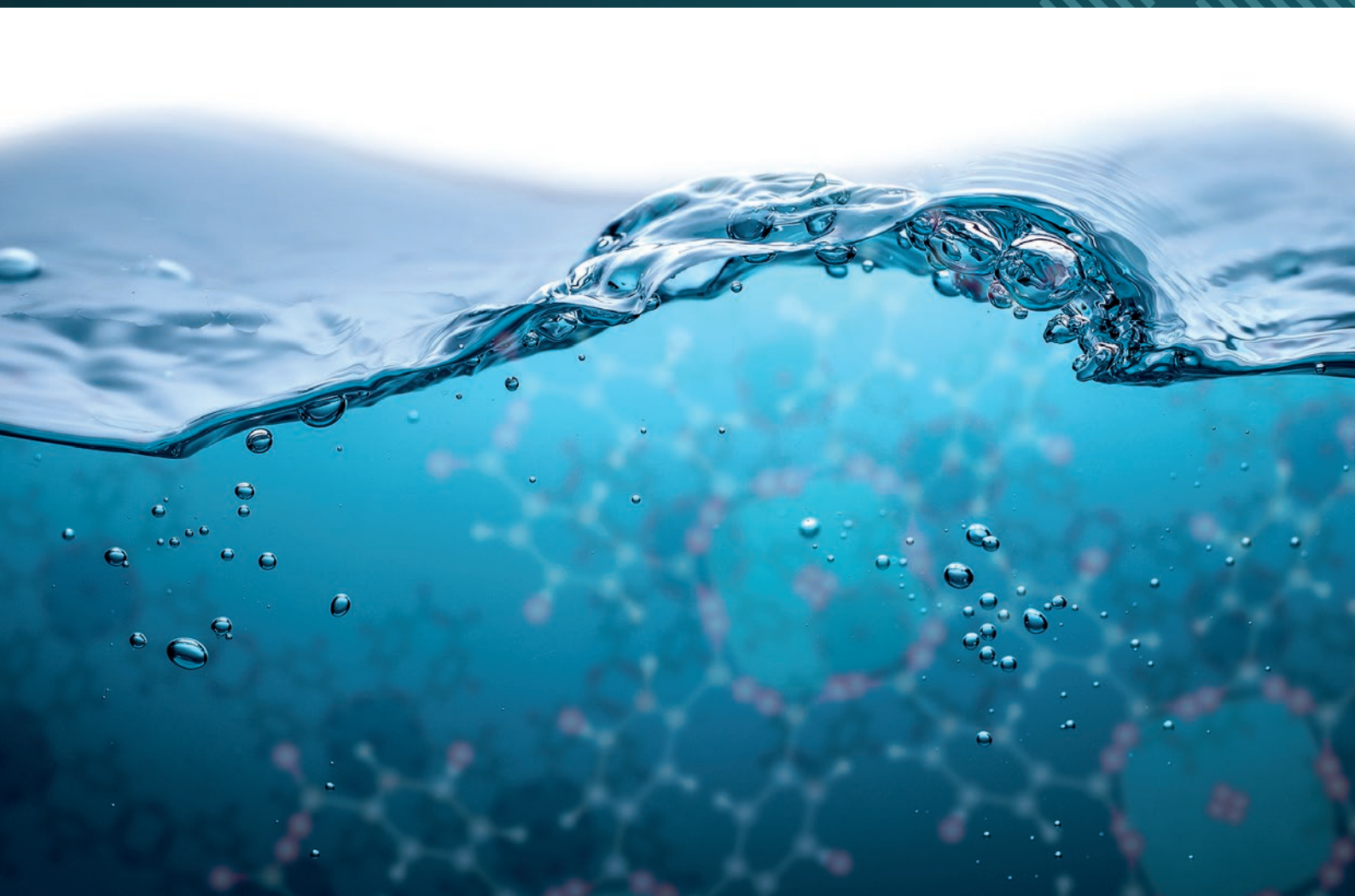




**Malvern  
Panalytical**  
a spectris company

# Analytical solutions for Metal Organic Framework materials

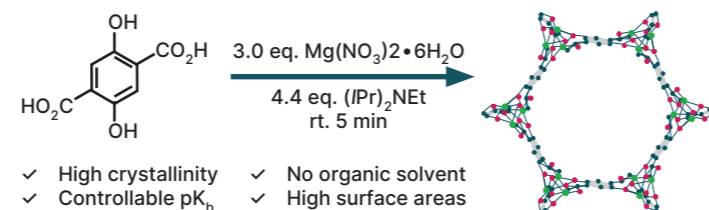
Grow your knowledge. Build robust structures.



# The world of Metal Organic Frameworks

## What are MOFs?

Metal Organic Frameworks (MOFs) are hybrid crystals created by self-assembling organic and inorganic molecules. MOFs contain nano-scale interconnected voids – giving them unrivaled potential for trapping, storing and catalyzing ions and molecules. Adjusting their local environment (such as the electric field, temperature, pressure or chemistry) can enable the adsorption or release of many molecules or ions.



### The hybrid structure of MOFs

Organic molecules can form large, interpenetrating lattices made of a framework of large pores. But, on their own, they are relatively fragile. Reacting these molecules with metal ions creates MOFs: strong covalent structures that are more durable than organic crystals. These materials can be synthesized on surfaces as thin films, obtained from solution as dry powders, or used as nanoparticles in liquids.

## What are they used for?

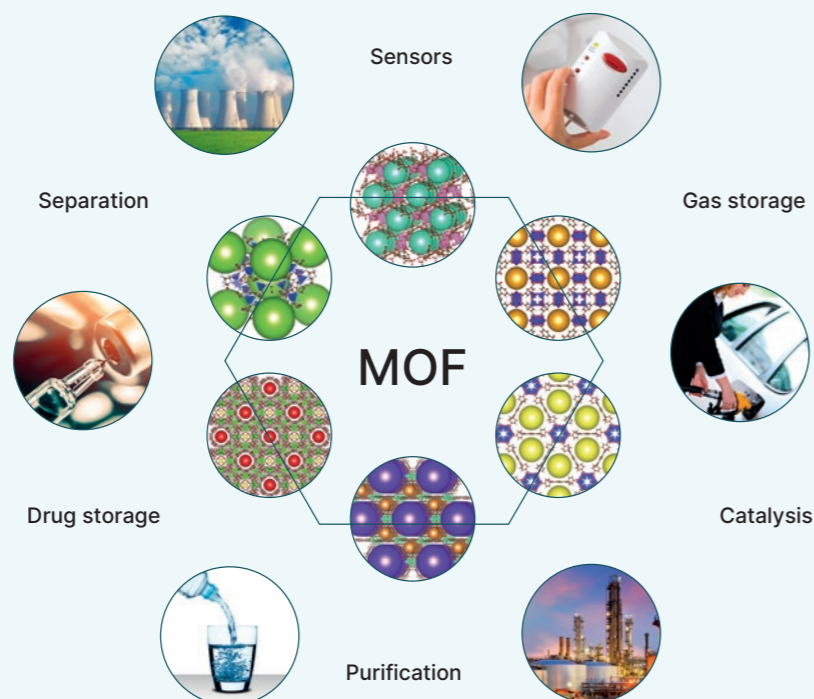
Thanks to their collection abilities, MOFs are ideal for several applications, including gas storage and separation, liquid separation and purification, electrochemical energy storage, catalysis, or drug delivery. They are also useful for gas sensor devices that measure and respond to this capture process.

