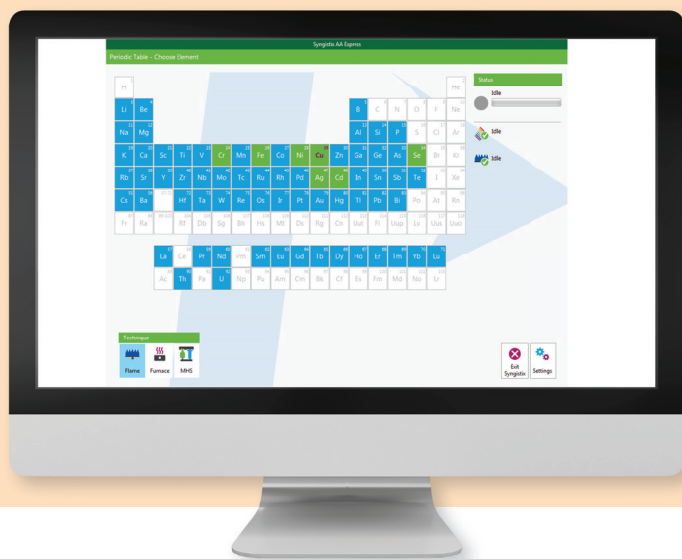


Atomic Absorption

Key Features:

- Streamlined workflow
- Optimized user interface
- Simplified reporting
- “At Your Service” function for efficient troubleshooting



Syngistix for AA Express Software – Only Three Steps to Reliable Results

Syngistix™ for AA Express, the latest addition to PerkinElmer’s Syngistix atomic spectroscopy software platform, is the newest instrument control software option for the PinAAcle 500, 900F and 900H AA spectrometers. Leveraging the icon-based design of

the Syngistix platform, Syngistix for AA Express delivers an even more streamlined workflow with only three steps to results, so novice and experienced users alike can truly enjoy quick hands-on and seamless operation. Syngistix for AA Express reduces complexity, speeds analysis time, and provides an unmatched user experience.

Syngistix for AA Express software supports PerkinElmer’s high-performance PinAAcle 500 and 900F flame systems, as well as the PinAAcle 900H flame/furnace system. These proven systems come standard with continuum source background correction and the flexibility of flame only or flame/furnace combination, delivering intuitive, highly efficient solutions that simplify your journey from sample to results—even with challenging matrices. They feature cutting-edge fiber optics for better detection limits and superior reliability. Ideal for laboratories running routine applications that require flame or both flame and furnace AA, the PinAAcle AA spectrometers and Syngistix for AA Express combine to deliver extremely easy-to-operate yet stable AA systems.

Optimized Interface, Faster Results

Whether you are an experienced or first time user of AA, Syngistix for AA Express will take you from setup to analysis seamlessly, in just three easy steps:

1. Choose your element
2. Set up your lamp
3. And analyze your samples

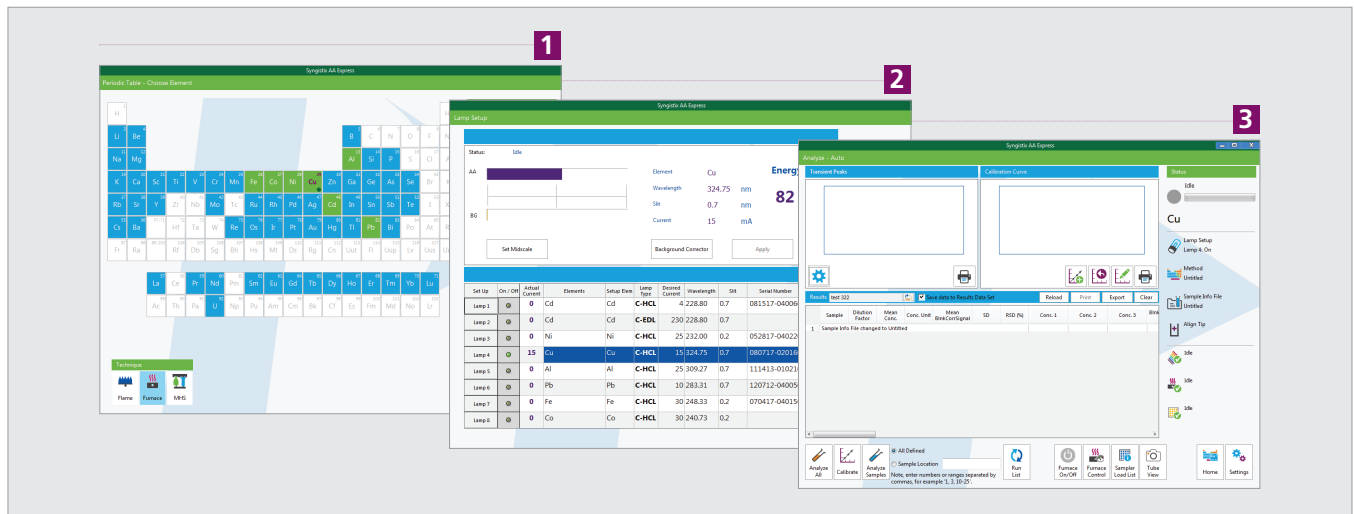


Figure 1. Example captures of the three steps with Syngistix for AA Express software.

