

25TH ORE AND MINERALS ANALYSIS (OMA) VIRTUAL EVENT

November 24th - November 25th, 2020

Register now on our website

WHAT'S THE FUTURE OF TECHNOLOGY IN MINING?

At Malvern Panalytical, we know that reducing costs and improving efficiency in the mining sector is an ongoing challenge. In previous years, we've helped mining professionals learn about the latest technologies through our highly popular series of Ore and Minerals Analysis (OMA) workshops. But this year, Covid-19 restrictions mean that we can't meet in person, so we're moving the events online, with a series of two-day Virtual OMAs with two hours of presentations on each day.

The 25th OMA virtual bonus edition is a great way of staying up-to-date with the latest advances in the field of mining and minerals. So join us to keep yourself informed on topics including:

Day 1: In the Field and Mine:

Advanced solutions for exploration and grade monitoring.

Day 2: In process and laboratory:

Adding efficiency and automation to processing and quality control.

This event brings together experts from the industry as well as our own experts to share knowledge for exploration to process optimization and environmental monitoring by improving the analysis of raw materials and process streams.

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 will be presented.

There is no fee for this event – register today and attend one or both days.

Summary

Date: November 24th - November 25th, 2020
Time: 10:00 - 12:30 Western European Time [West Europe]
Location: VIRTUAL
Language: English

www.malvernpanalytical.com/virtual-OMA



Malvern Panalytical
a spectris company

Day 1 (Exploration & Grade Monitoring)		
Time	Topic	Presenter
10.00 AM	Welcome, Value of analytical monitoring in mining	Dr. Uwe König , Malvern Panalytical, Business Development Mining & Metals (Netherlands)
	Importance of reconciled quantitative chemical and mineralogical information in mining projects	Prof. Pertti Lamberg PhD , Anglo American, CEO of AA Sakatti Mining Oy, (Finland)
	Taking your analysis to the field with portable NIR and remote sensing	Robert Cocciardi , Malvern Panalytical, Sales Representative and Channel Manager LATAM - NIR, (Canada)
	Data modelling for better mine site mapping and planning. Digitising your mine.	Dr. Sasha Pontual , IMDEX (New Zealand)
	Q&A	
	Growing demand for ore and competitive prices: Improve your grade control and optimise every dollar with on-line elemental, particle size, mineralogical analysis	Marek Gnyppek , Malvern Panalytical, Business Manager Eastern Europe (Poland)
	Geometallurgy characterization potential to be linked to process performance	Prof. Simon Michaux , GTK, Associate Professor Geometallurgy (Finland)
	Geometallurgy in Action - On-line gold mineralogy for process control	Dr. Jussi Liipo , Metso Outotec, Director, Geometallurgy & Mineralogy (Finland)
	Asset management - Realize maximum value during process monitoring	Martijn Fransen PhD , Malvern Panalytical, Director Digital Solutions (Netherlands)
12.30 PM	Q&A and closing	

Day 2 (Ore processing, QC, Environment)		
Time	Topic	Presenter
10.00 AM	Efficient minerals processing - what's particle size monitoring got to do with it?	Ulf Willén , Sector Sales Manager - Primary Materials - Western Europe, Malvern Panalytical (Sweden)
	Mineralogical monitoring applications - Case studies on how companies are using X-ray diffraction (XRD)	Dr. Sabine Verryn , XRD Analytical & Consulting (South Africa)
	Determination of lithium in different ores determined by X-ray diffraction (XRD) and statistical methods	Prof. Dr. Dr. Herbert Pöllmann , Martin-Luther-University Halle-Wittenberg (Germany)
	Keeping up with increased demand and working in Covid-19 safe distancing measures: why you should re-think robo-lab automation and on-line analysis for liquor monitoring	Michel Choua , Malvern Panalytical (Netherlands)
	Q&A	
	Advances in elemental analysis using XRF - increased throughput with combined ED and WDXRF technologies, cluster analysis, small spot analysis	Armand Jonkers , Malvern Panalytical, Application Specialist (Netherlands)
	Analysis of Ferro Chrome production samples using ED-XRF spectrometry	Hendrika Bosch , Glencore - Rustenburg smelter, Works Chemist (South Africa)
	The importance of accurate analysis for the monitoring of mining and beneficiation processes in Iron Ore and other industries	Italian Mashego , Anglo American, Section Manager, Laboratory Sishen Mine (South Africa)
	How optimizing your sample preparation leads to better control of your process	Dr. José Eduardo F.C. Gardolinski , Malvern Panalytical, Laboratory Manager (Netherlands)
12.30 PM	Q&A and closing	

MALVERN PANALYTICAL

Groveswood 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

www.malvernpanalytical.com/virtual-OMA