



Advanced Materials characterization toolbox

product – measurement – technology – applications – features



- Laser diffraction
- Light scattering
- Imaging
- Near Infra-red
- X-ray fluorescence
- X-ray diffraction
- Sample preparation

Particle size distribution, shape and concentration



Product	Mastersizer	Hydro Insight	Spraytec	Insittec
What does it measure?	Particle size	Particle size, shape, appearance	Particle size	Particle size
Technologies used	Laser Diffraction	Dynamic imaging	Laser Diffraction	Laser Diffraction
What is it used for?	<ul style="list-style-type: none"> Measuring the size distribution of suspensions, emulsions and dry powders from 10nm to 3500µm Controlling powder properties such as wettability, bulk density, powder flow and solubility Optimizing suspension and emulsion rheology 	<ul style="list-style-type: none"> Sits alongside Mastersizer 3000 collecting images of particles Providing quantitative information about particle size and shape covering the size range from 1µm to 800µm Helps in method design and simplifies troubleshooting 	<ul style="list-style-type: none"> Measuring the size distribution of sprays and aerosols from 0.1µm to 2000µm Defining the deposition pattern and bioavailability of drug materials delivered using pump sprays and inhalers Understanding the environmental impact of spraying in coating Consumer or agrochemical applications Resolve the fluctuations in droplet size during the rapid firing of automotive fuel injector systems 	<ul style="list-style-type: none"> Online continuous particle size analysis needed for efficient, cost-effective monitoring and control of industrial processes Suitable for the widest variety of process streams from dry powders to hot sticky slurries, sprays and emulsions, whether milligrams of material or hundreds of tonnes per hour Insittec systems measure particles in the size range 0.1 micron to 2.5 mm
What is special about this product?	<ul style="list-style-type: none"> World's most popular particle sizing instrument Class-leading particle sizing performance in a compact footprint Intuitive software with built-in expertise to ease your workload Flexible reporting to display your data the way you want Fast smart swap between wet and dry units Rapid and effective wet dispersion for dispersions and emulsions Fast, reliable particle size measurement of fragile and cohesive dry powders Mastersizer know-how throughout 	<ul style="list-style-type: none"> High-speed, high resolution dynamic imaging technology 127-frames-per-second digital camera with up to 5 megapixel resolution Imaging of individual particles and liquid particle dispersions Thumbnail images saved for post-run viewing Quantitative data on particle size and shape, including particle width and elongation data 	<ul style="list-style-type: none"> Measurement across a wide size range without requiring constant optics changes Resolve rapid changes in droplet size over time, by measuring up to 10,000 measurements a second Deliver accurate, concentration-independent results using a patented multiple scattering analysis Characterize wide spray plumes without risking optical contamination Reveal dynamic changes in spray particle size through the unique size history analysis software 	<ul style="list-style-type: none"> Industrially robust technologically proven measures particles in the size range 0.1 to 2500 µm delivers real-time monitoring and control Base model hardware manufactured to GAMP5 standards and compatible with CIP/SIP requirements to meet specific manufacturing specifications Easy to use software and fully automated operation to minimize training requirements and release operator time Integration with existing control platforms to simplify development of automated control protocols High reliability of >95% with little downtime, minimal maintenance and maximum ROI



Particle size, shape, chemistry, concentration, molecular weight, Formulation stability and intrinsic viscosity



