

# Small spot mapping

X-ray fluorescence as a process control tool for chemical characterization of macro-segregation occurring during continuous casting in steel

## Zetium

### Introduction

The continuous casting of slabs is aimed to produce a product with a certain chemical composition, geometry and surface quality, without any or a minimum acceptable level of external and internal defects. One of the most unpredictable defects of the slabs is centerline segregation, which has undesired effects on further processing of the slabs and hence on the possible uses of the final product.

Segregation can occur at two different scales: micro-segregation, leading to variations in composition in the micrometer range and macro-segregation, involving chemical variations over the length of the slab (Figure 1 and Figure 2). Micro-segregation can be removed by homogenization heat treatments, but it is practically impossible to remove macro-segregations. It is of critical importance to identify macro-segregation to be able to predict its effect on physical and chemical properties.

X-ray fluorescence systems equipped for small spot analysis and elemental distribution mapping can be employed to effectively control segregation during the production process. This application note demonstrates the capability of practical, fast elemental distribution mapping of two steel samples showing chemical segregation, analyzed for quality control purposes.



### Zetium wavelength dispersive XRF spectrometer

The Zetium spectrometer is a fully integrated XRF analyzer, consisting of an advanced analytical heart, integrated X-Y sample handler and software.

Sensitive optics and outstanding stability provide accuracy and precision.

The SuperQ<sup>®</sup> analysis software, which is delivered with the system, has an intuitive user interface and makes accurate quantitative analysis easy to perform.

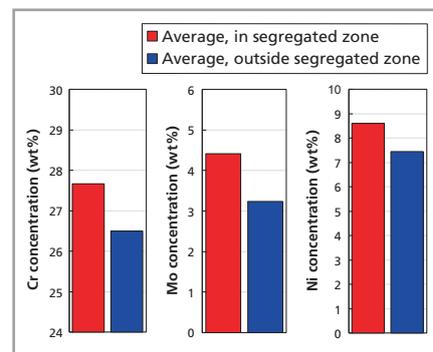


Figure 1. Variations in Cr, Mo and Ni concentrations between bulk slab and segregated zones occurred at casting

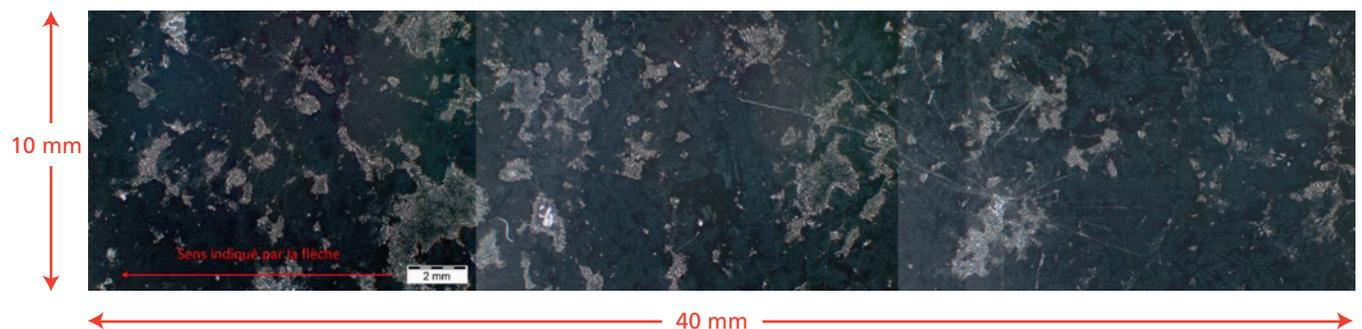


Figure 2. Scanning electron microscopy (SEM) micrograph, highlighting the surface of a slab sample after etching. White regions indicate segregation of Cr, Ni and Mo from the bulk.

### Instrumentation and software

Measurements were performed using a Zetium XRF spectrometer configured with a 4 kW Rh-anode SST R-MAX X-ray tube and the WD core for WDXRF analysis. To enable the spot analysis and mapping capability, the spectrometer was equipped with an ED core, high-precision translation mechanics for sample positioning and a high-resolution camera for sample imaging. Using the ED core offers the advantage of performing simultaneous analysis for all elements present, whereas using the WD core would imply already knowing which elements are present and repeating each spot analysis for the number of elements needed to be mapped.

