both amide I and amide II bands are sensitive to the secondary structure composition of a protein,^{3,4} although the amide II band is widely viewed as a less useful predictor for quantifying the secondary structure of proteins.

The shifts in the amide I band are often small compared to the intrinsic width of the band, resulting in one broad peak instead of a series of resolved peaks for each type of the secondary structure. Mathematical procedures such as Fourier self-deconvolution and second derivatives can be used to resolve the overlapping bands for the quantitative analysis of protein secondary structure.³ Table 1 shows the secondary structure band assignments for proteins in water. Note that all assignments are depicted as a range, as the exact position of each peak varies from protein to protein due to the differences in hydrogen bonding interactions and the environment of the proteins.

Secondary structure	Band assignment in water
α -Helix	1,648–1,657 cm^{-1}
β -Sheet (high-frequency component)	1,623–1,641 cm^{-1} 1,674–1,695 cm^{-1}
Random	1,642–1,657 cm^{-1}
Coils	1,662–1,686 cm^{-1}

Table 1. Secondary structure band assignments for protein in water.²

With a range of sampling techniques, including transmission, ATR, and infrared reflection absorption spectroscopy (IRRAS), FTIR is particularly advantageous in terms of its versatility and general applicability compared to other analytical techniques for protein secondary structure analysis. Protein sample forms suitable for FTIR analysis include lyophilized powders, water solution, and colloids, to name a few. We report herein two examples of protein secondary structure determination using transmission FTIR and ATR, respectively. Both methods are fast, consume a minute amount of sample, and require minimal sample preparation.

Experiment

All proteins were procured from Sigma-Aldrich (MO, USA) and used as received. For the transmission studies, a BioCell™ Calcium Fluoride Cell (Biotools, Jupiter, FL) was used, and all measurements were carried out at ambient temperature. A 10 μL protein solution was placed at the center of the window,

and the protein solution was sandwiched between the two CaF_2 windows, and placed in the holder. The concentration of protein tested was between 6 and 12 mg/mL. A 6 μm path length was created by sandwiching the two CaF_2 windows. CaF_2 windows are suited for water-based sample analysis. As water has a significant absorption peak at 1,645 cm^{-1} region, a small path length of 6 μm can effectively avoid saturated water peaks.

A purged Nicolet iS10 FTIR Spectrometer, equipped with a DTGS detector, was used for transmission analysis. The scan parameters used were 256 scans with a resolution of 4 cm^{-1} . The Thermo Scientific Smart OMNI-Transmission™ Accessory allows for a quick purge of the chamber, eliminating the need for water vapor subtraction in most analyses. Secondary structure analysis of the buffer-subtracted spectra was carried out using the built-in feature of the PROTA-3S™ FT-IR Protein Structure Analysis Software. Secondary structure calculation in PROTA-3S software is based on a database of 47 secondary structures (for more information visit www.btools.com).

For ATR analysis, a ConcentratIR2™ Multiple Reflection ATR Accessory (Harrick Scientific Products, Inc. Pleasantville, NY) with diamond crystal was used in a Nicolet iS50 FTIR spectrometer equipped with a mercuric cadmium telluride (MCT) detector. The diamond ATR has ten internal reflections with a nominal angle of incidence of 45 degrees. A 10 μL protein solution in phosphate buffer was dried on the surface of the ATR crystal under a stream of nitrogen. Scan parameters used were 256 scans and a resolution of 4 cm^{-1} . Secondary structure determination was carried out using the peak resolve feature of the OMNIC™ Software.

Results and discussion

Transmission-FTIR with Bio Cell

Figure 1 shows the overlay of three FTIR spectra: phosphate buffer, cytochrome C at 6 mg/mL and 12 mg/mL in phosphate buffer, respectively. At first glance, the spectra are predominantly water bands. The three spectra show little difference, even at a high protein concentration of 12 mg/mL.