## How can we help?

At Malvern Panalytical, our global network of over 170 scientists is continuously working to support customers with advanced materials like MOFs.

Our solutions deliver in-depth materials insights, from nanoscale to macroscale. Whatever the project, product or process, we'll help you to optimize your material.

On the following pages, we hope you'll find solutions that could benefit you. If you want to dive deeper into their specific benefits, don't hesitate to get in touch!

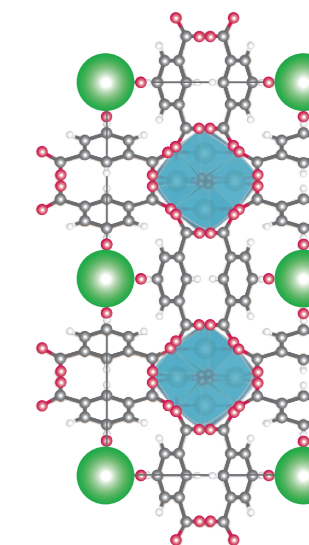


# Where materials analysis comes in

## Crystal structures: Key to collection

What exactly makes MOFs so good at collection? Their complex **crystal structures!** These structures contain voids, and the interconnectivity between these voids makes MOFs nanoporous. The exact structure depends on the MOF's constituent molecules and how it is synthesized. By tailoring this, you can optimize size, connectivity and **binding affinity** to attract specific molecules or ions.

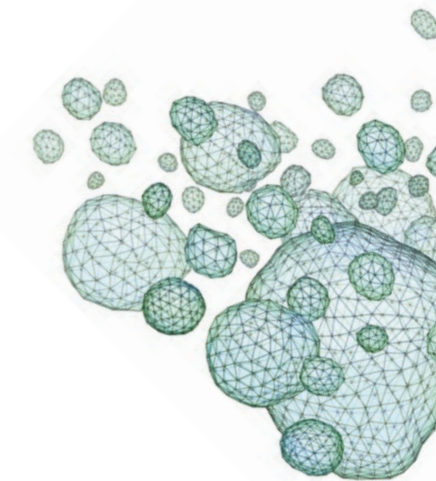
That means you need a deep understanding of the crystalline lattice structure and how it interacts with other molecular and ionic species. And for this, you need a strong understanding of the **structural changes** that result from variations in the host environment.



## Particle morphology for maximum performance

How do you maximize the performance of a MOF powder to ensure robust, consistently processable products? By optimizing the physical properties of the particles!

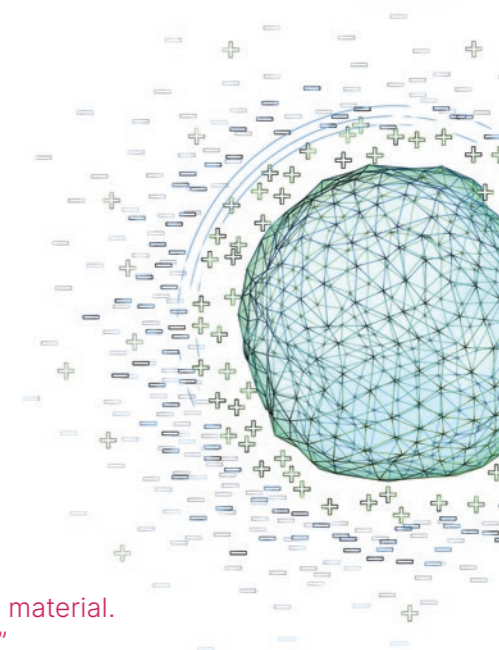
Specifically, the particles' total surface area determines how effectively gas or liquid can flow through and how quickly the MOF can trap or release molecules or ions. So, for the best MOF product, particles must remain a **constant size, intact** and not **agglomerated**. The powder's flow and packing also have to be controlled. And for this, you need to understand the **size and shape** of your MOF particles.



## Nanoparticles in liquid: hydrodynamic performance

One way to measure how effectively a MOF can harvest or release ions is to measure the **elemental concentration** of ions in residual concentrations in its carrier liquid.

When formulated as **nanoparticles** in suspension, MOFs can carry a whole range of active and targeted pharmaceutical drugs. That's because **surface modification** enables these nanoparticles to access cells. Understanding the **hydrodynamic behavior** of these surface-activated MOF nanoparticles is essential for all stages of the process.



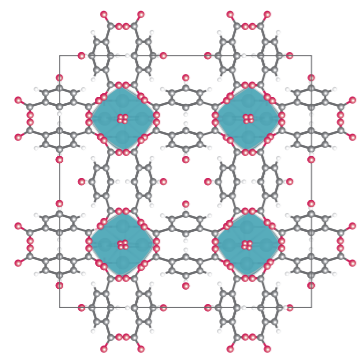
"These are really exciting times – I have never seen such capabilities in a material. The selectivity, the robustness, the scale of response: it's mind-blowing."

Dr Ana Morfesis –  
Principal Scientist, Malvern Panalytical

# Crystal structure

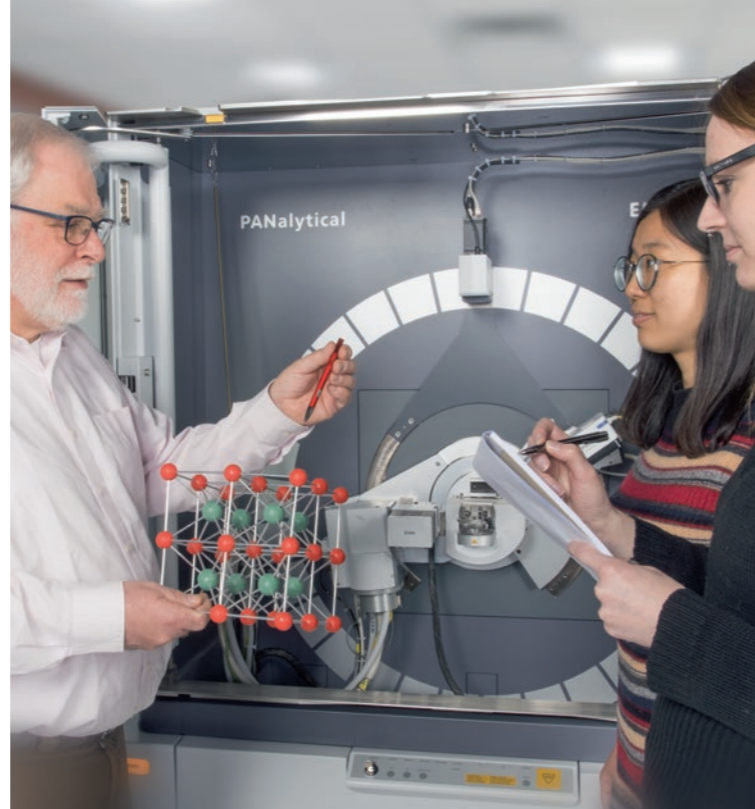
## Why is crystal structure analysis important?

