



# HIAC 9703+ Liquid Particle Counting System

Easily configurable for maximum flexibility



Be Right™

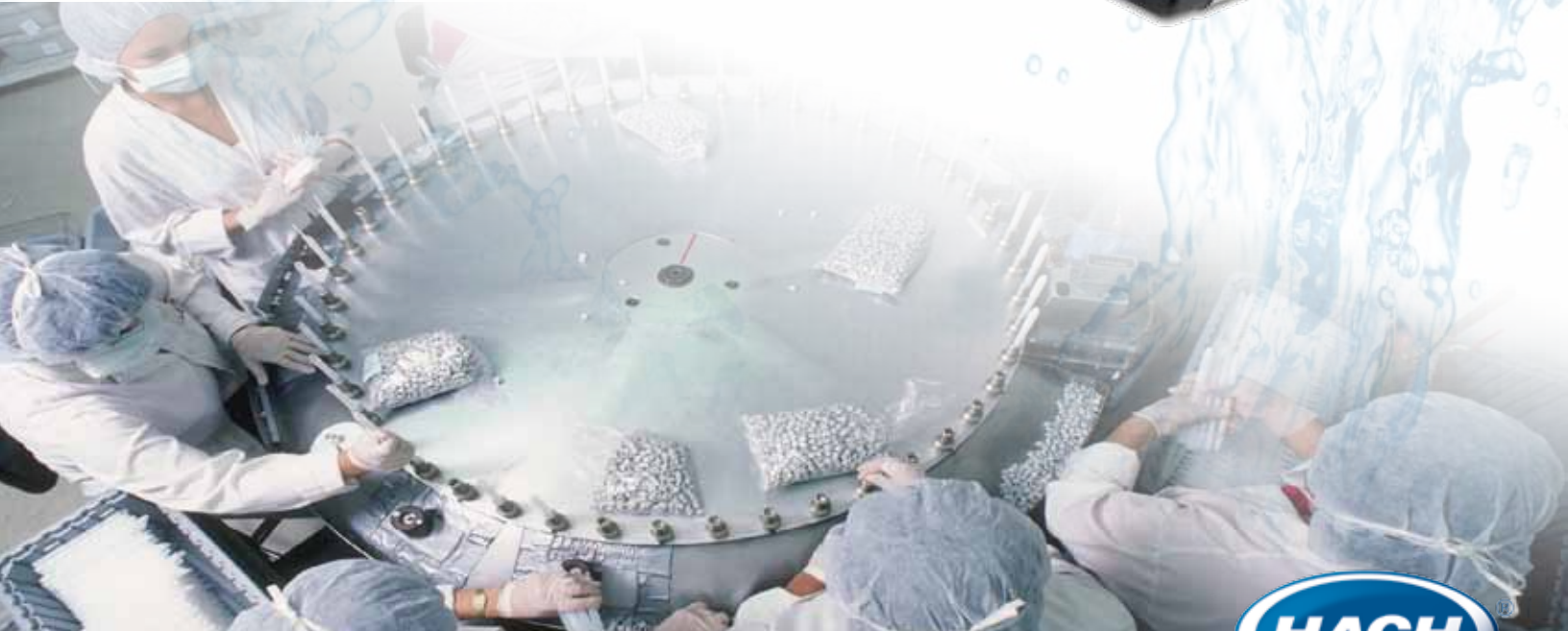
***Efficient Compliance***

***Highly Flexible***

***Advanced Diagnostics***

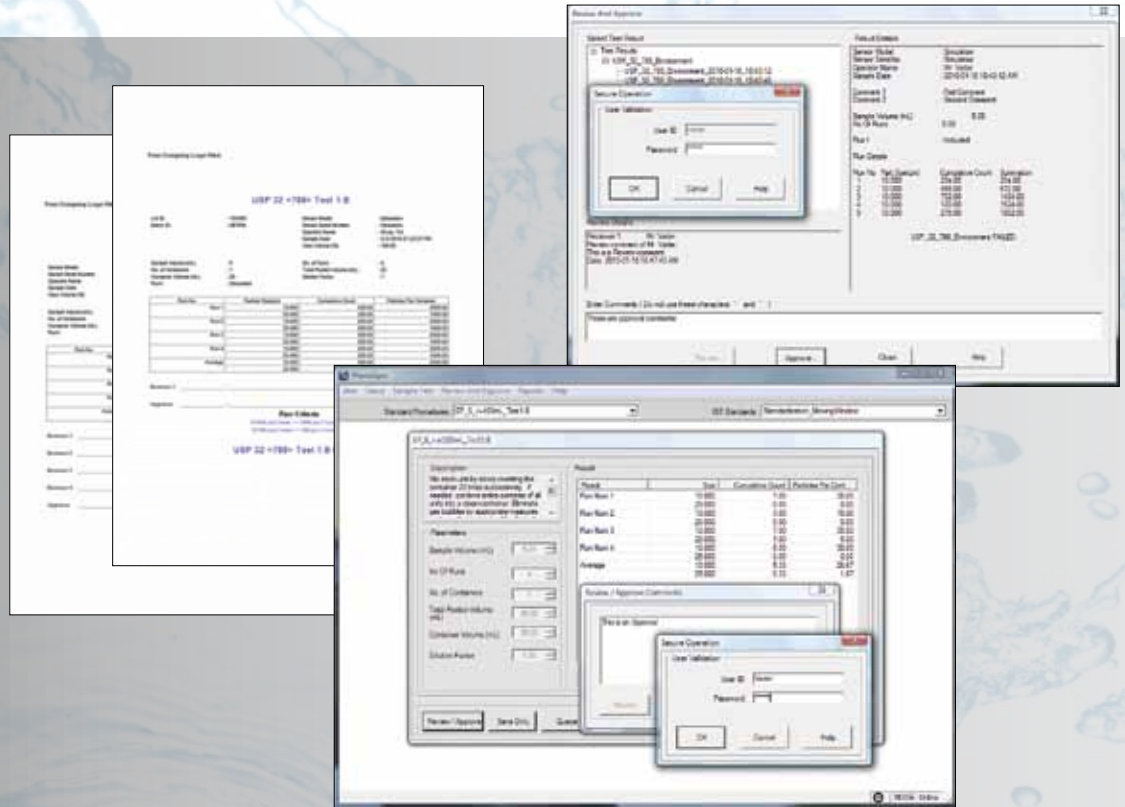
***Time Saving, Faster and Easier***

***Protects Your Investment***



***Be Right™***

# HIAC just got easier!



***Use less sample, save valuable time – protect your investment***

- Industry's smallest tare volume
- Small vial clamp ensures that your sample does not spill during testing

***Enjoy more walk away time***

- Automated flushing and cleaning routines



# PharmSpec

## ***Customized test routines and reporting***

- Customized particle size, sample volume and run test recipes for your unique application needs
- User configurable reports

## ***Added flexibility to configure security access***

- Allows administration to define multi-level security access
- Meets requirements for 21 CFR Part 11 compliance

