

ZETASIZER NANOSAMPLER

Automated size measurement for the Zetasizer Nano

 PARTICLE SIZE

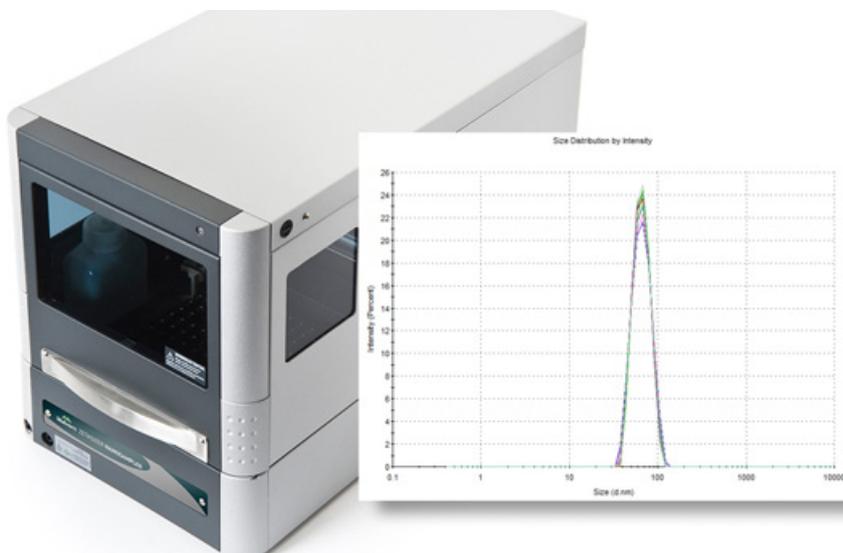
 MOLECULAR SIZE

Key Benefits of the NanoSampler

- Automated measurements reduces operator bias and improves laboratory efficiency
- Unattended multivariate analysis simplifies exploration of the effect of key formulation parameters
- Automates reproducibility studies
- Improved sample throughput
- Reduced operator time
- Maximises the return on your Zetasizer investment
- Simple change over from automated to manual measurements

Introducing the Zetasizer NanoSampler - Providing automated sample delivery for the Zetasizer Nano.

The Zetasizer Nano is the market-leading DLS system with unrivalled performance, made accessible to users of all abilities. Innovations such as the unique NIBS technology push the boundaries of detection limits while simultaneously removing user-influenced variables, thereby achieving better quality, more reproducible data with simplified operation. With the introduction of the new Zetasizer NanoSampler, these benefits can now be applied to an even larger sample set.



30 replicate sample loadings of a 60nm polystyrene latex standard show a standard deviation of only 0.287.

The NanoSampler adds automation and unattended operation in a versatile, compact package, maximising the productivity of your Zetasizer Nano. This is all achieved while maintaining highly precise and reproducible results (as shown in the graph), with confidence in quality of data assured by a carryover of less than 0.05%. The system can be used with a wide range of viscosities and has excellent solvent compatibility.

While the benefits of automated operation are useful for all users, the NanoSampler is ideally suited to laboratories where a large number of measurements are required and reproducibility or multi-variate studies are important.

The NanoSampler can accommodate up to 96 sample vials, accurately delivering up to 3 sequential aliquots of each sample into the Zetasizer Nano for automated particle or molecular size measurements. All operation remains within the easy-to-use Zetasizer software using the simple Standard Operating Procedure (SOP) interface.

The NanoSampler is a powerful addition to the Zetasizer Nano range, reducing the time to achieve the results you require from large sample sets by maximising instrument throughput. The NanoSampler is delivered complete with the software, analysis cell and fittings to attach to the Zetasizer Nano instrument, and, of course, is covered by the industry leading world-wide Malvern applications, support and service.

Parameter	Specification
Sample capacity	96 x 1.5mL vials in 2 plates
Injections per vial	1-3 programmable
Carryover	<0.05% *
Minimum sample volume for single loading	750µL
Maximum sample viscosity	6.44cP
Dimensions and weight (w x d x h)	300mm x 510mm x 360mm; 19Kg
Communication port	USB 2.0
Power requirements and consumption	100 - 240Vac +/- 10%, 50 - 60Hz; 200VA

***Sample dependent**

CONSUMABLES

The NanoSampler has been designed with your time, efficiency and reproducibility in mind. Malvern consumables will ensure that every measurement is made effectively when you load your samples into the NanoSampler. Vials made to exacting specifications avoid variation due to poor seating, and high quality tubing will eliminate failure points from your sample measurement process. In addition, ensuring that your instrument tubing is clean and free of any bends or twists will limit any contamination as well as deliver exactly the same volume with every aliquot. Speak with your local Malvern representative or call our support team to discuss choosing the right Malvern consumables for your needs.

Part Number	Description	Qty
BOT0051	2mL Vial and Snap cap	100
ZEN4070	Cell and Tubing for NanoSampler connection	1
ZEN4071	Tubing for NanoSampler connection	1

SUPPORT AND VALIDATION

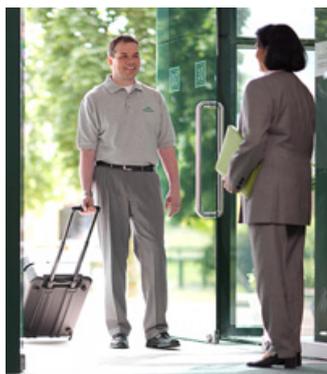
Malvern Applications and Systems Support Worldwide

Malvern instruments quality doesn't end with our stringent Quality Control testing; we provide extensive instrument and application support through our industry leading, 'time proven' Global Support Network.

Our highly trained Malvern Technical experts are just a phone call away from assessing and often resolving your issue. Our field based engineers are also on hand; ready to be dispatched to address those more challenging issues that require that professional, hands-on attention.

Choose from our carefully structured service plans that range from a total support, all inclusive offering to a simple annual service visit providing peace of mind with your instruments performance, condition, and reliability.

Making your annual financial support planning that job you no longer dread, your instrument will be kept in perfect working order and perform as well as the day it was installed. Speak with your local Malvern representative or call our support team to discuss options and how best we can support you.



Malvern Instruments Limited
Groewood Road, Malvern,
Worcestershire, UK, WR14 1XZ

Tel +44 1684 892456
Fax +44 1684 892789

www.malvern.com

Malvern Instruments Worldwide
Sales and service centers in over
65 countries; for details visit:
www.malvern.com/contact

© Malvern Instruments Ltd 2014

Malvern Instruments is part of Spectris plc, the Precision Instrumentation and Controls Company.
Spectris and the Spectris logo are Trade Marks of Spectris plc.



All information supplied within is correct at time of publication.

Malvern Instruments pursues a policy of continual improvement due to technical development. We therefore reserve the right to deviate from information, descriptions, and specifications in this publication without notice. Malvern Instruments shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance or use of this material.

Malvern and the 'hills' logo are International Trade Marks owned by Malvern Instruments Ltd.