To understand your MOF structures in detail and check that your target MOF has been synthesized correctly, you need to be able to identify any changes to the crystal structure – quickly and easily.



## What causes structural changes?

Structural variations can happen when raw materials are changed or if the material synthesis route is different. Long-term cycling or environmental conditions may also cause degradation in the lattice. Measurements can show whether your MOF's pores are empty or partially filled.

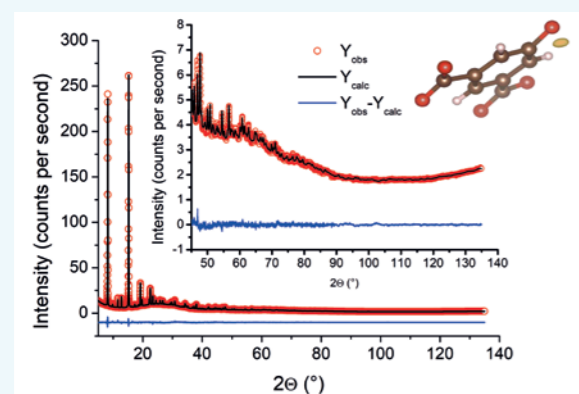


## X-ray diffraction: A certified solution

X-ray powder diffraction is a well-established way to measure and verify crystal structure. Our Empyrean and Aeris X-ray diffractometers can provide the highest-quality powder diffraction data. Count on their resolution and sensitivity to pin down even the finest structural details.

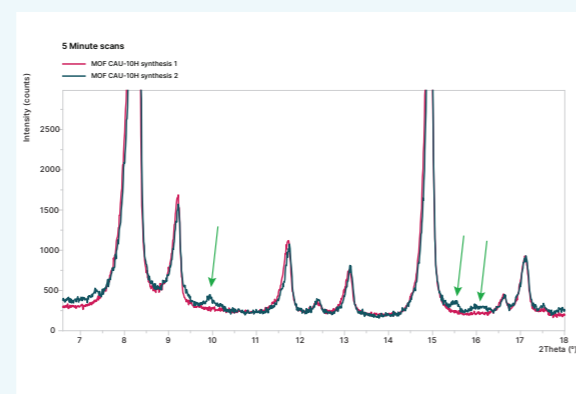
## Empyrean: Structural solutions

The excellent fit of this high resolution and low-background transmission data demonstrates that great data quality leads to an accurately determined structural model. The result even locates the hydrogen atom of the MOF's phenolic group.



## Aeris: Rapid fingerprinting

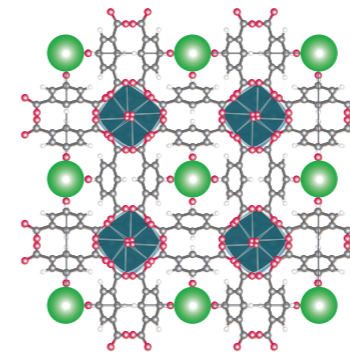
The powder diffraction scan below was obtained in a 5-minute scan on the Aeris diffractometer. It shows two MOF samples with nominally the same constituents, synthesized by two different routes. The green arrows show extra reflections, indicating that Sample 2 is not pure enough.



# Binding affinity

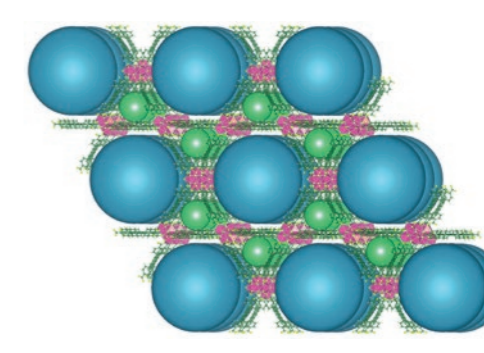
## Why analyze crystal structure during occupancy cycling?

By studying the MOF's crystal structure during occupancy cycling, you can understand how morphology, packing, and use environment affect its performance – and design your use environment accordingly.



## Binding affinity at different sites

What are the benefits of understanding binding affinity at specific binding sites in your MOF pores? Better MOF structures for specific adsorption species, and long-term robustness in various use environments.



## Integrated non-ambient measurements

Our X-ray diffractometers let you study the crystal structure changes in your MOFs – whatever the operating conditions. Choose from several integrated non-ambient chambers and reaction cells, and rapidly identify points of change with HighScore software, incorporating 2D and 3D graphics and cluster analysis.

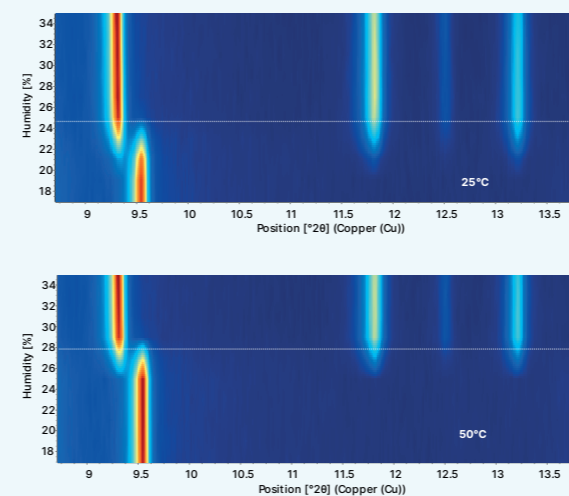
## A complete thermodynamic profile

By measuring heat transfer during binding, you can accurately determine binding constants ( $K_D$ ), reaction stoichiometry (N), enthalpy ( $\Delta H$ ), and entropy ( $\Delta S$ ). Our ultrasensitive Microcal PEAQ-ITC microcalorimeter makes measuring these thermodynamic changes easy – giving you a complete thermodynamic profile of adsorption and desorption.

