

21st Ore and Mineral Analysis – Virtual event

21st / 22nd July 2020



Ore and Mineral Analysis (OMA)

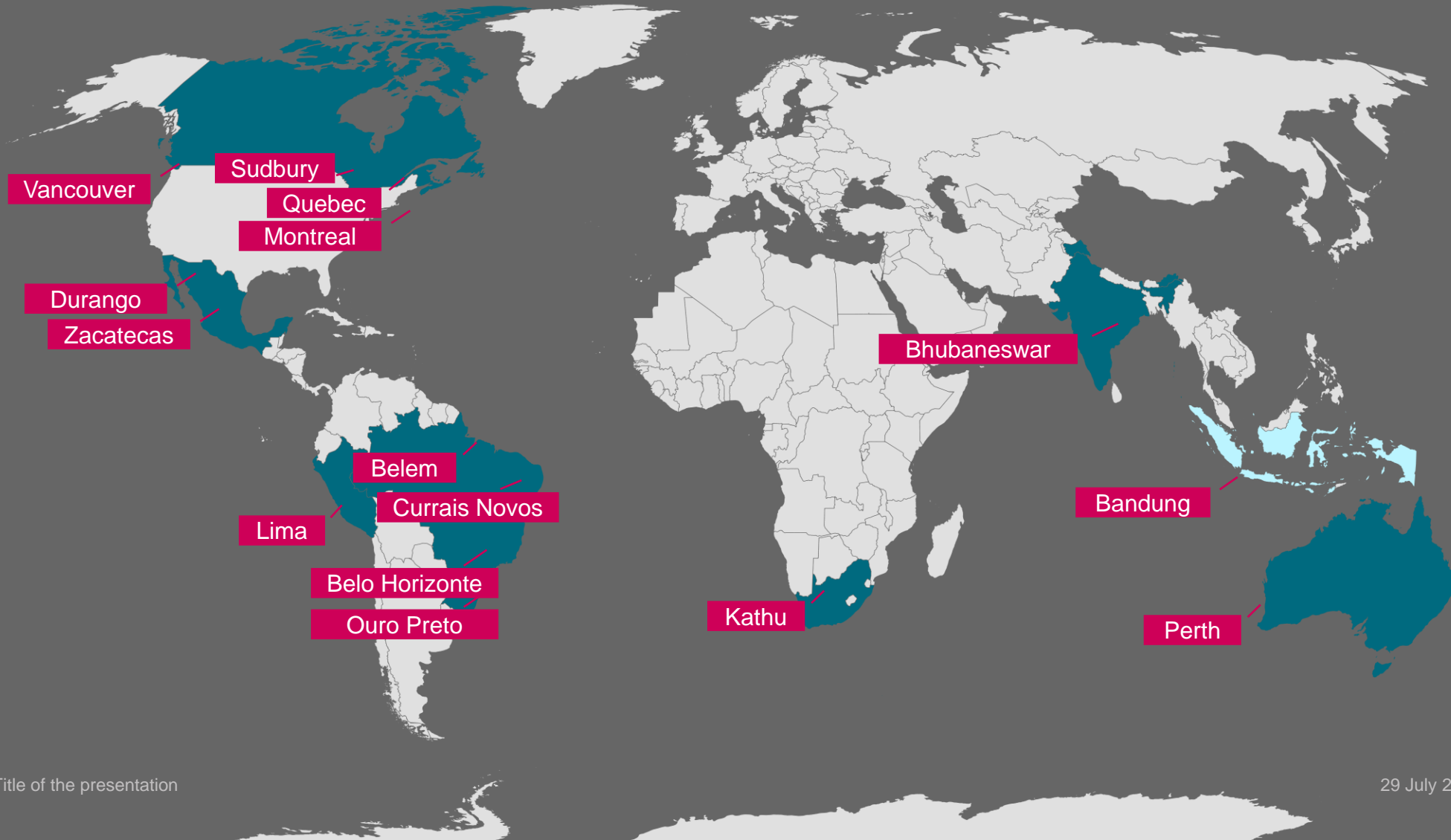
Workshop's



1st OMA 2015 Brazil, Belo Horizonte In cooperation with: Universidade Federal Minas Gerais 	2nd OMA 2016 Brazil, Nova Lima In cooperation with: Vale 	3rd OMA 2016 Peru, Lima In cooperation with: Enviroequip 	4th OMA 2017 Mexico, Zacatecas In cooperation with: Universidad Autonoma de Zacatecas 	5th OMA 2017 Brazil, Belo Horizonte In cooperation with: SGS Geosol 	6th OMA 2017 Canada, Sudbury In cooperation with: Laurentian Univ. / Goodman School of mining 	7th OMA 2017 Canada, Quebec In cooperation with: Claisse 	8th OMA 2017 Brazil, Belem In cooperation with: Universidade Federal do Para
