

THE SMOOTH OPERATION OF THE EFFICIENT LUBRICANT TESTING LAB

Avio® 500 ICP-OES Oils

- Vertical plasma torch ensures fast, stable operation, dependable results, and low cost per analysis
- Patented Flat Plate™ plasma technology uses half the argon of load-coil systems with the same robust plasma
- Simultaneously collects data on more than 73 elements (UV and visible wavelengths) in seconds for exceptional sample throughput
- Complies with: ASTM D5185, D4951 and D6130

LPC 500™ Liquid Particle Counter

- Inflow sensor measuring particle size and count as the sample makes its way to the Avio 500
- No additional sample preparation required
- Combined sample-to-sample analysis of 45 sec
- Smallest automated particle counter footprint in the industry

OilPrep™ 4/8 Oil Diluters

- High speed dilution of oil samples with a solvent
- Preparation for particle counting by SPOS and elemental analysis by ICP-OES
- Sample throughput between 200 and 500 samples per hour

Clarus® 590/690 GC

- Fast fuel analysis with unique PreVent backflush technology
- Robust, integrated liquid autosampler provides reliability with small sample sizes
- Deeper, more specific insights into the state of your oil
- Detection of potentially damaging compounds
- Precise results on fuel or coolant contamination levels
- More precise and effective scheduling of preventative maintenance
- Complies with ASTM D7593

PerkinElmer Consumables

- Engineered for your instruments
- Tested and validated to ensure accurate, repeatable results, on-time, every time
- Designed to deliver reliable performance, control operating costs and maximize uptime

Power Conditioner

- Ensures a stable power supply
- Recommended for most instrumentation
- Provides clean power at 120, 208, or 240 VAC

Teledyne CETAC Oils 7400 Autosampler with Dedicated Sample Introduction Systems

- Oils 7400 Autosampler for easy switching between oil and coolant testing
 - Features mixing paddle to stir sample prior to analysis
- ASXpress Plus for higher throughput: 2 samples/min for 23-element analysis
- Particle Xpress for LPC 500 coupling to Avio 500, offering particle and elemental analysis in the same run
 - Features switching valve with syringe assist

OilExpress™ 4 Oil Condition Monitoring System

- Automated FT-IR analysis of oils
- Detection of degradation and contamination properties of lubricating fluids
- Provides undiluted samples of used oil to the Spectrum Two FT-IR
- Sample throughput of up to 100 samples per hour
- Complies with: JOAP, ASTM E2414 and D7418

Spectrum Two™ FT-IR Lubricant Analysis System

- User-friendly design with dedicated in-service lubricants analysis protocols
- Clear, high-quality results, regardless of the level of expertise
- Designed to handle demanding quantitative applications
- Easy setup and reporting to get your lab up and running faster

TurboMatrix™ Headspace Sampler for GC

- Determines glycol contamination in in-service oils
- Complies with ASTM D7922

www.perkinelmer.com/lubricants