## Empyrean: Capture and release of water

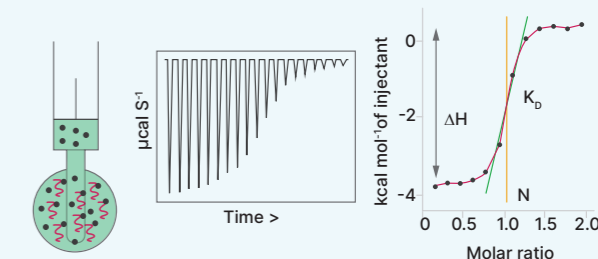
Because the MOF CAU-10-H can store water molecules ( $H_2O$ ), it has potential applications in water harvesting.

These isoline plots show variable *in situ* relative humidity data collected at two temperatures, demonstrating the effect of temperature on the adsorption and desorption of water within the sample.



## Basics of ITC experiment

Universal technique based on heat detection



Integration of heats are used to extract affinity (KD), stoichiometry (N) and binding enthalpy ( $\Delta H$ ) using appropriate binding model

## ITC: Example report

The MOF NU-1000 can effectively extract uremic toxins<sup>1</sup> – both when free in aqueous media and when bound to human serum albumin (HSA). These authors measured the interactions of the toxins at different binding sites in the MOF using ITC, revealing both entropically and enthalpically driven adsorption at two different pore sites.

<sup>1</sup> Cell Reports Physical Science, Volume 1, Issue 1, 22 January 2020, 100006: <https://doi.org/10.1016/j.xcrp.2019.100006>

# Particle size

## More optimized product formulation

MOFs are often produced in powder form, as they precipitate from a reaction in a solvent. The physical form must remain reproducible and stable. This means the powder particles must stay a constant size, intact, and not agglomerated throughout their use. Fast particle size distribution measurements can help you quickly understand these process parameters to deliver optimum product size.

## The right size for the right application

Each MOF has an optimum particle size – and size distribution – for its intended application:

- Bulk MOF powder particles designed to interact with gases can be 1 to 100s of microns large
- For batteries, smaller particles (20 nm to 1 micron) reduce ionic diffusion lengths. They must be distributed evenly in thin multi-layered structures.
- For drug delivery in biological systems, MOF nanocrystals (<100nm) in suspension are used. They can be small enough even to penetrate the blood brain barrier.

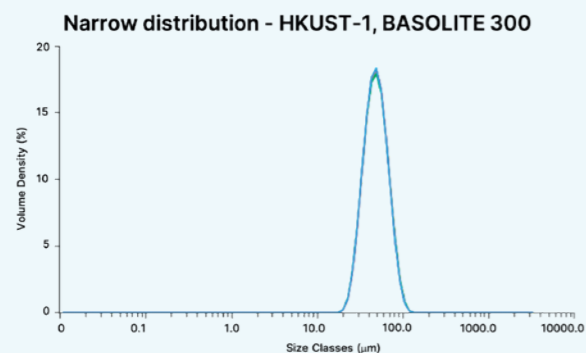
## Smarter particle sizing: What's the best option?

But which particle sizing method is right for you? Each instrument in the table has different benefits – and they can also measure additional attributes. Ultimately, your research, development, or production requirements will determine the best instrument and method. And with this many options, we're sure to find your match!

Partical Size Range	0.1nm	1nm	10nm	100nm	1µm	10µm	100µm	1mm	10mm
Dynamic Light Scattering <1nm to >1µm			Zetasizer						
Nanoparticle Tracking Analysis <100nm to >1mm			NanoSight						
Laser Diffraction <100nm to >3.5mm			Mastersizer						
Automated Imaging <1µm to >1.3mm					Morphology				
SAXS <1nm to >1µm		Empyrean							

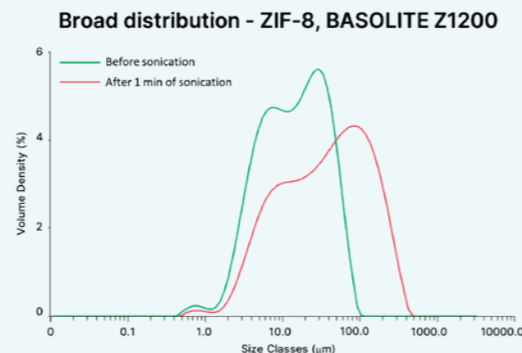
## Mastersizer 3000: Particle size distribution

This figure shows the results of a laser diffraction measurement using the Mastersizer 3000. The sample data from 5 measurements reveal a narrow and reproducible particle size distribution peaking at 48µm.



## Mastersizer 3000: Particle agglomeration

This figure shows the results of a laser diffraction measurement using the Mastersizer 3000. The sample was measured in a dispersion revealing a broad distribution of fine and agglomerated particles. The mean particle size reduces after sonication, showing a reduction in agglomeration.



# Particle shape

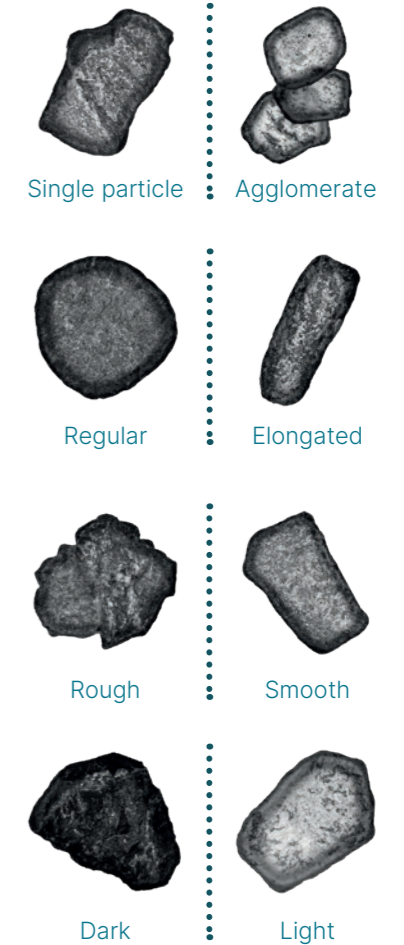
## The importance of particle shape

In powder, particle shape affects packing, flow, and the chance of agglomeration. In dispersion, particle shape can influence the stability of a formulation. Imaging is a great way to measure particle size and shape to understand or monitor performance.

## Tracking particle degradation

Repeated cycling, along with the influx and outflux of ions and molecules, can cause particle degradation or incomplete emptying of the MOF, leaving a residual occupancy that increases over time.

You can track this particle degradation using imaging with Raman spectroscopy chemical analysis, giving vital clues about your products' shelf-life. Raman can help you see what proportion of your particles have an adequate loading of, for instance, active pharmaceutical ingredient (API).



