



# ICP-MS RESEARCH CONTEST

## Could a FREE Multi-Quadrupole ICP-MS System Bring Your Analysis to the Next Level?

We know the challenges – and aspirations – of today's research labs. And that's the whole idea behind our **ICP-MS Research Contest**. It's our chance to find worthy inorganic research projects to sponsor. And it's *your* chance to take your trace-level analysis research to new heights with the use of the multi-award winning NexION® 5000 Multi-Quadrupole ICP-MS.

Simply submit your research project plan that shows how a NexION® 5000 ICP-MS could help in your research. Your project could be chosen to receive a **FREE\* PerkinElmer NexION® 5000 Multi-Quadrupole ICP-MS system** for an entire year.

\*Terms and conditions apply.



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)

### 3 Easy Steps to Apply

1. Visit [perkinelmer.com/AcademicContestUSD2021](http://perkinelmer.com/AcademicContestUSD2021) to explore our ICP-MS technology and learn more about the **research contest** program.
2. **Submit your project plan** and show how a NexION® 5000 ICP-MS could help in your research.
3. **Reap the rewards!**  
A single grand prize winner will receive a FREE NexION® 5000 ICP-MS system for one year to complete their research. A single runner up will be awarded a 50% discount on a new ICP-MS instrument. All eligible participants will receive a 35% discount on a new ICP-MS instrument.

*Entries are due October 1, 2021. Learn more details and apply today!*

**APPLY NOW**



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