Portability and precision: from handheld to lab analyzers, we have you covered

The iron and steel manufacturing processing

Stage 1 **Raw Materials Processing**

Weight verification

Elemental analysis using OES, XRF, XRD

Handheld XRF analysis

Online elemental analysis of coal

Density, flow and level measurement

Online gas and moisture analysis

Radiation measurement and protection

Particulate monitoring

Laboratory informatics

Stage 2 **Steel Making and Casting**

Elemental analysis using OES, XRF, XRD

Handheld XRF analysis

Elemental analysis including carbon content with LIBS

Flow measurement

Online gas analysis

Radiation measurement and protection

Particulate monitoring

Laboratory informatics

X-ray source/flux stability

OES Inclusion monitoring

Stage 3

Billet/Ingot/Final Product inspection

Certificate of analysis using OES

Elemental analysis including carbon content with LIBS

Handheld XRF analysis for product identification

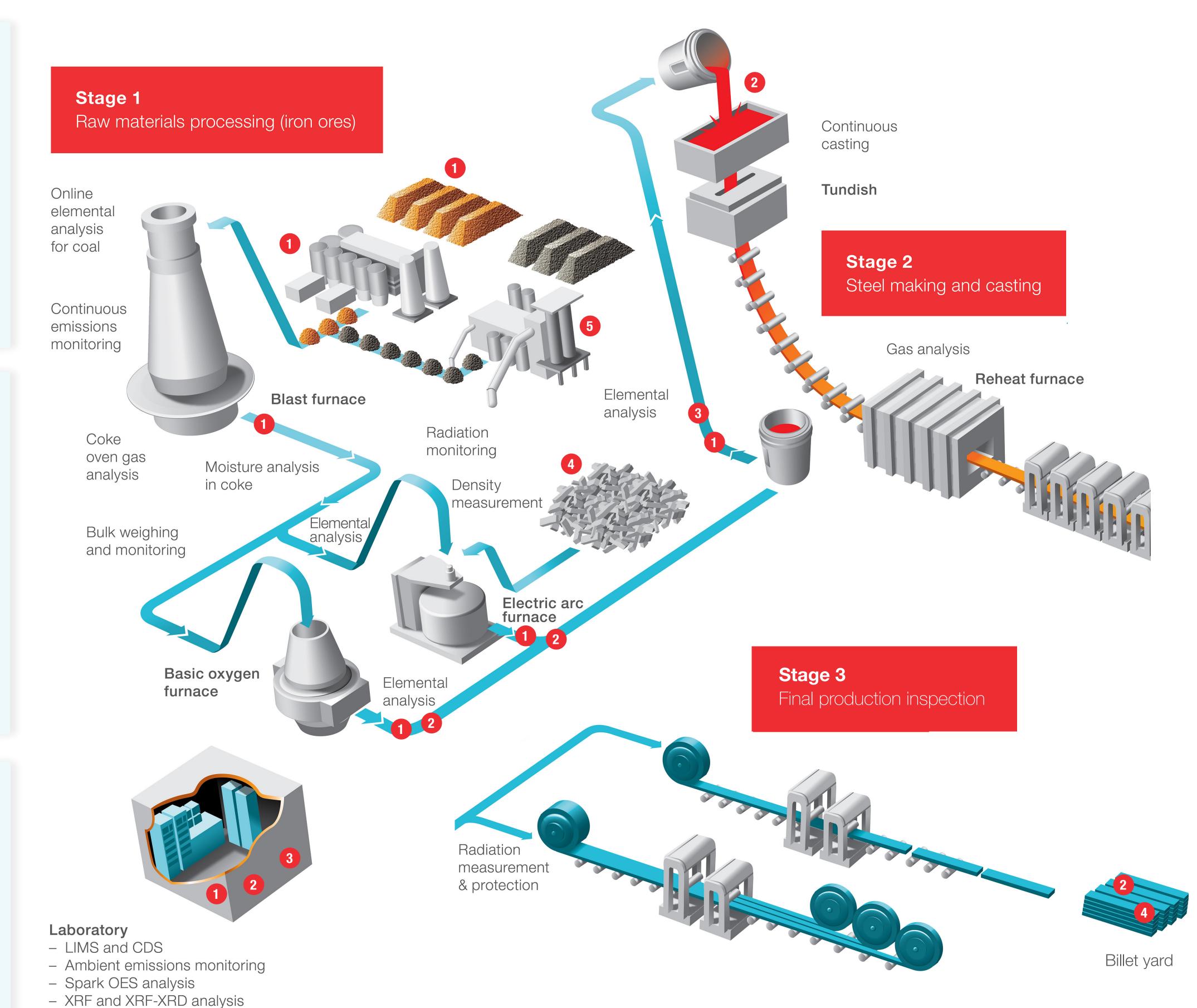
Laboratory informatics

Gas analysis

Radiation measurement and protection

Data acquisition & management

Particulate monitoring



Thermo Fisher SCIENTIFIC

Thermo Fisher products are useful at this stage



1 ARL 9900 XRF Analyzer





2 ARL iSpark Plus and ARL easySpark OES Analyzers



3 ARL SMS Automated Analyzers





4 Niton XL5 Plus and Niton Apollo Handheld Analyzers



6 CB Omni Agile Analyzer

Learn more at thermofisher.com/niton-arl

Automated sample preparation