9th OMA 2018 Canada, Vancouver In cooperation with: University of British Columbia 	10th OMA 2018 Brazil, Currais Novos In cooperation with: IFRN 	11th OMA 2018 Canada, Quebec In cooperation with: Claisse 	12th OMA 2018 South Africa, Kathu In cooperation with: Industrial Analytical (Pty) Ltd 	13th OMA 2019 Mexico, Durango In cooperation with: University of Durango 	14th OMA 2019 India, Bhubaneswar In cooperation with: IMMT 		
15th OMA 2019 Brazil, Ouro Preto In cooperation with: Universidade Federal do Ouro Preto 	16th OMA 2019 Australia, Perth In combination with: Opening new application laboratory 	17th OMA 2019 Indonesia, Bandung In combination with: Center for Mineral Research - tekMIRA 	18th OMA 2019 Canada, Montreal In combination with: Mc Gill University 	19th OMA 2019 Canada, Sudbury In combination with: Laurentian University 	20th OMA 2019 Brazil, São Paulo In combination with: LCT / USP 		

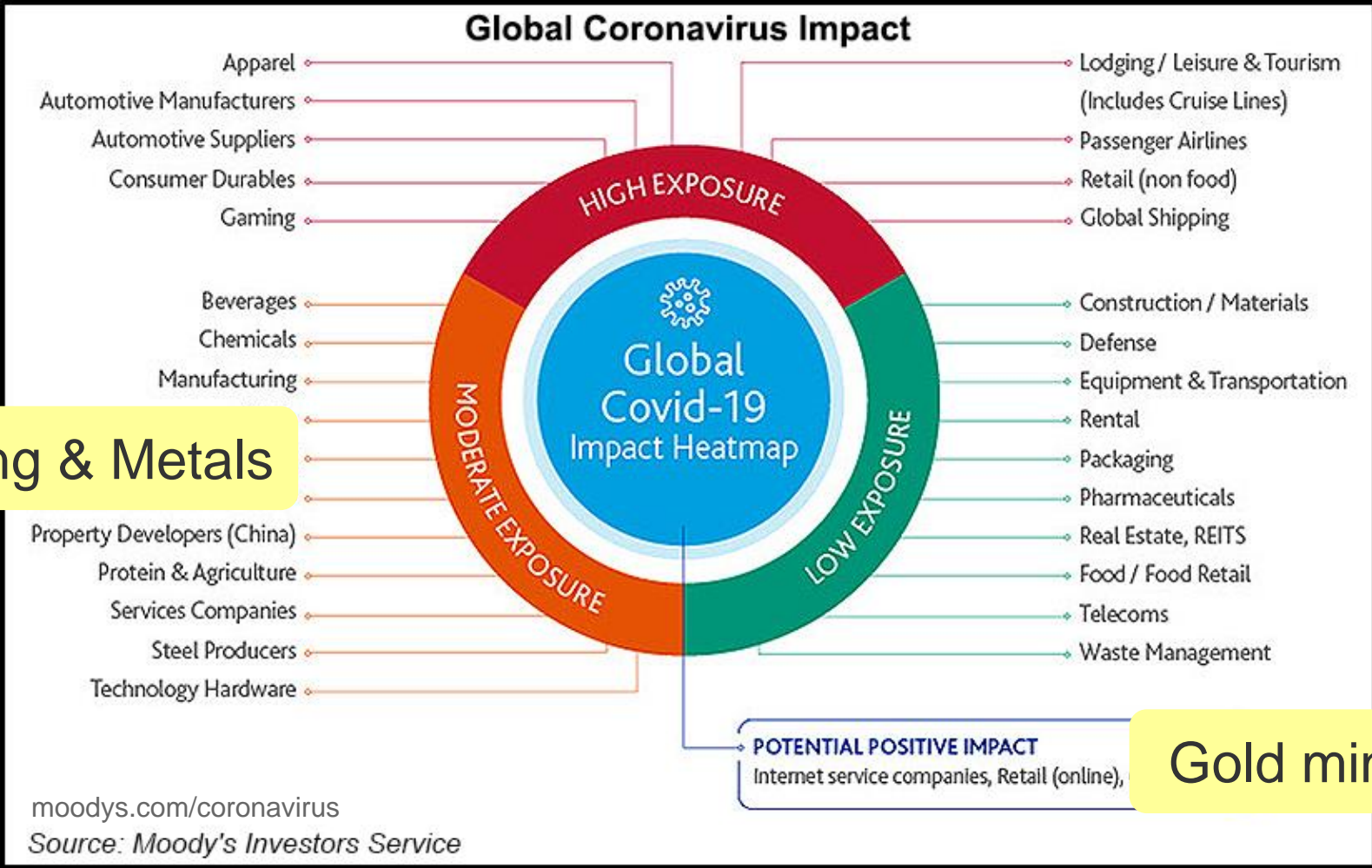
Ore and Mineral Analysis (OMA)

