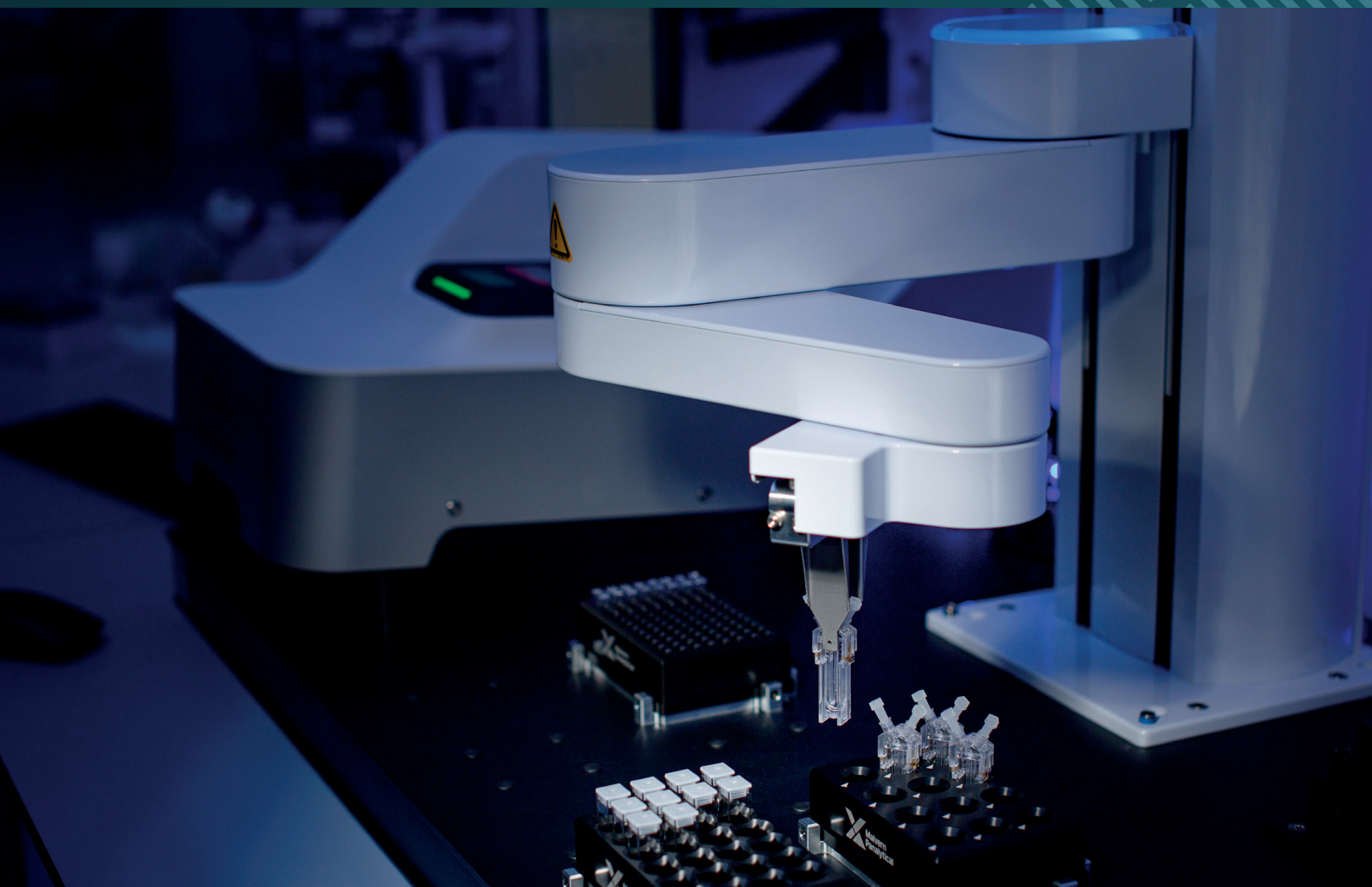




**Malvern
Panalytical**
a spectris company

Zetasizer sample assistant

Empowering scientists, one automated test at a time



Zetasizer Sample Assistant

Significantly improve your system utilization with the Sample Assistant – a new accessory for every Zetasizer Advance system. It frees the operator from manually changing samples.

The result? The same great data quality and versatility you expect from your Zetasizer Advance with even greater efficiency.