## Automated imaging for advanced particle characterization

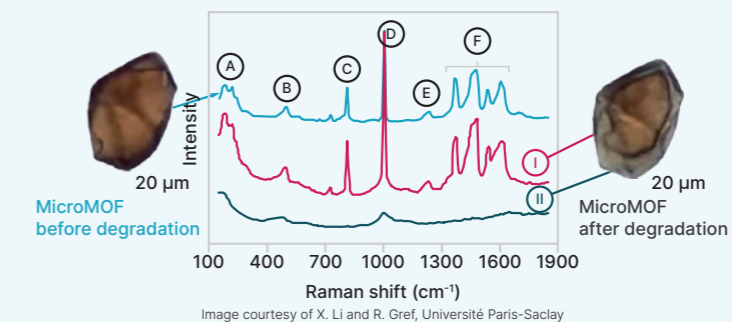
Automated imaging gives you both instant and long-term information on the physical robustness and performance of your individual MOF particles. With Morphologi 4, you can scan 10,000 – 500,000 particles and get a full set of size and shape results for each particle. All in just 30 minutes!

What's more, you can opt for your results to show only what's most relevant to you. And, with Morphologi 4-ID, you can return to any individual particle for more detailed analysis, including Raman chemical profiling.

## Morphologi 4-ID: Raman spectroscopy

This figure shows a published result of some imaging and Raman results demonstrating the degradation of biocompatible MOF crystals in a biological medium. The MOF crystals are being studied for their potential in drug delivery applications. It is important that the particles are dissolved in the body during drug release. The authors studied new particles before degradation

and then particles in biological medium (I) which, over time, developed denuded regions around the edges (II). Raman spectroscopy revealed that the degraded region was losing the organic ligand in the MOF<sup>2</sup>. Further studies showed the eventual collapse of the inorganic matrix as it dissolved in the biological medium.

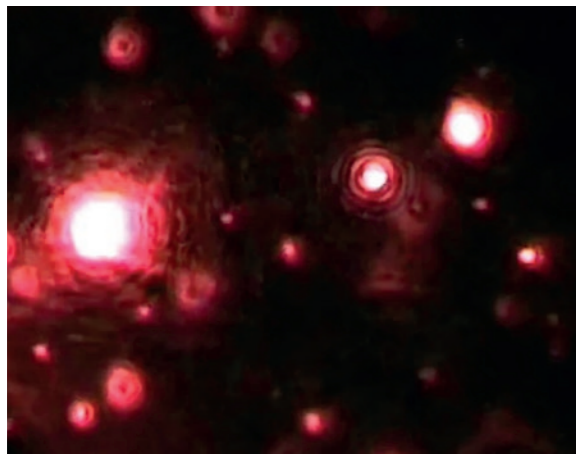


<sup>2</sup> Int. J. Pharm. 2017531(2): 424-432 <https://doi.org/10.1016/j.ijpharm.2017.05.056>

# Nanoparticles in liquids

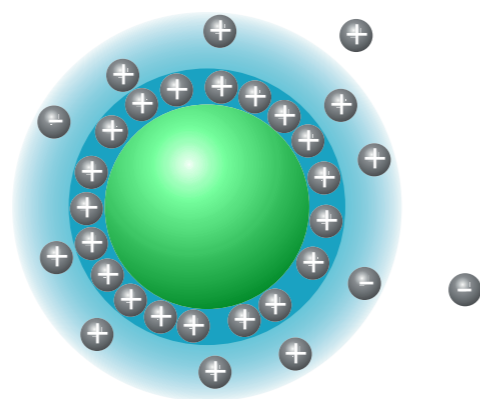
## Why study nanoparticle stability?

Understanding the size and surface charge of MOF nanoparticles in solution is critical to controlling mobility and aggregation. Controlling particle size in bioactive drugs is particularly important, as it ensures that drug carriers can pass through biological barriers and reach their intended destination.



## Modifying nanoparticle surfaces

Modifying the surface of MOF nanoparticles can help optimize their behavior in liquids. In drug delivery, for instance, the right surface enables access to cells. In water treatment, adjusting surface charge can optimize the absorption of toxic pollutants<sup>3</sup>.

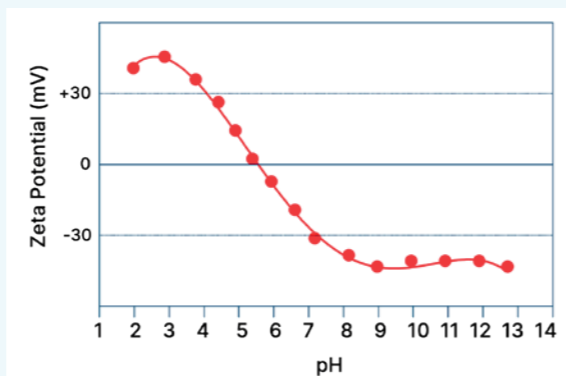


## Light scattering for in-depth nanoparticle analysis

You can measure the size of nanoparticles in liquids using Dynamic Light Scattering (DLS) on the Zetasizer Advance. Or, to analyze particle drift for Zeta potential, use Electrophoretic Light Scattering (ELS). And Nanosight NS300 complements these measurements, providing real-time imaging of particle movement, with tracking analysis for both sizing and concentration.

## Example applications

- For battery and electronic applications, MOF particles must remain structurally stable under environmental changes. Zeta potential is a useful measure of this stability<sup>4,5</sup> when pH values change.
- MOFs can be used as catalysts to degrade pharmaceuticals and cosmetics and prevent them from contaminating waste. Dynamic light scattering can qualify and control the synthesis of suitable catalytic MOF particles<sup>6</sup>.



<sup>3</sup> Journal of Environmental Sciences, Volume 80, June 2019, Pages 197-207: <https://doi.org/10.1016/j.jes.2018.12.013>

<sup>4</sup> Front. Chem., 04 August 2020: <https://doi.org/10.3389/fchem.2020.00617>

<sup>5</sup> Nano Energy, Volume 65, November 2019, 104032: <https://doi.org/10.1016/j.nanoen.2019.104032>

<sup>6</sup> Chemosphere, Volume 196, April 2018, Pages 105-114: <https://doi.org/10.1016/j.chemosphere.2017.12.164>

# Nanoparticles in liquids

## Know your purity levels

Knowing the purity of your starting materials and intermediate precursors can be important in developing your MOF synthesis and product processing.



## Monitoring impurities from catalyst residuals

MOFs can be used as catalysts – for example, in the production of carbon nanotubes.<sup>7</sup> Catalyst residues can cause unwanted impurities in nanomaterials, affecting their properties and the final application. XRF analysis can monitor these impurities.



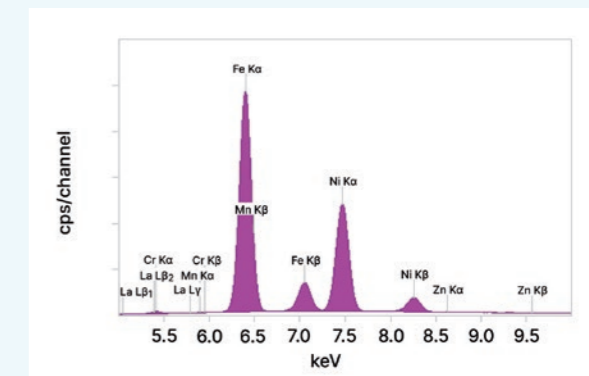
## Assessing ionic uptake and expulsion

You can measure how effectively your MOF particles trap and release ions from solutions by measuring ion concentrations – for example in filtered solutions, or (for batteries) in separators and electrolytes.



