

Great Technology Makes All the Difference

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

The **All Matrix Solution (AMS)** system lets you run samples with high total dissolved solids (TDS) and without manual dilution. And the solid-state free-running (RF) generator delivers superior plasma control and stability to handle these samples.

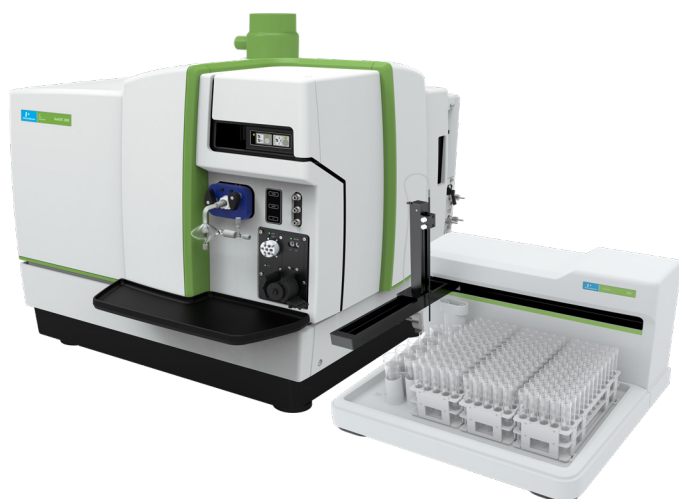
LumiCoil™ is a revolutionary RF coil that's guaranteed for life and requires no water or gas cooling. Coupled with an innovative, patented plasma generator, it enhances analytical performance as well as reliability while lowering operating costs.

The combination of **Triple Cone Interface and Quadrupole Ion Deflector** ensures that unionized material never reaches the cell or detector. This combination eliminates the need for maintenance behind the cones without sacrificing stability or accuracy.

Universal Cell Technology™ delivers best-in-class interference removal – and real confidence in your data. Its three modes of operation (standard, collision, and reaction) combine collision-cell simplicity with reaction-cell efficiency, all in a single instrument.

The exclusive **Extended Dynamic Range** functionality allows you to measure elements with low and high concentrations in the same sample in the same run. This extends your linear dynamic range by up to twelve orders of magnitude, optimizing your productivity and preserving the lifetime of the detector.

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. Plus, it features preloaded testing methods, eliminating the need for method development in various applications.



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

For more information about our NexION ICP-MS high-throughput solution, visit www.perkinelmer.com/NexIONHTS

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.

75973

PKI

DRAFT