



XPS Evaluation of Wear Resistant Coatings

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

Thermo Scientific™ K-Alpha™ X-ray Photoelectron Spectrometer (XPS) System, was used to correlate the behavior of wear resistant coatings with their elemental and chemical composition.

Introduction

Additives in lubricating oils play an important role in forming friction modifying layers on metal surfaces. Overbased calcium sulfonate additives are used to deposit wear resistant layers on steel surfaces, for example, and zinc dialkyldiphosphonate (ZDDP) has been observed to form a protective, glassy phosphonate coating on surfaces under tribological load.¹ Layers formed on surfaces under load are unlikely to have a uniform distribution across the sample. Optical images of worn surfaces (Figure 1) show darker and brighter tracks (50–100 μm wide) caused by non-uniform processes during tribological testing.



Figure 1: Optical images taken by the instrument of the analyzed surfaces

The samples used in tribological testing may have a large surface area (several cm^2) compared to the size of a typical high quality X-ray spot (30–400 μm). An experimental protocol which includes both micro and macro analyses is therefore required to characterize the wear tracks. Small spot X-ray analysis can be used to detect and distinguish different chemical bonding states within narrow interesting features and large area mapping can be used to investigate the distribution of these states across the surface.

Thermo Scientific K-Alpha is the ideal analytical tool for investigating the elemental and chemical composition of these surfaces, combining imaging capability with surface sensitivity and chemical selectivity. The design philosophy of K-Alpha means that even when the anti-wear coatings are electrically insulating or have magnetizable substrates, such as steel, straightforward analysis with high sensitivity is still possible even for small spot analysis.

Experimental and Results

Three steel samples were tested for friction stability. One of the samples, labelled as GOOD_{OLD}, had already been under consistent tribological load throughout its lifetime, whereas the other two samples (GOOD_{NEW} and BAD_{NEW}) had not been worn prior to the test. The GOOD_{OLD} sample was tested to see whether aging under load had detrimentally affected its friction stability behavior. It was found that the GOOD_{NEW} sample behaved very well under tribological test. One of the samples (BAD_{NEW}), however, was found to have abnormal

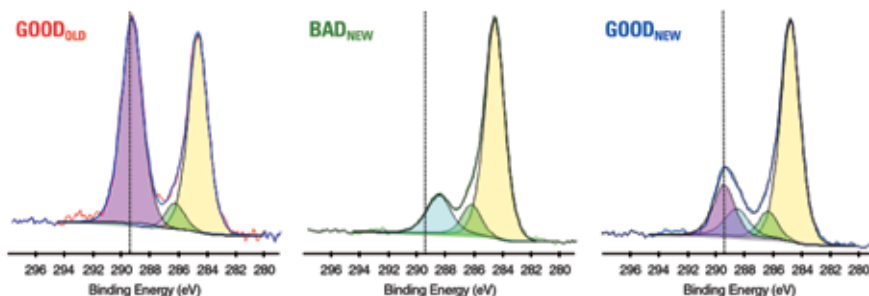


Figure 3: Chemometric analysis of carbon



Thermo Scientific K-Alpha XPS

friction properties. K-Alpha XPS was used to correlate surface composition and chemistry to these different tribological characteristics. Survey spectra (Figure 2) were acquired over wide areas on each sample to identify and quantify the elements present on the surfaces. Although the oil contained a high

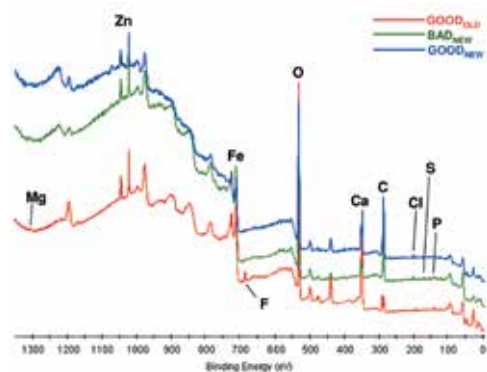


Figure 2: Elemental identification of the elements on the steel surfaces

level of ZDDP additive, the actual amount of zinc on the samples after tribological testing is quite small. (See quantification in Table 1.) In contrast, the GOOD_{OLD} and GOOD_{NEW} samples have a significant quantity of calcium. The source of calcium was the overbased detergents that had been mixed, in differing ratios, with the ZDDP into the oil. The GOOD_{OLD} sample was found to have significantly more calcium on the surface than the BAD_{NEW} and GOOD_{NEW} samples, but significantly less total carbon. The carbon chemistry on each sample was therefore investigated more thoroughly.