Workshop's



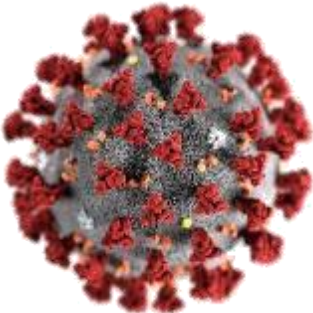
Mining industry

Covid-19 impact



Mining & Metals

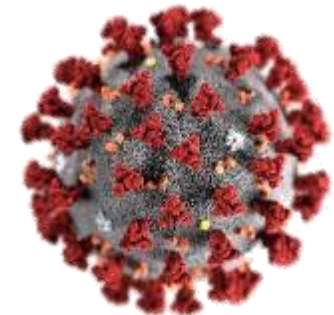
Gold mining



Mining industry

Covid-19 impact

- Isolated outbreaks and government mandated shutdowns
 - E.g. South Africa and Peru
- Demand for commodities like Cu, Fe, Zn remain low
 - Exception gold, benefiting from higher levels of uncertainty
- Lower energy prices impact mining industry positive
 - Energy typically constituting 20-25% of direct operating costs
- Companies likely to use crisis as a catalyst to rethink:
 - How and where work is done,
 - How to improve the ability to collaborate remotely,
 - How to accelerate adoption of automation and digital capabilities and
 - How to operate safe with a minimal workforce or no human intervention.



Mining Industry

Trends



GEOPOLITICAL

Global

- 2016 turning point for mining sector
- Challenges from technological disruption, CO₂ restrictions and environmental restrictions, health and safety



Rebound expected beginning 2021

Increased need for digitalization / automation

Mining Industry

Trends



GEOPOLITICAL

Global

- 2016 turning point for mining sector
- Challenges from technological disruption, CO₂ restrictions and environmental restrictions, health and safety



Rebound expected beginning 2021

Increased need for digitalization / automation

GEOLOGICAL

Mining industry

- Decreasing grades require more frequent and accurate monitoring, remote locations
- Reduce costs - Stabilize margins – Increase efficiency
- Increase of secondary and recycled materials in supply chain

Mining Industry

Trends



GEOPOLITICAL

Global

- 2016 turning point for mining sector
- Challenges from technological disruption, CO₂ restrictions and environmental restrictions, health and safety

Rebound expected beginning 2021

Increased need for digitalization / automation

GEOLOGICAL

Mining industry

- **Decreasing grades require more frequent and accurate monitoring**, remote locations
- Reduce costs - Stabilize margins – Increase efficiency
- Increase of secondary and recycled materials in supply chain

