

CUBIX³

3rd generation of successful industrial XRD applications

Cement
Iron
Minerals
Pharma
Potflux



CUBIX³

Ultimate speed for high-throughput production control

Innovation technology



The CubiX³ builds the next step in PANalytical's excellent reputation for fast, push-button mineralogical analysis.

Equipped with PANalytical's X'Celerator the CubiX³ is the fastest and most accurate diffractometer for production control. It yields an enormous gain in speed – up to 150 times faster than conventional detection technology for a complete diffraction pattern.

With the R&D 100 Award winning X'Celerator detector you achieve the ultimate speed, resolution, data quality and ease of use.

Smart process control

X-ray diffraction is the most efficient method of directly obtaining mineralogical information of different materials. Direct information about the mineralogical composition of your materials enables you to realize a constant quality and maximize the efficiency of your production processes. Analysis results can be obtained within a few minutes, thanks to solid-state X-ray detection technology and modern software algorithms.

The whole process of sample preparation, measurement, analysis and reporting can be fully automated and set up for push-button operation, 24 hours per day, 7 days per week. The reliable and rugged construction of the hardware ensures maximum uptime. The exceptionally high repeatability and accuracy of the measurements gives precise and reliable results throughout the system's life time.

CubiX³ advantages

- **Fast and precise production control**
 - High-speed data collection
 - Fast sample throughput
 - Rapid data analysis
- **Easy integration in automation environments**
- **Easy sample preparation**
- **No chemicals required**
- **Push-button operation**
- **Traceability in controlled environments, meets 21 CFR Part 11 requirements**
- **Fully automated measurement, analysis and reporting. No specialist XRD knowledge required**

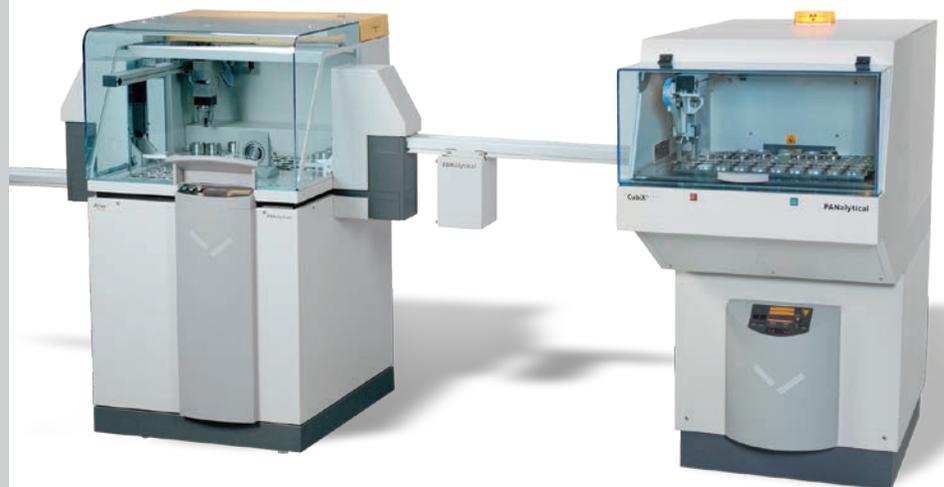


CUBIX³

TWIN solution

PANalytical TWIN solutions combine a PANalytical XRF and a PANalytical XRD system. It uses the single interface of the SuperVisor software to announce all samples just once, to route the sample to one or both systems and to control the actual measurements. The results can be merged before transmission into a joint analysis report. They are stored, printed or transmitted separately in different formats on demand.

- One user interface for both XRD and XRF
- Robot or conveyor connections possible
- Two complete systems for simultaneous use
- One sample for both methods, one result
- Support of 40 mm and 51 mm sample holder rings



Intelligent automation

PANalytical offers all types of full laboratory automation and integration solutions. We also work together with major suppliers of laboratory automation. A completely integrated solution combines the analysis systems with the sample preparation and automatically takes care of incoming sample materials. It also cleans and recycles sample holders, achieving the shortest throughput time.