Carbon Bonding States Quantification

Chemical State	Relative Amounts of Carbon Bonding States		
	GOOD _{OLD}	BAD _{NEW}	GOOD _{NEW}
C-C	44.03	74.07	65.97
C-O	5.52	9.81	7.68
C=O	1.13	15.61	10.04
CO ₃ ²⁻	49.32	0.51	16.31

The chemical bonding states of carbon on each surface were analyzed by acquiring high energy resolution carbon spectra (Figure 3). Four chemical states were observed: organically derived C-C, C-O and C=O, and inorganic carbonate. A detailed analysis of the carbon chemistry is possible by fitting the data with a series of synthetic Lorentzian-Gaussian peakshapes. This allows a full quantification of the chemical states on the sample surfaces. The GOOD_{OLD} sample has significant amounts of both carbonate and organic carbon. Peak fitting confirms that the GOOD_{NEW} sample has a mixture of carbonate and C=O but that the BAD_{NEW} sample has virtually no carbonate.

Atomic Concentration			
Element	GOOD _{OLD}	BAD _{NEW}	GOOD _{NEW}
P	0.29	0.28	0.21
S	0.29	0.59	–
Cl	0.22	0.80	0.83
C	15.96	39.86	40.10
Ca	14.12	3.79	8.31
O	57.73	45.12	44.71
F	1.50	–	0.28
Fe	6.74	7.67	4.28
Zn	3.03	1.63	1.28
Na	–	0.26	–
Mg	0.12	(tr)	–

Table 1: Elemental surface quantification of samples

XPS imaging of the GOOD_{OLD} sample revealed well defined wear tracks, with alternating regions of high and low carbonate (Figure 4). These regions correlate with high and low calcium concentrations, indicating that the calcium carbonate is formed in the tracks. The thickness of the carbonate film at each point in the map was measured simultaneously with the composition. In the carbonate tracks, the film can be as thick as 87 Å. XPS imaging of the GOOD_{NEW} sample also revealed thick calcium carbonate tracks. These tracks were slightly thicker than for the GOOD_{OLD} sample.

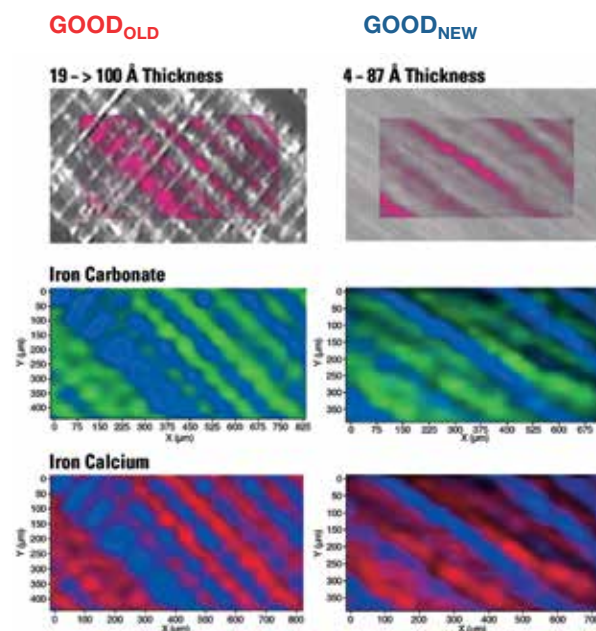


Figure 4: XPS imaging measurements of carbonate film thickness and surface composition

Summary

Thermo Scientific K-Alpha was used to comprehensively characterize the surfaces of three coated steel samples. Two of the samples, labelled GOOD_{OLD} and GOOD_{NEW}, were known to have good friction stability properties and XPS showed the presence of calcium carbonate tracks on these samples. A third sample, which was known to have poorer friction stability properties, labelled BAD_{NEW}, had virtually no calcium carbonate on its surface.

K-Alpha XPS analysis showed that the correct ratio of ZDDP to calcium detergent will result in the formation of calcium carbonate during tribological load, and it is this carbonate that confers good friction stability properties. Calcium carbonate, however, did not form on the BAD_{NEW} sample, indicating an inappropriate ratio of ZDDP to detergent in the oil formulation.

