

# Rare Earth Elements Turnkey Solution

Contact us for more information

## Overview

Malvern Panalytical offers a comprehensive turnkey solution for the elemental analysis of all Rare Earth Elements (REE) in ores and concentrates at highest accuracies in a vast range of materials.

## Components

### 1. Standards

- Synthetic calibration materials tailored for REE analysis
- All 17 REE covered, in total 40 elements included
  - REE analysis in a wide variety of materials



### 2. Fusion

- Optimized Sample preparation recipes
- Enhances the precision and reliability of the analysis



### 3. XRF Method

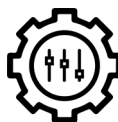
- Universal calibration over wide concentration ranges
- Optimized measurement conditions for REE
- Method validation

**XRF**  
Method

## Benefits



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



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



- **Wide Coverage:** Covers 40 elements including all 17 REEs, across wide concentration ranges.



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

## Applications

Analysis of rare earth elements in various matrices. Suitable for industries requiring accurate and reliable REE measurements.

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

Compound	Max conc.% *)
BaO	40
La2O3	40
CeO2	60
Pr6O11	20
Nd2O3	25
Sm2O3	15
Eu2O3	3
Gd2O3	5.1
Tb4O7	1.5
Dy2O3	7.5
Ho2O3	2.1
Er2O3	7.5
Tm2O3	1.5
Yb2O3	7.5
Lu2O3	1
Hf2O2	10
Ta2O5	3
PbO	10
ThO2	5
U3O8	1.5

\*) minimum concentrations start at the LOQ, typically ~0,01 %



### SCAN QR CODE

For more information on how our REE Canned Solution can benefit your operations, please contact our sales team.



**Malvern Panalytical**  
a spectris company