## ***Save time utilizing pre-configured compendial test recipes***

### **Eliminate manual calculations – Pass/Fail limits**

#### **Injectable Drugs**

- USP <788>
- EP <2.9.19>
- JP <11>, <24>
- KP <52>

#### **Ophthalmic**

- USP <789>

#### **Inhalants**

- USP<788>/Levalbuterol Inhalation

## ***Manage all of your application needs***

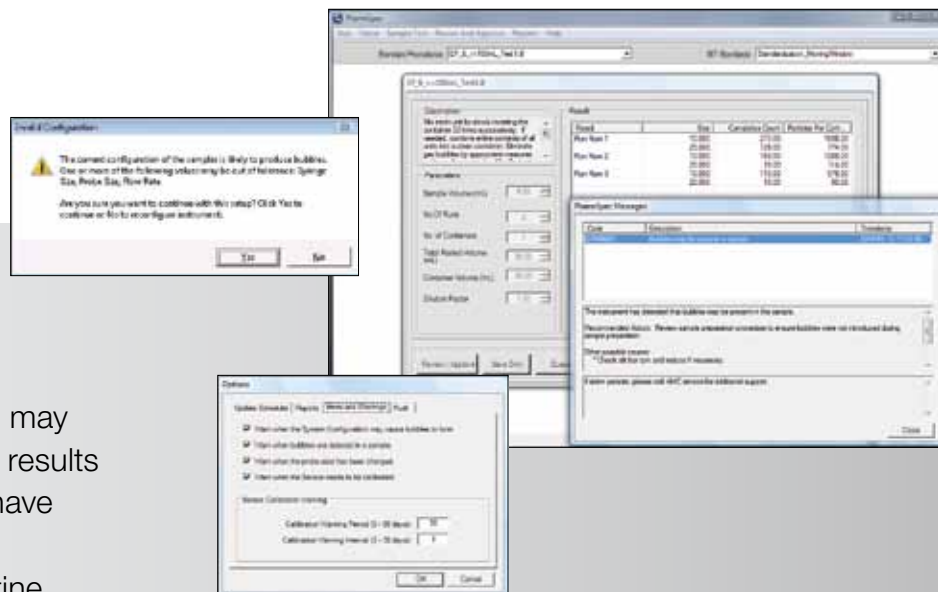
### **Designed to handle the full range from very small (1 ml) to very large (>1000 ml) volumes**

- Small volume injectable (SVI)
- Large volume injectable (LVI)
- Research and Development
- Troubleshooting
- Stability testing

### **Easily configurable for maximum flexibility**

- Interchangeable sensors from 0.5  $\mu\text{m}$  to 600  $\mu\text{m}$
- Interchangeable sampling probes
- Interchangeable syringes – based on sampling size
- Configurable flow rates from 10 ml/min to 100 ml/min





## Eliminate uncertainty from data anomalies

- Alarm notification when bubbles may have impacted particle counting results
- Operator alerted if the sensors have been contaminated
- Advanced notification when routine service or calibration is needed
- Operator notification when regulatory updates become available
- All alarms provide recommended actions

## Eliminate anxiety of data loss

- Data loss recovery feature protects test data in the event of a power interruption or loss
- PharmSpec allows the operator to schedule data base back up to a secure network location

## Turn-key validation and after sale service

Comprehensive local and on-site HIAC service programs with multiple field service offices and authorized distributors are available throughout the world

### Avoid costly down time

- On-site service significantly reduces down time – no shipping of your unit required
- Notification with a phone call when your unit is in need of routine service

### Instrument Validation Services

- SOP assistance and IQ/OQ validation services
- Utilize our expertise to get your system out-of-the-box and ready-to-use
- Ensure the accuracy of your instrument with documented validation and ISO 21501-3 compliant calibrations

## Specifications

### 9703+ Sampler

<i>Temperature Range</i>	5 to 40°C (40 to 104°F) 0 to 80% relative humidity, non-condensing
<i>Temperature Range of Sample</i>	5 to 40°C (40 to 104°F)
<i>Viscosity Limit</i>	<50 cP
<i>Power Requirement</i>	100/230 VAC, 50/60 Hz
<i>Maximum Power</i>	67 W at 115/230 VAC
<i>Dimensions</i>	343 W X 337 D X 482 H mm (13.5 X 13.3 X 19.0 inches)
<i>Weight</i>	10.7 kg (23.6 lbs)
<i>Sample Bottle Clearance</i>	153 mm (6.02 inches)
<i>Volume Accuracy</i>	> 95%
<i>Flow Rate Accuracy</i>	> 95%
<i>Sample Flow Rate</i>	10 to 100 ml/min; Actual flow rate for the system is determined by the sensor's flow rate with step-by-step instructions and Pass/Fail results.
<i>Tare Volumes for Probes</i>	0.091 ml for 81 L X 1.2 ID mm (3.2 X 0.047 inches) 0.172 ml for 154 L X 1.2 ID mm (6.1 X 0.047 inches) 1.57 ml for 140 L X 6.35 ID mm (5.5 X 0.25 inches) - large bore

