

Base Metal Sulfides Turnkey Solution

Contact us for more information

Overview

Malvern Panalytical offers a comprehensive turnkey solution for the elemental analysis of base metal sulfides, designed to achieve the highest accuracies in a wide range of materials such as ores, concentrates and tailings.

Components

1. Standards

- Synthetic reference materials tailored for base metal sulfides analysis.
- Covers a variety of base metal sulfides (Cu, Zn, Pb, Mo), ensuring accurate and consistent measurements.



2. Fusion

- Optimized sample preparation recipes.
- Enhances the precision and reliability of the analysis.
- Protects Pt ware during aggressive oxidation process.



3. XRF Method

- retains sulfur in the bead
- Universal calibration over wide concentration ranges.
- Optimized measurement conditions for base metal sulfides.
- Method validation.

XRF
Method



SCAN QR CODE

For more information on how our Base Metal Sulfides Expert Analytical Kit can benefit your operations, please contact our sales team.

Benefits



- **Expertise:** Leverages Malvern Panalytical's extensive expertise in base metal sulfides analysis.



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



- **Wide Coverage:** Covers a variety of base metal sulfides, including Cu, Zn, Mo, Pb, and more.



- **Validated Solution:** Tested on external Certified Reference Materials (CRMs) to prove "fit for purpose" and analytical performance, for defensible analysis.

Applications

- Analysis of base metal sulfides in various matrices.
- Suitable for industries requiring accurate and reliable base metal sulfides measurements.

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



Malvern Panalytical
a spectris company