

LiPPS Analysis of Chemical Bonding States in an Ionic Liquid

Paul Mack, Thermo Fisher Scientific, East Grinstead, West Sussex, UK

Key Words

Carbon Chemistry, Elemental Composition, Surface Analysis, Tribology, XPS

The Thermo Scientific™ K-Alpha™ XPS was used to analyze the elemental and chemical composition of a worn steel surface, coated with an ionic liquid. Different points on the worn surface were found to have different carbon chemistry.

Introduction

Ionic liquids have recently been identified as potential lubricants for metal/metal contacts. Tribological experiments are used to study the wear and friction behavior of the liquids, but interpreting the data from such experiments requires a detailed understanding of the surface chemistry. Liquid Phase Photoelectron Spectroscopy (LiPPS), a variant of X-ray photoelectron spectroscopy (XPS), is the ideal analytical technique for this purpose, combining surface sensitivity with chemical selectivity.

To fully characterize the surfaces formed during tribological experiments it is necessary to detect and distinguish different chemical bonding states within small, localized areas of the sample. Additionally, the ionic liquid may be deposited on a magnetic substrate, such as steel, which further increases the demands of the analysis. This requires an XPS tool which combines small spot capability with high sensitivity and excellent energy resolution, but which can achieve this on magnetic samples.

The Thermo Scientific K-Alpha was used to investigate the wear surface formed by the tribological action of a brass rod on a steel disc. The ionic liquid 1-hexyl-3-methylimidazolium tris(pentafluoroethyl)trifluorophosphate, was used as the lubricant. The chemical structure of the ionic liquid is shown in Figure 1.

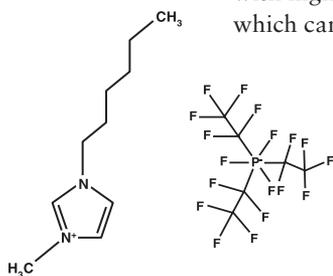


Figure 1: Chemical structure of 1-hexyl-3-methylimidazolium tris(pentafluoroethyl)trifluorophosphate



Experimental

An optical image of the worn steel surface is shown in Figure 2. The image was acquired using the live Reflex optics system, unique to K-Alpha. With this system, the user can confidently select the analysis points for large

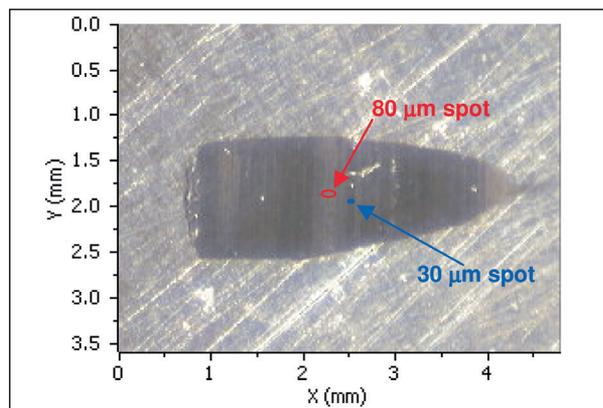


Figure 2: K-Alpha live optical image of tribologically induced wear track. The ellipses indicate the positions and sizes of the analyzed areas.

or small area XPS analysis. The XPS probe is accurately replicated on the live optical view by a graphical marker, allowing the user to easily select the most appropriate probe size for the feature of interest.

The wear track on the steel surface has dark and light stripes of differing widths, which are easily observed in the K-Alpha optical image. Using the graphical probe marker, it was straightforward to select 30 μm and 80 μm spots respectively for analyzing the differently-sized dark and light stripes (see Figure 2).

Results

Elemental Composition of Worn Surface Lubricated with Ionic Liquid

LiPPS survey spectra (Figure 3) were acquired from the areas marked in the optical image. Survey spectra allow the elemental composition of the worn surface to be quantified at the two different analysis points (see Table 1).

Atomic Concentration		
Element	Light Stripe	Dark Stripe
P	3.20	3.05
C	43.05	45.05
N	4.47	4.48
O	–	3.13
F	49.27	44.29

Table 1: Elemental quantification of worn steel surface at two analysis points

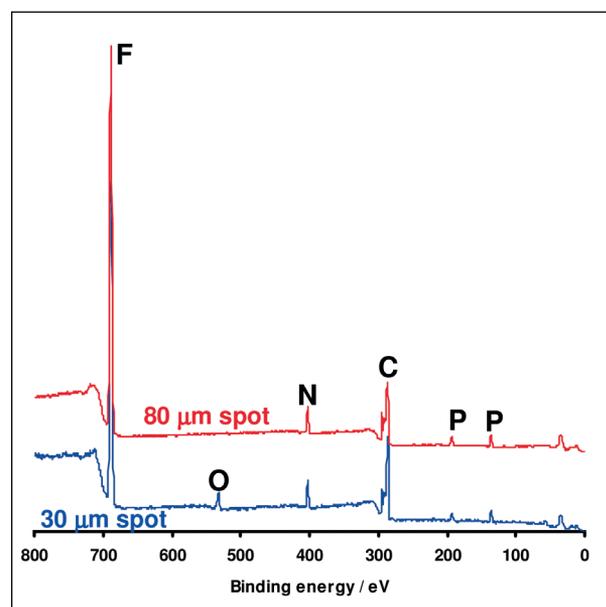


Figure 3: XPS spectra acquired from darker (blue) and lighter (red) stripes in the wear track of the steel disc. Spectra have been normalized and offset for clarity

No spectral features associated with the steel substrate were observed, indicating that the ionic liquid layer at each point was sufficiently thick to completely attenuate photoelectrons from the substrate. It can be concluded, therefore, that the ionic liquid layer is at least 10 nm thick.

