

OilPrep Family of Oil Diluters

Infrared Spectroscopy



Introduction

The OilPrep™ family of oil diluters automatically dilutes oil samples with kerosene (V-Solv™), in preparation for wear metals analysis by inductively coupled plasma (ICP), such as with the PerkinElmer® Optima™ ICP-OES. The OilPrep oil diluter configuration is capable of diluting up to 300 samples/hour using multiple (4 or 8) disposable tips. The Varispan® arm delivers the flexibility to access original sample containers, whereas ultrasonic liquid-level detection provides unattended dilution of samples of variable volume. OilPrep 4 can be upgraded to the OilExpress™ 4 Oil Condition Monitoring System to provide FT-IR analysis and oil dilution in a single platform. The PerkinElmer OilPrep oil diluters provide a turnkey solution for improving productivity in wear metals testing labs while decreasing labor and material costs.

Specifications

	OilPrep 8	OilPrep 4	OilPrep 4 XL
Weight (kg)	125	97	118
Dimensions (mm) excluding PC, bottles and cabling	D=810 W=1580 H=810	D=800 W=1000 H=850	D=800 W=1450 H=850
Maximum capacity:			
17 mm tubes	96	192	384
4 oz bottles	96	32	64
Throughput samples/hr (oil 120 cSt)	300	260	260
Oil Viscosity Range (cSt at 40 °C)	5-500		

Specifications

	OilPrep 8	OilPrep 4	OilPrep 4 XL
Possible upgrades	No	Yes • OilExpress 4 • Liquid-level sensor arm	Yes • OilExpress 4 XL • Liquid-level sensor arm
Tips model	1000 µL disposable tip	1000 µL, wide bore disposable tip	
Ultrasonic liquid sensor arm for variable volume measurement	Included	Optional. Part numbers available for 17 mm tubes and 4 oz bottles	
Maximum bottle height, all racks	100 mm		
Maximum bottle diameter for 32-position rack	50 mm O.D.		
Maximum bottle diameter for 96-position rack	17 mm O.D.		
Tubing material	Tubing FEP (Fluorinated Ethylene Propylene)		
Solvent for dilution	Kerosene (V-Solv™)		

General features

	OilPrep 8	OilPrep 4	OilPrep 4 XL
Software	Oil dilution: OilPrep (WinPREP™) Oil Dilution Software		
Operating ambient temperatures	15-35 °C		
Operation relative humidity	80% at 30 °C non-condensing		
Safety and certifications	<ul style="list-style-type: none"> • Emergency Stop button on diluter • Clear Safety front-panel • CE and CSA certifications 		
Power	100/120 and 200/230 VAC; universal voltage power supply		