

SIMPLIFIED IN-SERVICE LUBRICANT ANALYSIS

What Support Does Your In-Service Lubricants Lab Need?

Your in-service lubricant testing lab demands high productivity and reliable results. Does your lab measure up to today's standards with:

- Effortless high-throughput oil analysis
- Reliability and cost efficiency operations for highly automated, rapid, reliable oil condition monitoring results
- Recognized industry standard protocols such as ASTM® and JOAP

With a full suite of instrumentation like Spectrum Two™ FT-IR and OilExpress 4 systems, consumables, training, and service, our combined oil condition monitoring package is designed to support every key aspect of your oil analyses.

Need Infrared Analysis for 100 Samples Per Hour?

The OilExpress 4 system adapts to your laboratory's needs, from dozens to thousands of samples per day. The system minimizes your operational costs by increasing throughput to reduce cost per sample and offers significant savings from decreased solvent waste.

Upgrade your lab today and experience the following benefits:

- Modular solution adapts to changing sampling and throughput needs – simply upgrade the OilExpress 4 system rather than purchasing additional instruments
- Lower solvent consumption – up to 80% reduction in solvent and waste volumes compared with other systems
- Increased throughput – up to 100 samples analyzed per hour with OilExpress 4 DUO. Up to 96 4 oz bottles, or up to 384 17 mm tubes in a single unattended batch on the OilExpress 4 XL system



Spectrum Two™



OilExpress 4™

What Consumables Do I Need for My Lubricant Analyses?

You invest great efforts into your in-service oil monitoring program – and we do the same with our consumables and accessories, tested and validated to fit your lubricant analysis needs. That's why we offer a full range of quality consumables designed only for lubricant analysis. Browse our consumables portfolio for products that offer reliable performance, control of operating costs, and maximized uptime of your instruments. Like our trusted instruments, our consumables offer the best performance over and over.

Confidently select your OilExpress 4 consumables and accessories here:

- ▶ [FT-IR Accessories](#)
- ▶ [Pipettes and Disposable Tips](#)
- ▶ [OilExpress Supplies](#)
- ▶ [Sampling and Starter Kits](#)
- ▶ [Materials Characterization Reference Materials](#)



Complete your solution with OneSource® Asset Management Services offers a single point of contact and custom service plans built to suit your needs. Some services include:

- Improved Service Levels: Proactive monitoring and management of service providers
- Reduced Administration Burden: One contract and PO to manage, one contact to request service
- Global Coverage: 1600 extensively trained and experienced team members located worldwide

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com

For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

23623

PKI

Some Lubricants Applications to Suit Your Lab

[Lubricant Testing Solutions Interactive Brochure](#)

[On-Site Oil Condition
Monitoring Technical Note](#)

[Determination of Benzene in
Gasoline Application Note](#)