The main elemental difference between the two points is the presence of oxygen in the light stripe. The quantities of the other elements identified are similar in each stripe. There is an observable difference in the shape of the $C1s$ peak, suggesting the possibility of different carbon chemistry in the darker and lighter stripes.

Chemical Bonding States in Worn Surface Lubricated with Ionic Liquid

The carbon chemistry at the lighter and darker stripes was investigated using the high energy resolution of K-Alpha. Carbon spectra acquired at the two analysis points allow the user to distinguish between the different bonding states on the worn surface (Figure 4).

The carbon chemistry at the lighter and darker stripes was investigated using the high energy resolution of K-Alpha. Carbon spectra acquired at the two analysis points allow the user to distinguish between the different bonding states on the worn surface (Figure 4).

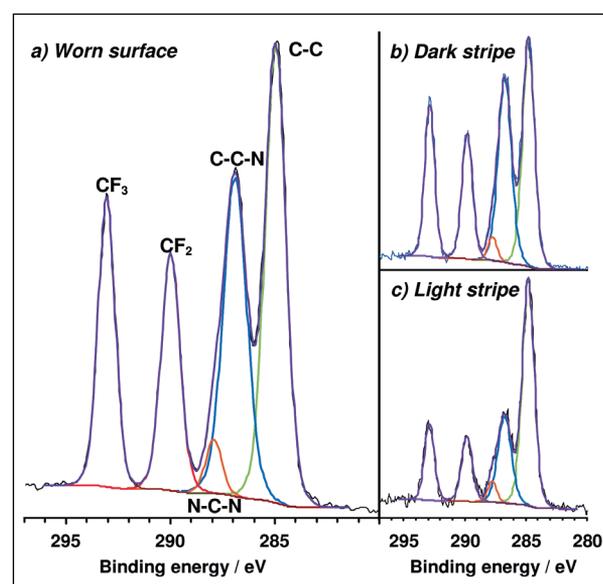


Figure 4: High energy resolution $C1s$ spectra acquired from a) worn surface using 400 μm X-ray spot b) darker stripe and c) lighter stripe on worn surface

The peaks labelled $C-C$ and $C-C-N$ or $N-C-N$ are respectively due to the alkyl chain and imidazolium ring in the cationic component of the ionic liquid.¹ (A small amount of the $C-C$ peak is also due to surface contamination.) The energy resolution of the K-Alpha is sufficient to distinguish between the two different chemical states of carbon in the imidazolium ring, which consists of carbon bonded to single nitrogen atom and carbon bonded to two nitrogen atoms.

The CF_3 and CF_2 peaks are due to the two different carbon bonding states in the anionic fluorophosphate component of the ionic liquid. The main difference between the carbon chemistry of the pale and dark stripes is the increased amount of aliphatic carbon in the latter (indicated by the larger $C-C$ component in Figure 4).

Summary

LiPPs was used to analyze the elemental composition and carbon chemistry of a worn steel surface formed during a tribological experiment. The surface was coated with the ionic liquid 1-hexyl-3-methylimidazolium tris(pentafluoroethyl)trifluorophosphate. Different points on the worn surface were found to have different carbon chemistry, and the thickness of the ionic liquid layer was determined to be at least 10 nm.

Acknowledgement

Thermo Fisher Scientific wishes to thank the Austrian Center of Competence for Tribology (www.ac2t.at) for the ionic liquid sample and helpful discussions.

Reference

1. S. Caporali *et al.*, *Journal of Electron Spectroscopy and Related Phenomena* 151 (2006) 4-8.

www.thermoscientific.com

©2011-2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Africa +43 1 333 50 34 0	Denmark +45 70 23 62 60	India +91 22 6742 9494	New Zealand +64 9 980 6700
Australia +61 3 9757 4300	Europe-Other +43 1 333 50 34 0	Italy +39 02 950 591	Russia/CIS +43 1 333 50 34 0
Austria +43 810 282 206	Finland/Norway/Sweden	Japan +81 45 453 9100	Spain +34 914 845 965
Belgium +32 53 73 42 41	+46 8 556 468 00	Latin America +1 561 688 8700	Switzerland +41 61 716 77 00
Canada +1 800 530 8447	France +33 1 60 92 48 00	Middle East +43 1 333 50 34 0	UK +44 1442 233555
China +86 21 6865 4588	Germany +49 6103 408 1014	Netherlands +31 76 579 55 55	USA +1 800 532 4752