All measurements were conducted using the state-of-the-art SuperQ software package.

### Sample preparation

Two steel samples named Brame 1 and Brame 2 are the object of this study (Figure 3). Samples were placed into a special mapping sample holder and imaged using a high-resolution camera. The samples were then loaded into the measurement position using a special turret mechanism and analyzed using SuperQ software.

### Measurement conditions

In order to map the elemental distribution of a small area, a conventional calibration on a small spot was set up using 11 standards, including six NiFeCo setup samples and two standards from the ECRM series (298/1 and 289-1). The two samples were measured against the calibration using a step size of 250 microns and respectively analyzed in 1269 spots for sample Brame 1 and 1932 spots for sample Brame 2.

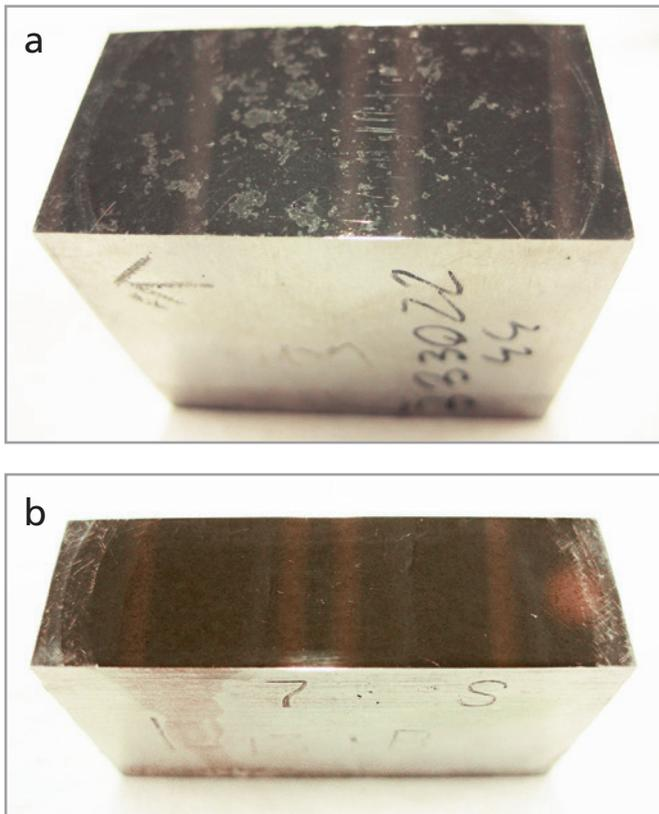


Figure 3. a) Sample Brame 1 analyzed for a 6.5 mm x 11.5 mm mapped area; b) sample Brame 2 analyzed for 10 mm x 11 mm mapped area

### Results

Figure 4 shows calibration lines for Fe, Ni, Cr and Mo with RMS and K values as a result of a NiFeCo calibration performed on small spot mapping (0.5 mm).

