

# Improved Measurement Accuracy of Wear Metals in Lubricant Oils by ICP-OES

Following ASTM D5185-18 method using an Agilent 5800 RV ICP-OES and IntelliQuant

## Author

Alejandro Amorin  
Agilent Technologies, Inc.

## Get the right answer the first time

The Agilent 5800 Radial View (RV) ICP-OES incorporates smart capabilities, including IntelliQuant, Outlier Conditional Formatting (OCF), and Early Maintenance Feedback (EMF), that deliver deeper insight into methods, samples, and operational status. This insight enables new proactive approaches to reduce remeasurement and downtime, bringing you greater confidence in your results.

Identifying wear metals in lubricant oils can prevent damage to important equipment. Many labs use ASTM D5185-18 standard method for the rapid determination of 22 elements in lubricating oils. Although the ICP-OES method is well-established, labs sometimes need to remeasure samples. In one case, an oils analysis lab had problems with carbon forming in the torch during analysis by ICP-OES. Replacing the torch with an Agilent Easy-fit fully demountable torch for oil analysis reduced the carbon build-up.

In another example, a small contracting lab specializing in the analysis of lubricating oils occasionally received specialty oils containing a high Mo content. Colloidal Mo compounds are often used as antifriction additives in lubricant oils and greases. At high concentrations, Mo will interfere with other elements at the typical emission lines specified in ASTM D5185-18. IntelliQuant is a software function that can accurately identify potential interferences on analytical wavelengths. The function will suggest the most appropriate wavelength based on the elements present in the sample, without remeasuring the samples.

## Automatic flagging of outlying results

The OCF feature in the ICP Expert software makes it quick and easy to find potentially problematic results in a large data set. To simulate an oil sample containing a high concentration of Mo, engine oil samples were spiked at 250 mg/kg and analyzed using IntelliQuant. As shown in Figure 1, OCF highlighted inconsistencies in the results between three Si wavelengths, indicating a potential interference on one of the wavelengths.

Rack:Tube	Solution Label	Outlier Summary	Si 251.432 nm ppm	Si 251.611 nm ppm	Si 288.158 nm ppm	V 309.310 nm ppm	V 311.837 nm ppm
3.24	EngineOil-A21_Spkd-1	F	2.9625	2.9717	2.9796	2.3414	2.3338
3.24	EngineOil-A21_Spkd-2	F	2.9865	2.9878	3.0141	2.3536	2.3380
3.24	EngineOil-A21_Spkd-3	F	3.0333	3.0252	3.0538	2.3683	2.3503
3.25	EngineOil-A21_Mo10_Spkd-1	F	2.9581	3.0367	3.0053	2.3503	2.3314
3.25	EngineOil-A21_Mo10_Spkd-2	F	2.9561	3.0136	2.9878	2.3138	2.2879
3.25	EngineOil-A21_Mo10_Spkd-3	F	2.9801	3.0195	2.9881	2.3302	2.3055
3.26	EngineOil-A21_Mo100_Spkd-1	F	2.9539	3.5467	2.9931	2.3196	2.2878
3.26	EngineOil-A21_Mo100_Spkd-2	F	2.9382	3.5247	2.9709	2.3009	2.2674
3.26	EngineOil-A21_Mo100_Spkd-3	F	2.9704	3.5642	2.9952	2.2982	2.2720
3.27	EngineOil-A21_Mo250_Spkd-1	F	2.9277	4.3820	2.9956	2.2628	2.2340
3.27	EngineOil-A21_Mo250_Spkd-2	F	2.9004	4.3732	3.0015	2.2350	2.2110
3.27	EngineOil-A21_Mo250_Spkd-3	F	2.9986	4.4219	3.0785	2.2438	2.2083

Figure 1. Flagging of Mo-spiked oil sample results informs the user of a difference in the Si concentrations determined using the different wavelengths.

IntelliQuant uses data analytics to rank different emission wavelengths for the same element and identifies likely interferences. By hovering over the "?" symbol next to three Si lines, IntelliQuant identified an interference on Si 251.611 from Mo 251.609, and Si 184.685 and Si 185.005 were considered concentration outliers (Figure 2). IntelliQuant suggested 251.432 nm as the best wavelength to use for the quantitative analysis of Si, indicated by the highest confidence ranking (five stars).

## Oil SRM recoveries

The 5800 RV ICP-OES fitted with an integrated AVS 7 valve system was used for the fast analysis of lubricating oil samples following ASTM D5185-18. The sample introduction system included the Easy-fit fully demountable RV torch fitted with a removable 1.4 mm i.d. quartz injector. NIST 1085c Wear Metals in Lubricant Oil SRM was used to validate the method. As shown in Table 1, recoveries for all elements were within  $\pm 10\%$  of the certified or reference values. The data shows the effectiveness of the 5800 RV ICP-OES for the analysis of wear metals and additives in lubricating oils according to the ASTM method.

## Minimizing downtime

To keep the 5800 running over long periods, smart EMF diagnostics and counters alert the analyst when maintenance is needed. Ensuring peak instrument performance minimizes QC failures and reduces the need for sample reruns.

[www.agilent.com/chem](http://www.agilent.com/chem)

DE.7688425926

This information is subject to change without notice.

© Agilent Technologies, Inc. 2020  
Printed in the USA, April 6, 2020  
5994-1646EN

Element	Used	Flags	Wavelength	Rating	Concentration	Intensity	Background
Si			251.611	* ?	2.86	12124.5	1353.8
			288.158	***	1.96	7084.7	1576.1
			185.005	* ?	2.27	2420.2	716.3
	✓		251.432	*****	1.92	2718.5	1368.6
			184.685	* ?	1.75	1110.4	783.3
			252.411	***	1.99	2088.5	1121.1

Analyte: Si(251.611)  
Confidence: very weak  
Interference: Mo(251.609)  
Confidence: very strong

Analyte: Si(185.005)  
Confidence: very weak  
Concentration outlier

Analyte: Si(184.685)  
Confidence: very weak  
Concentration outlier

Figure 2. IntelliQuant ranks different emission wavelengths for the same element using a star-ranking system. Hovering over the "?" symbol displays reasons for the poor rating on a wavelength.

Table 1. Recoveries of elements in oil SRM.

Element and Wavelength (nm)	Certified Concentration (mg/kg)	Measured Concentration (mg/kg)	Recovery %
Ag 328.068	298	295	99
Al 167.019	292	309	106
B 249.678	304	280	92
Ba 233.527	306	289	94
Ca 422.673	299	305	102
Cd 214.439	301	311	103
Cr 267.716	302	293	97
Cu 324.754	298	307	103
Fe 259.940	301	299	99
K 766.491	295	303	103
Mg 280.270	300	294	98
Mn 257.610	299	300	100
Mo 202.032	299	296	99
Na 588.995	300	299	100
Ni 231.604	306	289	94
P 177.434	304	286	94
Pb 220.353	303	293	97
Si 251.432	293	286	98
Sn 189.925	298	289	97
Ti 334.941	300	302	101
V 311.837	285*	288	101
Zn 213.857	285	275	97

\*Reference value

Full details of this application can be found in Agilent publication [5994-1671EN](#).