





Mining QC Strategy Checklist XRF vs. XRD


Quick Reference for Lab & Field Use

 Purpose of the analysis	XRF	XRD
Elemental anomalies (e.g. presence of Co)	✓	
Elemental composition (e.g. % Ni, ppm REE)	✓	
Mineral identification (e.g. spodumene or petalite)		✓
Mineralogical composition (e.g. % chalcopyrite)		✓
Trace element detection	✓	
Detect oxidation stages (e.g. Fe ²⁺ vs Fe ³⁺)		✓
Distinguish crystalline vs amorphous		✓
Clay identification (kaolinite vs illite)		✓

 Sample requirements	XRF	XRD
Loose powder (dry & fine, < 75 μm)	✓	✓
Pressed pellets	✓	✓
Fused beads	✓	
Liquids	✓	
Sample homogeneity critical	✓	✓

 Operating context	XRF	XRD
In-field (portable)	✓	
Laboratory	✓	✓
Process (real-time)	✓	
Automation	✓	✓

 Stage in the mining process	XRF	XRD
Exploration phase	✓✓✓	✓✓
Extraction operation	✓	
Process control (e.g. blending, geometallurgy)	✓✓✓	✓✓✓
Research	✓	✓✓✓
Environmental screening	✓✓✓	✓

 Tips and Pitfalls
!! Calibrate the instrument regularly using CRMs similar in composition to your ore.
!! XRD quantifies amorphous content but cannot identify its composition.
✓ Use XRF and XRD together for optimal geometallurgical or process analysis.