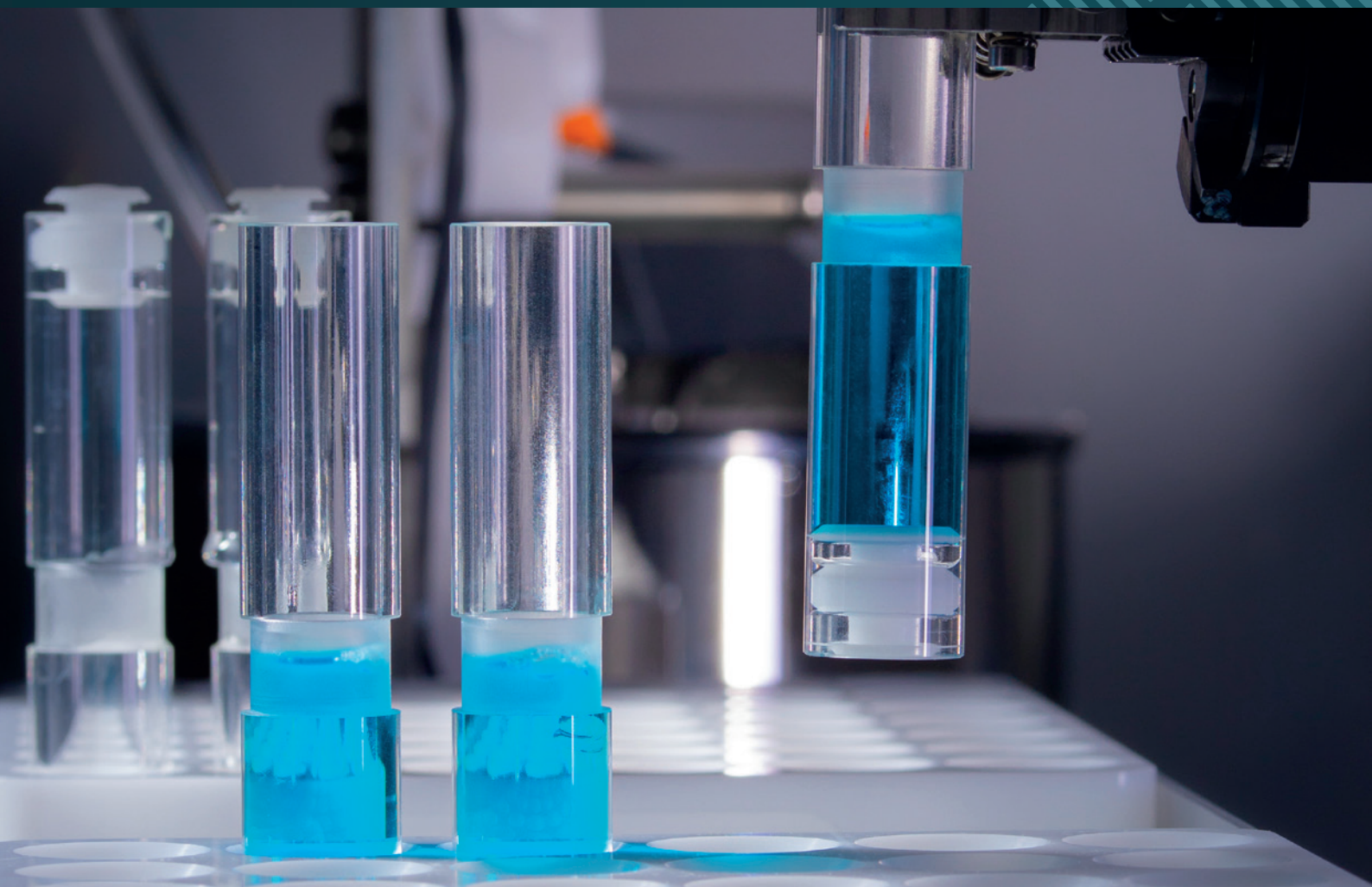




**Malvern
Panalytical**
a spectris company

Mastersizer Auto-Lab

Measuring your samples so you don't have to



Your trusted Mastersizer... now automated

The Mastersizer range of laser diffraction particle size analysers set the standard for delivering rapid, accurate particle size distributions for both wet and dry dispersions. But deep understanding of your samples takes time. The Mastersizer Auto-Lab enables you to minimize manual operating time to add value where you need it most.

Big on size, small on time

Now the Mastersizer can be automated with the Mastersizer Auto-Lab giving you the same great measurement flexibility but without the need for dedicating hours testing samples. Meaning you can dedicate your time where you can add the most value to your business.

