



March 30th, 2024
Crystal Girod, Biomek Product Manager
Biomek FX_P and Biomek NX_P End of Service and Support

Dear Valued Biomek Automated Workstation Customer,

Since 1935, Beckman Coulter Life Sciences has had a relentless commitment to innovation. Our teams are always on a journey of continuous improvement to bring you the best products and service possible.

This letter is to inform you of the End of Service and Support date effective December 31st, 2028 for the following:

- Biomek FX_P Single Arm System and Single Multichannel Pipettor (A31842)
- Biomek FX_P Single Arm System with Span-8 Pipettor (A31843)
- Biomek FX_P Dual Arm System, Multichannel Pipettor and Span-8 Pipettors (A31844)
- Biomek FX_P Dual Arm System with Dual Multichannel Pipettors (A31845)
- Biomek NX_P Multichannel Laboratory Automation Workstation with Integrated Gripper (A31841)
- Biomek NX_P Automated Workstation, Span-8 With Gripper (A31840)
- Biomek NX_P Span-8 Without Gripper (A31839)

As part of our customer support commitment, Beckman Coulter will make its best effort to support the instruments through the End of Service and Support date. Preventive maintenance services will continue to be offered until December 31st, 2028.

As we will not be able to offer service and support after December 31st, 2028 for instruments listed above, we realize you may have questions about this change. We encourage you to contact your local Beckman Coulter Life Sciences service or sales representative to learn more about alternative solutions.

Kind Regards,

Crystal Girod
Biomek Product Manager, Beckman Coulter Life Sciences
cgrod01@beckman.com