## Elemental insights with X-ray fluorescence

Our X-ray fluorescence (XRF) instruments can help determine your sample's chemical composition – whether it's solid, liquid, slurry, or loose powder. You can also use our sample preparation equipment to make peroxide solutions for ICP elemental analysis.



XRF spectrum of carbon nanotube sample showing residual impurities

<sup>7</sup> Reference: Inorg. Chem. 2019, 58, 5, 3227–3236 : <https://pubs.acs.org/doi/10.1021/acs.inorgchem.8b03318>

# Crystal structure



## Research

You need a MOF crystal that fulfils your application requirements – without fail:

- A robust crystal structure that adsorbs the correct molecules or ions
- The right adsorption capacity – obtained through the right physical morphology
- Understanding of the process, with matching fluid flow and composition (liquid or gas)
- Characterization and fine-tuning of adsorption and release parameters



## Development

You need an upscaled process that achieves:

- The correct crystal structure
- The required physical morphology
- A robustly contained material
- Reproducible adsorption and release parameters



## Production

And you need product quality, every time:

- Testing for chemical impurities
- Fast quality assurance of crystal structure
- Fast quality assurance of sample morphology
- Effective in-line measurements
- Automated measurement and analysis

Let us help you

# Our solutions

## Crystal structure

Let our **Empyrean** and **Aeris** X-ray diffractometers take your crystal structure analysis to the next level.



## Binding affinity

The ultrasensitive microcalorimetry of our **Microcal PEAQ-ITC** gives you direct measurements of binding affinity ( $K_D$ ), stoichiometry ( $n$ ), enthalpy ( $\Delta H$ ), and entropy ( $\Delta S$ ).



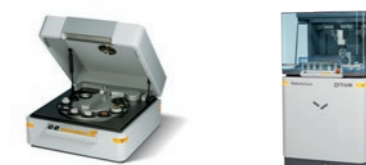
## Particle size and shape

Use our **Mastersizer 3000** for particle size distribution with laser diffraction, **Morphologi 4** for particle size and shape analysis, and **Morphologi 4-ID** to analyze particulate chemistry with automated imaging and morphologically directed Raman spectroscopy (MDRS).



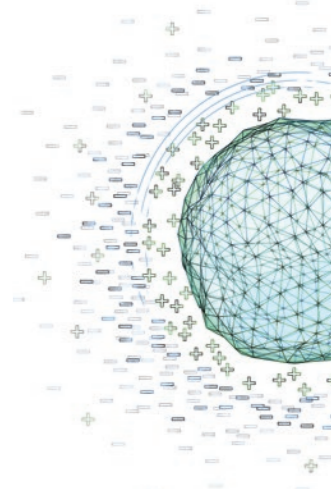
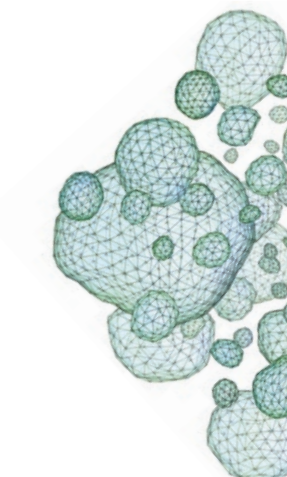
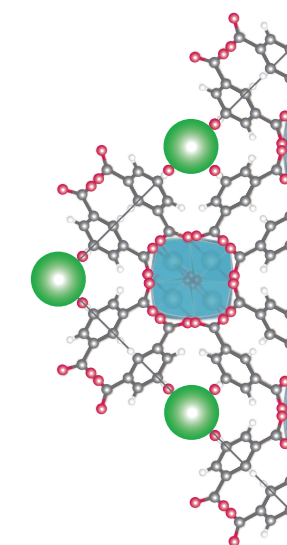
## Elemental analysis

The X-ray fluorescence (XRF) spectroscopy of **Epsilon** and **Zetium** makes elemental analysis of MOFs and their fluid environments easier than ever.



## Nanoparticles

Need to analyze MOF nanoparticles in suspension? **Zetasizer Advance** has you covered with dynamic light scattering (DLS) and electrophoretic light scattering (ELS). For particle size and concentration analysis, try **NanoSight NS300** using Nano Tracking Analysis (NTA).



These are just some of the measurements and analyses you can perform with Empyrean – get in touch to discover the full range of possibilities.

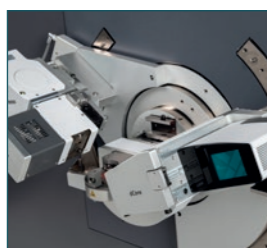


## Empyrean

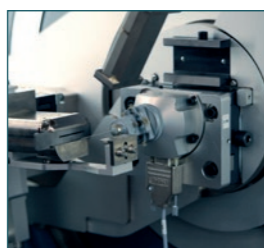
### Empyrean is for you if you want:

- The world's most accurate **high-resolution goniometer** for state-of-the-art diffraction and scattering experiments.
- A full choice of Universal PreFIX **optics and stages** for fast, accurate component exchange.
- **iCore and dCore** optics for multiple measurements with batch automation.
- 0D, 1D and 2D **solid-state detectors** – fully integrated, low-noise, and high-resolution.
- **Advanced configurations** for specialized measurements (capillary transmission, SAXS, GISAXS, HR-XRD, PDF and CT).
- The widest range of **sample holders and sample stages** for manual and automated sample exchange.
- Fully integrated **non-ambient sample chambers**.
- Bespoke **design and integration** for sample mounting and *in situ* experiments.
- **HighScore and HighScore Plus**: user-friendly software for phase analysis, Rietveld refinement, fingerprinting, cluster analysis, non-ambient studies – and more!
- Fully **automated analysis routines** – for individual measurements or large data sets.

### And for MOFS take a look at:



Low-background, high-resolution **iCore and dCore** detectors reveal even the smallest peaks in automated measurements.