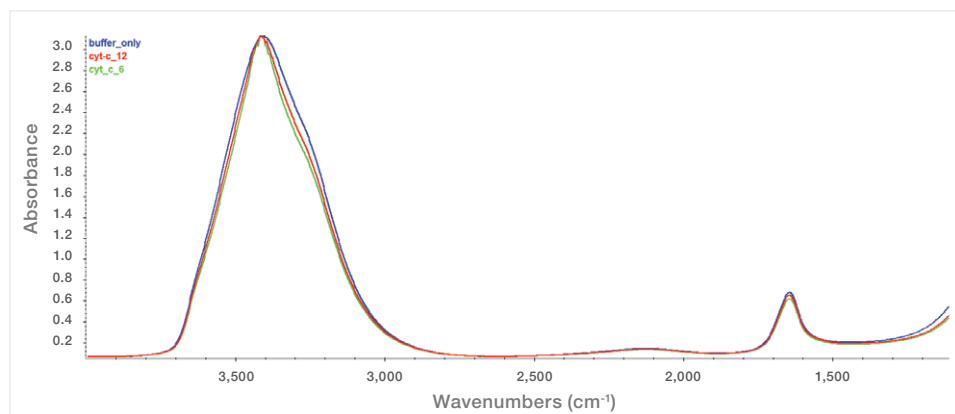


Figure 1. Transmission-FTIR spectra for cytochrome C in phosphate buffer (cytc_12) at 12 mg/mL and 6 mg/mL (cytc_6), and phosphate buffer blank.

Next, the buffer spectrum was subtracted from the raw protein spectra using the PROTA-3S software, and the results are shown in Figures 2A (cytochrome C) and 2B (concanavalin). The amide I and II peaks are clearly discernible in both spectra. The amide I peak position for cytochrome C spectra is $1,654\text{ cm}^{-1}$, suggesting an α -helix dominant secondary structure. For concanavalin A, the amide I peak centers at $1,633\text{ cm}^{-1}$, and there is also a noticeable shoulder peak at $1,690\text{ cm}^{-1}$ (red circle), indicative of the β -sheet component and its associated high-frequency component.²

Table 2 summarizes the secondary structure prediction using the PROTA-3S software. The cytochrome C has 45% α -helix and 5% β -sheet, whereas concanavalin A has 42% β -sheet and 4% α -helix. Differences in secondary structure composition between X-ray and FTIR data are likely due to the physicochemical state of the protein samples such as crystalline versus solution, temperature, pH, buffer conditions, etc. Furthermore, different prediction algorithms could have slightly varying outputs.⁷ Notwithstanding the differences in analytical technique, sample state, and prediction algorithms, the secondary structure elucidation by FTIR using PROTA-3S software is largely in line with that from X-ray. Transmission-FTIR measurements combined with PROTA-3S software offer a facile and fast means to analyze the secondary structure of proteins in solution^{2,3} with minimal sample prep.

ATR-FTIR with ConcentratIR2 Accessory

When the quantity and concentration of protein are limited, FTIR measurements with the ConcentratIR2 Multiple Reflection ATR offer a better alternative than transmission-FTIR spectroscopy. The unique design of this ATR accessory allows for the direct measurement of protein powders, gels, solutions as well as proteins dried on the ATR surface. When concentrating proteins on the crystal surface, caution should be exercised in buffer selection since buffer will also concentrate on the surface of the crystal.

Only those buffers with minimum or no peaks in the amide I and II region should be selected. Figure 3 shows the ATR-FTIR spectra of BSA in phosphate buffer, dried on the crystal from a 1 mg/mL solution. In addition to the amide I and II bands, there are spectral features of the side chain, such as $1,515\text{ cm}^{-1}$ from tyrosine and $1,498\text{ cm}^{-1}$ from aspartic acid. Side chain peaks are critical for the elucidation of protonation and de-protonation states of various amino acids.²

Protein	α -Helix (%)		β -Sheet (%)		Random (%)	
	FTIR	X-ray	FTIR	X-ray	FTIR	X-ray
Cytochrome C	45	41	5	0	50	59
Concanavalin A	4	0	42	48	54	52

Table 2. Comparison of secondary structure calculation from FTIR (PROTA-3S) and X-ray data.

