Application Note: 52080

Analysis of a Wound Dressing

Paul Mack, John Wolstenholme, Andy Wright, Thermo Fisher Scientific, East Grinstead, UK

Key Words

- ESCALAB 250Xi
- Spectroscopic Imaging
- Surface Analysis
- Wound Dressing
- XPS

The Thermo Scientific ESCALAB 250Xi has been used to measure the silver distribution in a wound dressing. This analysis requires the ability to produce images over a large area and at high resolution over a small area. The quantity of silver present in these dressings is small so high sensitivity is essential.

Introduction

Silver containing wound dressings are now widely used to minimize the occurrence of bacterial infection in certain types of wound. The usefulness of these dressings depends upon the chemical and physical nature of the silver, which includes the way in which the silver is distributed over the dressing.

The spectroscopic imaging capability in the ESCALAB 250Xi makes this an ideal instrument for assessing the quality of this type of material. It can be used, for example, to determine the distribution of the silver within the dressing on both a large and a small scale. Large scale spectroscopic imaging can be accomplished by stage mapping and small-scale, high-resolution imaging makes use of the parallel imaging capability of the ESCALAB 250Xi. The concentration of silver in these materials is very low, requiring the high sensitivity available from the ESCALAB 250Xi.

Experimental

Figure 1 shows the optical view of the sample of dressing in the analysis position. The woven nature of the sample is clear in this image.



Figure 1: Optical view of the wound dressing

A set of large-area spectrum images was acquired from this sample. This was accomplished by setting the monochromator spot size to $200 \ \mu\text{m}$ and scanning the sample under the monochromator beam. A spectrum was acquired for Ag 3d and C 1s at each pixel.

Using the large area images, a position was defined for small-area, high-resolution, parallel imaging. This method ensures that the position for the high-resolution map can be selected quickly and accurately. This is a powerful capability provided by the Thermo Scientific Avantage data system.

For the high-resolution measurements, the parallel imaging was combined with spectroscopic imaging. This ensures that, for each element, there was a spectrum acquired at each pixel of the image, enabling accurate quantification of the images.

Results

Figure 2 shows the quantified images from the large area mapping measurement from the C 1s peaks associated with C-C and C-O. Figure 3 is an overlay of quantified images of O 1s and Ag 3d from the same area as the images in Figure 2. Comparison of these images shows that the silver is associated with the areas of the dressing having a low concentration of oxygen.



Figure 2: Quantified chemical state images from the wound dressing. The image on the left shows the distribution of C-C bonds and the image on the right shows the distribution of the C-O bonds. The intensity scale is calibrated in atomic percent.





Figure 3: An overlay of the quantified Ag 3d and 0 1s images from the same area as that shown in Figure 2 $\,$



Figure 4: An overlay of the high-resolution images from C 1s and Ag 3d acquired using the parallel imaging mode These large area maps can then be used to define the analysis position for the small-area, high-resolution, parallel imaging. Figure 4 shows an overlay of C 1s and Ag 3d parallel images clearly showing the distribution of silver in the dressing. This shows that the distribution of the silver is well-defined on a microscopic scale.

Summary

ESCALAB 250Xi has been successfully used to map quantitatively the low concentration of silver on a wound dressing. The results reveal that the distribution of silver is well defined on both a millimeter and a micrometer scale.

The ability to select the position of the high resolution images directly from the large area image is a rapid and accurate method for producing these quantitative images. In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

Africa-Other +27 11 570 1840 Australia +61 3 9757 4300 Austria **Belgium** +32 53 73 42 41 **Canada** +1 800 530 8447 **China** +86 10 8419 3588 **Denmark** +45 70 23 62 60 **Europe-Other** +43 1 333 50 34 0 Finland/Norway/ Sweden +46 8 556 468 00 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 +1 561 688 8700 Middle East Netherlands New Zealand +64 9 980 6700 Russia/CIS **South Africa** +27 11 570 1840 **Spain** +34 914 845 965 Switzerland **UK** +44 1442 233555 **USA** +1 800 532 4752



G Systems Ltd. Trading as ermo Fisher Scientific, East instead, UK is ISO Certified.

AN52080_E 04/11N



www.thermoscientific.com

©2011 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.