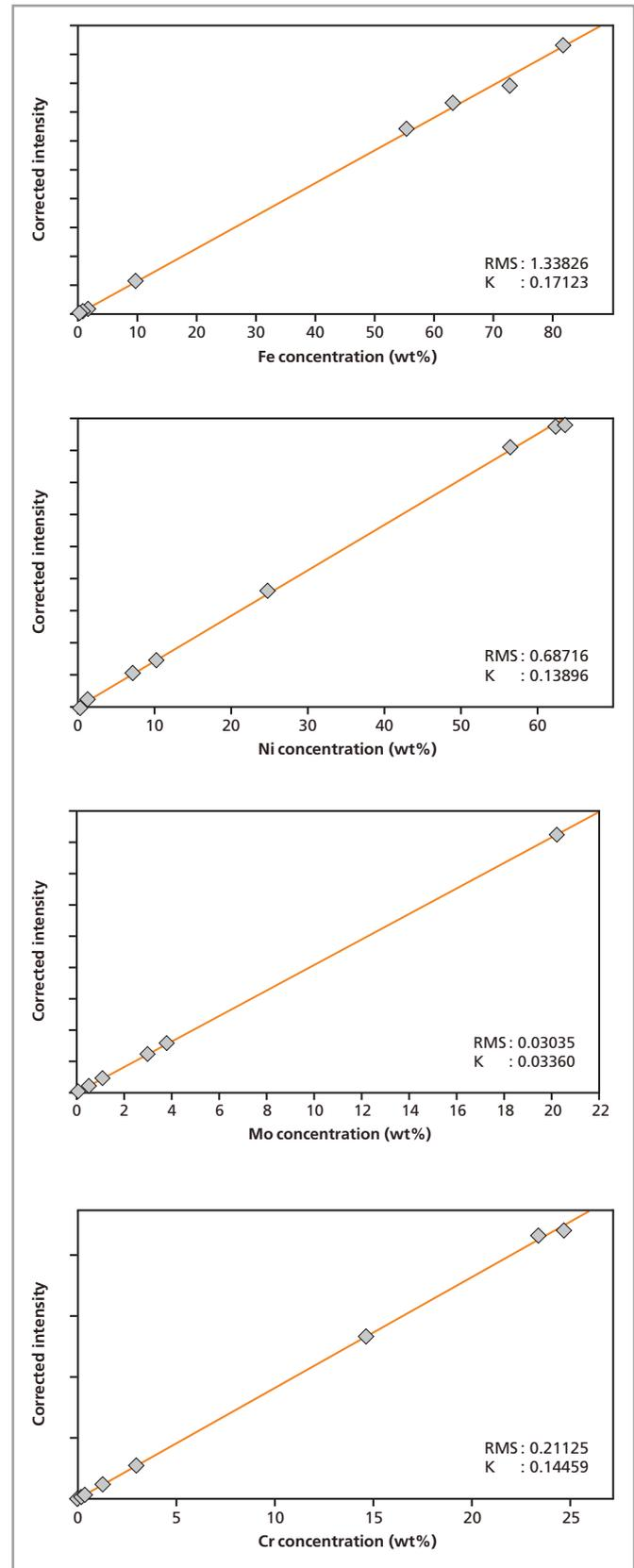


Figure 4. Calibration lines on 0.5 mm spot size for Fe, Ni, Mo and Cr

Compositional images (Figure 5) show the variations in elemental distribution for the two samples in the mapped areas, clearly more evident for sample Brame 1. Compositional variations measured for the mapped areas of both samples are reported in Table 1.

The elemental distribution mapping can help to identify and localize the chemical segregation, which can ultimately alter chemical and physical properties of the final product and therefore cause failure to meet product specifications.

a

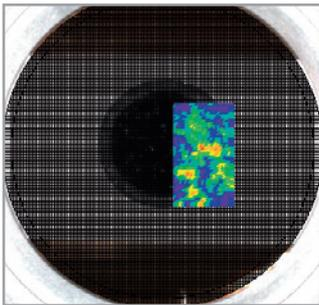
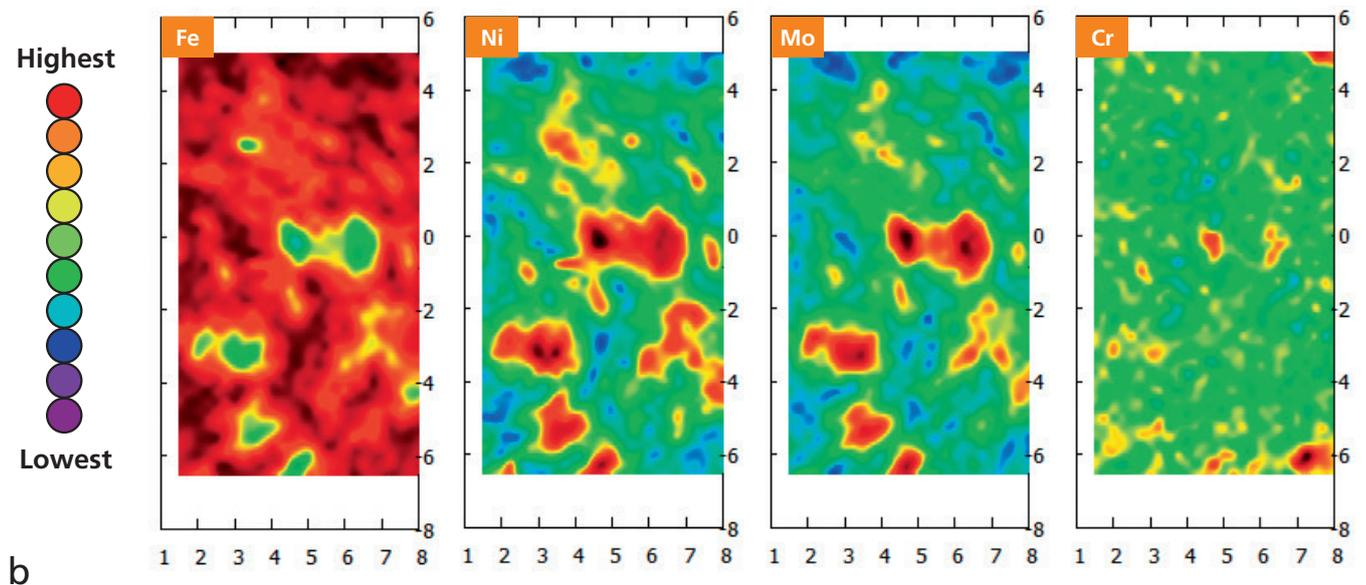


