

Iron Ore and DRI Expert Analytical Kit

Contact us for more information

Overview

Malvern Panalytical offers a comprehensive turnkey solution for the elemental analysis of iron ore concentrates and direct reduced iron (DRI), designed to achieve the highest accuracies.

Components

1. Standards

A dedicated suite of standards will be assembled to specifically address the customer's list of elements and concentrations, utilizing:

- The WROXI CRM solution module,
- External Certified Reference Materials (CRMs),
- Synthetic standards specifically made for the customer
- Or a combination of the above



2. Fusion

- Optimized sample preparation recipes.
- Enhances the precision and reliability of the analysis.



3. XRF Method

- Calibration fully matched to the customers materials and compositions.
- Optimized measurement conditions for iron ore, concentrates and DRI.
- Method validation.

XRF
Method



SCAN QR CODE

For more information on how our Iron Ore and DRI Expert Analytical Kit can benefit your operations, please contact our sales team.

Benefits



- **Expertise:** Leverages Malvern Panalytical's extensive expertise in iron ore and DRI analysis.



- **Customization:** Includes customization options to meet specific customer needs.



- **Wide Coverage:** Tailored to cover the specific elements required by the customer.



- **Validated Solution:** Reduces production costs and improves price determination accuracy.

Applications

Analysis of iron ores/concentrates and DRI/HBI. Suitable for industries requiring accurate and reliable iron ore and DRI measurements.

<https://www2.malvernpanalytical.com/contactus>



Malvern Panalytical
a spectris company