

## LABStartUp Program

# GET 15% OFF

## SAMPLE PREPARATION SYSTEMS

### Apply Now!

Simplify your sample preparation process and save.

Offer valid through December 31, 2022

### LABStartUp Savings – Apply Now and Get 15% Off!



OMNI International, the newest member of the PerkinElmer family, provides a uniquely diverse line of laboratory homogenizers. OMNI streamlines your sample preparation process with easy to use solutions designed for grinding, lysing, and homogenizing samples prior to downstream analysis. Fast, accurate, repeatable results every time.

To apply, complete the online joining form for the LABStartUp program, and our sales representatives will contact you.

[JOIN LABSTARTUP – APPLY NOW](#)



**15% Off Homogenizers  
and Automated Sample  
Prep Instruments**

## Laboratory Essentials



### Rotor-Stator Homogenizers

Fast. Powerful. Affordable.

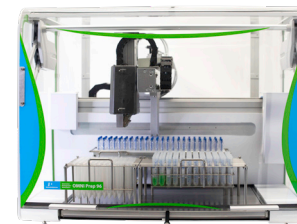
## High-Throughput



### Bead Mill Homogenizers

Versatile. Accurate. Repeatable Results.

## Automation



### Homogenizer Workstations

Set and Walkaway. Time Saver.



**How to Design Your Experiment.**

**Shop online today at [www.omni-inc.com](http://www.omni-inc.com)**

**\*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.