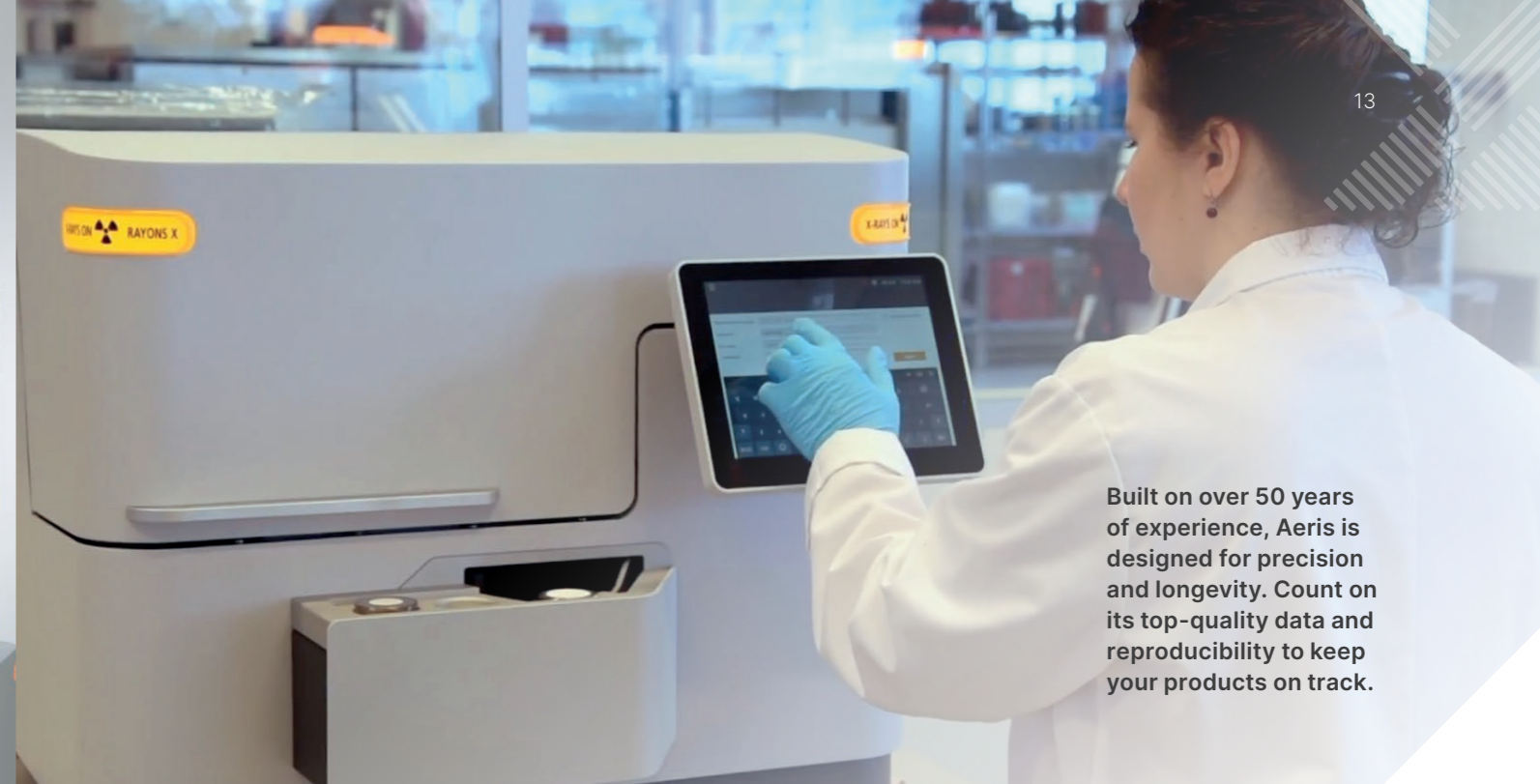
Your choice of **optics and configurations** to minimize texture – and the best low-angle peak shapes for Rietveld fitting.



Integrated **non-ambient stages** for controlled, synchronized measurements – every time.



All the **sample holder types** you could want – and up to 48 samples loaded for automated batch analysis.



Built on over 50 years of experience, Aeris is designed for precision and longevity. Count on its top-quality data and reproducibility to keep your products on track.

## Aeris

### Aeris is for you if you want:

- An easy-to-operate instrument combining **one-step sample loading** with **simple push-button operation**.
- The quality of a floor-standing diffractometer in **compact** form.
- Unrivaled **low detection limits** for small powder diffraction peaks and material variations.
- High-resolution 1D and 2D **solid-state detectors**.
- **Powder diffraction** measurements in reflection, transmission and grazing incidence.
- Safe, precise **external sample loading** anytime. No need to disturb ongoing measurements!
- The widest range of **sample holders and sample stages** – with various options for manual and automated sample exchange.
- An uncoupled **theta-2theta goniometer** to keep your samples safe and your instrument clean.
- Rapid instant results with **RoboRiet automated analysis**. Wave goodbye to waiting times!
- Integrated **non-ambient temperature chamber**.
- **Full integration** into an automated lab or processing plant.

### And for MOFS take a look at:



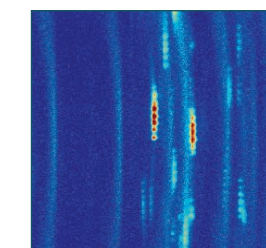
With **push-button powder diffraction**, you don't need to be an expert to manage quality control.



**Automation** streamlines your processes and cuts your production overheads. No more unnecessary expenses!



**Grazing-incidence diffraction** improves detection limits: Get ready to analyze even thinner films and coatings!



**2D capability** lets you check your crystallites' processing quality – large or small, random or preferred orientation.

A rugged, easy-to-use design with intelligent software and high sensitivity gives you all the thermodynamic parameters you'll need.



## MicroCal PEAQ-ITC

### MicroCal PEAQ-ITC is for you if you want:

- An **easy-to-operate**, ergonomically designed instrument. Say hello to a smoother experience!
- **High-quality, reproducible results** with minimal experimental artifacts.
- High **signal-to-noise** – for greater confidence in your data's quality and relevance.
- An integrated **washing module** with quick disconnect fittings and liquid sensors.
- **Full cleaning** of both sample cell and titration syringe – even at elevated temperatures.
- Robust injection syringe **filling and cleaning** system.
- **Hastelloy® sample cell**: suitable for a wide range of compounds.
- **Instrument control software** with guided workflows, step-by-step tutorials, and maintenance alerts.
- Analysis software with **experimental design simulation, batch evaluation** of large data sets, **automated assessment** of data quality, and a **streamlined user interface**.
- **Presentation-quality graphics**: sharing your findings just got easier.

### And for MOFS take a look at:



Robust injection syringe **filling and cleaning** system.



Unique **pipette and titration syringe**.



Integrated **washing module** with liquid sensors.



MicroCal PEAQ-ITC **Automated**.

The Mastersizer 3000's versatility means you can create the ideal setup for consistent results – again and again.



