

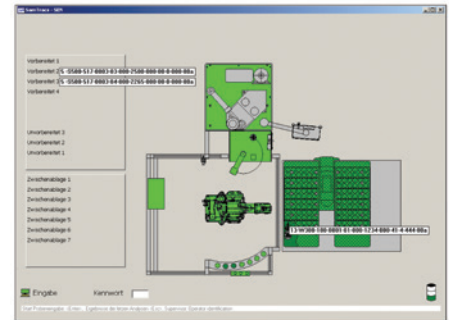


# The automated laboratory

## Charge correction

In this automation the PANalytical Charge Correction software is integrated with the process flow. From customized sample announcement to full metals accounting the complete production process is integrated in SamTracs. After sample announcement the samples are milled and analyzed on a PANalytical Magix FAST. On completion of the analysis charge correction is started. Charge correction determines based on the furnace parameters, actual furnace weight and the alloy whether or not correction is needed (and possible). The operator still has the possibility to fine tune the proposal. All data is stored and various reports can be printed based on production codes or over periods of times. As the complete process flow is integrated also the metal accounting is based on the data collected by SamTracs. Also a connection to the plants ERP (SAP) implementation is part of the project.

- Small setup < 10 m<sup>2</sup>
- Charge correction integrated
- Completely adapted to the end-user requirements
- Customer specific sample announcement screens



## The concept

Within an automated laboratory, the complete sample transportation preparation-and analysis process can be automated. Our automation projects can cover all steps involved in the process control and quality control, summarized as follows:

- Sample treatment before sending
- Sample transportation to and within the laboratory
- Preparation of the sample material
- Sample analysis and results distribution
- Control of the automated laboratory
- Container

The automated laboratory is a multi disciplinary strategy to increase the productivity and reduce lab process cycle times of our customers by making the best use of technology. Based on customer needs, the automated laboratory is designed for the customer and with the customer.

PANalytical is not only the world's leader supplier for X-ray instruments, but recognize also the added value of X-ray systems in an integrated automated laboratory. Since 1994 we have built automation projects, surpassing 90 installations worldwide.

## 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