



**Malvern
Panalytical**

EPSILON 4 FOOD & ENVIRONMENT



TRUST YOUR PRODUCT QUALITY

Analyze accurately and frequently

Growing populations increase the need for a better handling of our earth's resources. The Epsilon 4, an energy dispersive X-ray fluorescence (EDXRF) spectrometer, is a powerful analytical tool that can obtain useful elemental information to maximize crop growth and ensure product and environmental safety.

The latest advances in excitation and detection technology in Epsilon 4 opens possibilities for challenging applications that are traditionally performed by ICP and AAS. Switching from ICP to EDXRF significantly reduces the use of consumables, utilities and time . Discover how elemental analysis with Epsilon 4 adds value to the food production process.

Epsilon 4's value for Food & Environment

- Suitable for wide range of sample types: air filters, liquids, powders
- Low cost of ownership
- Low infrastructural requirements, ideal for at-line analysis
- Ready for any sample using standardless solution Omnian
- Simple, fast and safe sample preparation
- Non-destructive analysis
- Analysis from Carbon to Americium, from sub-ppm's to 100 wt% concentrations



Animal Feed

Control nutrients in animal feed accurately to keep cattle healthy and minimize manure greenhouse gasses.

Air Quality

Monitor the elemental composition of particle matter in ambient air, while complying to international norms

EPA IO 3.3, EN14902, HJ829-2017

Nutrient quality analysis

Frequent and accurate analysis during production processes, for example the addition of nutrients in milk powder, ensures product quality and consistency

Soil & Fertilizer analysis

Analyze the soil and match fertilizers to optimize growth conditions for crops. Also assure there are no toxic elements affecting the crops.

ISO 18227, ISO 15309 and ASTM C1255

Total salt concentration analysis (Na, Cl, K)

Derive NaCl content in food in one measurement.

Toxic element screening

Use Epsilon 4's high sensitivity to screen for toxic elements in raw materials and food products.



ACCURATE AND NORM COMPLIANT TRACE ANALYSIS AIR FILTERS AND SOILS

Epsilon 4 can accurately quantify many elements down to trace level concentrations according to stringent international test methods and norms, like EPA IO-3.3 for particles on air filters. Below are two application examples that demonstrate the analytical capability of Epsilon 4.

Elemental analysis of air filters according to US EPA method IO-3.3

Detection limits are an important measure of an instrument's performance. The detection limits (LLD) for this application were calculated from 20 replicate measurements of a blank sample and are based on 1 sigma (as specified in method IO-3.3).

Figure 1 shows a comparison between the LLD values reported in the EPA method and the LLD values obtained by Epsilon 4, in 45 minutes. All LLD values are smaller than the EPA limits.

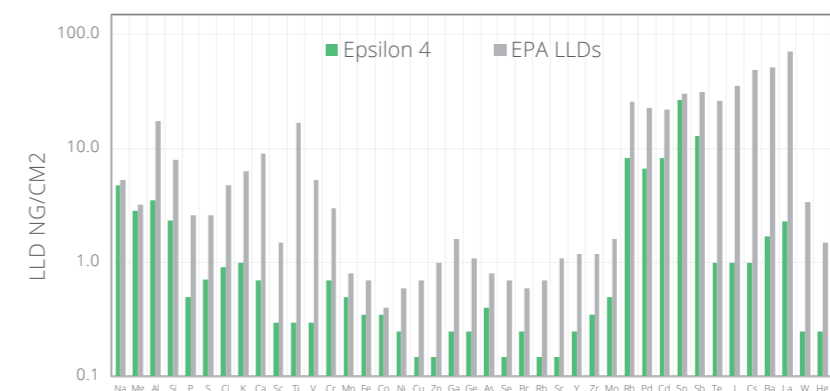


Figure 1. Comparison between the LLDs obtained using the Epsilon 4 and the required LLDs of the EPA IO-3.3 method

Elemental analysis of nutrients and toxic elements in soils

Twenty-six certified soil samples, measured as pressed pellets, were used for the application set-up of Epsilon 4. To test the accuracy and precision of the method, one Chinese soil standard GSS8 was measured five times consecutively. The certified and average measured concentrations, together with the standard deviations, are presented in **Table 1**, demonstrating excellent accuracy and precision, in 30 minutes measuring time.