Plus, the user interface has been optimized for seamless operation, further streamlining the process:

- The side bar (A) offers shortcuts to the different steps of the analysis.
- The instrument status bar (B) provides real-time information of the instrument during the analysis.
- The icon bar (C) at the bottom displays most of the functions to avoid switching back and forth between different interfaces.

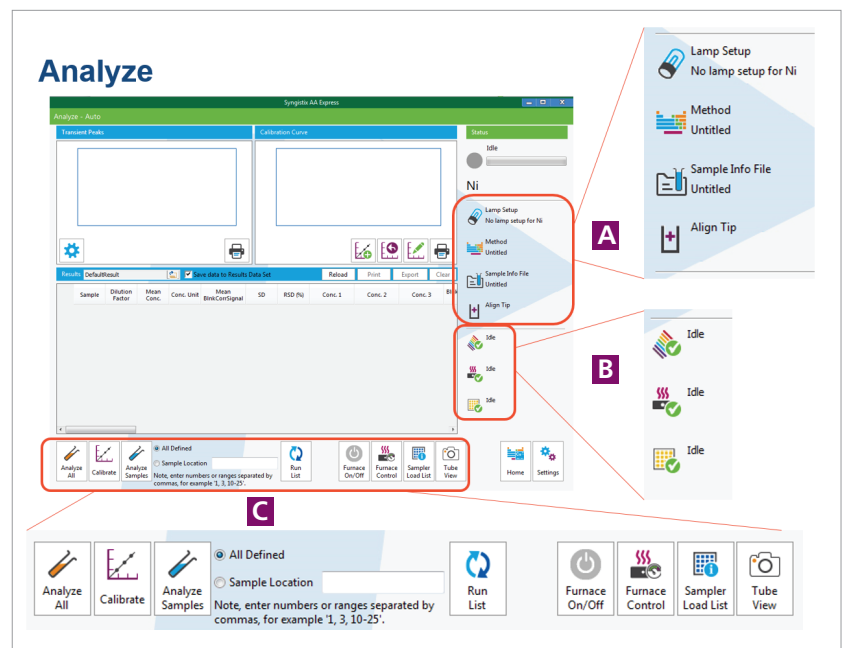


Figure 2. Analyze screen in Syngistix for AA Express software, displaying the easy navigation.

Simplified One-Page Report

At the end of a run, printing out a comprehensive report with information from different windows can be time consuming and sometimes inefficient. With Syngistix for AA Express, reporting also is more streamlined – when your run is complete, the software will automatically generate a simplified report including Method, Calibration, and Results, all on a single page.

“At Your Service” Function for Easy and Efficient Troubleshooting

In those cases where you need assistance in order to address a technical issue, Syngistix for AA Express software features a unique “At Your Service” function that allows the user to collect, in just one click, all of the information needed to facilitate system diagnostics and maintenance:

- Log information of base modules
- Result logs
- Lamp information, including wavelength, current, hours of usage, etc.
- Current active method and sample information
- Instrument information and logs
- Error logs
- Latest snapshots of furnace camera when the tip is in the tube, if furnace technique is in use

The software automatically packages all of the necessary files. At this point, you can choose what to do with the package – you may wish to forward it to your local Service Support Engineer, so that your situation is addressed efficiently, and you are able to return to testing quickly.