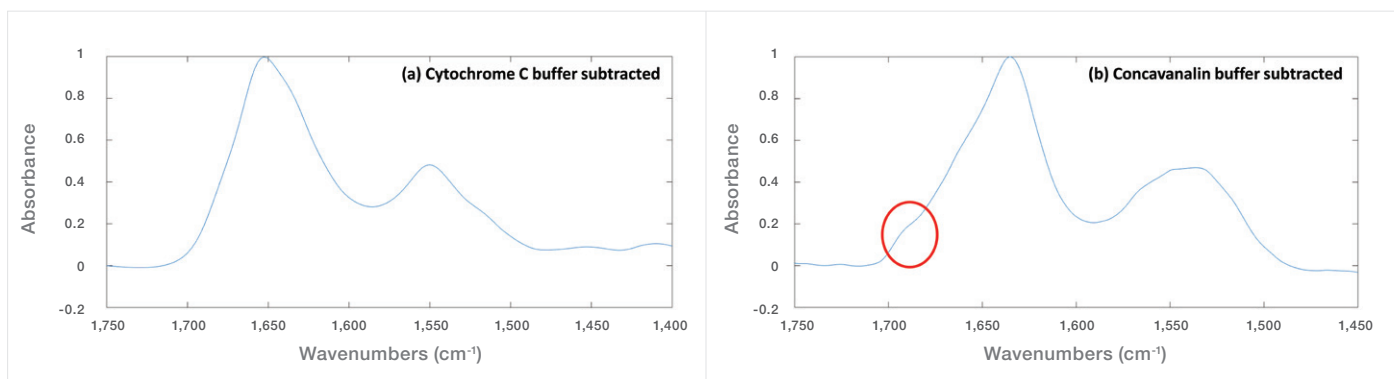


Figure 2. FTIR spectra of (a) cytochrome C and (b) concanavalin A after the buffer spectrum was subtracted using PROTA-3S software.

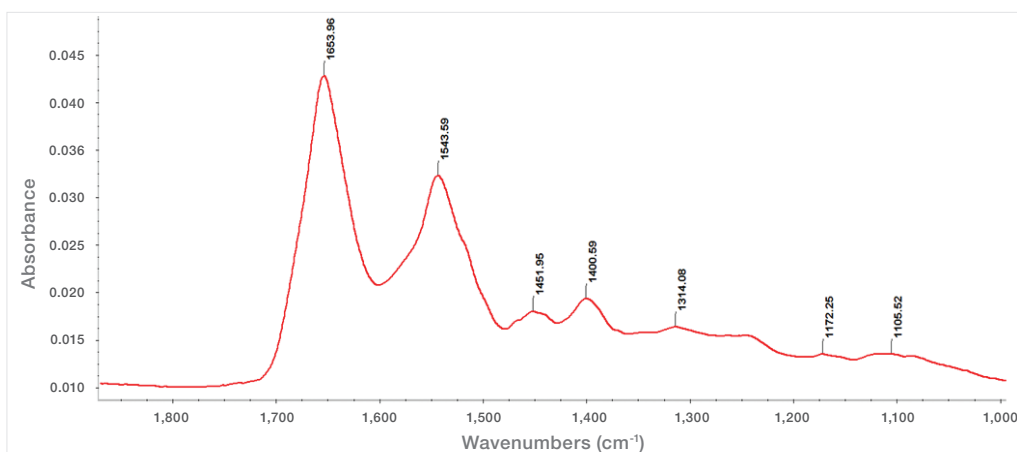


Figure 3. Amide I and II for 1 mg/mL BSA analyzed using ConcentratIR2 ATR on the Nicolet iS50 FTIR Spectrometer equipped with an MCT detector.

Peak deconvolution of the amide I peak (Figure 4) of BSA was carried out using the OMNIC software. It is important to note that second derivative analysis is often performed prior to deconvolution to clearly identify the peaks required for peak fitting.² In the current study, the second derivative peaks obtained (results not shown) are well correlated to the secondary structure peak assignments in Table 1. In order to obtain a good peak shape for peak fitting, a baseline correction on the amide I region was also performed. Baseline correction also effectively excluded the contributions from the amide II region. The deconvolution of amide I resulted in 5 peaks, and the area under each peak was then evaluated against the total area. Amide I peak deconvolution shows a secondary structure composition of 47% α -helix, 3% β -sheet, 24% coils, and 26% random, which is to published FTIR5 and X-ray data.

Conclusion

In this note, we have demonstrated two examples of protein secondary structure elucidation using FTIR spectroscopy. Transmission-FTIR measurements combined with PROTA-3S software provides a facile means to analyze secondary structure of proteins in solution with minimal sample preparation. When the quantity and concentration of protein are limited, ATR-FTIR offers a better alternative by drying the proteins in ATR crystals directly. The data were collected using an older model, the Nicolet iS10 Spectrometer. An improved model, the Nicolet iS20 Spectrometer, offers superior speed and performance over this predecessor model.