TECHNOLOGICAL

Industry 4.0

- Big data, predictive analytics
- **Multiple sensors, combined technologies**
- On-line analysis of ores (ore sorting), liquids and slurry
- Automated sample preparation and analysis (standardization, cost saving)

Analytical environments

Bring the laboratory to the sample



Field

- Geologists exploring new or existing deposits



Mine

- Mine operators, geologist monitoring ore grades and blending



Laboratory

- Chemists, mineralogists, quality managers analyzing routine samples (feed, concentrates, tailings)



Remote

- Operators analyzing on-site at remote areas or underground mines

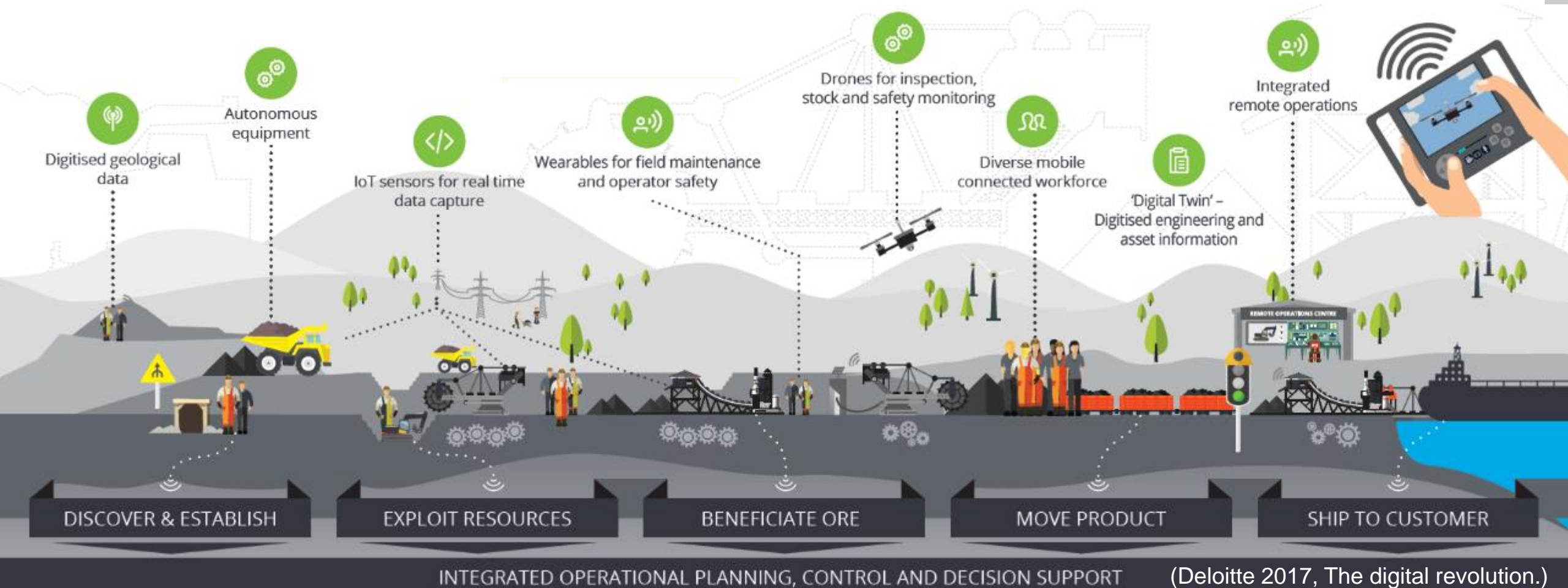


Automation

- Quality managers close to processing and beneficiation plants

Digital mine - future

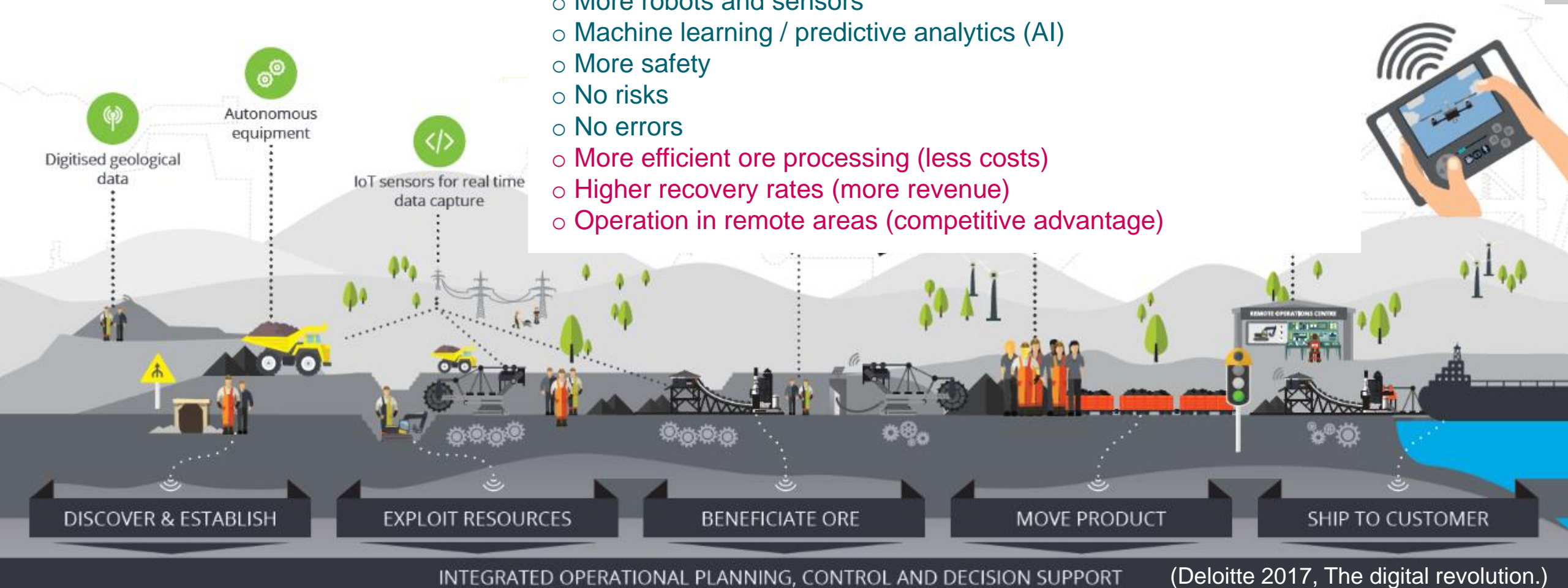
From ore discovery towards metal shipment



Digital mine - future

From ore discovery towards metal shipment

- Autonomous mining process
 - Less / no people
 - More robots and sensors
 - Machine learning / predictive analytics (AI)
 - More safety
 - No risks
 - No errors
 - More efficient ore processing (less costs)
 - Higher recovery rates (more revenue)
 - Operation in remote areas (competitive advantage)



Value predictive monitoring

Mining and metals industry - Examples