Table 1. Variation in elemental concentration in the mapped areas for samples Brame 1 and Brame 2.

Concentration (%)	Fe	Ni	Mo	Cr
Brame 1				
Min.	57	6.5	3.6	24.6
Max.	63	9.5	4.8	26.4
Brame 2				
Min.	67.8	9.7	1.85	15.9
Max.	70	10.7	2.15	16.5



b

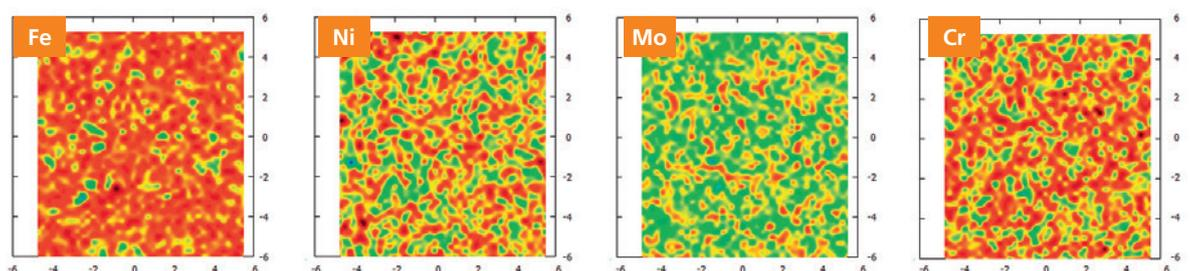
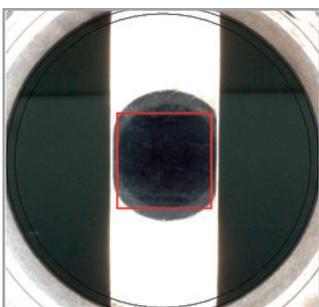


Figure 5. 2D contour plots for Fe, Ni, Mo and Cr in samples Brame 1(a) and Brame 2 (b) presenting and macro-segregation and a high-resolution image of the area mapped

## NiFeCo-FP specifications

### 1. Setup samples: elements and concentration ranges (wt %)

Al	0.01	–	6
Si	0.01	–	2.4
P	0.01	–	0.3
S**	0.01	–	0.1
Ti	0.01	–	4
V	0.01	–	3
Cr	0.01	–	24
Mn	0.01	–	15
Fe	0.01	–	78
Co	0.01	–	63
Ni	0.01	–	64
Cu	0.01	–	30
Y*	0.01	–	0.3
Zr	0.01	–	0.4
Nb	0.01	–	6.3
Mo	0.01	–	21
Hf*	0.01	–	1.3
Ta*	0.01	–	7.2
W	0.01	–	15
Re*	0.01	–	5.4
Pt*	0.01	–	0.3

Elements indicated with

\* were calibrated using in-house reference materials.