References

1. Elliott, A., Ambrose, E. J. Structure of synthetic polypeptides, *Nature* (1950) 165, 921-922.
2. Jackson, M., Mantsch, H.H. The use and misuse of FTIR spectroscopy in the determination of protein structure, *Crit. Rev. Biochem. Mol. Biol.* (1995) 30, 95-120.
3. Barth, A. Infrared spectroscopy of proteins, *Biochim. Biophys. Acta* (2007) 1767, 1073-1101.
4. Byler, D.M., Susi, H. Examination of the secondary structure of proteins by deconvolved FTIR spectra, *Biopolymers* (1986) 25, 469-487.
5. Surewicz, W.K., Mantsch, H.H. New insight into protein secondary structure from resolution-enhanced infrared spectra, *Biochim. Biophys. Acta* (1988) 952, 115-130.
6. Sukumaran, S., Hauser, K., Maier, E., Benz, R., Mantele, W. Tracking the unfolding and refolding pathways of outer membrane protein porin from *Paracoccus denitrificans*, *Biochemistry* (2006) 45, 3972-3980.
7. Klose, D., Janes R.W. 2Struc – the protein secondary structure analysis server, *Biophysical Journal* (2010) 98, 454-455.

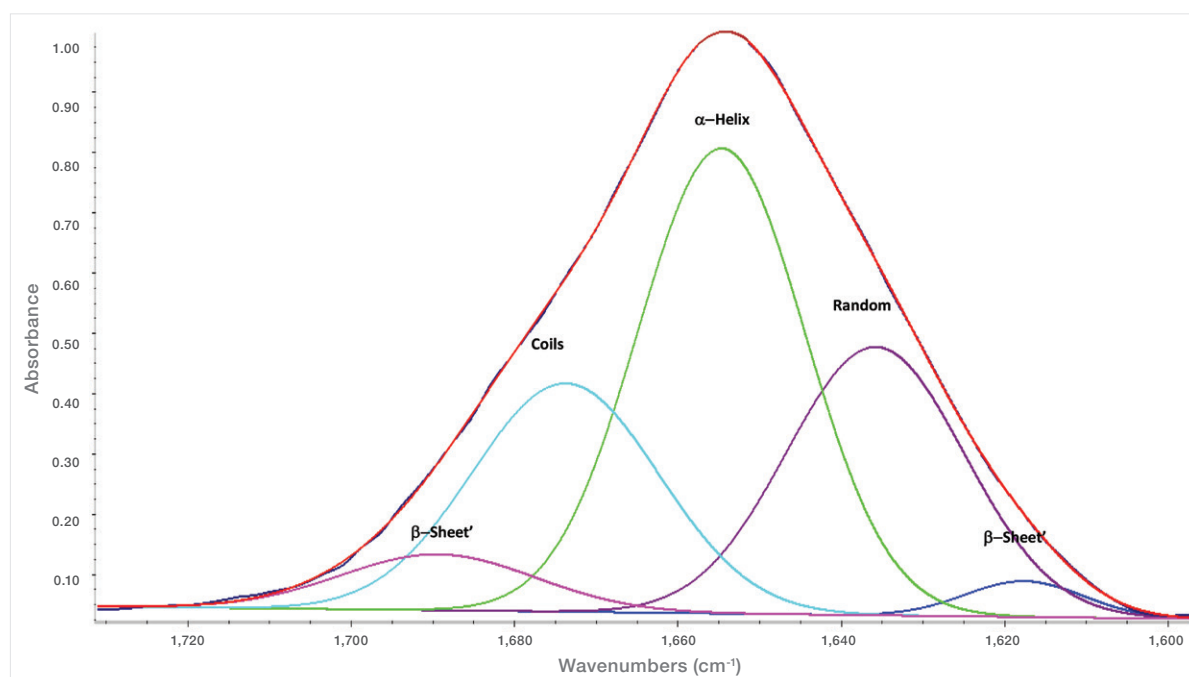


Figure 4: Peak deconvolution of amide I peak of BSA using Peak Resolve function of OMNIC software.

Learn more at thermofisher.com/brighteroutcomes