

# DIRECT FEEDBACK AND FAST COUNTERACTIONS

From bauxite to alumina to aluminium processing

Contact us for more information

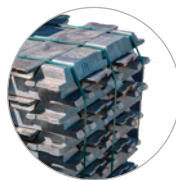
## REDUCE COSTS FOR REAGENTS AND ENERGY DURING MINING, REFINING AND SMELTING

Malvern Panalytical delivers innovative and unique analysis solutions to increase the efficiency of your bauxite mining operation, alumina refinery and aluminium smelters. Constant and accurate monitoring will enable significant savings in each step of the process.



### Bauxite mining

**Monitor** available alumina and reactive silica in the mine.  
**Enable** fast counteractions on changing bauxite composition.  
**Guarantee** ore quality to avoid penalties.



### Aluminium smelting

**Avoid** bath solidification with accurate and high frequent monitoring.  
**Minimize** energy consumption.  
**Automate** analysis for fast, safe and robust feedback.



### Alumina refining

**Reduce** consumption of caustic soda and energy.  
**Control** alumina particle size.  
**Optimize** red mud settling rates and flocculant consumption.

---

**Remote sensing (NIR) - On-line monitoring (CNA / NIR) - Elemental testing (XRF)**  
**Particle size and shape characterization (PSD)**  
**- Mineralogical analysis (XRD / NIR) - Sample preparation (Fusion)**

---

[www.malvernpanalytical.com/mining](http://www.malvernpanalytical.com/mining)



**Malvern Panalytical**  
a spectris company

## ANALYTICAL SOLUTIONS TO INCREASE PROFIT AND EFFICIENCY OF YOUR ALUMINIUM PRODUCTION

### DETECT AVAILABLE ALUMINA AND REACTIVE SILICA IN THE MINE

#### QualitySpec7000 cross-belt analyzer

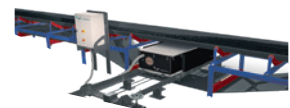
- Allows on-line mineralogical monitoring.
- Enables real-time analysis of available alumina and reactive silica.
- Operates under completely safe conditions.



### ENABLE OPTIMAL BLENDING OF THE ALUMINA REFINERY RAW MATERIAL FEED

#### CNA cross-belt analyzer

- Provides the elemental composition of your entire ore in real-time.
- Delivers reliable results independent from your belt width, load, variation or particle size.
- Operates under completely safe conditions.



### CONTROL MINERALOGY AND PROCESS PARAMETERS FAST AND EASY

#### Aeris

- Facilitates intuitive and accurate analysis of the mineral content of bauxite, alumina modifications ( $\alpha$ ,  $\gamma$ ) and bath composition.
- Operates in rough industrial environments and remote destinations.
- Automatable.



### ENHANCE LABORATORY PERFORMANCE WITH AUTOMATED ANALYSIS

#### Automation

- Guarantees safe sample taking, transportation, preparation and analysis.
- Turn-key laboratory solutions.



### ANALYZE IMPURITY LEVELS AND GUARANTEE QUALITY

#### Zetium

- Allows fast, accurate and simultaneous analysis of light and heavy elements.
- Tracks impurities to lowest levels.
- Designed for high-throughput environments.



### INCREASE LIFETIME OF MILLS AND ENSURE OPTIMAL REFINING AND SMELTING CONDITIONS

#### Mastersizer 3000

- Monitors green and accurate the particle size of bauxite and alumina.
- Enables wet and dry analysis
- Automatable.



## 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](http://www.malvernpanalytical.com)

[www.malvernpanalytical.com/mining](http://www.malvernpanalytical.com/mining)