

2830 ZT Wafer analyzer

Non-contact determination of layer thickness and composition



Supporting the semiconductor, data storage and saw device industry - now and in the future

Semiconductor and data storage technology has advanced rapidly over recent years, and the rate of change continues to gather momentum. This places considerable pressure on the metrology techniques used to control production developments and meet ever more rigorous demands.

Deposited thin films and multi-layers continue to decrease in thickness while increasing in complexity. As new technologies emerge at an escalating rate, yield improvement (during process ramp up) must evolve accordingly. Flexible and cost-effective process monitoring has become increasingly critical to obtain optimal process conditions and maximum yield.

X-ray fluorescence (XRF) spectroscopy meets the key requirements of semiconductor and hard disk manufacturers. The technique is inherently suited to thin film analysis, offering exceptional precision and sensitivity for the simultaneous determination of layer thickness and composition. Moreover, XRF is compatible with a broad range of applications, facilitates straightforward stack analysis and is synonymous with high uptime and low running costs.





XRF in surface acoustic wave (SAW) device applications

SAW devices are used in electronic components to provide a number of different functions such as RF filters, oscillators, and transformers. It is widely used in telecom networks, cellular base stations, satellite navigation systems, mobile phones, automotive, and TV set-top boxes.

As smartphones are reaching Gigabit LTE speeds in 5G technology, the number of cellular bands within the phones is also rapidly increasing to provide support (current smartphones already have ~40 cellular bands). Advanced filtering technology provided by these SAW devices is needed to support carrier aggregation across a wide variety of frequency bands while managing interference issues and delivering excellent radio performance.

The thickness control of the thin films in these SAW devices is of prime importance because a very small change in the thickness can already change the operating frequency of the device. In this application space, the 2830 ZT wavelength dispersive X-ray wafer analyzer is an excellent and proven solution to meet the most stringent requirements of top SAW device manufacturers.

The 2830 ZT is designed for maximum uptime and capability. It features the latest X-ray technology, including:

- Zeta technology, eliminating X-ray tube aging, greatly enhancing operational efficiency.
- 4 kW SST-mAX tube for operation at 160 mA current to give the highest sensitivities and lowest detection limits.
- FP Multi software allowing simultaneous analysis of film thickness and composition of stacks up to 16 layers.



Ultimate capability for measuring films thickness and composition

The 2830 ZT wavelength dispersive X-ray fluorescence (WDXRF) Wafer Analyzer is the successor to Malvern Panalytical's highly acclaimed PW2830. Designed specifically for the industry, it enables the determination of layer composition, thickness, dopant levels and surface uniformity for a wide range of wafers up to 300 mm.

The 2830 ZT Wafer Analyzer is equipped with Malvern Panalytical's advanced 4 kW SST-mAX X-ray tube, featuring groundbreaking ZETA technology which eliminates the effects of X-ray tube aging – by far the largest contributor to instrument drift. As a result, the 2830 ZT attains the highest levels of productivity and sensitivity, with excellent light element performance. What's more, thanks to ZETA technology, the 2830 ZT maintains these qualities throughout its entire lifetime.

Benefits

Unrivalled productivity

- Continuous capability and speed.
- Maximized uptime.
- Superior light element capability.

Fast and cost-effective multi-layer analysis

- Powerful Fundamental Parameter software.
- Analysis of stacks up to 16 layers.
- Simultaneous measurement of up to 24 elements.
- Intuitive operation.

Seamless fab integration

- Efficient, compact system design
- From manual to fully automated wafer loading.
- Compliance to all relevant standards.

A complete solution

- Excellent tool matching.
- Complete support.