Product	Morphologi 4/4-ID	Zetasizer	Nanosight	Product
What does it measure?	Chemical identification, Particle shape, Particle size	Zeta potential, Particle size, Particle concentration, Molecular size, Molecular weight, Protein mobility	Particle concentration, Particle size	What does it measure?
Technologies used	Image Analysis Raman Spectroscopy	Dynamic Light Scattering, Electrophoretic Light Scattering, Static Light Scattering	Nanoparticle Tracking Analysis	Technologies used
What is it used for?	<ul style="list-style-type: none"> Size measurement of non-spherical particles such as needle shaped crystals from 0.5µm to 1000µm Measurement of shape differences where particle size alone does not allow differentiation Detection and enumeration of agglomerates, oversized particles and contaminant particles Automation of manual methods such as microscopy Physical characterization of individual components within a mixture Cross-validation of particle size measurements such as laser diffraction 	<ul style="list-style-type: none"> Measuring the size of colloids, nanoparticles and molecules in solution from 0.6nm to 15µm Determining the molecular weight of polymers and proteins Optimizing and predicting dispersion stability of colloids and biomolecules Monitoring molecular aggregation and particle flocculation processes 	<ul style="list-style-type: none"> Detection and visualization of nanoparticle populations on a particle-by-particle basis from 10nm to 1µm Measurement of particle size and particle concentration Tracking of aggregation and dissolution processes in real time Applications include nanotoxicology, biomarker detection and drug delivery research Fluorescence mode to provide differentiation of labeled or naturally fluorescing particles 	What is it used for?
What is special about this product?	<ul style="list-style-type: none"> Measures particle size, shape and chemical identity in one platform Integrated dry powder dispersion unit automates sample preparation for consistent measurements Versatile sample presentation accessories for measuring suspended and filtered samples Simple SOP operation from sample dispersion through to size, shape and chemical analysis Automatic selection, targeting and chemical classification of thousands of individual particles Powerful and intuitive software interface making both visual and statistical interpretation of your data easier than ever 	<ul style="list-style-type: none"> Technologies include NIBS, M3-PALS, DLS, ELS Simplicity of operation means minimal training and robust results High sensitivity for nanoparticles, proteins, and macromolecules High optical quality and temperature control ensures accuracy and repeatability MPT-3 Autotitrator option Adaptive Correlation to enhance repeatability MADLS for angle independent size analysis with improved precision and resolution. Also used for accurate particle concentration measurements M3-PAL and constant current mode for Zeta potential to reduce errors 	<ul style="list-style-type: none"> Simultaneous measurement of multiple characteristics Visual validation of results gives extra confidence User friendly software with easy set-up of SOPs for routine use Minimal sample preparation Automated multiple sample analysis when used with a syringe pump or autosampler Minimal consumables reduce running costs on a day-to-day basis High resolution particle sizing technique, ideal for polydisperse systems 	What is special about this product?



Focus on semiconductors



Product	Automated X'Pert ³ MRD (XL)	2830ZT Wafer analyzer
What does it measure?	Wafer orientation and quality Thin film strain, composition Multilayer thickness Epilayer integrity	Elemental composition layer thickness
Technologies used	High-resolution X-ray diffraction (HR-XRD)	X-ray fluorescence
What is it used for?	<ul style="list-style-type: none"> QC and R&D of thin film layers and substrates Semiconductor single-crystal epitaxial layers Multicrystalline electro ceramics PPM-resolution lattice parameter and lattice parameter strain measurements Monolayer precision thickness measurements epitaxial layers Phase purity and orientation distribution in sputter deposited films 	<ul style="list-style-type: none"> Elemental QC of thin films (including Si and electro ceramics) Alloy composition and film thickness High precision high throughput measurement Elemental composition (for Elements heavier than Be)
What is special about this product?	<ul style="list-style-type: none"> Can be used in both lab and fab versions (upgradable from lab version) Can access coplanar, skew and in-plane geometry 200mm wafer mapping capability GEN series compatible waferloading and automation Clean room ready AMASS analysis software, with 'any orientation - any spacegroup' capability The only platform that has a 180 deg Chi cradle The only platform that can reach extremely high incident angles on 8" wafers Motorized divergence and antiscatter slits for monochromators can automatically vary the X-ray beam height Microbeam collimators can reduce the X-ray beam size to as small as 100 x 100 microns 4 detection modes fully automated 	<ul style="list-style-type: none"> This tool can measure up to 300mm wafer size up to 25 wafers per hour Several channels of the same element can be equipped to boost the element sensitivity



