



# RESULTS IN 3...2...1

## Meet Syngistix™ for ICP software for the Avio® Max ICP-OES series

For most labs conducting routine elemental analysis, time constraints and complex systems are the biggest challenges to overcome. Simplified and streamlined workflows are critical to success.

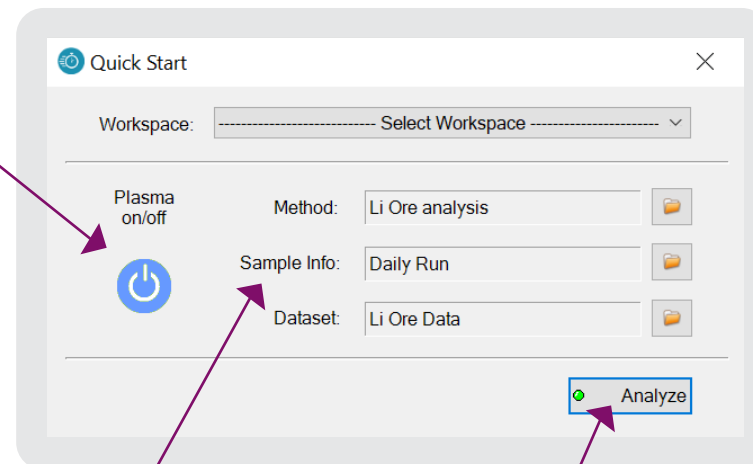
Syngistix for ICP software for the Avio Max ICP-OES series is engineered with an assortment of smart features, streamlining analysis with only **three clicks** from sample to results!

- Click 1: Ignite the plasma
- Click 2: Load the sample sequence
- Click 3: Analyze

**Simple – easy – fast!**

For more information, visit [www.perkinelmer.com/SyngistixICP](http://www.perkinelmer.com/SyngistixICP)

CLICK 1



CLICK 2

CLICK 3

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

  
**PerkinElmer**  
For the Better

For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2022, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.