

COMPLETE FLEXIBILITY IN OIL DILUTION

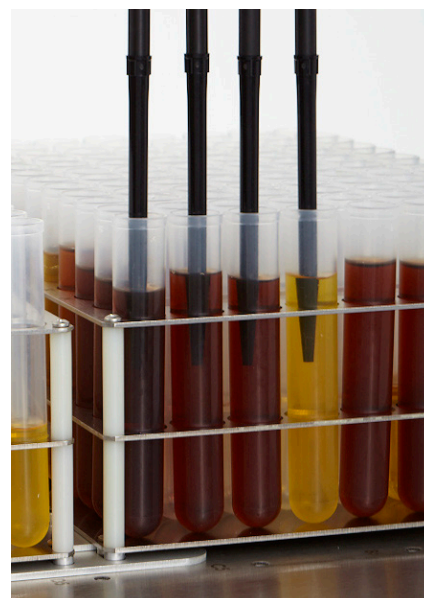
OilPrep 4 Oil Diluter

Need flexible sample preparation for wear metals and particle counting/sizing analyses?

The OilPrep™ dedicated four-tip oil diluter automatically dilutes oil samples with kerosene, in preparation for wear metals and particle counting/sizing analyses by inductively coupled plasma (ICP) and single particle optical sizing (SPOS), such as with the Avio® ICP-OES and LPC 500™ Liquid Particle Counter.

- Pipettes four samples simultaneously, improving laboratory productivity and efficiency
- Four-tip Varispan® arm allows the flexibility to sample from either tubes or bottles
- Ultrasonic Liquid-Level Sensing technology provides unattended dilution of samples of variable volume, saving time
- Disposable pipette tips minimize cross contamination and eliminate tip washing
- Easy-to-use software features predefined pipetting templates for optimized workflows
- Can be upgraded to OilExpress™ 4 Oil Condition Monitoring System to provide FT-IR analysis and oil dilution in a single platform

The PerkinElmer OilPrep 4 oil diluter provides a turnkey solution for anyone needing to increase their wear metals and particle counting/sizing testing throughput, while decreasing labor and material costs.



Why choose PerkinElmer?

The OilPrep 4 oil diluter delivers significant benefits:

- **High throughput** – Significantly improves productivity and efficiency versus manual or single-tip systems. Diluting up to 250 samples per hour, one OilPrep 4 oil diluter can prepare approximately 2000 samples in 8 hours.
- **Flexibility** – OilPrep 4-tip Varispan technology provides automatic computer-controlled variable sample probe spacing for multi-tipped processing of original sample containers. An OilPrep 8 configuration (eight-tip capacity) is also available.
- **Reduced solvent waste** – Use of disposable tips with OilPrep 4 oil diluter eliminates time-consuming tipwash steps required with fixed-tip systems, accelerating throughput and significantly reducing the volume of waste solvent.
- **Minimize cross contamination** – Use of disposable pipette tips also minimizes cross contamination reducing false-positive results that require significant time and expense to retest.
- **Optimized workflows** – WinPREP® software allows user flexibility in choosing vessel types (vials, bottles, tubes), assay setup, and provides potential for fully automated (analytical) integrations.
- **Time savings** – Unlike alternate technologies, PerkinElmer's Ultrasonic Liquid Level Sensing is not impacted by individual sample viscosities, color, or external lighting.



OilPrep 4 XL
Oil Diluter

Move away from manual or single-tip sample preparation methods and leverage complete flexibility with the OilPrep 4 oil diluter!

To find out more, please visit the web at www.perkinelmer.com/lubricants

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 ©2012-2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.