

Is Your Lab Equipped for Lubricant Analysis?

Avio® 500 ICP-OES Oils

- High throughput sample analysis, measuring concentrations of metals:
 - Lubricants - new and in-service
 - Fuels
 - Biofuels
 - Coolants
- Complies with:
 - ASTM D5185
 - ASTM D4951
 - ASTM D6130

OilPrep™ 4

- High speed dilution of oil samples with a solvent
- Preparation for metals analysis by Inductively Coupled Plasma (ICP)
- Sample throughput can be between 200 to 500 samples per hour

Clarus® 590 GC

- Automated GC system
- Analysis of fuel contamination in in-service oils, including:
 - Gasoline
 - Diesel
 - Biodiesel
- Complies with ASTM D7593

Large Screen Monitors

- Easy viewing of instrument software
- Shown: OilExpress 4 Software and TotalChrom Software

OilExpress™ 4

- Automated systems for FT-IR analysis of oils
- Determine key degradation and contamination properties of lubricating fluids
- Autosampler delivers undiluted samples of used oil to the Spectrum Two FT-IR
- Sample throughput can be up to 100 samples per hour
- Complies with:
 - JOAP
 - ASTM E2412
 - ASTM D7418

STA 6000

- Lubricant stability studies
- Complies with:
 - ASTM D6375
 - ASTM D2595
 - ASTM E1131
 - ASTM E1858

TurboMatrix™ Headspace Sampler for GC

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

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 ASXpress Plus Rapid Sample Introduction System

- 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
 - Features switching valve and vacuum assist

www.perkinelmer.com/lubricants