Precise and accurate results you can trust

We ensure every particle is measured by adding all the contents vial and using a plunger and rinse. This ensures no sampling errors, bias or chance of cross contamination. By utilising the same Standard Operating Procedures (SOPs) as the Manual Mastersizer you can be confident that our automated analysis will deliver precise and accurate results you can trust.



Up to 45 samples



Highest accuracy



Fast results



Robust and durable

Flexible to changing plans

What if I have urgent samples?

We know things don't always go to plan, so the solution supports walk up sample addition to the end of the run and priority analysis with 3 priority positions available. You don't need to wait for the sequence to finish or have to discard your samples to get the rapid insights you need to control your process.

Instrument Qualification and performance

Comprehensive installation acceptance documents are available to aid validation of the solution, ensuring you can rely on the results. Using the certified reference and calibration material standards for Mastersizer, we can confirm correct installation and track the ongoing performance of our analytical systems.

Our history of automation solutions

The Mastersizer Auto-Lab is developed by our Process and Automation Solutions (PAS) team, a mélange of Panalytical and Malvern engineers with a proven track record of producing automation solutions.

Our automation solutions are already established on products like Insitac inline particle size analysis, or Epsilon, Zetium, Axios, Aeris, and Empryan for XRD and XRF providing you with the capability to customise and integrate with your process, and workflow.

The Mastersizer Auto-Lab benefits you with:

- Fully enclosed and portable Mastersizer autosampler
- Run Mastersizer SOPs
- Measures whole vial to ensure no bias
- Removes chance of cross contamination
- Increase sample throughput
- Position your technicians where they can add more value to your business
- Spend less time worrying about operator error and mix-ups
- Conforms to <ISO13320>

Technical specifications

Description	
Number of samples	42 plus 3 in a priority tray
Sample size range	0.01 µm to 1500 µm (MV) 2100 µm (LV)
Time per measurement	5 minutes (SOP dependent)
Sample volume range	0.5 ml to 10 ml
Unit size (H x W x L)	1696 × 995 × 1370 (Including portable trolley stand)
Unit weight	428 kg (Including portable trolley stand)
Supply voltage	100-240 Vac +/- 10%

How much time can I save?

Want to automate your process with the exceptional capabilities of the Mastersizer particle sizer with autosampler, contact your local Malvern Panalytical representative to discuss how our automation solution could deliver for your process.



About Malvern Panalytical

We draw on the power of our analytical instruments and services to make the invisible visible and the impossible possible.

Through the chemical, physical and structural analysis of materials, our high precision analytical systems and top-notch services support our customers in creating a better world. We help them improve everything from the energies that power us and the materials we build with, to the medicines that cure us and the foods we enjoy.

We partner with many of the world's biggest companies, universities and research organizations. They value us not only for the power of our solutions, but also for the depth of our expertise, collaboration and integrity.

We are committed to Net Zero in our own operations by 2030 and in our total value chain by 2040. This is woven into the fabric of our business, and we help our employees and customers think about their part in creating a healthier, cleaner, and more productive world.

With over 2300 employees, we serve the world, and we are part of Spectris plc, the world-leading precision measurement group.

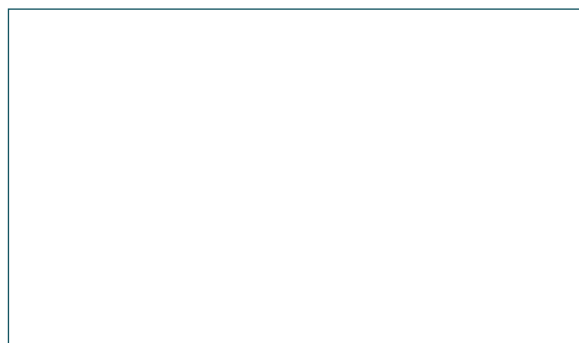
Malvern Panalytical. We're BIG on small™

Service & Support

Malvern Panalytical provides the global training, service and support you need to continuously drive your analytical processes at the highest level. We help you increase the return on your investment with us, and ensure that as your laboratory and analytical needs grow, we are there to support you.

Our worldwide team of specialists adds value to your business processes by ensuring applications expertise, rapid response and maximum instrument uptime.

- Local and remote support
- Full and flexible range of support agreements
- Compliance and validation support
- Onsite or classroom-based training courses
- e-Learning training courses and web seminars
- Sample and application consultancy



Malvern Panalytical

Groveswood Road, Malvern,
Worcestershire, WR14 1XZ,
United Kingdom

Tel. +44 1684 892456
Fax. +44 1684 892789

Lelyweg 1,
7602 EA Almelo,
The Netherlands

Tel. +31 546 534 444
Fax. +31 546 534 598

www.malvernpanalytical.com