



FLUID POWER CONTAMINATION MONITORING

INTUITIVE - IMMEDIATE - RELIABLE





**TAKE CONTROL OF YOUR
FLUID CLEANLINESS WITH
HIAC INSTRUMENTS**



Portable Liquid Particle Counter

HIAC PODS+

The HIAC PODS+ provides instant, proactive and predictive maintenance information. Sample your low-viscosity or high-viscosity fluids and report the data to ISO, NAS or SAE standards. The PODS+ has multi-liquid testing capabilities for hydraulics, oil, fuels, glycol and water.

Applications:

- Hydraulic fluids
- Fuels
- Oil
- Water
- Glycol



Key Benefits:

- Save time with custom sample recipes and sample test runs under 60 seconds
- Easily move from one sample point to another
- 1-button sampling, no training or instrument expertise required
- Onsite instant reports that eliminate lab wait times and lab fees
- Multi-liquid sampling capability (fuels, petroleum, water & glycols)
- Ability to report out to multiple standards from the same sample
- Detects moisture in petroleum-based fluids
- Reporting standards: ISO, SAE, NAS, ASTM, GOST, DEF STAN, NAVAIR, User defined

Laboratory Liquid Particle Counter

HIAC 8011+

The HIAC 8011+ is the culmination of over 30 years of particle counting expertise. Designed for ease of use, one-button sampling yields results in under 60 seconds. The sample management system ensures consistent, accurate data and is capable of testing fluids from 1cSt to 425cSt without dilution.

HIAC - the benchmark in industrial liquid particle counting, the brand you know and trust.

Applications:

- Lab sampling
- Large volume samples defined
- Corrosives
- Hydraulic fluids
- Fuel oils
- Water
- Alcohol
- Solvents
- Glycols



Key Benefits:

- Viscosity range from 1 - 425 cSt at room temperature
- 1-button sampling with results in under 60 seconds
- Standards: ISO, NAS, SAE, GOST, DOD and ASTM, User defined
- Size analysis range 0.5 μm to 600 μm (sensor dependent)
- Interchangeable sensors

Online Liquid Particle Counter

HIAC ROC

HIAC ROC particle counters are constructed for harsh environments. The HIAC ROC excels in high pressure and excels in high-pressure and high-temperature applications and offers carefree maintenance. It supports a wide range of industrial and mobile applications, multipoint system monitoring as well as point-of-use applications.






Applications:

- Multipoint system monitoring
- Hydraulic presses and machines
- Filter carts
- Fluid fill stations
- Hydraulic power units
- Reclamation station
- Component test stands

Key Benefits:

- Suitable for use on systems up to 500 bar 7500 PSI
- Long-life laser diode, required for 24-hour continuous on-line operations
- Local display offers ISO codes
- 24/7 trend monitoring

Take Control of Your Fluid Cleanliness with HIAC

	Product	Applications	Reporting Standards	Fluid Viscosity	Features
	HIAC PODS+ Portable Liquid Particle Counter	<ul style="list-style-type: none"> Hydraulic fluids Fuels Oil Water Glycol 	ISO 4406 SAE AS4059 NAS 1638 ASTM D7619-12 GOST 17216 DEF STAN 91-91 NAVAIR 01-1A User defined	1 to 425 cSt with shop air pressure at 100 psig 1 to 150 cSt with internal pump	ISO-MTD: 4, 6, 10, 14, 21, 25, 30, 38, and 70 µm
	HIAC 8011+ Laboratory Liquid Particle Counter	<ul style="list-style-type: none"> Lab sampling Large volume samples defined Corrosives Hydraulic fluids Fuel oils Water Alcohol Solvents Glycols 	ISO 4406 NAS 1638 SAE AS4059 GOST 17216 NAVAIR 01-1A DEFSTAN 91-91 ASTM D7619 User defined Raw counts	Capable of testing fluids from 1 cSt to 425 cSt without dilution	0.5 µm to 600 µm (sensor dependent) Interchangeable sensors
	HIAC ROC Online Liquid Particle Counter	<ul style="list-style-type: none"> Multipoint system monitoring Hydraulic presses and machines Filter carts Fluid fill stations Hydraulic power units Reclamation station Component test stands 	ISO 4406 NAS 1638 SAE AS4059	2 to 424 cSt Suitable for use on systems up to 500 bar 7500 PSI	24/7 trend monitoring

Discover more at beckman.com/HIAC



©2020 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are the trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com

PART-6467APP03.20