



**Malvern
Panalytical**
a spectris company

Zetium

Cement edition





Need to improve process efficiency and margins at your cement plant? The Cement edition of the Zetium spectrometer is your instrument. With robust, accurate, fast analysis of raw materials, fuels, raw mix, hot meal, clinker, and cement, it takes elemental composition analysis to the next level. Say hello to stable, optimized processing!

Elemental innovation

Continuous development, improved customer experience

Benefit-driven innovation on the Zetium Cement edition has made it the most powerful multipurpose tool for the analysis of cement and associated raw materials.

Elemental intelligence

Advanced analytical software for advanced analytical hardware

Now, our SuperQ software is enabling even more technology combinations and analytical possibilities for the Zetium. Its Virtual Analyst also makes setting up and operating the system simpler.

Elemental technology

60 years of experience: The ideal starting point

The Zetium Cement edition is the next generation in a long line of WDXRF platforms, including Axios, MagiX, and PW2400. Over the years, we've fine-tuned this proven technology – providing a strong foundation for the Zetium platform.

Elemental support

Reliable support – whatever the location

From service and training to laboratory analysis, we support you all the way. With a worldwide network of experienced engineers and application specialists, and full remote support, we'll make sure your spectrometer works at its best – wherever you are in the world.

Superior accuracy, simplicity, and reliability

By seamlessly combining our core and innovative technologies, the Zetium Cement edition delivers unrivaled analytical performance, speed, and robustness.

Virtual Analyst: Wave goodbye to guesswork

Analysis is a complex task with many critical choices and variables. SuperQ's Virtual Analyst software can help you make those choices – just like having an application specialist available 24/7. The software actively calculates your ideal measurement conditions, taking into account spectrometer configuration, sample information, line overlaps, required analytical range, detection limits, and precision. No more guesswork!

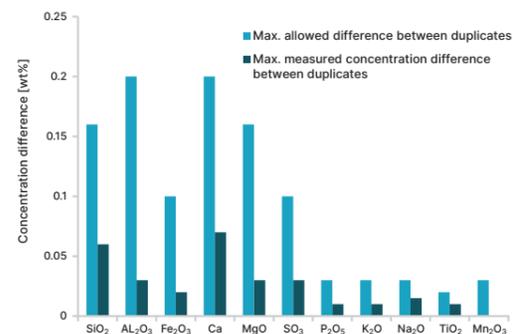
Maximum uptime, stability and precision

Instrument drift and contamination can affect the accuracy of your results. Luckily, the Zetium Cement edition's unique drift-free SST R X-ray tubes eliminate the

largest source of instrument drift, while its direct optical position sensing (DOPS) technology maximizes precision. And we've minimized sample contamination through a range of advanced technologies, including automatic sample-type recognition, turret-style sample loading, preloading dust removal, and CHI-BLUE tube window protection.

Stay ahead of ASTM and ISO compliance

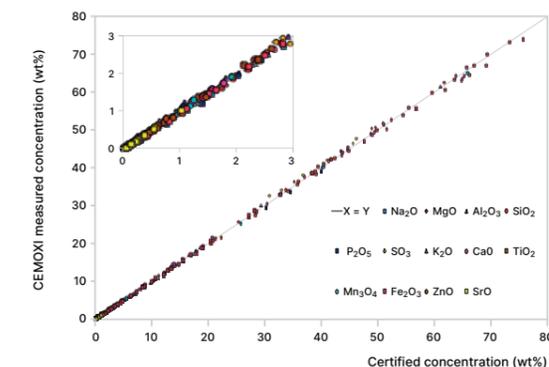
We know how important compliance is – so rest assured that the Zetium Cement edition satisfies ASTM C114-13 and ISO 29581-2 requirements for cement analysis. You can be fully confident in your analysis and product quality!



Cement Extension for WROXI-CRM

Preparing enough standards for reliable calibrations can be tricky – not to mention expensive. Cement Extension of WROXI Certified Reference Materials: a complete analytical solution for the analysis of 13 oxides in cement and related materials.

The Cement Extension contains nine synthetic Certified Reference Materials contains nine synthetic standards, reliable fusion recipes, an application template, and monitor samples. In other words, it's an out-of-the-box solution for the analysis of major and minor elements in raw materials, raw mix, clinker and cement.





Seamless integration in automation

Want to analyze the mineralogy of raw materials, raw meal, hot meal, clinker, additives, or cement, as well as their elemental composition? Just link the Zetium to our Aeris X-Ray diffractometer, and you'll be all set with reliable mineralogical feedback alongside your elemental data.

The power of combining technologies

You can easily integrate the Zetium with other instruments and sample preparation equipment via a belt connection. The result? An automated lab that supports both mineralogical and elemental analysis – which means higher productivity and lower process cycle times.

A process tailored to your needs

We make sample loading for these automated systems easy. You can access it from either side or from the back of the Zetium Cement edition, and choose from 51.5mm or 40mm sample holders. We can also include an optional sample inverter to make sure your sample is oriented correctly.

Interested in a turnkey automation solution?
Contact us today for a tailored quotation!



SMART MANAGER, SMART ZETIUM

The distractions and demands of quality control management can make it difficult to get the most out of your instruments. To make it easier, we've connected the Zetium to our Smart Manager platform – making it a Smart Zetium.

Meet your new trusted team member

This cloud-based 'control room' connects all your Zetium instruments, giving you a clear picture of their performance – wherever they are in the world. So you can keep on top of performance, optimize usage, reduce downtime, and unleash the potential of your data. Just like having a new, trusted member of your team!

Maximum data security

And rest assured: all your data remains yours, and is only visible to Malvern Panalytical. Smart Manager uses the latest Microsoft Azure cloud technology, ensuring that your data is safe and secure at all times. We don't collect data from your own samples unless you explicitly request it.

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.

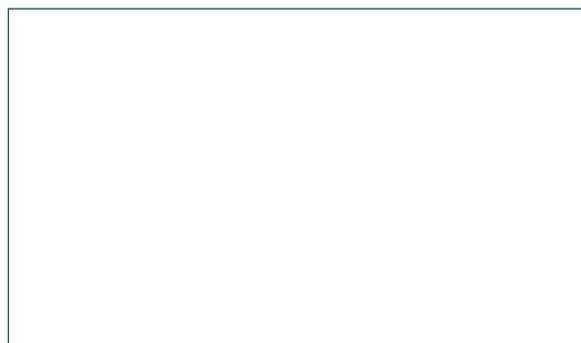
Malvern Panalytical. We're BIG on small™

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