2830 ZT application examples:

- Dielectrics: BPSG, PSG, BSG, ASG
- Doped polysilicon
- Barrier films and stacks: Ta, TaN, WC_xN_y , TiN, Si_3N_4
- Silicides/salicides: TiSi_x, CoSi_x, WSi_x, NiSi_x
- Metallization films and stacks: Cu, AlCu, Ti, W, Au, Pt, AuGe, Ag, Sn, etc
- Low-k dielectrics: FSG, SiOF
- High-k dielectrics: HfO₂, HfAlO_x, Ta₂O₅, etc
- TMR and GMR related films and stacks: CoFe(B), AIO_x, Ru, NiFe, IrMn, CrPtMn, etc
- PRAM and FeRAM films: PZT, SBT, BLT, GeSbTe
- SAW and BAW films and stacks: Al, Cu, AlCu, Ti, Ta, Pt
- SiGe
- Phase change materials: GeSbTe, InSbTe, SeAgGe, GeAsT

Unrivalled productivity

Continuous capability and speed

SST-mAX features groundbreaking ZETA technology which eliminates the effects of X-ray tube aging. 'New tube' performance is maintained throughout the tube's lifetime. Together with high sensitivity, ZETA technology ensures that rapid analysis and short measurement times are maintained across the lifetime of the tube. ZETA technology strongly reduces the need for drift correction and recalibration which increases productivity and uptime of the instrument.



Maximized uptime

Conventional X-ray tubes suffer tungsten evaporation, which causes deposits on the inside of the tube's beryllium window. Instrumentation using such tubes requires regular drift correction to compensate for decreasing intensity, especially for light elements. Implementation of the SST-mAX in the 2830 ZT solves this drift problem, thereby maximizing uptime and maintaining instrument precision over time.



Superior light element performance

The key features to maximizing sensitvity and stability of 2830 ZT for light elements include:

- 4 kW output SST-mAX, operating at a high current of 160 mA.
- Dedicated high performance channels for light elements in the range from boron to magnesium.
- Dry (oil-free) pumps maintaining a clean, sub-Pascal vacuum within the measurement chamber.



Save on costly downtime hours

Save time on instrument monitor corrections with the ultimate long-term stability of SST-mAX tube. Light elements such as boron and carbon often monitored on a daily basis can now be monitored weekly or bi-monthly, minimizing instrument downtime and thus, saving money.

Consistent and stable measurement results

The 2830 ZT is designed for maximum uptime and capability. It features the latest X-ray technology, including:

A range of critical features maximize reproducibility and minimize measurement times. For example, the laserassisted wafer height positioning system guarantees consistent and stable measurement results. This feature also makes it possible for 2830 ZT to analyze films on a large range of substrate materials in addition to Si wafers, including piezo- and pyroelectric, glass and quartz substrates.





Zeta technology enables long term stability and limits tube replacements. Save additionally on tube cost and downtime by upgrading to SST-mAX.





Fast and cost-effective multi-layer analysis

Powerful Fundamental Parameters

Analysis of stacks

FP modeling allows reliable thin film analysis. Malvern Panalytical's best-in-class algorithms model X-ray fluorescent behavior of the elements to calculate precise layer composition and thickness. This reduces the time and cost involved in setting up thin film and multi-layer analyses. System calibration requires only a minimum number of reference samples to analyze complex stacks of up to 16 layers.

Simultaneous measurement of up to 24 elements

The 2830 ZT can be fitted with up to 24 fixed measurement channels, from boron onwards. These high performance channels allow for multi-layer analysis according to specific user requirements. Furthermore, as new channels can be added at any time, the 2830 ZT offers powerful flexibility, allowing the user to keep up with the latest technologies.

Intuitive operation

The 2830 ZT utilizes Malvern Panalytical's proven SuperQ Thin Film software platform. This powerful, easy-touse thin film data collection and analysis package is renowned for its highly intuitive interface and menu. It enables simple setup of standard measurement protocols and individually customized programs. User-assigned function keys or large on-screen 'speed buttons' permit rapid initiation of measurement for single wafers or complete batches.

Keep track of your stack





Malvern Panalytical's SuperQ software makes it simple to set up spot patterns for uniformity measurement over the full surface area of wafers up to 300 mm.



Malvern Panalytical's Fundamental Parameter (FP) software package, FP Multi, is fully integrated into the SuperQ thin film platform of the 2830 ZT. It enables the direct calculation of chemical composition and thickness of layered materials and is ideal for analyzing both single layer thin films and complex multi-layer film stacks.

Uniformity patterns and wafer maps are easily defined with user-friendly spot location selection. Click-anddrag commands allow straightforward comparison with international wafer configuration standards.