Time

EXAMPLE

**IN-SITU measurements
in oil well drill cuttings**

Lab reports on a side wall core suite cost about \$40k – \$70k and information is **untimely**. In-situ analysis will reduce lab costs plus result in **more efficient use** of the rig.

Save (per drilling rig)
>100k USD/y

Downtime

EXAMPLE

**MINERLOGY
for ore processing**

A copper mine suffered from stopping ore processing due to NOT checking mineralogy

- Ration **oxides/sulfides**
- Presence of hard minerals during **milling**
- Missing awareness of **clay minerals**

Loss
>70m USD/y

OPEX

EXAMPLE

**ACID consumption
for ore processing**

Reduction of **1kg H2SO4 per ton ore** due to accurately and online monitoring of **clay minerals** results of huge savings.

Save
>3m USD/y

Penalties

EXAMPLE

**MOISTURE
for shipment of Fe-ore**

1% moisture deviation from the agreed-upon moisture content results in a **\$0.8 per ton penalty**.
2% lower target for moisture content of iron ores will cost less profit.

Loss
>69m USD/y

Energy

EXAMPLE

**COKE consumption
for sinter production**

Saving in **coke consumption** (1kg coke / ton sinter) due to frequent and online monitoring of sinter raw mix and sinter represents a significant saving in energy.

Save
>1.4m €/y

www.malvernpanalytical.com