

MAKE YOUR LAB BUDGET GO FURTHER

LABStartUp

With PerkinElmer's 0% Lease Financing Options You Can:



Accelerate Productivity

Whatever your budget, when your lab needs it.



Stretch Your Budget

With no upfront costs, you can pay monthly for up to 3 years.



Keep Innovating

Don't let budget constraints stand between you and science.

How It Works:

Order Today, Pay Over Time

Choose from select PerkinElmer Analytical Instruments.

Flexible Leasing

Finance 100% of your purchase with a 0% interest rate. Choose from 12, 18, 24 or 36 month contract options.

Finance Eligibility

Complete our credit application and you will receive a quick credit decision.

** Certain restrictions apply. The information contained herein is not a commitment. All financing is subject to credit approval, completion of all required documentation, terms and conditions and availability, each at the sole discretion of PerkinElmer. PerkinElmer does not provide legal, tax or accounting advice. You must obtain and rely on such advice from your own accountants, auditors, attorneys or other professional advisors. 12, 18, 24, and 36 month financing validity on quotes with validity through Dec 31, 2022. Only PerkinElmer Analytical instruments will be considered for this offer. Offer may not be applied to existing, pending, or prior orders. 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). 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 or changed without notice at any time. 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."*

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.

311611 (83411) PKI