Here's how Zetasizer Sample Assistant makes a difference:

- Improves system utilization by minimizing lost time between samples
- More efficient sample handling without compromising data quality
- Sample Assistant works together with your existing methods to simplify method transfer
Need a new method? You can create them in the Method Builder as normal
- Use multiple methods, even on the same sample tray
- Straightforward workflow guides users through maintenance with ease.

Bespoke grippers

Nimble and precise: allows the Assistant to hold your samples securely.



Co-bot

Works safely around people. Changes samples for you with smooth and precise movements.

Space for your Zetasizer

Holds your Zetasizer in the correct position for the co-bot, allowing walk-up use when needed.

Baseboard

Everything has its place on the baseboard: the instrument, co-bot, and storage.

Sample trays

Three types of trays to support Standard and Universal sizing cells and Zeta cells.

Tray bays

Hold any tray in a configuration that works for you. Trays can be held in up to six different positions.

Tool holder

Securely stores datum tools and cell adapters; ready for when the Assistant requires them.

Providing the flexibility you need

The great advantage of the Zetasizer is that the entire system is designed holistically. From the instrument itself to the sample cells, everything about the Zetasizer is designed to ensure optimal data quality. So what's the missing piece of the puzzle?

Until now, you had to be present to operate the instrument and achieve that unparalleled data quality. From now on, thanks to the Sample Assistant users can enjoy the same, high quality data without the need to attend the instrument, having the chance to do other tasks while the Assistant is dealing with the samples.

The Sample Assistant solves both problems: high-quality sample data that inspires confidence, without the need to be present. It is also easy to upgrade, with no instrument modifications, no time lost when cleaning between samples, and no risk of cross-contamination.

Improve Utilization

Minimizes the time lost between samples. Sample Assistant is always ready.

Extends the workday or week. Your Zetasizer with Sample Assistant can run unattended at any time.

Reduces time lost from cleaning flow cells between samples.



Standard Tray

Our standard tray holds samples in presented in 10x10mm cuvettes, such as DTS0012 disposable cuvettes, holds up to 24 cuvettes.



Zeta Cell Tray

Our zeta cell tray is designed to hold up to 12 Zeta capillary cells, such as DTS1070, in the correct orientation.



Universal Cell Tray

Our universal cell tray is designed to hold up to 96 of our Universal disposable sizing cells and it honours the plan geometry of ANSI/SLAS 96-well plates to provide improved compatibility with multi-pipettes and fluid handling systems.

Operates Safely

Works around you: detects obstructions and stops safely.

Makes it easy to recover from a system failure: your data is safe and the last known state is preserved.

Descriptions	Specification
Weight (without Zetasizer)	45kg
Weight (with Zetasizer)	65kg
Dimension (l x w x h)	600 x 900 x 734mm
Power	110 – 240V AC 50/60Hz
Communications	Via Ethernet (USB – Ethernet converter supplied)

Main System	Description
Zetasizer Sample Assistant (base configuration)	Includes: baseboard, stands, 2 x tray holders, 1 x standard tray, tool holder with datum tool, tray datum tool and all fixings
Optional Accessories	Description
Universal Sizing Cell kit	1 x tray holder, 1 x Universal Cell Tray, 2 x Universal Cells, Cell Sealing Jig, 1 x pack cell sealing film (25 sheets)
Zeta cell tray	Cell tray to hold DTS1070 cells
Standard Cell tray	Cell tray to hold standard 10x10mm cuvettes
Universal Cell Tray	96-position cell tray
Tray holder	Additional tray holder with fixings



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.

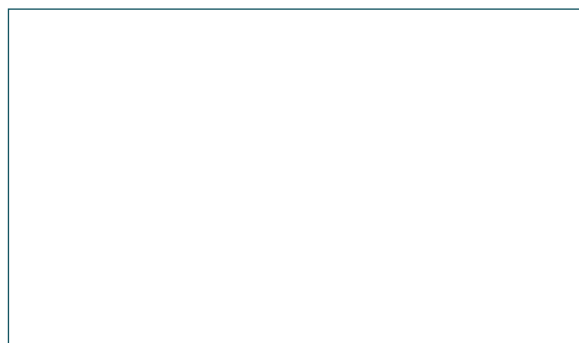
Malvern Panalytical. We're BIG on small™

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