



**Malvern
Panalytical**
a spectris company

**REDUCE ANALYSIS TIMES
BY UP TO 80% FOR INCLUSIONS
AND ELEMENTAL MAPPING**

www.malvernpanalytical.com/HiPer

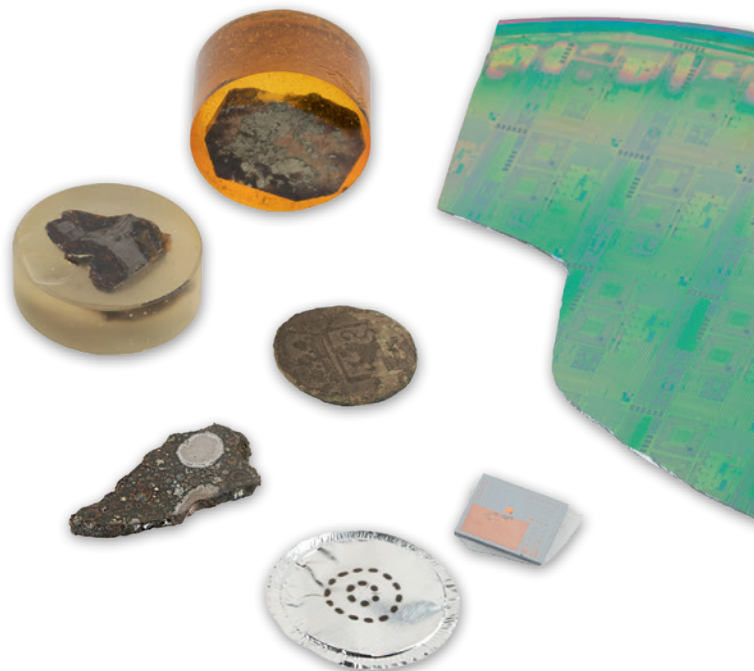
HiPer SMALL SPOT MAPPING

HiPer Small Spot Mapping is the latest addition to the Zetium platform and its SumXcore technology – an integration of WDXRF, EDXRF and XRD.

With the unique X-ray optics of HiPer Small Spot Mapping, measurement times can be reduced by up to 80%. This puts the Zetium in a class of its own with respect to analytical power, speed and task flexibility in a wide range of environments.

- Small spot mapping available on your spectrometer for production control
- Inclusion analysis in just minutes or map over 100 mm² in only a few hours
- Cover Na to Am and screen for unexpected elements
- Schedule and run small spot mapping on sample batches

www.malvernpanalytical.com/HiPer

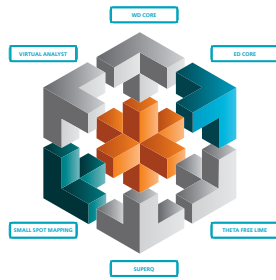


ZETIUM - ELEMENTAL EXCELLENCE

Elemental innovation

Continuous development, improved customer experience

The Zetium platform embodies SumXcore technology - an integration of WDXRF, EDXRF and XRD, creating a revolutionary step in materials analysis.



Elemental technology

60 years of experience and heritage - the ideal starting point

Zetium is designed to set new standards in terms of analytical power, usability and sustainability, to users in process optimization, quality control and research.

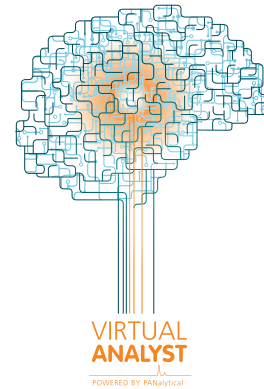
- Flexible sample handling
- Intelligent sample introduction
- Unrivalled accuracy & reproducibility



Elemental intelligence

SuperQ, evolutionary software for a revolutionary platform

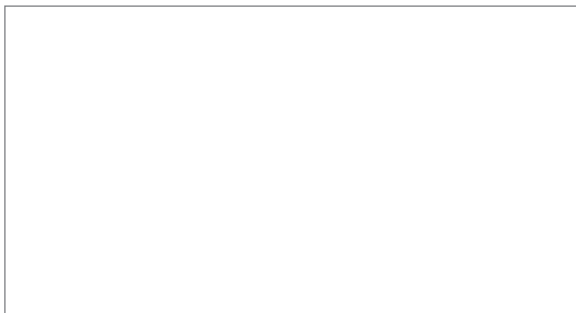
With the integrated intelligence of the Virtual Analyst, SuperQ software provides guidance on the optimal setup for each experiment.



Elemental support

Transparent and reliable

- Service
- Expertise
- Training & education
- Analysis & standards preparation



MALVERN PANALYTICAL

Groewood 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

info@malvernpanalytical.com
www.malvernpanalytical.com