



# POWER UP YOUR LAB

## Maximize Your Grant Funding

### Take advantage of our LABStartUp Grant Match

**LAB**StartUp

At PerkinElmer we are committed to supporting the discovery and advancements in science programs and understand the demand for reliable systems in budget conscious laboratories. Whether you are teaching the next generation of scientists or doing applied or fundamental research and development, we are offering you the opportunity to maximize the return on your investment with our **Grant Match Program** that could help you fund your laboratory.

Whether you are interested in atomic spectroscopy, chromatography or molecular spectroscopy, you will have the opportunity to build your teaching lab and leverage your grant funding to create a state-of-the-art lab environment for your students and faculty.

- 50% grant match = 33.33% instrument discount
- 60% grant match = 37.50% instrument discount
- 70% grant match = 41.20% instrument discount
- 80% grant match = 44.40% instrument discount
- 100% grant match = 50.00% instrument discount

*Qualifying Products must be purchased at list price only, and all products must be included on a single purchase order. Offer applies to qualifying products only.*

### Take advantage of this fantastic opportunity to start your lab with best-in-class technology

#### Step 1: Select the number of instruments your lab requires and PerkinElmer will match your grant.\*

- 1 instrument – 50% grant match\*
- 2 instruments – 60% grant match\*
- 3 instruments – 70% grant match\*
- 4 instruments – 80% grant match\*
- 5 instruments – 100% grant match\*

#### Step 2: Select the instruments your lab requires to receive your grant match percentage from Step 1.\*\*

- NexION® 1000, 2000, 5000 ICP-MS
- Avio® 220 Max, 550 Max, 560 Max ICP-OES
- PinAAcle™ 500, 900 AA
- TGA 4000, 8000
- DSC 4000, 6000, 8000, 8500
- STA 6000, 8000
- CHN 2400 Organic Elemental Analyzer
- LAMBDA™ 365, 465, 850+ UV/Vis Spectrophotometer
- LAMBDA™ 1050+ UV/Vis Spectrophotometer
- Spectrum™ Two FT-IR Spectrometer
- Spectrum™ 3 MIR/NIR Spectrometer
- Spotlight™ 150, 200, & 400 FT-IR
- FL 6500, 8500 Fluorescence Spectrophotometer
- Clarus® 590, 690 GC
- Torion® T-9 Portable GC/MS
- TurboMatrix™ System 350 ATD & Head Space
- QSight® Triple Quad 120, 210, 220, 420 LC/MS/MS
- Flexar & LC 300 HPLC/UHPLC

\* Certain restrictions apply. Offer is valid only with Qualifying Products orders. Applicable to quotes with validity through December 31, 2022 and may not be applied to existing, pending, or prior orders. Offer cannot be combined with any other promotions, discounts, volume purchase agreements, or other existing contracts. Used, refurbished or demonstration instruments do not qualify. Service parts, service contracts, training, software, accessories, consumables, and other products do not qualify for this offer. Other restrictions may apply. All sales are subject to PerkinElmer's general terms and conditions of sale (terms and conditions of sale may vary by country, please refer to the quote provided for the applicable terms and conditions of sale). Taxes and shipping are not applicable toward qualifying purchase amounts. Shipping charges apply to all purchases. Healthcare professionals, which include without limitation, and "Health Care Practitioner" under the Massachusetts Marketing Code of Conduct Law (105 C.M.R. §907.004) are not eligible for and may not participate in this promotion. This promotion/incentive may be discontinued, suspended, or amended by PerkinElmer in its sole discretion at any time without notice. Offer void where prohibited, licensed, or restricted by country, federal, state, provincial, or local laws or regulation or agency/institutional policy. Offer is expressly not valid in any country, state, province, or locality that requires prior registration of such offer. Other restrictions may apply. For details, please contact your sales representative. No cash or cash equivalent.

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)



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.