



a | 5350 Lakeview Parkway South Dr. | Indianapolis, IN 46268
o | 317.808.4200 or 800.352.4975
f | 317.808.4395
w | beckman.com

PRODUCT OBSOLESCENCE NOTIFICATION

June 2, 2022

Dear Valued Customer,

This letter is to inform you that Beckman Coulter Life Sciences will obsolete the Biomek 4000 Laboratory Automation Workstation product line, effective **December 30, 2022**.

This obsolescence applies to all hardware configurations and accessories, including:

- Biomek 4000 Automated Workstation (A99749)
- Biomek 4000 Automated PCR Set-up Package (B22637)
- Biomek 4000 Automated PCR Clean-up Package (B22638)
- Biomek 4000 Genomic Workstation Package (B22640)
- Biomek 4000 Liquid Handling Package (B22867)
- Biomek 4000 Workstation Package for CE SDS Sample Prep (B37830)
- Biomek 4000 NGS Workstation (B42929)
- Biomek 4000 HLA Prep Workstation (C23350)

We plan to support the Biomek 4000 instruments for at least five years from the date of obsolescence. A formal notification will be sent one year prior to the discontinuation of support. **Please contact your local service representative with any questions.**

Since the introduction of the Biomek 1000 instruments over 30 years ago, Beckman Coulter has continued to innovate new automation solutions to accelerate scientific discoveries. The Biomek i-Series instruments with improved capacity, flexibility, and usability—represents the latest step in this evolution.

While we're proud of the continued advancement of our automated platforms, we understand that this obsolescence may impact our current customers. Please be assured that most existing Biomek instruments' methods can be adapted to the new Biomek i-Series Workstations, and we're confident you'll benefit from the new platform enhancements if you choose to upgrade to an Biomek i-Series workstation in the future. Please go to Beckman.com or contact your local sales representative for more information.

Thank you for your continued partnership and support of Beckman Coulter Life Sciences automation.

Sincerely,

Crystal Girod,
Commercial Product Manager, Liquid Handling

© 2022 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the Stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.