Measurement programs are quickly selected via clearly labeled "speed buttons" on the built-in touch screen monitor.

Seamless fab integration

Efficient, compact system design

8

With a footprint comparable to that of 200 mm XRF wafer analyzers, the 2830 ZT occupies a minimum amount of valuable cleanroom space. A very narrow face print contributes further to space savings. The 2830 ZT can even be flush-mounted into a cleanroom wall, leaving only the loading port panel apparent within the workspace.

In accordance with normal fab procedures, the 2830 ZT measures wafers in a surface-up orientation. A transfer robot and notch/flat alignment unit position samples on the x-y wafer stage. Vacuum pumps can be installed remotely for convenient maintenance.

The system integrates readily into plant-wide process control and quality assurance networks, using SECS/ GEM protocols to communicate with the host computer.

From manual wafer loading to full automation

2830 ZT accepts 100 mm to 300 mm wafers, and smaller wafers or fragments can be measured using adaptors. To ensure integration into any fabrication environment, various wafer-handling options are available:

- Manual loading
- Open cassette loading with the robot in the cleanroom atmosphere
- Single or dual FOUP loading with robot and alignment unit in an ISO class 1 mini-environment
- Single or dual SMIF loading with robot and alignment unit in an ISO class 1 mini-environment
- Any combination of SMIF, FOUP and open cassette loading

Single load port FOUP and SMIF systems can be equipped with an internal reference position which can harbor a cassette with frequently used reference wafers. The reference position accommodates three wafer sizes in a range of 100 to 300 mm.



Improving safety in semiconductor manufacturing

As the semiconductor industry expands and gets more complex, the safety requirements in the manufacturing environment becomes even stricter. More and more semiconductor fabs require full compliancy to the SEMI-S2 standard.

What is the SEMI-S2 standard?

SEMI-S2 is in fact the most well-known standard in semiconductor manufacturing equipment for EHS. It covers 22 specific EHS categories, including regulatory requirements; electrical, mechanical, fire, chemical, radiation, noise, and ergonomics hazards; emergency shutdown, ventilation, and exhaust specifications; hazard warnings and more.

2830 ZT SEMI-S2 compliance

The 2830 ZT Wafer Analyzer has been certified as compliant to the latest SEMI-S2 standard. This means that the 2830 ZT instrument can be trusted as being a completely safe instrument, helping you to ensure and improve safety in the semiconductor manufacturing environment.



Compliance to all relevant standards

Malvern Panalytical's PW2830 was the first X-ray metrology system to comply with all 300 mm standards. This led to its rapid establishment as the system of choice for automated wafer analysis. Now, the 2830 ZT builds on this reputation, and is designed to accommodate all relevant industry norms and standards.

Importantly, the 2830 ZT is fully compliant with SECS/GEM (Semiconductor Equipment Communication Standards/Generic Equipment Model), GEM300, and AMHS (Automatic Material Handling System).



9

Completing the solution

Excellent tool matching



Metrology accuracy, reproducibility and repeatability are essential for accurate results within a fab. However, many organizations operate fabs at multiple locations. With different fabs making the same product, consistency of results is essential. 2830 ZT can be placed at different fab locations and matched to a 'gold standard' reference tool using a universal set of reference wafers.

This provides easy multi-production line quality control of thin film deposition and other processes. Excellent tool matching is the result of the high quality optical components, in combination with robust software algorithms.

Complete support

Malvern Panalytical understands the importance of continuity in production control. The 2830 ZT is based on proven concepts, established throughout the company's XRF thin film metrology product line.

However, technology is only part of a successful process control solution. To make sure that users maximize the benefit of the 2830 ZT, Malvern Panalytical provides excellent technical, software and application support.





2830 ZT users can take advantage of a large knowledge base when creating new applications, training new users, or solving technical issues.

Malvern Panalytical takes great pride in offering outstanding customer service, and values the trust users repeatedly place in the organization. A range of service contracts, including 24/7 coverage in most areas, is available, delivered by the largest X-ray support and service force in the market.



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



correctness or completeness

accuracy,

Malvern Panalytical

Grovewood 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