



LC 300 HPLC & UHPLC RESEARCH *CONTEST*

Could a Free HPLC or UHPLC 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 **HPLC & UHPLC Research Contest**. It's our chance to find worthy research projects to sponsor. And it's your change to take your chromatography analyses to new heights with the use of the new LC 300 HPLC or LC 300 UHPLC System. Simply submit your research project plan telling us how a LC 300 HPLC or LC 300 UHPLC could help in your research. Your project could be chosen to win the Grand Prize of a **FREE* PerkinElmer LC 300 HPLC or LC 300 UHPLC system** for a year or win big discounts in our runner up prizes.



Applying Is Easy as 1, 2, 3

1. Visit www.perkinelmer.com/AcademicContestEUR2021 to explore our LC solution technology and learn more about the research contest program.
2. **Submit your project plan** and tell us how a LC300 HPLC or LC 300 UHPLC could help in your research.
3. **Reap the rewards!** Two Grand Price Winners will receive either an LC 300 HPLC or LC 300 UHPLC system for one year to support their research. A runner up will be awarded a 50% discount on any new PerkinElmer HPLC or UHPLC system. All eligible participants will receive a voucher for a 35% discount on any new PerkinElmer HPLC or UHPLC system.

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

*Terms and conditions apply. Participating countries: Austria, Belgium, Croatia, Czech Republic, Denmark, Estonia, Finland, France, Germany, Greece, Hungary, Ireland, Italy, Latvia, Liechtenstein, Lithuania, Luxembourg, Monaco, Netherlands, Norway, Poland, Portugal, Romania, Serbia, Slovakia, Slovenia, Spain, Sweden, Switzerland, Turkey, UK (England, Scotland, Wales & Northern Ireland)

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.