

HIAC GlyCount Glycol Liquid Particle Counter

Features

- With Transaqua HT, Oceanic HW540, Oceanic HW443 and Pelagic 100
- Efficient and intuitive to use
- Immediate laboratory-quality on-site results
- Reports SAE and ISO cleanliness classifications
- Standard bottle and on-line modes
- Multiple language support
- Reduce cycle time for filtration processes
- Monitor real-time system operations
- Track on-line system cleanliness at the point-of-use

Applications

- Glycol hydraulic fluid contamination monitoring
- Cleanliness certification for subsea petroleum production equipment

Rapid and accurate results in liquid analysis

The new HIAC GlyCount provides rapid test results that are as accurate as a conventional laboratory microscope or liquid particle counter analyses. The portable HIAC GlyCount delivers cleanliness results at the point-of-use under actual operating conditions.

With a concentration limit of 30,000 particles/mL, the HIAC GlyCount monitors the dirtiest glycol fluids. Superior optics and design provide eight channels for reporting particle counting data. Temperature measurement is also available for assessing fluid conditions.

The HIAC GlyCount features a wide array of reporting formats, including NAS 1638, SAE AS 4059 and ISO 4406. The HIAC GlyCount can be calibrated with PSL spheres in water or ISO-MTD in glycol.

The instrument contains a buffer for storing 500 analysis results. A rugged carrying case is also included for your portable offshore needs.



Hach, the makers of HIAC liquid particle counters, introduces a new development in glycol cleanliness monitoring. The portable HIAC GlyCount™ measures, stores, and reports glycol fluid cleanliness in a fraction of the time needed for traditional analysis methods.

This new portable HIAC instrument analyzes glycol fluids and coolants in bottle sampling or on-line modes.



Product Specifications

<i>Number of Channels</i>	8
<i>Flow Rate</i>	50 mL/min standard
<i>Calibration</i>	PSL Calibration and ISO-MTD in glycol
<i>Counting Efficiency</i>	Meets JIS B9925:1997
<i>Concentration Limit</i>	20,000 particles/mL at 5% coincidence loss (per ISO 11171) 30,000 particles/mL at 10% coincidence
<i>Fluid Temperature Range</i>	0 to 65°C at 25°C ambient (32 to 150°F at 77°F ambient)
<i>Measured Fluid Temperature</i>	0 to 65°C, ±0.5°C (32 to 150°F, ±0.9°F)
<i>Viscosity Range</i>	2 to 50 cSt
<i>Wetted Materials</i>	Brass, stainless steel, sapphire, PTFE, and Aflas®
<i>Cleanliness Classification</i>	SAE AS 4059, NAS 1638, ISO 4406-1991, ISO 4406.2-1999
<i>Data Storage</i>	500 Sample Records
<i>Dimensions</i>	17.8 D x 33.0 W x 35.6 H cm (7 x 12.5 x 14 inches)
<i>Weight</i>	9.5 kg (21 lbs)
<i>Serial Communication</i>	RS-232
<i>Bottle Operation</i>	Sample Volume: 3 runs (averaged) of 5, 10 or 20 mL/run, programmable Purge Volume: 15 to 30 mL, programmable Pressure Cartridge: CO ₂ , replaceable, rechargeable Operating Time: 60 samples per cartridge (120 mL sample bottle) Shop Air: 90 to 110 psi (6.2 to 7.6 bar) clean, dry
<i>On-line Operation</i>	Sampler Volume: 3 runs (averaged) of 5, 10, 20, 50 or 100 mL/run, programmable Purge Volume: 25 to 999 mL, programmable Fluid Pressure: 100 to 3000 psi (7 to 207 bar)
<i>Power</i>	+24 VDC @ 2A maximum
<i>AC/Battery Adapter</i>	Universal 100 to 240 VAC, 50 to 60 Hz, 60 W
<i>Battery</i>	Nickel-Metal Hydride
<i>Battery Operating Time</i>	100 samples or 4 hours continuous
<i>Battery Recharge Time</i>	2.5 hours
<i>Ambient Operating Conditions</i>	0 to 50°C (32 to 122°F), 20 to 85% relative humidity, non-condensing
<i>GlyCount Storage Conditions</i>	-40 to 70°C (-40 to 158°F), up to 98% relative humidity, non-condensing
<i>Accessories Included</i>	Ultrasonic Bath, Carrying Case, High Pressure Hose Adapter, CO ₂ Bottles, Hand Pump, Sample Bottles, PODS Control Software
<i>Optional Accessories</i>	Additional Sample Bottles (glass) Additional CO ₂ Bottles

Beckman Coulter Life Sciences
250 S Kraemer Blvd
Brea, CA 97821 USA
Telephone: 800-866-7889
E-mail: insidesalesgp@beckman.com
www.particle.com

CE CLASS 1 LASER PRODUCT
Contact manufacturer for complete compliance details

Lit No. 5222
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.