\*\* The high sulfur concentration sample (0.1 wt%) is the glass monitor.

### 2. Presentation:

- Robust aluminium carrying case containing setup samples plus user information and concentrations on CD-ROM
- 6 calibration setup samples, calibrated against more than 120 CRMs
- 2 drift monitors

### 3. Materials that can be analyzed:

- Stainless steel
- High-speed steel
- Low-alloy steel
- Tool steel
- Mild steel
- Nimonic steels
- Inconel
- High-Mn steel
- Ni alloy

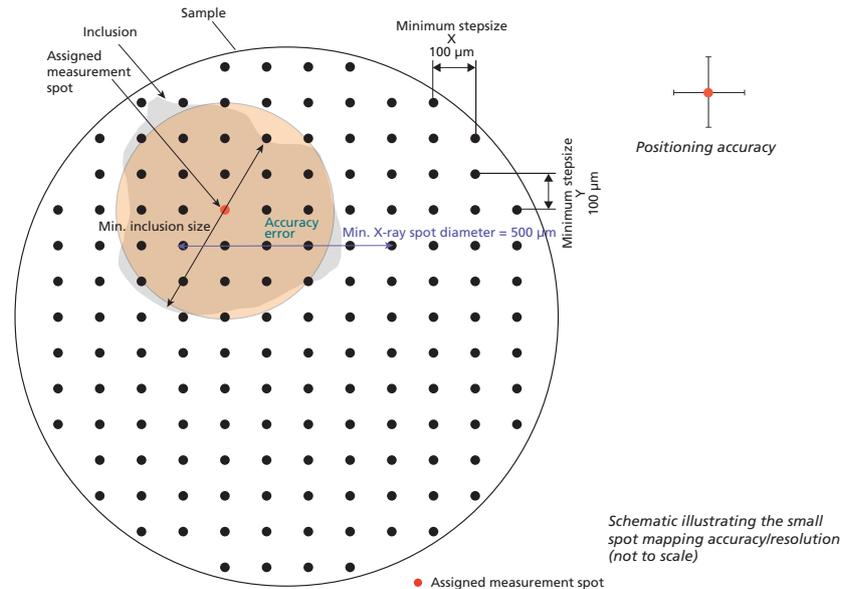


Note: NiFeCo-FP is not intended for the analysis of slow-cooled cast alloys.

### Setup of small spot analysis and mapping

The small spot analysis and mapping module can accommodate samples with diameters up to 35 mm. The measurement spots have a diameter of 500 µm (FWHM) and a minimum intervening step size of < 100 µm (see schematic drawing), resulting in a maximum of 24,000 spots per sample.

Sample preparation is simple and samples are mounted in a dedicated holder, specially designed for irregularly shaped samples of varying size. This holder incorporates a sample clamping device, ensuring that small samples can be placed into the instrument without compromising positioning accuracy or damaging the sample.



### Conclusions

The results clearly demonstrate that the integration of EDXRF (equipped to perform small spot analysis and mapping) in the Zetium spectrometer delivers fast and accurate analysis of a steel sample which experienced macro-segregation during casting. This analysis performed as routine test may be helpful to improve process control.

The possibility to perform small spot analysis and mapping using the ED core enables fast simultaneous analysis of all elements present without compromising the analytical performance of the WDXRF optics. Furthermore it is considerably faster than using a conventional WDXRF, where each element present needs to be measured sequentially. The inclusion of the ED core has other advantages, such as fast sample screening, with or without Omnian, and identification and flagging of unexpected elements in process control.

**Acknowledgement:** We thank Aperam, Isbergues in France for providing samples used in the present application

### Global and near



**PANalytical B.V.**  
Lelyweg 1, 7602 EA Almelo  
P.O. Box 13, 7600 AA Almelo  
The Netherlands  
T +31 (0) 546 534 444  
F +31 (0) 546 534 598  
info@panalytical.com  
www.panalytical.com

**Regional sales offices**  
Americas  
T +1 508 647 1100  
F +1 508 647 1115

Europe, Middle East, Africa  
T +31 (0) 546 834 444  
F +31 (0) 546 834 499

Asia Pacific  
T +65 6741 2868  
F +65 6741 2166