



a | 5350 Lakeview Parkway South Dr. | Indianapolis, IN 46268  
o | 317.808.4200 or 800.352.4975  
f | 317.808.4395  
w | [beckman.com](http://beckman.com)

## **PRODUCT OBSOLESCENCE NOTIFICATION**

November 20<sup>th</sup>, 2025

Dear Valued Customer,

This letter is to inform you that Beckman Coulter Life Sciences will obsolete the Biomek NGenius Next Generation Library Prep System, effective December 31<sup>st</sup>, 2025.

This obsolescence applies to all hardware configurations listed below:

- Biomek NGenius Next Generation Library Prep System (C62703)
- Biomek NGenius Next Generation Library Prep System (D03220)

We plan to support the Biomek NGenius instrument for at least five (5) years from the effective date of obsolescence. Please contact your local service representative with any questions regarding continued support.

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 many of the existing Biomek NGenius methods can be adapted to the current Biomek i-Series Workstations (Biomek i3, i5 and i7). We are confident you will benefit from the new platform enhancements should you choose to upgrade to a Biomek i-Series workstation in the future. For more information, please visit [Beckman.com](http://Beckman.com) or contact your local sales representative.

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

Sincerely,

*Kyle Curry*

Kyle Curry,  
Senior Product Manager, Liquid Handling

© 2025 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.