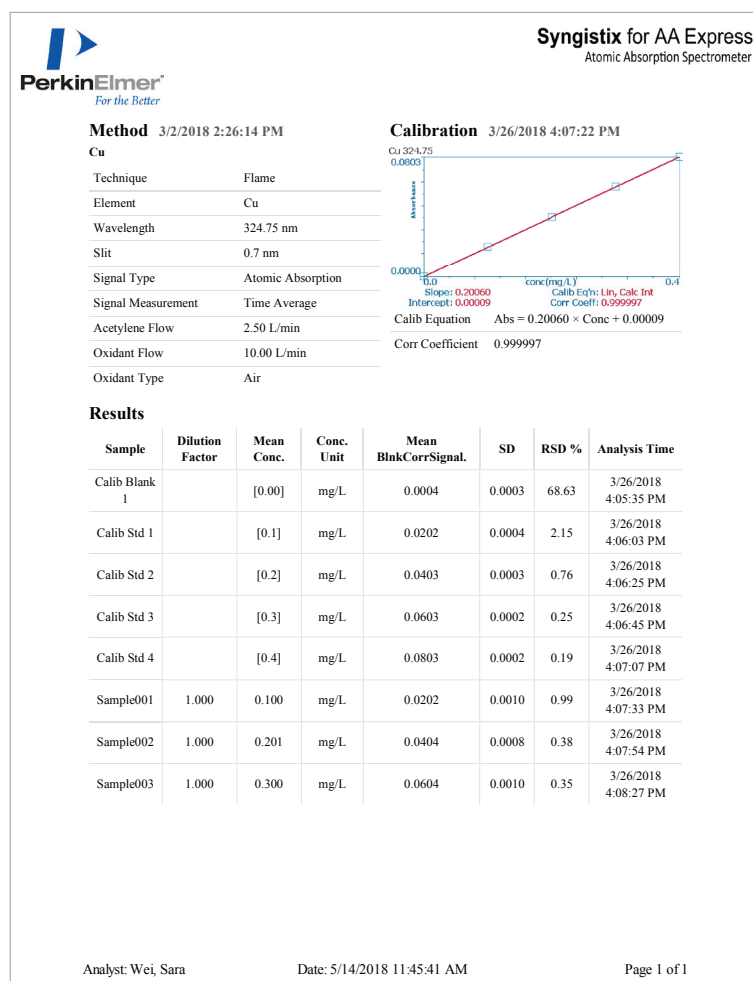


Figure 3. Example single-page report generated by Syngistix for AA Express software.

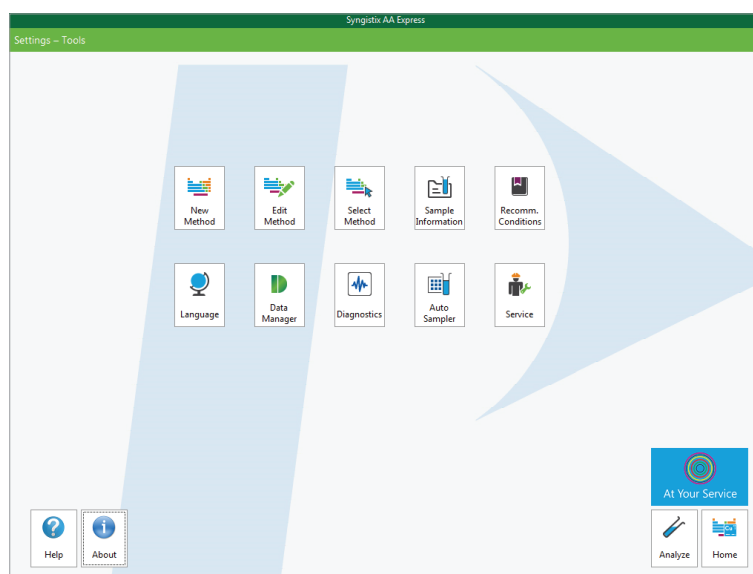


Figure 4. “At Your Service” button in Syngistix for AA Express software.

Powerful Software Options for Diverse Needs

We understand that different labs have different requirements. For the PinAAcle 500, 900F and 900H AA spectrometers, you can select from a full suite of powerful software tools, starting with the innovative and easy-to-use Syngistix for AA Express, the more comprehensive Syngistix for AA standard software, or you can add on the special Syngistix Enhanced Security™ option that fulfills the special needs of highly regulated laboratories, such as those operating under the rules of 21 CFR Part 11.

Feature	Syngistix for AA Standard	Syngistix for AA Express
Flame and HGA Control Capability	Yes	Yes
TubeView Furnace Camera	Yes	Yes
Database Manager	Yes	Yes
Method Editor	Yes	Simplified
FIMS/FIAS/FAST Flame Support	Yes	
Data Reprocessing	Yes	
Automatic QC	Yes	
Multi-element Methods and Analysis	Yes	
Streamlined UI Workflow for First-Time User		Yes
Simplified Report Format		Yes
Supports Syngistix Enhanced Security Add-On	Yes	

Ordering Information

Model	Part Number
PinAAcle 500 Flame AA Spectrometer with Touchscreen	N3180011
PinAAcle 900F Flame AA Spectrometer	N3200051
PinAAcle 900H Combined Flame/ Furnace AA Spectrometer	N3200073
Syngistix for AA Express Software	N1010308
Syngistix for AA Standard Software	N1010302
Syngistix for AA Enhanced Security Software	N1010314

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2021, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

297956 (014126_01)

PKI