Molecular and crystalline structure, composition, crystalline quality

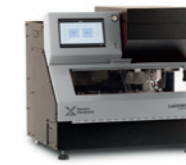


Product	OMNISEC	Empyrean	X'Pert ³ MRD (XL)	Aeris	FieldSpec - LabSpec	Product
What does it measure?	Absolute molecular weight (Mw), molecular size, intrinsic viscosity (IV), branching and other polymer parameters	Crystalline phase crystallographic structure materials microstructure Thin film structure Epitaxial strain	Thin film structure, epitaxial strain, materials microstructure	Crystalline phase	Materials properties, molecular structure, multiple constituents	What does it measure?
Technologies used	Gel Permeation Chromatography, Size Exclusion Chromatography, Static Light Scattering	X-ray diffraction, reflectometry, SAXS, CT	X-ray diffraction, reflectometry	X-ray diffraction, reflectometry	Near-infrared spectroscopy (NIR)	Technologies used
What is it used for?	<ul style="list-style-type: none"> Determining absolute and relative Mw and Mw distribution of synthetic and natural polymers Measuring IV to investigate molecular structure and branching Assessing polymer degradation Controlling mechanical properties of plastics Characterizing and quantifying components in blends Measuring sample concentration e.g., polymer additives in fuel 	<ul style="list-style-type: none"> Phase Identification Phase quantification Thin Film characterization Epitaxial layer analysis Bulk material characterization Non-Ambient phase change Battery in-operando study In-situ phase change Stress and residual strain Crystallographic texture 3D flaw detection Nanoparticle analysis 	<ul style="list-style-type: none"> Measuring Epitaxial film composition and thickness from rocking curves and reciprocal space maps Analyzing full-sized semiconductor wafers Speciality thin films, strain, preferred orientation, stress, texture, requiring large 4-circle goniometer All standard methods to ultra-high resolution 	<ul style="list-style-type: none"> Phase identification and quantification in reflection, transmission or grazing incidence Thin film analysis Stress measurement 	<ul style="list-style-type: none"> Non-destructive analysis of a wide range of materials in the laboratory or in the field Quality control of any type of product and requires no sample preparation Uses statistical analysis to correlate spectra to any number of materials properties 	What is it used for?
What is special about this product?	<ul style="list-style-type: none"> OMNISEC REVEAL – a fully integrated temperature controlled multidetector module with minimized inter-detector volumes for maximum stability and sensitivity Market leading light scattering detector for absolute Mw High sensitivity RI detector, for low concentrations Self-balancing viscometer for accuracy in IV Wide UV/VIS wavelength range for all applications Maximum sensitivity detectors with MALS capability Intuitive, advanced and automated software OMNISEC RESOLVE - an integrated pump, degasser, autosampler and column oven in a single unit 	<ul style="list-style-type: none"> Most versatile and productive XRD system with the highest data and product quality on the market Highest angular resolution of any laboratory powder diffractometer and delivers data closest to synchrotron quality Widest range of non-ambient and in-situ environments All relevant diffraction geometries (reflection, transmission, capillary, microdiffraction, Debye-Scherrer) with batch automation Wide selection of components to match every customer Cost-effective options for SAXS and CT configurations Customizable and special solutions Full selection of analysis software packages including HighScore+ 	<ul style="list-style-type: none"> New high resolution goniometer using Heidenhain encoders Improved accuracy and faster positioning Rapid tool-free exchange of tube position from point to line focus Pneumatic shutters and beam attenuators Longer lifetime of incident beam components with CRISP* including a lead-free tube tower. Second generation of PreFIX for even more accurate optics positioning Full selection of analysis software packages including AMASS 	<ul style="list-style-type: none"> Compact X-ray diffractometer External loading intuitive operation, accessible to non-experts Touch screen user interface lets you proceed effortlessly through the measurement process of your samples Low cost of ownership, limited infrastructure requirements no need for compressed air or external cooling, lowest power consumption X-ray tube has a virtually unlimited lifetime Automation capabilities HighScore software for ease of analysis Upgradable for power, grazing incidence, transmission, non-ambience 	<ul style="list-style-type: none"> Can be calibrated for a large range of materials' properties Highly advanced statistical analysis software to correlate spectra to any number of materials properties Spectral resolution 3nm - 10nm Wavelength range 350nm - 2500nm Non-destructive method requiring little or no sample preparation Wide spectral range can measure multiple properties simultaneously 	What is special about this product?

Ask about specials and automation!