Table 1. Accuracy and repeatability results obtained by measuring a soil standard GSS8 five times, consecutively.

Elements & Compounds	Certified values (mg/kg)	Results \pm std dev (mg/kg)	Elements & Compounds	Certified values (mg/kg)	Results \pm std dev (mg/kg)	Elements & Compounds	Certified values (mg/kg)	Results \pm std dev (mg/kg)
Na ₂ O (%)	1.72	1.33 \pm 0.02	Fe ₂ O (%)	4.48	4.35 \pm 0.01	Nb	15	14.2 \pm 0.1
MgO (%)	2.38	2.28 \pm 0.01	Co	12.7	12.2 \pm 0.1	Mo	1.16	1.06 \pm 0.08
Al ₂ O ₃ (%)	11.92	12.10 \pm 0.02	Ni	31.5	31.1 \pm 0.4	Cd	0.13	< 1
SiO ₂ (%)	58.60	54.84 \pm 0.05	Cu	24.3	22.0 \pm 0.3	Sn	2.8	2.0 \pm 0.3
P ₂ O ₅ (%)	0.18	0.18 \pm 0.01	Zn	68	59.6 \pm 0.3	Sb	1.04	1.1 \pm 0.4
SO ₃ (%)	0.03	0.05 \pm 0.01	Ga	14.8	16.9 \pm 0.2	Ba	480	499 \pm 2
K ₂ O (%)	2.42	2.35 \pm 0.01	As	12.7	12.6 \pm 0.2	Ce	66	60 \pm 6
CaO (%)	8.27	8.30 \pm 0.01	Br	2.6	2.6 \pm 0.1	W	1.7	1.9 \pm 0.3
Ti	3837	3625 \pm 4	Rb	96	95.5 \pm 0.1	Pb	21	19.1 \pm 0.2
V	81.4	83.1 \pm 1.4	Sr	236	227.3 \pm 0.2	Bi	0.3	1.1 \pm 0.1
Cr	68	72.1 \pm 0.8	Y	26	24.6 \pm 0.1	Th	11.8	11.7 \pm 0.4
MnO (%)	0.08	0.08 \pm 0.01	Zr	229	227.7 \pm 0.5	U	2.7	11.5 \pm 0.4

FAST, REPRODUCIBLE ANALYTICAL METHOD

Compared to other analytical techniques XRF requires little or no sample preparation

XRF is an ideal means of determining the chemical composition of all kinds of materials. Measurements in Epsilon 4 are carried out directly on the solid material (or liquid) with little to no sample preparation. There is no need for any dilution or digestion and therefore no disposal of chemical waste.

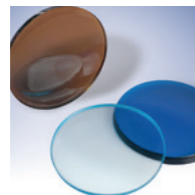
Epsilon 4 spectrometers can handle a large variety of sample types weighing from a few milligrams to larger bulk samples. Samples can be measured as:

- Solids
- Pressed powders
- Loose powders
- Liquids
- Fused beads
- Slurries
- Granules
- Filters
- Films and coatings

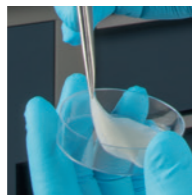
LIQUIDS



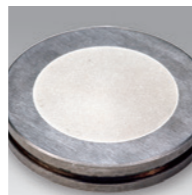
SOLIDS



AIR FILTERS



POWDERS



TAILORED SOLUTIONS THROUGH EXPERTISE

Experienced Malvern Panalytical staff work in close cooperation with you to provide not only training but also tailored analytical programs and procedures, balancing throughput and accuracy while minimizing set-up and running costs.

Access to the right calibration samples is key in XRF. Malvern Panalytical helps in obtaining or creating the standards you need. We provide total solutions including standards for several key applications. We can also generate suites of in-house standards by certifying your materials through our ISO 17025 certified laboratory.

Sample preparation, although typically straightforward for XRF, is an important factor in the overall analytical precision and accuracy. Sample preparation needs to be quick, robust and reproducible, and the choice of sample preparation technique starts with your requirements and materials.

Combine XRF with the analytical solutions Malvern Panalytical offers, like near infrared spectroscopy, particle size analysis and X-ray diffraction. Our experts can advise you which approach suits best given your material types and analytical requirements.

