

Ferro Alloys Turnkey Solution

Contact us for more information

Overview

Malvern Panalytical offers comprehensive turnkey solutions for the elemental analysis of ferro alloys, designed to achieve the highest accuracies.

Components

1. Standards

- Always customer specific: mix of international Certified Reference Materials and synthetic standards, tailored for your specific need.
- Covers a variety of ferro alloys, ensuring accurate and consistent measurements.



2. Fusion

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



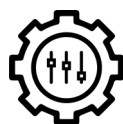
3. XRF Method

- Optimized measurement conditions for ferro alloys.
- Method validation.

XRF
Method



Benefits



- **Expertise:** Leverages Malvern Panalytical's extensive expertise in ferro alloys analysis.

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



- **Wide Coverage:** Covers a variety of ferro alloys, including FeMn, FeSi, SiMn, FeMo, FeNi, FeNb, FeP, FeTi, FeW, FeV, and more.



- **Validated Solution:** Tested on CRMs to prove "fit for purpose" and analytical performance, for defensible analysis.



SCAN QR CODE

For more information on how our Ferro Alloys Expert Analytical Kit can benefit your operations, please contact our sales team.

Applications

- Analysis of a variety of ferro alloys
- Suitable for industries requiring accurate and reliable ferro alloys measurements.

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



**Malvern
Panalytical**
a spectris company