- Universal Automation Interface
- Compatible with all laboratory automation systems
- Fits into any automation environment
- Extensible concept
- Results ready for transfer to a LIMS system



Making the most of your investment

The purchase of an XRD system will add value to a company for many years. To make even more out of your investment we provide a range of EXPERTISE programs that are designed to ensure that the capabilities of an X-ray system can be fully realized. They provide priority access to the company's unrivalled application knowledge and experience in analytical X-ray methodologies.



EXPERTISE programs

Gain priority access to PANalytical's unrivalled expertise in:

- Sample preparation methodology
- Applications development
- Good Laboratory Practice (GLP) consultation
- Optimization of calibration maintenance
- Tailored training packages

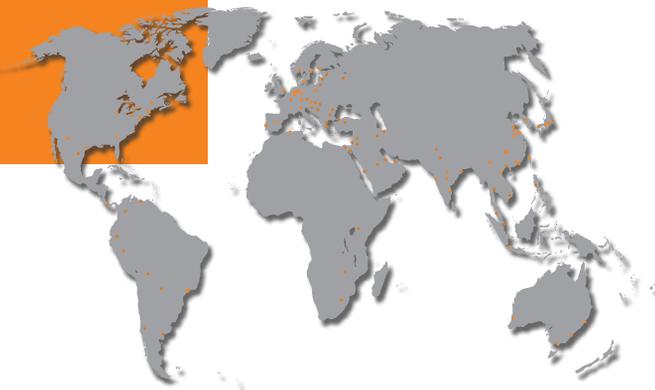
Customer support and application laboratory

PANalytical instruments play a key role in industry and research. High uptime, consistency and efficient use are essential. We place a high priority on customer support, and are proud of our extensive global network of sales and service centers.

Support facilities include:

- Preventive maintenance and repair
- Remote diagnostics
- Installation and commissioning
- Application courses
- Application support

Global and near



PANalytical is one of the world's leading suppliers of analytical instrumentation and software for X-ray diffraction (XRD) and X-ray fluorescence spectrometry (XRF). The materials characterization equipment is used for scientific research and development, for industrial process control applications and for semiconductor metrology.

During the last decade PANalytical has added a variety of other analysis techniques to their product portfolio. Optical emission spectrometry (OBLF GmbH, Germany), pulsed fast thermal neutron activation (Sodern, France) and near-infrared (ASD Inc.) capabilities together with XRD and XRF can provide customers with tailor-made analytical solutions for the characterization of a wide range of products such as cement, metals, nanomaterials, polymers and many more.

PANalytical's headquarters are in Almelo, the Netherlands. Fully equipped application laboratories are established in Japan, China, the USA, and the Netherlands. PANalytical's research activities are based in Almelo (NL) and on the campus of the University of Sussex in Brighton (UK). Supply and competence centers are located on two sites in the Netherlands: Almelo (development and production of X-ray instruments) and Eindhoven (development and production of X-ray tubes) and in Boulder, USA (development and production of near-infrared instruments). A sales and service network in more than 60 countries ensures unrivalled levels of customer support.

The company is certified in accordance with ISO 9001 and ISO 14001.

Visit our website at www.panalytical.com for more information about our activities.

PANalytical is part of Spectris plc, the productivity-enhancing instrumentation and controls company.

PANalytical B.V.

Lelyweg 1, 7602 EA Almelo
 P.O. Box 13, 7600 AA Almelo
 The Netherlands
 T +31 (0) 546 534 444
 F +31 (0) 546 534 598
info@panalytical.com
www.panalytical.com

Regional sales offices

Americas

T +1 508 647 1100
 F +1 508 647 1115

Europe, Middle East, Africa

T +31 (0) 546 834 444
 F +31 (0) 546 834 969

Asia Pacific

T +65 6741 2868
 F +65 6741 2166