

Great Technology Makes All the Difference

What's more, the Avio series ICP-OES leverages a host of other unique and proprietary technologies that deliver seamless operation, so you can focus on what matters most – your results.

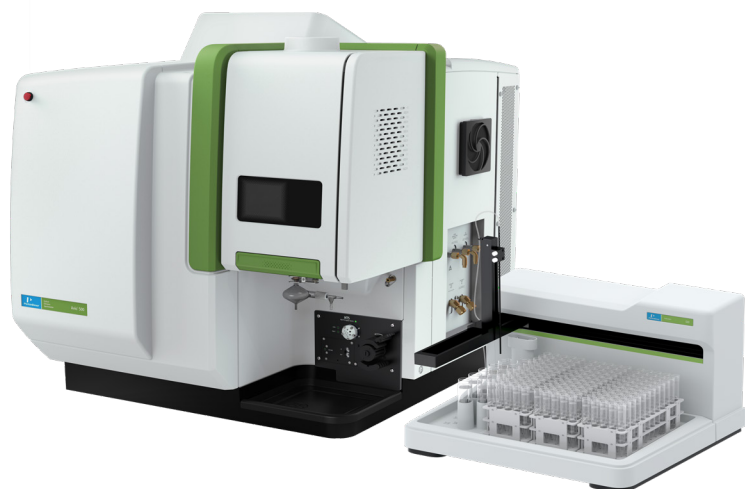
Our patented **Flat Plate™** Plasma Technology delivers a more robust, stable plasma, and uses half the argon of load-coil ICP systems (at least 50% less than other systems).

Vertical Plasma Torch Design delivers unrivaled matrix flexibility and reduced maintenance. Its design lets you make adjustments quickly and easily, even while the system is running. The removable injector is independent of the torch, simplifying maintenance and reducing the potential for breakage.

Dual View delivers full wavelength coverage, with no loss of light or sensitivity. Even elements at high (>500 nm) or low (<200 nm) wavelengths can be measured with complete confidence in the same run.

Color PlasmaCam™ offers continuous viewing of the plasma, simplifying method development and enabling remote diagnostic capabilities.

And our **Syngistix™** software provides an intuitive interface that mirrors your workflow, with left-to-right icon-based navigation that guides you through your analysis – and because it's a cross-platform solution, it eases the transition from technique to technique. What's more, it features preloaded testing methods, eliminating the need for method development in many applications.



HTS and the S20 Series Autosamplers integrate with the Avio ICP-OES, Syngistix Software delivering seamless operation and faster turnaround.

For more information about our Avio ICP-OES high-throughput solution, visit www.perkinelmer.com/AvioHTS

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 © 2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

75984

PKI

DRAFT