### PharmSpec 3 Software

<i>Processor</i>	Pentium® 1.0 GHz
<i>Operating System</i>	Windows® XP with Service Pack 3 or later Windows Vista® with Service Pack 2 or later Windows 7 platforms
<i>Memory</i>	512 MB of RAM
<i>Hard Disk</i>	2 GB of available hard-disk space
<i>Drives</i>	CD drive, DVD drive, USB port or network connection is required for installation
<i>Display</i>	Super VGA video adapter and monitor, 1024 x 768 or higher resolution
<i>Power Supply</i>	Uninterrupted power supply (UPS)
<i>Removable Drive</i>	Removable drive for backing up data, such as a recordable CD or a network server that can back up the computer hard drive
<i>Printer</i>	Color printer capable of printing at 300 dpi
<i>COM Port</i>	USB port to connect to the particle counting sensor
<i>Software</i>	Internet Explorer 6.0 or higher and Microsoft® Excel®, Word® or Adobe® Acrobat® (to access PharmSpec reports)

### Typical System Configurations - other options available

<i>Part Number</i>	<i>Description</i>
2089105-01	9703+ Sampler
2089250-150	HRLD-150 Sensor
700060GP	PharmSpec 3 Software
DOC026.53.80099	Validation Documents - English (other languages available)
2089007-025	25 ml Calibration Flow Rate (other flow rates available)

<i>Standard Sensors</i>	<i>Range</i>
HRLD-150	1.3 µm to 150 µm
HRLD-150JA	1.3 µm to 150 µm (chemical compatibility)
HRLD-400	2 µm to 400 µm
HRLD-600JS	2 µm to 600 µm (chemical compatibility)

<i>Compliance Certifications</i>	CE Compliant per EMC and Low Voltage Directives Class 1 Laser Product Complies with IEC/EN 60825-1 and 21 CFR 1040.10 pursuant to Laser Notice 50 **
----------------------------------	---

\*\* Contact manufacturer for complete compliance details

Lit. No. 5223

F10 Printed in USA

© Hach Company 2010. All rights reserved.

In the interest of improving and updating its equipment, Hach Company reserves the right to alter specifications to equipment at any time.

**At Hach, it's about learning from our customers and providing the right answers.**

**Keep it pure.**

**Make it simple.**

**Be right.**

For current price information, technical support, and ordering assistance, contact the Hach office or distributor serving your area.

*In the United States, contact:*

HACH COMPANY World Headquarters  
PO Box 608  
Loveland, Colorado 80539 USA  
Telephone: 800-866-7889  
Fax: 970-461-3914  
E-mail: customersupport@hach.com  
**www.particle.com**

*U.S. exporters and customers in Canada, Latin America, sub-Saharan Africa, Asia, and Australia/ New Zealand, contact:*

HACH COMPANY World Headquarters  
PO Box 608  
Loveland, Colorado 80539 USA  
Telephone: 970-663-9760  
Fax: 970-461-3914  
E-mail: intl@hach.com  
**www.particle.com**

*In Europe, the Middle East, and Mediterranean Africa, contact:*

HACH LANGE GmbH  
Willstätterstraße 11  
D-40549 Düsseldorf, GERMANY  
Tel: +49 (0) 211 5288-0  
Fax: +49 (0) 211 5288-143  
E-mail: info@hach-lange.de  
**www.particle.com**



**Be Right™**