

ORE AND MINERALS ANALYSIS (OMA) WORLDWIDE WORKSHOPS

Reflect – Refresh – Recover

Please join one of
our OMA workshops

Throughout the year Malvern Panalytical organizes multiple OMA workshops all over the world. During these free-of-charge seminars, dedicated to the mining industry, we discuss how to achieve better recovery rates, reduce downtime and increase productivity.

The one-day workshops bring together experts from the mining industry, research institutions and Malvern Panalytical, as supplier of the newest technologies, to create value for mining and metal production.

News and advances in X-ray Fluorescence, X-Ray Diffraction, On-line monitoring (CNA), Particle Size and morphology analysis, Near Infrared (NIR) characterization and sample preparation are topics to be presented.

“Delegates appreciate the close collaboration of the host organization, usually a research institution or a mining company with our specialists, who address challenges the mining industry faces daily”.

Uwe König,
Global Mining Segment Manager,
Malvern Panalytical

Visit www.malvernpanalytical.com/OMA to find a workshop in your area.

Costs

The events have no costs; however, it is limited to a maximum of attendees. So book now.

Background OMA workshops

The first Ore & Minerals Analysis (OMA) workshop was held in August 2015 at the Physics Department of the Federal University of Minas Gerais (UFMG) in Belo Horizonte, Brazil. In November 2018 the first OMA workshop held outside of the American continents was hosted in Kathu, the iron ore capital of the Northern Cape province of South Africa. After then more countries were added to host the OMA workshop. The 20th OMA workshop will go back to its origin and we will celebrate in Brazil.

www.malvernpanalytical.com/OMA



**Malvern
Panalytical**
a spectris company