

HIAC™ ROC Remote On-line Counter

Simple, trouble-free, and affordable: The HIAC ROC allows you to spend your time preventing problems and less time fixing them.

Benefits

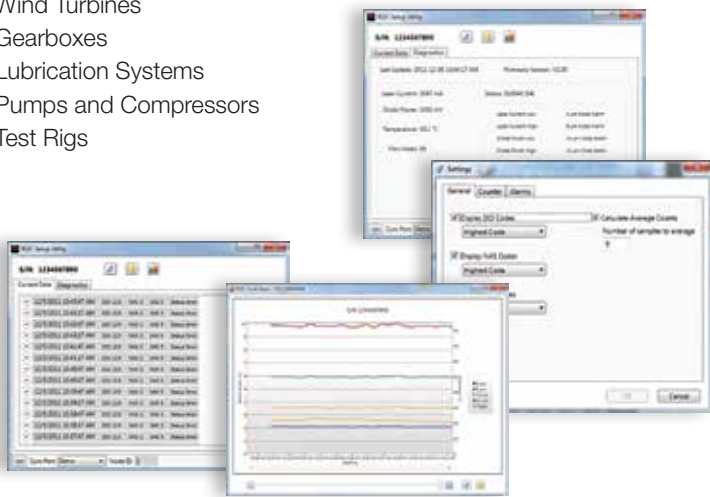
- Can fit into any application, 2-424 cSt
- Data available on highly visual local display in either ISO, NAS or SAE Reporting Codes
- Easily adapt to filter carts with auto stop when oil is clean
- High temperature and pressure capabilities for harsh environments
- Designed for continuous online, maintenance free operation

Software

- Easily configure the HIAC ROC with your PC to fit your specific application
- Log and export data for customized reporting
- Trend data in real time for proactive maintenance
- Connect to a serial port from your Windows® PC using included utility program

Applications / Industries

- Flushing Stands
- Earth Moving Machinery
- Wind Turbines
- Gearboxes
- Lubrication Systems
- Pumps and Compressors
- Test Rigs



Extend machine life, reduce downtime

Constructed for harsh environments, the HIAC ROC excels in high pressure and high temperature applications and offers carefree maintenance. The large flow path minimizes blockages during operation; the lack of moving parts makes maintenance trouble-free. The HIAC ROC on-line particle monitor serves a wide range of industrial and mobile applications: multipoint system monitoring as well as point-of-use applications including hydraulic presses and machines, filter carts, fluid fill stations, hydraulic power units, reclamation stations, and component test stands.



Particle contamination levels provide easy analysis of a machine's condition. With hydraulics and high-speed rotational machinery, particle contamination leads to machine failure, downtime, and maintenance costs. Detecting failure mechanisms such as early detection of oil contaminants allows maintenance personnel to increase machine life and reliability.



Product Specifications

Sizes	4, 6, 14, and 21 μm (ISO MTD)
Light Source	Laser diode
Performance Verification	Optional validation certificate available (ISO MTD at 2.8 mg/L concentration)
Accuracy	0.5 ISO code (min. concentration ISO MTD 2.8 mg/L , max. ISO code = 24)
Display	Optional local display presents ISO, NAS or SAE codes, alarms, fluid temperature and status
Power	9 to 33 VDC, 150 mA (power must be supplied to instrument for operation)
Output	RS-232, RS-485, ground on alarm output
Reports	ISO 4406 NAS 1638, SAE 4059
Fitting Connections	SAE -4 and -8
Sensor Flow Rate	50 to 500 mL/min (0.01 to 0.1 gal/min) through view area
Fluid Compatibility	Hydraulic and lubrication oils, mineral, synthetic (Phosphate Ester option)
Environment	
Operating Temperature	-10 to 60°C (-14 to 140°F)
Storage Temperature	-40 to 85°C (-40 to 185°F) 97% relative humidity, non-condensing
IP Rating	IP66
Wetted Path	Anodized aluminum (standard black), 300 series stainless steel, Sapphire, Brass, Steel, Aflas
Viscosity	2 cSt to 424 cSt viscosities tested at ambient temperature: 25°C +/- 2°C (Model dependant)
Sample Pressure	20-7250 psi (Model dependant)

Beckman Coulter Life Sciences
250 S Kraemer Blvd
Brea, CA 92821 USA
Telephone: 800-866-7889
E-mail: insidesalesgp@beckman.com
www.particle.com



CE CLASS 1 LASER PRODUCT

Lit5224
G9 Printed in USA
© Beckman Coulter Life Sciences 2013. All rights reserved.
In the interest of improving and updating its equipment, Beckman Coulter reserves the right to alter specifications to equipment at any time.

**BECKMAN
COULTER**
Life Sciences