## Mastersizer 3000

### Mastersizer 3000 is for you if you want:

- Robust, rapid **particle size measurements** spanning 10nm – 3500µm in a single optical path.
- Intuitive software with **built-in expertise** – for faster analysis than ever!
- Measurement of **dry powders, suspensions and emulsions**, with multiple plug-and-play dispersion units.
- Easy, reproducible operating procedures via the **Measurement Manager** interface.
- **Live feedback** on dispersion and sonication, for optimum particle conditions.
- A full **real-time view** of instrument controls, live data, and trends.
- **Automatic cleaning** and **easy switching** from liquid to dry dispersions. No additional stress necessary!
- Flexible **dry dispersion** – for robust or fragile materials, and whatever the sample volume.
- **Fully editable reporting** – so the benefits continue after analysis.

### And for MOFS take a look at:



**Aero funnel sample feeder** for Aero dry powder dispersion units. Enjoy rapid quality control for your bulk powder samples!



**Hydro EV** – a unique dip-in wet sample dispersion unit, used with standard laboratory glassware for a wide range of dispersant volumes.



**Hydro Insight** - real-time images of liquid particle dispersions and individual particles, with quantitative data on particle shape.



Ask us about **in-line** and **at-line** particle sizers and **bespoke auto-sampling solutions!**



Automated particle imaging lets you discover the physical properties of each one of your MOF particles – now and over time.

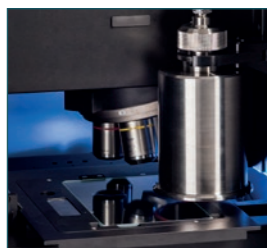


## Morphologi 4-ID

### Morphologi 4 and Morphologi 4-ID are for you if you want:

- To scan **10,000 – 500,000** particles from **0.5 to 1300µm** – with full results of up to 20 parameters in **30 minutes**.
- Results that present only what's **relevant** to you.
- To analyze individual particles in more **detail** with Raman chemical profiling.
- A **high-resolution microscope** for the highest-quality particle images.
- Automated **'sharp-edge' analysis** to detect low-contrast particles.
- An integrated **dry powder dispersion unit** for reproducible sample dispersion.
- Advanced **data exploration tools** – for maximum sample knowledge.

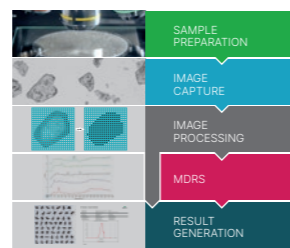
### And for MOFs take a look at:



Integrated **dry powder dispersion unit**, making reproducible sample preparation easy.



**Wet dispersion cell** – so you can analyze 2-6 mL of your sample.



**Imaging workflow and SOP control** – from sample dispersion to data analysis.



A great choice of **sample holders**, including **slide holders** for traditional microscopy.

With the Zetasizer Advance's wealth of measurement options, you know you're getting the best results for your unique samples.



## Zetasizer Advance

### Zetasizer Advance is for you if you want:

- A **wider range** of sample concentration and particle sizes – thanks to **Non-Invasive Back Scatter (NIBS) technology**.
- The **highest-resolution sizing data** from combined **Multi-Angle Dynamic Light Scattering (MADLS®) technology**.
- Accurate size measurements of **1nm-10µm particles**.
- A confident understanding of your sample's **stability** and **aggregation** potential.
- **Constant current mode** – say goodbye to data errors from electrode polarization.
- **Repeatability** with minimal sample preparation, thanks to the **Adaptive Correlation algorithm**.
- A **sample-centric workflow** to automate even the most complex analyses.
- **Instant automatic feedback** and **actionable advice**.

### And for MOFs take a look at:



The **ZSU1002 cell** for sample volumes as low as 3µL and particles up to 10µm. Say hello to a wider dynamic concentration range!



Our **completely disposable cell** – no more cross-contamination between measurements!



The **MPT-3 titrator** automates your titrations for Zeta or size vs. pH.



Our **dip cell** lets you measure zeta potential in both **aqueous and non-aqueous** dispersants.

With Nanosight NS300, you can understand your nanoparticle suspensions in more depth than ever.



## NanoSight

### NanoSight NS300 is for you if you want:

- Particle size images and measurements from **10-1 $\mu$ m**, at up to **1  $\times$  10<sup>9</sup> particles/mL**.
- **Particle-by-particle, high-resolution** particle size data, plus **concentration measurements** for colloidal suspensions or nanoparticle solutions.
- **Quantification** of primary sample components, aggregates, and fluorescently tagged particles.
- High-resolution information on **monomodal and polydisperse systems** – across all kinds of materials.

### And for MOFS take a look at:



Comprehensive **real-time data viewing** and analysis.



NanoSight **syringe pump** lets new particles flow continuously into the sample chamber.



NanoSight **Sample Assistant**: enjoy precise, reproducible, automated sample loading from a 96-well plate.



**Low-volume flow cell** – delivering data from as little as **250 $\mu$ L**.

Whatever your requirements, Epsilon has something for you – from the simplest monitoring to the most demanding measurements.



## Epsilon range

### Epsilon 1 or Epsilon 4 is for you if you want:

- **Non-destructive detection** of elements from C-Am, with detection limits from 1 ppm – 100%.
- A **unique dynamic range** and **high elemental resolution** – whatever the sample matrix.
- **Easy management** of powders, solids, films, filters, and solutions.
- **Fast, accurate** analysis, at-line or in-line.
- Advanced **data treatment algorithms**.
- Optional **standardless analysis** with Omnic software.
- A pass-fail **fingerprint program** for quick quality control.
- A **full workflow package**: from auto-sampling and sample preparation to measurement and analysis.
- Automation options.

### And for MOFS take a look at:



Epsilon 4's **10-position carousel, 50-100 position sample changer**, and unlimited **belt automation**. The possibilities are endless!



**Xflow automated elemental analysis** of liquids – whether it's on-line or at-line.



**Fusion** sample preparation for XRF and ICP analysis.



Choose from a wide range of our **certified reference materials (CRMs)**. We'll even create bespoke CRMs for your unique applications.



## Why choose us?

**When you make the invisible visible, the impossible is possible.**

Our analytical systems and services help our customers to create a better world. Through chemical, physical and structural analysis of materials, they 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.

With over 2200 employees, we serve the world, and we are part of Spectris plc, the world-leading precision measurements group.

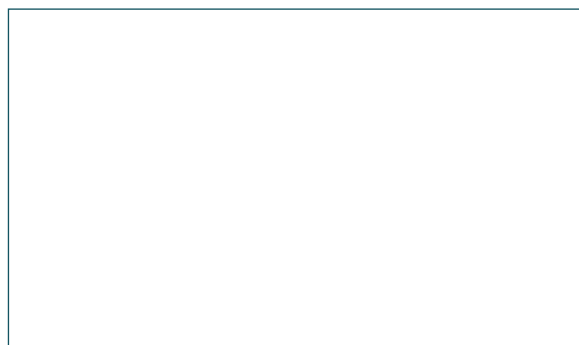
**Malvern Analytical. We're BIG on small™**

## Service & Support

Malvern Analytical 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 Analytical

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.malvernanalytical.com](http://www.malvernanalytical.com)