

# ORDERING ONLINE: IT'S THAT EASY



Sign Up and Get Started Today at [www.perkinelmer.com/register](http://www.perkinelmer.com/register)

## Quick and Efficient Three Steps - that's how easy it is to buy online

1. Visit [www.perkinelmer.com](http://www.perkinelmer.com)
2. Click on "Login/Register" at the top of the web page to register.  
All you need is your PerkinElmer account number.
3. Search for the product you need and add to your cart

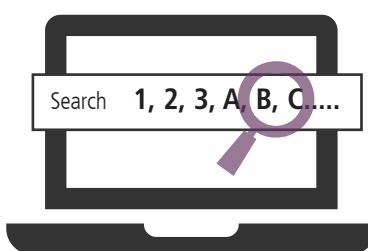
### Benefits of ordering online:

- Real time product availability
- Immediate order processing
- Order confirmations
- Shipment notifications
- Quick re-order
- Save items to favorites list
- Tracking information and updates
- All orders available in "my account" order history

Step 1: Quick registration.

Step 2: Shop by product number, alphabetical listing, or by technology.

Step 3: Checkout via your shopping cart.



The power and convenience of real-time purchasing information are now available online.  
When will the product ship? Shipment date of each item is visible in the shopping cart.

Shop Online today!  
<https://www.perkinelmer.com/buy>

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 ©2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.