Tap into our knowledge network through our global Expertise Centers to optimize your complete analytical process, including sample preparation methods and equipment.



Our aim is to make Epsilon 4 an essential part of your elemental analysis

The added value for you is what drives us:

- The largest support network in the industry
- Training programs customized to your needs
- Reference materials
 - *Certified reference materials (CRMs)*
 - *Synthetic reference materials tailored to your requirements*
- Analytical services
 - *Certify your samples through our ISO 17025 certified laboratory*
- Consultancy
 - *Norm compliance*
 - *Laboratory information management*
 - *Process automation*
 - *standard operating procedures*
 - *Interlaboratory standardization*



MEASURE IT IN YOUR OWN LANGUAGE

1. Load your sample
2. Select required method
3. Enter relevant sample information
4. Just click Measure
 - 測量
 - 測定
 - Mesurer
 - Messung
 - Mesure
 - Zmierzyć
 - Medida
 - Измерить
 - Médir



ENHANCE YOUR ANALYSIS THROUGH SOFTWARE OPTIONS

Five industry software options are available to further enhance the capabilities of Epsilon 4: Omnian, Stratos, Oil-Trace, Enhanced Data Security and FingerPrint.

These dedicated options add new functional dimensions to benchtop spectrometry and take the hard work out of regulatory compliance.

Elemental screening OMNIAN



Our powerful Omnian software is ideal when there is no conventional calibration established for materials that require analysis. When faced with non-routine samples or materials for which there are no certified reference materials, Omnian provides excellent insight into the elemental composition. Designed to provide fast and reliable quantification, Omnian's advanced fundamental parameters (FP) algorithm automatically deals with the analytical challenges posed by samples of widely differing types.



Enhanced data security AUDIT TRAIL SOFTWARE



The enhanced data security software option is designed for GMP and GLP environments, and enables the user to comply with FDA 21CFR Part 11. The software includes every feature required to satisfy the strict environmental protocols, like multiple security levels, log in with user identification, reporting of date and time, results storing, extensive audit trails and LIMS integration

Pass/Fail analysis FINGERPRINT



FingerPrint is a material type confirmation routine that uses a rapid statistical analysis of the spectrum for a simple PASS/FAIL answer. Spectra used for the FingerPrint routine can also be used for conventional compositional determination for a more complete diagnostic analysis.



WHY CHOOSE MALVERN PANALYTICAL?

We are global leaders in materials characterization, creating superior, customer-focused solutions and services which supply tangible economic impact through chemical, physical and structural analysis.

Our aim is to help you develop better quality products and get them to market faster. Our solutions support excellence in research, and help maximize productivity and process efficiency.

Malvern Panalytical is part of Spectris, the productivity-enhancing instruments and controls company.

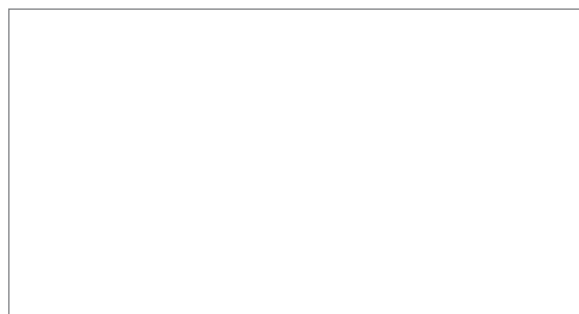
www.spectris.com

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

Groewood Road, Malvern,
Worcestershire, WR14 1XZ,
United Kingdom

Tel. +44 1684 892456
Fax. +44 1684 892789

Lelyweg 1, 7602 EA Almelo,
Netherlands

Tel. +31 546 534 444
Fax. +31 54 534 598

info@malvernpanalytical.com
www.malvernpanalytical.com

Disclaimer: Although diligent care has been used to ensure that the information in this material is accurate, nothing herein can be construed to imply any representation or warranty as to the accuracy, correctness or completeness of this information and we shall not be liable for errors contained herein or for damages in connection with the use of this material. Malvern Panalytical reserves the right to change the content in this material at any time without notice.
Copyright: © 2018 Malvern Panalytical. This publication or any portion thereof may not be copied or transmitted without our express written permission. 9498 707 6611 PNI 1453

www.panalytical.com/epsilon4