Elemental composition, film thickness

Sample preparation for XRF and ICP



Product	Zetium	Epsilon 4	Epsilon 1	Epsilon Xflow	LeNeo	FORJ	Eagon2	LeDoser 12	Product
What does it measure?	Elemental composition Film thickness	Elemental composition Film thickness	Elemental composition Film thickness	Elemental composition QC of liquids	<ul style="list-style-type: none"> Automated fused bead machine, for sample preparation that provides a factor x10 improvement in accuracy for XRF measurements Automated preparation of peroxide and borate solutions for ICP. Decrease the sample preparation time, achieves complete dissolution, increase users' safety 			Automated sample weighing and borate flux dispensing for sample preparation by fusion	What does it provide?
Technologies used	X-ray fluorescence (XRF)	X-ray fluorescence (XRF)	X-ray fluorescence (XRF)	X-ray fluorescence (XRF)	Automated borate fusion			High precision autonomous weighing	Technologies used
What is it used for?	<ul style="list-style-type: none"> Quantification of elements Be-Am with superior lowest limits of detection for most elements All materials in solid, powder or liquid form Detection and measurement of elements in thin films including film thickness measurement Small spot mapping for all elements across the range 	<ul style="list-style-type: none"> Quantification of elements B-Am with lowest limits of detection for most elements All materials in solid, powder or liquid form Detection and measurement of elements in thin films including film thickness measurement 	<ul style="list-style-type: none"> Quantification of elements from Na to Am with lowest limits of detection for most elements Online elemental analysis of solids, powders and liquids 	<ul style="list-style-type: none"> Online elemental analysis of liquids 	<ul style="list-style-type: none"> Preparing glass disks for elemental analysis using XRF Preparing peroxide or borate solutions for elemental analysis using ICP 	<ul style="list-style-type: none"> Preparing glass disks for elemental analysis using XRF Preparing peroxide or borate solutions for elemental analysis using ICP 	<ul style="list-style-type: none"> Preparing glass discs for elemental analysis using XRF 	<ul style="list-style-type: none"> Eliminates repetition of operators' tasks, reduces fatigue and errors. Increases productivity 	What is it used for?
What is special about this product?	<ul style="list-style-type: none"> Combination of ED and WD reduce measurement time up to 50% Batch automation Simple & Intuitive SuperQ software with the Virtual Analyst FastScan Omnia program standardless analysis Largest range of bespoke and in-house certified calibration standards Dust removal device minimizes contamination and maximizes instrument uptime SST R-mAX tube with CHI-BLUE window coating for increased X-ray tube durability and less drifting Small-volume airlock design for rapid cycling of samples into vacuum, or low He consumption for liquids analysis Supported by expertise and CRMs for all materials types 	<ul style="list-style-type: none"> Can handle a large variety of sample sizes, from less than a gram up to larger bulk samples, including irregularly shaped objects Small footprint allows it to be sited near to, or even next to, the production line for process control Automatable Close coupling of Tube-sample-detector for optimized sensitivity Automatic and builtin drift monitor for best accuracy Unique combination of 10 position sample changer with spinner Creating unlimited applications Automatic Program Selection (APS) for easy operation Supported by expertise and CRMs for all materials types 	<ul style="list-style-type: none"> All measurements in air. No need for helium or vacuum pump Highest analytical performance in its class Completely X-ray safe operation Built-in drift monitor for best accuracy and easy operation Creating unlimited applications Automatic Program Selection (APS) for easy operation Supported by expertise and CRMs for all materials types 	<ul style="list-style-type: none"> Instantaneous results for feedback into control systems Supported by expertise and CRMs for all materials types Ask about bespoke in-line solutions! 	<ul style="list-style-type: none"> 3 preparation modes in 1 instrument 1 fusion position Small and compact, fits in limited space Ready to use right out of the box Self-installed Absolute safety for the operator Quick and easy replacement of internal refractory plates 	<ul style="list-style-type: none"> 6 positions for fusion Resistance-based heating system Completely closed furnace Cold to cold operation without risks of injuries for the operators A patented handling mechanism transfers the crucibles and molds into the furnace, improving fusion cycle time and robustness Heat-drop and heat-ramp functions to enhance oxidation conditions while boosting production rate Sample agitation by swirling to increased sample homogeneity, dissolution speed and fusion success rate Optimized chamber design ensuring heat homogeneity across the different fusion positions Pre-heat and shut-off timer to program heating periods and ensuring power saving Supercapacitor for safe closure of furnace door and continuation of fusion cycle in case of brief (< 30 seconds) power cut 	<ul style="list-style-type: none"> Absolute safety for the operator 2 fusion positions Optional exhaust adapter for minimum infrastructure requirements Casting dish sensors: no possible damage of the instrument related to pouring without the platinum mold Non-wetting agent pills injection for optimized fusion method efficiency Pause and inspection function to visualize the fusion process during the fusion cycle 	<ul style="list-style-type: none"> Synchronizes the sample preparation process and enables the samples to be ready just in time for the subsequent fusion and analysis steps 12 positions Saves 90% of labor time related to the weighing step Fully adaptable to your standard operating procedure (SOP) Eliminates data transfer errors because it is LIMS ready and has sample tracking option Self-installed 	What is special about this product?

Ask about smart manager!

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



www.malvernpanalytical.com

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